1 October 2024

Structural basis for saxitoxin congener binding and neutralization by anuran saxiphilins

Sandra Zakrzewska¹, Samantha A. Nixon¹, Zhou Chen^{1#}, Holly S. Hajare², Elizabeth R. Park², John V. Mulcahy², Kandis M. Arlinghaus³, Eduard Neu⁴, Kirill Konovalov⁴, Davide Provasi⁴, Tod A. Leighfield³, Marta Filizola⁴, J. Du Bois², and Daniel L. Minor, Jr.^{1, 5-8*}

¹Cardiovascular Research Institute

⁵Departments of Biochemistry and Biophysics, and Cellular and Molecular Pharmacology

⁶California Institute for Quantitative Biomedical Research

⁷Kavli Institute for Fundamental Neuroscience

University of California, San Francisco, CA 94158-9001 USA

⁸Molecular Biophysics and Integrated Bio-imaging Division Lawrence Berkeley National Laboratory, Berkeley, CA 94720 USA

²Department of Chemistry Stanford University, Stanford, CA 94305 USA ³National Oceanic and Atmospheric Administration National Centers for Coastal Ocean Science Charleston, SC 29412 USA ⁴Department of Pharmacological Sciences

Ichan School of Medicine at Mount Sinai New York, NY 10029 USA

*Correspondence to: daniel.minor@ucsf.edu

[#]Present address: Department of Anatomy and Physiology Shanghai Jiao Tong University School of Medicine Shanghai, 200025, China

Keywords: Saxitoxin, Toxin resistance, Saxiphilin, Gonyautoxins, Neosaxitoxin, Paralytic Shellfish Poisoning (PSP)

Short Title: STX congener binding to saxiphilins

1 October 2024

Abstract

Dinoflagellates and cyanobacteria in harmful 'red tide' algal blooms produce saxitoxin (STX) and \sim 50 congeners that block voltage-gated sodium channel (Na_v) function and disrupt bioelectrical signals¹⁻⁴. Consuming seafood carrying these lethal toxins causes paralytic shellfish poisoning (PSP), a growing public health hazard due to climate change⁵⁻ ⁷ that motivates efforts to detect these toxins and counteract their noxious effects. Although structural studies of Navs^{8,9} and anuran soluble STX binding proteins known as saxiphilins (Sxphs)^{10,11} revealed convergent binding modes for the bis-guanidinium STX core^{10,11}, the structural basis for STX congener recognition is unknown. Here, we show that American bullfrog (Rana catesbeiana) RcSxph^{10,11} and High Himalaya frog (Nanorana parkeri) NpSxph¹⁰ use a pre-organized pocket to sequester STX congeners through a binding mode shared with STX. This 'lock and key' recognition vields a tradeoff between a relatively rigid high-affinity toxin binding site in which bound waters are crucial and the ability of Sxphs to accommodate STX congener modifications. Importantly, functional studies show that Sxphs act as 'toxin sponges' that reverse Nav block by multiple STX congeners and can detect these bis-quanidinium toxins in a radioligand receptor binding assay (RBA) for PSP toxin environmental testing^{12,13}. Our findings establish how Sxphs sequester diverse neurotoxins and reveal structural factors underlying STX congener binding differences between Sxphs and Navs that are rooted in the distinct toxin binding orientations on these two targets. These insights expand the molecular foundation required for understanding toxin sponge action and for guiding development of new means to monitor PSTs and mitigate their harmful effects.

1 October 2024

Introduction

Saxitoxin (STX) is the archetype of a family of lethal, paralytic, small-molecule guanidinium neurotoxins that block voltage-gated sodium channels (Na_vs) and their bioelectrical signals in nerve and muscle^{1,5,14} and is a chemical weapon^{5,14}. Diverse species of dinoflagellates and cyanobacteria in marine and fresh water harmful algal blooms produce this toxin class as cocktails of STX and STX congeners, some of which match or exceed STX toxicity¹⁻⁴. Accumulation of STX and related toxins in seafood causes paralytic shellfish poisoning (PSP) and poses an increasing threat to commercial fishing and public health from climate change driven harmful algal blooms⁵⁻⁷. Consequently, there are extensive worldwide efforts to detect contamination of food and water supplies by PSP toxins^{15,16}. Structural studies of STX bound to Na_vs^{8,9} and frog saxiphilins (Sxphs)^{10,11}, a class of soluble high affinity STX binding proteins^{10,17}, reveal a common STX recognition mode^{10,11} by which these two structurally and functionally unrelated proteins use a relatively rigid binding site to coordinate the di-cationic, bis-guanidinium core of the toxin through a similar set of ionic and cation- π interactions^{10,11}. How STX congeners bind to these targets and whether differences in binding affinity arise from changes to the STX binding mode or by some other means is unknown.

Natural modifications to STX yield a family of ~50 STX congeners, collectively known as paralytic shellfish toxins (PSTs)^{1,3,5}, that affect toxin affinity for Na_Vs^{5,14} and frog plasma containing Sxph activity¹⁷. STX is commonly modified at three positions (Fig. 1A): R1, the carbamoyl site; R2, the six membered guanidinium ring N1 position; and R3, the C11 carbon. R1 modifications include acylation (acetate, mono- and bis-hydroxybenzoate), decarbamoylation (e.g. decarbamoyl STX, dcSTX) and sulfation (e.g. gonyautoxin 5, GTX5, also known as B1). R2 hydroxylation creates neosaxitoxin, neoSTX, a more effective Na_V blocker¹⁸ and more potent poison^{1,3} than STX. R3 can be modified by carboxymethylation (e.g. 11-saxitoxin ethanoic acid) and sulfation (e.g. gonyautoxin 2/3, GTX2/3), creating two interconverting epimers^{1,5,14}. Combinations of these alterations further diversify the STX family. The chemical complexity and variable proportions of STX congeners in natural PST samples poses challenges for quantitative testing¹² and for developing PSP treatments.

Anuran Sxphs from diverse frogs and toads comprise a family of soluble, transferrin-like 'toxin sponge' proteins that contain a single high-affinity (Kd ~ nM) STX binding site^{10,11,17}. This property allows Sxphs to rescue Na_Vs from STX block^{10,19} and likely contributes to the ability of some frogs to resist STX poisoning^{17,19-21}. Competitive radioligand binding studies of bullfrog¹⁷ and cane toad plasma²² suggest that Sxphs can bind select STX congeners. However, there have been no direct

1 October 2024

studies of STX congener binding to purified Sxphs and no structural information exists for STX congener binding to either Sxphs^{10,11} or Na_vs^{8,9}, yielding a gap in understanding STX congener recognition.

To evaluate how STX congeners interact with their molecular targets, we investigated toxin binding to RcSxph from the American bullfrog (Rana catesbeiana)^{10,11} and NpSxph from the High Himalaya frog (*Nanorana parkeri*)¹⁰, two previously characterized Sxphs that share a common toxin binding pocket but that have different STX affinities¹⁰. Our studies using a suite of thermofluor (TF), fluorescence polarization (FP), and isothermal titration calorimetry (ITC) binding assays^{10,23} show that both Sxphs have similar selectivity profiles ranging from enhanced binding of GTX5 to extremely weak affinity for the neoSTX series. High resolution X-ray crystal structures of NpSxph:STX congener complexes together with molecular dynamics (MD) simulations reveal toxin binding poses identical to STX and identify a network of ordered water molecules involved in toxin recognition. These findings highlight the ability of the pre-organized, relatively rigid STX binding pocket to accommodate many STX modifications and uncover structural differences that affect the binding preferences of Sxphs and Na_vs for different STX congeners. Importantly, functional studies using two-electrode voltage clamp (TEVC) and planar patch clamp experiments establish that Sxphs act as 'toxin sponges'¹⁹ that can reverse Na_V block by various STX congeners. Further, we demonstrate that Sxphs can substitute for rat brain homogenates in a widely used radioligand receptor binding assay (RBA) for PSTs^{12,13}. Together, our findings establish a molecular foundation for STX congener binding to Sxphs that should inform studies of 'toxin sponge'-based resistance mechanisms^{19,24-26} and enable efforts to develop new biologics to detect and neutralize STX and related PSTs.

1 October 2024

Results

RcSxph and NpSxph bind diverse STX congeners but not neosaxitoxins

To profile *Rc*Sxph^{11,27} and *Np*Sxph¹⁰ binding to different STX congeners, we used a thermofluor (TF) assay^{10,23} in which toxin binding causes a concentration-dependent change in the apparent melting temperature (Tm) of the target protein (Figs. 1A-E). We found that decarbamyolation at R1 (dcSTX) and alterations such as carbamate sulfation (GTX5) at R1, C11 sulfation at R3 (GTX2/3), combined C11 sulfation and decarbamoylation (dcGTX2/3), and combined carbamate and C11 sulfation (C1/C2) (Fig. 1A) shift the apparent melting temperature (Δ Tm) (Table S1) of both Sxphs (Figs. 1B-E, Table S1). Among the various congeners, GTX5 behaved most like STX, having a similar Δ Tm for both Sxphs (Δ Tm = 3.7 ± 0.1 and 3.6 ± 0.2°C, and 3.3 ± 0.1 and $3.2 \pm 0.2^{\circ}$ C, for GTX5 and STX, and RcSxph and NpSxph, respectively). By contrast, dcSTX, GTX2/3, dcGTX2/3, and C1/C2 exhibited reduced Δ Tm changes, indicating that these congeners bind more weakly than STX (Δ Tm = 3.1 ± 0.1, 2.7 ± 0.1, 2.3 ± 0.1, and 2.6 ± 0.1°C, for *Rc*Sxph, respectively and $\Delta Tm = 2.2 \pm 0.1$, 2.2 ± 0.1 , 1.4 ± 0.1 , and $1.5 \pm 0.1^{\circ}C$ for NpSxph, respectively), (Figs. 1B-C, Table S1). RcSxph and NpSxph show similar Δ Tm concentration dependences for this toxin panel (Figs. 1D-E). Importantly, these data establish that the Sxph TF assay can discriminate among the various modified forms of STX and serve as a facile means to assess Sxph:STX congener interactions.

Hydroxylation of the six membered guanidinium ring at N1 (the R2 site) yields neoSTX^{3,5}, a toxin that is a more effective Na_V blocker¹⁸ and more potent poison^{1,3} than STX. To address whether *Rc*Sxph and *Np*Sxph bind N1 hydroxylated toxin congeners, we used the TF assay to profile neoSTX, dc-neoSTX, GTX6, and GTX1/4 (Figs. 1B-E and S1A-D, Table S1). In contrast to the results obtained with STX congeners, neither neoSTX nor its derivatives affected Δ Tm substantially (Figs. S1A-D, Table S1), indicating that N1 hydroxylation potently interferes with binding to *Rc*Sxph and *Np*Sxph. Although none of these toxins altered the *Np*Sxph Δ Tm, neoSTX, dc-neoSTX, and GTX6 showed some minor Δ Tm changes for *Rc*Sxph at the highest tested concentrations (>10 µM) (Fig. S1C). The *Rc*Sxph Δ Tm rank order for dcSTX, GTX5 (B1), C1, and neoSTX (Table S1) agrees with radioligand binding competition studies of *R. catesbeiana* plasma¹⁷ and reinforces prior conclusions that N1 (neoSTX) and C11 (C1) modifications have the largest impact on *Rc*Sxph binding¹⁷.

To measure the affinity of STX and neoSTX congeners to *Rc*Sxph and *Np*Sxph, we modified our toxin binding FP assay^{10,23} into a fluorescence polarization competition (FPc) assay²⁸ that

1 October 2024

measures the ability of unlabeled toxins to compete with a fluorescein-labeled STX derivative. F-STX¹⁰ (Figs. 2A-B). Control experiments with STX yielded apparent affinities for RcSxph and NpSxph (Kd = 8.9 ± 1.6 and 8.7 ± 1.7 nM, respectively) that correspond with direct FP measurements (Kd = 7.4 ± 2.6 and 0.5 ± 0.3 nM, respectively). Subsequent FPc profiling of STX congeners (Figs. 2A-B, Table S2) revealed trends that follow the TF results (Figs. 2C-D). As expected from its STX-like effects on Δ Tm, GTX5 had an affinity similar to STX (Kd = 9.6 ± 0.8 and 8.9 \pm 1.6 nM and 7.3 \pm 1.8 and 8.7 \pm 1.7 nM, for GTX5 and STX and RcSxph and NpSxph, respectively), whereas dcGTX2/3, the toxin having the smallest Δ Tm, had the lowest affinity (180 ± 12 and 568 ± 68 nM for RcSxph and NpSxph, respectively) (Table S2). neoSTX, dc-neoSTX, and GTX1/4 bound RcSxph and NpSxph weakly, having affinities >5 µM (Figs. S1E-F, Table S2). In this panel, GTX6 was notable for displaying a weak but measurable affinity compared to other neoSTX congeners (1516 ± 222 and 1239 ± 578 nM for RcSxph and NpSxph, respectively). However, this effect could be attributed to a ~0.5% contamination of the strong binder, GTX5, in the commercial toxin sample used in this assay (Table S2). These results highlight the sensitivity of our assay and the ability of Sxphs to select between toxins in mixtures. The excellent agreement between Δ Tm and the $\Delta\Delta$ G values (Figs. 2C-D) obtained by FPc aligns with TF-FP correlations observed in RcSxph STX binding pocket alanine scans¹⁰ and demonstrates that ΔTm provides an accurate rank order assessment of STX congener affinity differences. Further, there is a strong correlation in STX congener binding between RcSxph and NpSxph (Fig. 2E) that aligns with the similar STX binding pockets found in these two Sxphs¹⁰.

RcSxph and NpSxph discriminate among STX congeners in a receptor binding assay

Testing for PST food contamination often relies on a radioligand receptor binding assay (RBA) using tritiated STX ([³H]STX) and rat brain homogenate^{12,29}. As Sxphs exhibit high thermostability, having Tms ~>50°C¹⁰, we wanted to examine whether purified Sxphs could be used in place of brain homogenates. RBAs using *Rc*Sxph (Figs. 2F, S2A-C) and *Np*Sxph (Figs. 2G, S2D-F) show that both proteins detect STX and its congeners consistent with their performances in the TF and FPc assays. The data reflect clear differences among the various toxins, including those from the neoSTX series (Table S3). Importantly, none of the toxins bind to *Rc*Sxph E540A, a mutant having ~2000-fold lower affinity for STX than *Rc*Sxph¹⁰ (Fig. S2G). This result demonstrates that, as with the TF and FP assays¹⁰, toxin detection relies on an Sxph STX binding pocket competent for toxin binding. RBA for *Rc*Sxph and *Np*Sxph and rat brain homogenate display comparable binding affinities, spanning 3-4 orders of magnitude (Kds ~1 nM–1 µM and ~3 nM-10 µM for rat brain homogenate and Sxphs, respectively) (Figs. S2H-I). Rat brain homogenates have a similar

1 October 2024

dynamic range for both the STX series (STX, dcSTX, GTX5, GTX2/3, dcGTX2/3, and C1/C2) and the neoSTX series (neoSTX, dc-neoSTX, GTX6, and GTX1/4) (Figs. S2H-K), whereas Sxphs are better at discriminating between these two toxin groups (Figs. S2A-F, Table S3). The RBA binding affinities and $\Delta\Delta G$ changes are well-correlated with those measured by FPc (Figs. 2H-I, Tables S2-S3) providing further validation of the TF and FPc assay trends (Figs. 2C-D). Together, these data demonstrate that purified Sxphs can serve as viable replacements for rat brain homogenate in the RBA PST detection assay.

Thermodynamics of toxin binding reveal critical moieties for Sxph interaction

To examine the thermodynamics of STX congener binding, we conducted isothermal titration calorimetry (ITC) experiments to measure the binding of dcSTX, GTX2/3, dcGTX2/3, GTX5, and C1/C2 to *Np*Sxph directly (Table 1, Figs 3A-E). As expected from prior studies with STX^{10,11,17,30}, this STX congener set binds *Np*Sxph with a 1:1 stoichiometry (Table 1). GTX5 showed thermodynamic binding parameters similar to those of STX, in line with the comparable affinities of these two toxins, whereas the other toxins showed reduced Kds and concomitant Δ H and Δ S changes. The $\Delta\Delta$ G values determined by ITC follow those from the FPc assay, but show systematically tighter binding (Fig. 3F, Tables 1 and S2).

Comparison of the *Np*Sxph:STX congener binding free energy reveals interesting trends regarding the various STX modifications. Carbamate removal (dcSTX and dcGTX2/3) reduces binding by ~1 kcal mol⁻¹ consistent with this moiety interacting with key STX binding determinants lle559, Phe562, and Pro728 within the *Np*Sxph binding pocket¹⁰. C11 sulfation is uniformly detrimental, causing $\Delta\Delta$ G changes of ~2-3 kcal mol⁻¹ for GTX2/3 vs. STX, dcGTX2/3 vs. dcSTX, and C1/C2 vs. GTX5. This modification makes close contacts with the binding pocket Gly798-Val799 backbone. By contrast, R1 carbamate sulfation (Fig. 1A) is favorable compared to STX ($\Delta\Delta$ G = -0.2 kcal mol⁻¹ for GTX5 vs. STX) but reduces binding when C11 is also sulfated ($\Delta\Delta$ G = 0.7 kcal mol⁻¹ for C1/C2 vs. GTX2/3). Hence, each modification evokes specific changes to the Sxph:toxin interaction. Some are straightforward to interpret, such as a reduction in binding enthalpy for dcSTX versus STX consistent with the fewer dcSTX contacts to the Sxph binding site, whereas others, such as the C11 sulfated toxins, show more complex enthalpy-entropy compensation effects suggesting that changes in water solvation of the toxin and binding pocket are important^{31,32}. The *Np*Sxph TF, FPc, RBA, and ITC results are in excellent agreement, and together provide a versatile assay suite for characterizing Sxph:toxin interactions²³.

NpSxph:STX congener structures reveal similar interactions and a common water network

1 October 2024

To visualize Sxph:STX congener interactions, we co-crystallized and determined the X-ray crystal structures of *Np*Sxph:dcSTX, *Np*Sxph:GTX2, *Np*Sxph:dcGTX2, *Np*Sxph:GTX5, and *Np*Sxph:C1 at resolutions of 1.90Å, 1.95Å, 1.80Å, 1.90Å, and 1.90Å, respectively, using molecular replacement (Table S4). The maps show excellent quality densities that correspond to the unique shape of each bound toxin (Figs. 4A-E) and show that all STX congeners bind *Np*Sxph with a 1:1 stoichiometry matching the functional studies (Table 1). GTX2/3, dcGTX2/3, and C1/C2 are epimeric mixtures at the C11 sulfated position^{1,5,33}. Because it was not possible to discern the bound epimer from the density, we assigned the density to the most prevalent epimer in the sample (GTX2, dcGTX2, and C1) (Figs. 4B, D, and E).

Comparisons of the *Np*Sxph:STX congener structures with the *Np*Sxph:STX complex (PDB: 8D6M)¹⁰ show that there are no large-scale changes in the Sxph core (N- and C-lobes) and thyroglobulin domains (Thy1-1 and Thy1-2) (root mean square deviation of C α positions (RMSD_{C α}) = 0.229Å, 0.259Å, 0.223Å, 0.235Å, and 0.239Å for dcSTX, GTX2, dcGTX2, GTX5, and C1, respectively) (Fig. 4F). Superposition of STX and STX congeners highlights that all have a similar binding pose involving identical interactions of the tricyclic bis-guanidinium core with *Np*Sxph Glu541, Asp786, Asp795 and Tyr796 that follow the STX molecular recognition fingerprint¹⁰ (Figs. 4G, S5A-F). In general, there are only minor changes in the position of the toxin among the structures. The most variation occurs at the carbamate (Cb) (Fig. 4H), with the largest shifts observed for the sulfocarbamate toxins GTX5 and C1 (~0.9Å and ~1.8Å, respectively).

Despite the variety of STX core modifications, the *Np*Sxph STX binding pocket residues display minimal structural movement between the apo¹⁰ and STX congener bound forms (Movie S1). In all structures, Asp786 coordinates the five-membered ring as seen in the *Np*Sxph:STX complex¹⁰ (Figs. S4A-F) and is repositioned from the outward-facing rotamer of the apo structure (Fig. S4G)¹⁰. In the *Np*Sxph:GTX5 and *Np*Sxph:C1 complexes, Arg719 moves by ~5Å towards the Cb sulfate (Figs. 4G, S4C, and F, Movie S1) and together with Arg566 forms an electrostatically positive ridge near the Cb sulfate (Fig. 4I). Both residues are conserved among Sxph sequences¹⁰.

The high resolution structures also revealed a set of water-mediated networks that contribute to *Np*Sxph:toxin interactions. Two water molecules, Wat541 and Wat795, occupy the same positions in the *Np*Sxph:STX complex¹⁰ and all *Np*Sxph:STX congener structures and form interactions with key hot-spot residues¹⁰ (Figs. S3A-E, S4A-G, and S5A-F). Wat541 is deeply buried in the binding pocket, and bridges the Glu541 sidechain, the residue that contributes the most binding

1 October 2024

energy to STX core recognition¹⁰, with the backbone carbonyl of Glu784 (Fig. S4H). Notably, this water bridge is absent in the apo structure¹⁰ (Fig. S4G). Instead, there is an axial water (Wat541a) on the opposite side of Glu541 (Fig. S4I). The second key water molecule, Wat795, bridges the sidechain and backbone oxygens of Asp795 with STX OH-14 and STX N9 (Fig. S4J) and is also lacking in the apo structure (Fig. S4G). The Asp795 site contributes nearly as much binding energy to toxin binding as the Glu541 site (~4.5 kcal mol⁻¹)¹⁰. Thus, the two most important residues for binding the STX core, Glu541 and Asp795, rely on water-mediated interactions. All structures, except those with C1 and GTX2, revealed a third water molecule, Wat786, coordinated with the Asp786 sidechain and the N7 position on the five-membered toxin ring (Figs. S3A-C, S4J, and S5). This Asp786 residue is one of the few residues that move between the apo and bound forms^{10,11} but contributes little to toxin binding energetics¹⁰.

To examine the roles of these water molecules further, we conducted a set of MD simulations totaling 5 µs for NpSxph:STX and each of the NpSxph:STX congener complexes, as well as models of complexes with the C11 epimers GTX3, dcGTX3, and C2. The STX core (Figs. S6A-B) and STX binding pocket (Figs. S6C-D) show remarkably low mobility leading to very stable toxin:protein contacts that persist throughout the simulations, regardless of toxin type (Fig. S6E). The similar behavior of the GTX2/3, dcGTX2/3, and C1/C2 epimers (Figs. S6A-C) indicates that there is little discrimination between these forms as underscored by their similar contacts in all trajectories (Fig. S6E). The stability of the Sxph:toxin interactions extends to the key water molecules Wat541 and Wat795, that coordinate Glu541 and Asp795 and their interactions with the toxin (Figs. S6E-F). Notably, Wat795 shows lower occupancy in the C11 sulfated toxin complexes (Fig. S6F), suggesting that part of the affinity loss caused by the sulfate modification derives from its effect on water occupancy at this site. This Wat795 behavior agrees with the observation that C11 sulfated congeners show a more favorable binding entropy (Table 1) and supports the idea that toxin and binding pocket solvation changes contribute to toxin affinity. In line with the lesser energetic importance of the Asp786 position¹⁰, its bound water, Wat786, exhibits low occupancy (Fig. S6H). Further, the absence of Wat786 is frequently correlated with the Asp786 movement away from the toxin and hydration of the coordination site by bulk water. Taken together, the X-ray data and MD simulations highlight the structurally rigid, pre-organized nature of the Sxph binding pocket and identify a shared set of protein-toxin and water-mediated contacts that engage diverse STX congeners having affinities that vary by ~two orders of magnitude.

Sxphs rescue Navs from STX congener block

1 October 2024

*Rc*Sxph is a 'toxin sponge' that can reverse STX inhibition of Na_Vs by competing for STX^{10,19}. As this ability is linked to Sxph affinity for STX¹⁰ and *Np*Sxph binds STX better than *Rc*Sxph by ~15 fold¹⁰, we wanted to test whether *Np*Sxph could act as a more effective toxin sponge. To this end, we used a previously established two-electrode voltage clamp (TEVC) assay^{10,19} in which 100 nM STX yields 90% block (IC₉₀) of the Golden poison frog *Phyllobates terribilis* skeletal muscle channel Na_V1.4 (*Pt*Na_V1.4) expressed in *Xenopus laevis* oocytes to measure the effect of Sxph addition (Figs. 5A-B, S7A-B). Consistent with their shared low nanomolar STX affinity¹⁰, Sxph titration showed that 2:1 Sxph:toxin ratios were sufficient for both *Np*Sxph and *Rc*Sxph to reverse STX block completely (Figs. 5A-C), agreeing with prior *Rc*Sxph studies¹⁹. Notably, the Sxph ratio required for 50% rescue of the blocked current was better for *Np*Sxph (Effective Rescue Ratio 50, ERR₅₀ = 0.5 ± 0.1 for *Np*Sxph and 0.7 ± 0.2 for *Rc*Sxph) (Figs. 5A-C, S7A-B), in agreement with its higher STX affinity¹⁰.

As *Np*Sxph and *Rc*Sxph^{10,19} both act as effective anti-STX toxin sponges (Figs. 5A-C) and also have strong binding affinity for some STX congeners (Tables 1, S1-S3), we next asked whether these proteins could act more broadly to rescue Na_Vs from different STX family members. We first measured the *Pt*Na_V1.4 response to dcSTX, GTX2/3, and GTX5 (Figs. 5D-E) using TEVC. All of these toxins were weaker blockers relative to STX (IC₅₀ = 12.6 ± 1.4 nM¹⁹) having a rank order of STX<GTX2/3<dcSTX<<GTX5 (IC₅₀ = 27.5 ± 2.1, 144.4 ± 29.7, and 2290 ± 563 nM, respectively) (Figs. 5D-E, Table S5) similar to mammalian Na_V1.4³³⁻³⁶. We then tested whether a 3:1 *Np*Sxph:toxin ratio could neutralize GTX2/3 and dcSTX block of *Pt*Na_V1.4 at their respective IC₉₀s (200 nM and 800 nM, respectively). GTX5 was not tested because its weak IC₉₀ (>10 µM) required prohibitively large amounts of *Np*Sxph and toxin in this assay. These experiments revealed that *Np*Sxph reverses channel inhibition by both GTX2/3 (Fig. 5F) and dcSTX (Fig. 5G). Further, in line with its lower affinity for GTX2/3 relative to STX, *Np*Sxph had an increased ERR₅₀ for this toxin (ERR₅₀ = 2.1 ± 0.7 and 0.5 ± 0.1, respectively) (Figs. 5E and S7C), providing further support that rescue effectiveness depends on Sxph toxin affinity¹⁰.

Comparison of rescue kinetics revealed unexpected differences between dcSTX and GTX2/3. A 3:1 *Np*Sxph:toxin ratio caused rapid reversal of dcSTX block and complete current recovery within one minute (102.6 \pm 1.2%, n=6) (Figs. 5G, S7D). By contrast, reversal of GTX2/3 block using the same Sxph:toxin ratio was slower and biphasic showing an initial fast recovery phase yielding 87.4 \pm 6.5% (n=6) current recovery after 5 minutes (Figs. S7C and E) followed by a slower recovery phase that reached 91.8 \pm 2.8% (n=2) at 30 minutes (Figs. 5F and S7E). To test if the two phases resulted from contaminants in the naturally isolated GTX2/3 sample (GTX2/3_{nat}), we

1 October 2024

repeated the experiment using pure, synthetic GTX2/3 (GTX2/3_{syn}). These experiments showed that the 3:1 *Np*Sxph:GTX2/3_{syn} ratio had identical reversal behavior to GTX2/3_{nat}, displaying a biphasic response reaching 87.3 \pm 3.3% (n=5) current rescue at 5 minutes and a slower rescue phase that reached 94.4 \pm 3.3% (n=5) at 30 minutes (Figs. S7F-G). Having eliminated toxin contaminants as the source of the two phases, we next tested whether *Rc*Sxph would behave comparably given that it has a similar GTX2/3 binding affinity to *Np*Sxph (Tables S2-S3). Matching the effect of *Np*Sxph, a 3:1 *Rc*Sxph:GTX2/3_{syn} ratio was sufficient for complete rescue of channel block (Figs. S7H-I). However, *Rc*Sxph showed a faster current recovery during the first phase (97.2 \pm 4.5% (n=4) after 5 minutes) and complete rescue after 10 minutes (100.6 \pm 2.8%, n=4) (Fig. S7H). Structural comparison suggests that the GTX2/3 selectivity differences may originate from more electropositive surface near the C11 sulfate found in the *Rc*Sxph STX binding pocket compared to *Np*Sxph (Fig. S7J). Together, these results demonstrate that Sxphs rescue Na_V from STX congener block and uncover functional differences between Sxphs having similar STX binding sites^{10,11} (Tables S1-S3, and S5).

Sxph:Toxin binding neutralizes human Nav poisoning by diverse STX congeners

To overcome the sample limitations posed by the TEVC assay and enable investigation of a wider range of toxins, we established a planar whole cell patch-clamp assay using mammalian cells stably expressing *Homo sapiens* Na_V1.4 (*Hs*Na_V1.4) that required ~10,000-fold lower test sample volumes than the TEVC assay. Profiling the IC₅₀ values of STX and five congeners against *Hs*Na_V1.4 (Figs. 6A-B, Table S5) showed that low nanomolar concentrations of STX, GTX2/3, dcSTX, and dcGTX2/3 inhibited *Hs*Na_V1.4 (IC₅₀ = 3.0 ± 1.6, 6.8 ± 1.1, 14.7 ± 8.7, and 31.8 ± 10.6 nM respectively) consistent with prior measurements for *Hs*Na_V1.4⁴. These measurements also revealed modest changes in potency for decarbamoylation and C11 sulfation that match prior studies^{14,34}. By contrast, carbamate sulfation on GTX5 and C1/C2 reduced *Hs*Na_V1.4 block by ~2 orders of magnitude relative to STX (IC₅₀ = 335.3 ± 55.4 and 151.5 ± 37 nM, respectively) (Figs. 6A-B, Table S5) consistent with previous reports^{14,34}.

We then assessed the ability of *Np*Sxph to prevent *Hs*Na_V1.4 block by STX, dcSTX, GTX2/3, dcGTX2/3, GTX5, and C1/C2 at their respective IC₉₀ values (Table S5). Comparing the *Hs*Na_V1.4 sodium currents in the presence of a 5:1 *Np*Sxph:toxin mixture against those elicited in the absence and presence of each toxin at its IC₉₀ concentration revealed that *Np*Sxph neutralized all tested STX congeners (Figs. 6C-D). As with the TEVC rescue experiments¹⁰ (Fig. 5C), the degree of neutralization followed toxin binding affinity trends. We observed nearly complete neutralization of STX, GTX5, C1/C2, and dcSTX (93.7 \pm 6.2, n=5; 99.2 \pm 1.4, n=4; 90.4 \pm 5.8,

1 October 2024

n=6; and 88.3 \pm 5.5%, n=7; respectively) (Figs. 6C–D) and partial neutralization of GTX2/3 and dcGTX2/3 (53.2 \pm 9.5 , n=7; and 59.9 \pm 7.0%, n=7; respectively) (Fig. 6E). The lower fractional recovery for the C11 monosulfated toxins in the planar patch-clamp recordings highlights that rescue is governed by the relative affinities of a given toxin for the target channel and Sxph. Together with the TEVC studies (Figs. 5F-G), these data demonstrate the capacity of Sxphs to act as broad-spectrum 'toxin sponges' that can neutralize multiple, naturally-occurring STX congeners and reverse their inhibition of Na_vs.

Divergent STX congener affinities highlight Sxphs and Navs toxin binding factors

Structural and binding studies have uncovered a remarkable convergence in the way Sxphs and Na_vs recognize the STX core^{10,11}. Our studies here reveal that despite this common recognition framework, there are clear differences in how R1, R2, and R3 toxin modifications (Fig. 1A) affect interaction with each target. The sulfocarbamate (GTX5) has minimal effect on Sxph binding (Tables S2-S3) but is detrimental to Na_V1.4 binding (>3 kcal mol⁻¹) (Table S3). Conversely, C11 sulfation (GTX2/3) has modest effects on channel interactions but destabilizes binding to Sxphs by ~1 kcal mol⁻¹. Whereas N1 hydroxylation (neoSTX) increases toxin affinity (Table S3) and ability to block the channel¹⁸, this modification severely compromises *Rc*Sxph and *Np*Sxph binding (>3.7 kcal mol⁻¹) (Table S2-S3). There are no Na_vs:STX congener structures. Thus, we used the NpSxph:C1 structure, having a toxin bearing sulfation at both the carbamate and C11 to compare the toxin pose with that found in the human Na_V1.7:STX complex⁸ (Fig. 7A-B) to gain insight into the STX congener binding differences between the two targets. The position of the DI and DII residues that engage the STX core sets a lateral toxin orientation across the Nav pore ring that leads to sulfocarbamate clashes with the P1 and P2 helices and Nav selectivity filter of the DIII pore module and exposure of the C11 sulfate to the extracellular solvent facing side of the pore above the DIV the selectivity filter (Figs. 7A-B). The putative sulfocarbamate-DIII clash and accommodation of the C11 sulfate are consistent with the effects of these modifications on Na_V (Tables S3 and S5), common effects of DIII sequence variations on STX and GTX3 block³⁷, and energetic coupling between carbamate modification and DIII mutants³⁶. By contrast, the Sxph residues that coordinate the STX core sit at the bottom of a deep pocket (Fig. 7C-D). This positioning sets a toxin orientation that exposes both the carbamovl and C11 sulfate groups to the solvent-exposed opening of the binding site and explains why both toxin forms bind Sxphs. Notably, N1 hydroxylation appears to create a clash with the key binding position Glu541¹⁰ in the deepest part of the Sxph binding site (Figs. 7C-D and S8), offering a rationale for why neoSTX and its congeners are poor RcSxph and NpSxph binders (Tables 1 and S1-S3). Together, these comparisons show that the orientation of the STX core components with respect to other binding

1 October 2024

target elements has a strong influence on how STX targets interact with various STX core modifications and identify key structural features that differentiate how Na_Vs and Sxphs bind STX congeners.

1 October 2024

Discussion

Sxphs^{10,11} and Na_vs^{8,9,38} bind STX with high affinity by means of a rigid binding site that uses a convergent molecular recognition strategy to engage the STX bis-guanidinium core^{10,11} even though each presents the shared STX binding motif on different scaffolds. How STX congener modifications affect binding interactions with these two classes of PST binding proteins and whether Sxph^{10,11} and Nav^{8,9,38} high affinity STX sites can be reshaped to accommodate the various STX congener modifications have been unclear. Our data show that the most prevalent STX congener alterations (decarbamoylation and carbamate sulfation at R1, N1 hydroxylation at R2, and C11 sulfation at R3) (Fig. 1A) have profound effects on RcSxph and NpSxph toxin affinities that range over of ~3 orders of magnitude (Tables 1, S1-S3). Strikingly, despite these large affinity differences, our structural analysis shows that STX congeners bearing various combinations of decarbamovlation, carbamate sulfation, and C11 sulfation bind with poses matching STX (Figs. 4G-H, S6A-B). Further, similar to STX¹⁰, dcSTX, GTX2/3, dcGTX2/3, GTX5, and C1/C2 binding induces minimal conformational change in the Sxph toxin binding pocket, apart from the movement of an aspartate (NpSxph Asp786) to coordinate the five-membered guanidinium ring of the toxin (Fig.S4 Movie S1). The strongest binding perturbations are caused by N1 hydroxylation in the neoSTX series (Tables S1-S3) that creates an apparent clash in a deeply buried part of the binding pocket (Fig. S8). Thus, the 'lock and key' mode of toxin binding used by Sxph strongly influences toxin preferences.

Our structural studies (Figs. S3-S5) and MD simulations (Fig. S6) identified a set of common water molecules that are crucial to toxin binding. These ordered waters, absent from the apo-*Np*Sxph structure¹⁰ (Fig. S4G), buttress toxin interactions by stabilizing the residue that interacts with the N1 nitrogen of the six membered guanidinium ring, *Np*Sxph Glu541, and by bridging *Np*Sxph Asp795 and the toxin hydrated ketone (Figs. S4A-J). Notably, these conserved Sxph residues are most critical for STX binding affinity¹⁰, each contributing >4 kcal mol⁻¹ binding energy. Hence, maintaining 'lock and key' binding comes at the cost of reduced affinity for decarbamoylated, C11 sulfated, and N1 hydroxylated congeners (Tables 1, S1-S3) as each of these modifications affects interactions with various elements of the Sxph pocket. These affinity differences arise from loss of carbamate contacts¹⁰ (dcSTX and dcGTX2/3), destabilization of the Asp795 water-mediated bridge that affects binding entropy (Table 1) paired with close contacts to the binding pocket Gly798-Val799 backbone (C1/C2, GTX2/3, and dcGTX2/3) (Fig. S7J), and an apparent clash of R2 modified toxins with a key STX-binding residue, *Np*Sxph Glu541 (Fig. S8). By contrast, modifications that face the solvent-exposed mouth of binding pocket, such

1 October 2024

as carbamoylsulfation (GTX5) or linker addition¹⁰, are well tolerated. Together, these data highlight the tradeoff between the rigidity of the high-affinity the STX site^{10,11} in which water has a key role in mediating toxin binding and the ability of *Rc*Sxph and *Np*Sxph to accommodate STX congener modifications.

Sxphs and Navs have divergent rank orders of STX congener affinities that suggest binding site differences beyond the convergent framework used to recognize the STX bis-guanidinium core^{10,11}. The GTX5 sulfocarbamate is well-tolerated by Sxphs but not Na_vs, whereas Na_vs are better at accommodating the C11 sulfation of GTX2/3 than Sxphs (Tables S1-S3 and S5). Comparison of the Sxph and Nav toxin binding sites indicates that these differences are rooted in the divergent orientation of the toxins in the two binding sites that is set by the location of the core STX recognition elements. In Navs, the toxin sits laterally across the ring of the channel pore such that Domains I and II elements engage the bis-guanidinium core while DIII and DIV frame the carbamate and C11 sites (Figs. 7A-B). By contrast, Sxphs present a vase-like binding cavity in which the STX core binding residues are at the bottom of a deep pocket from which the carbamate and C11 sites project to the solvent exposed opening (Fig. 7C-D). The resulting orientation differences lead to sulfocarbamate clashes with Nay Domain III selectivity filter (Figs. 7A-B) that align with evidence for carbamate-DIII energetic coupling³⁶ and strong loss of binding (Table S3). Such clashes with carbamate modifications are absent in Sxphs (Figs. 7C-D). Indeed, Sxphs allow large modifications at this site with no consequence to affinity¹⁰. Conversely, in Na_vs, C11 sulfation appears to be positioned in a solvent exposed part of the channel Domain IV selectivity filter free from steric clashes (Fig. 7A-B), consistent with its modest reduction in Nav block (~2-fold, e.g. STX vs. GTX2/3 Table S5)³⁷. This situation differs from Sxph where C11 sulfation yields close contacts with a binding pocket wall (Figs. 7C-D) and affects the stability of the water molecule that bridges the Asp795 and the toxin (Fig. S6F). These comparisons underscore the importance of toxin orientation with respect to the surrounding features of its binding site and the crucial role of water in mediating receptor-toxin interactions. Both factors are likely to be important in designing Sxphs or other proteins capable of recognizing various STX congeners. Whether STX congeners bind to Navs in orientations that differ from STX and whether water-mediated interactions like those seen in Sxph:toxin complexes are critical for high affinity Nav binding remain key undressed questions.

Naturally occurring PST mixtures comprising STX and its congeners are potent Na_V blockers^{1-4,14}, making such STX variants the concern of public health efforts to mitigate the effects of PSP outbreaks. Deciphering how STX and its congeners interact with their molecular targets is

1 October 2024

important for developing new ways to detect and neutralize PSTs, particularly as much seafood monitoring largely relies on a ~90 year old mouse lethality assay^{39,40}. A major challenge is developing reagents that can capture and distinguish different STX congeners. In this regard, it is notable that Sxphs are very stable having Tms ~>50°C¹⁰, produce toxin binding curves of comparable quality to rat brain homogenate in the RBA^{12,13} (Fig. S2, Table S3), and show a broader dynamic range in this assay. Hence, our results demonstrate that Sxphs can serve as an attractive, thermostable alternative for PST detection that merits further study.

Sxphs are thought to act as 'toxin sponges' that protect some frogs from STX poisoning by competing with Na_vs for toxin binding^{17,19}. Our studies show that *Rc*Sxph and *Np*Sxph rescue both amphibian and human Na_v1.4 from STX congener block (Figs. 5-6). Similar to STX¹⁰, the ability of these proteins to sequester toxin congeners is linked to Sxph:toxin affinity, as weaker Sxph binders such as GTX2/3 require higher Sxph:toxin ratios to affect rescue (Figs. 6D-E). Further, RcSxph and NpSxph show distinct capacities to discriminate STX from dcSTX (Table S2) and different kinetic behaviors in GTX2/3 rescue assays (e.g. Figs. S7G-I). These two Sxphs vary at a key position that influences toxin binding affinity (Tyr558 and Ile559, respectively)¹⁰, providing support for the idea that natural Sxph STX binding pocket variations shape STX congener selectivity¹⁰. NeoSTX series members comprise some of the most potent STX congeners^{1,3,5,18} but do not have appreciable affinity for RcSxph or NpSxph. Given Sxph STX binding pocket sequence diversity¹⁰ and evidence for neoSTX binding activity in cane toad (*Rhinella marina*) plasma²², an organism that has a Sxph¹⁰, profiling the STX congener binding of diverse Sxphs could provide a path to identify Sxphs capable of neutralizing neoSTX series toxins. Together, our results establish the necessary foundation for understanding how Sxphs sequester diverse neurotoxins and highlight the key role of water in toxin binding. This framework is important for understanding toxin sponge resistance mechanisms^{17,19} and for the development of new means to detect and neutralize diverse types of PSTs.

1 October 2024

Acknowledgements

We thank M. Huynh and A. Bara for technical help and K. Brejc for comments on the manuscript. This work was supported by grants DoD HDTRA-1-19-1-0040, HDTRA-1-21-1-10011, and HDTRA-1-23-1-0026 to D.L.M., and NIH-NIGMS R01-GM117263-05 to J.D. E.N. was supported by a fellowship from the Studienstiftung des deutschen Volkes and acknowledges the use of computational resources provided in part by the Erlangen National High Performance Computing Center (NHR@FAU) through the b132dc grant. Additional computations utilized resources from Scientific Computing at the Icahn School of Medicine at Mount Sinai, supported by the Clinical and Translational Science Awards (CTSA) grant UL1TR004419 from the National Center for Advancing Translational Sciences, as well as the Office of Research Infrastructure at the National Institutes of Health under award numbers S100D026880 and S100D030463. This research was supported in part by an appointment to the NOAA Research Participation Program administered by the Oak Ridge Institute for Science and Education (ORISE) through an interagency agreement between the U.S. Department of Energy (DOE) and NOAA.

Author Contributions

S.Z., S.A.N., Z.C., J.D., and D.L.M. conceived the study and designed the experiments. S.Z. and Z.C. performed the TF assay and produced and purified Sxphs. S.Z. established the FPc assay, performed the ITC studies, and determined the crystal structures of *Np*Sxph complexes. K.M.A. and T.A.L. designed and conducted the RBA experiments. E.N. designed, ran, and analyzed the molecular dynamics simulations. K.K. and D.P. contributed additional analyses of the simulations. S.A.N. performed two-electrode voltage clamp electrophysiology experiments. S.Z. and S.A.N. performed whole-cell patch clamp toxin concentration-response experiments. S.Z. established and performed human Na_V neutralization experiments. H.S.H., E.R.P., and J.V.M. synthesized and quantified samples of F-STX, STX, and GTX2/3. S.Z., S.A.N., Z.C., K.M.A., E.N., K.K., D.P. M.F., and D.L.M. analyzed data. T.A.L., M.F., J.D., and D.L.M. provided guidance and support. S.Z., S.A.N., J.D., and D.L.M. wrote the paper.

Competing interests

J.D. is a cofounder and holds equity shares in SiteOne Therapeutics, Inc., a start-up company interested in developing subtype-selective modulators of sodium channels.

The scientific results and conclusions, as well as any views or opinions expressed herein, are those of the authors and do not necessarily reflect the policies and views of the Department of Commerce (DOC), NOAA, DOE, or ORAU/ORISE.

1 October 2024

The other authors declare no competing interests.

Data and materials availability

Coordinates and structure factors and for *Np*Sxph:dcSTX (PDB:8V68), *Np*Sxph:GTX2 (PDB:8V69), *Np*Sxph:dcGTX2 (PDB:8V65), *Np*Sxph:GTX5 (PDB:8V66), and *Np*Sxph:C1 (PDB:8V67) are deposited with the RCSB and will be released upon publication.

Requests for material should be sent to D.L.M.

1 October 2024

Materials and Methods

Sxph expression and purification

Rana catesbeiana saxiphilin (RcSxph) (GenBank: U05246.1) and Nanorana parkeri saxiphilin (NpSxph) (GenBank: XM 018555331.1) were produced by expression in insect cells using baculoviruses and were purified as previously described^{10,11,23}. In brief, Sxphs carrying a C-terminal 3C protease cleavage site, green fluorescent protein (GFP), and a His₁₀ tag in series were expressed in Spodoptera frugiperda (Sf9) cells using a baculovirus expression system²³. After addition of P2/P3 baculovirus to Sf9 cells at the dilution ratio of 1:50 (v/v), cells were incubated in a non-humidified New Brunswick Innova 44 incubator (Eppendorf, cat. no. M1282-0010) at 27°C, shaking at 130 rpm for 72h. Expressed Sxphs were secreted into the growth media. Cells were harvested by centrifugation. The supernatant was adjusted to pH 8.0 with a final concentration of 50 mM Tris-HCl and treated with 1 mM NiCl₂ and 5 mM CaCl₂ to precipitate contaminants. Precipitants were removed by centrifugation, and the clarified supernatant was incubated with antiGFP nanobody-conjugated Sepharose resin for 5 hours at room temperature (23±2°C). The resin was washed with 20 column volumes of a wash buffer containing 300 mM NaCl and 30 mM Tris-HCl (pH 7.4). After purification with antiGFP nanobody resin, protein samples were treated with 3C protease (0.2 mg mL⁻¹ in the wash buffer) overnight at 4°C to remove the GFP-His tag from Sxphs. The cleaved eluates were collected and purified by size exclusion chromatography (SEC) using a Superdex 200 10/300 GL column (Cytiva). For the TF, FP, RBA, ITC assays, and electrophysiology experiments, Sxphs were purified using a final SEC step in 150 mM NaCl, 10 mM HEPES (pH 7.4). Structure determination of NpSxph complexes with STX congeners, was done following the same methods, except for the final SEC buffer containing 30 mM NaCl, 10 mM HEPES (pH 7.4). Protein concentrations were determined by measuring UV absorbance at 280 nm using the following extinction coefficients calculated using the ExPASY server (https://web.expasy.org/protparam/): RcSxph and RcSxph E540A mutant 96,365 M⁻¹ cm⁻¹; *Np*Sxph 108,980 M⁻¹ cm^{-1 23}.

Toxin preparation

Saxitoxin (STX) and fluorescein-labeled saxitoxin (F-STX) were synthesized, purified, and validated as outlined in ^{10,23,41,42}. STX and F-STX powders were directly dissolved with MilliQ water to make 1 mM and 1 µM stocks, respectively. Decarbamoylsaxitoxin (dcSTX, cat. no. dcSTX-c), gonyautoxin2/3 (GTX2/3, cat. no. GTX2/3-d), decarbamoylgonyautoxin-2/3 (dcGTX2/3, cat. no. dcGTX2/3-d), gonyautoxin-5 (GTX5, cat. no. GTX5-d), N-sulfocarbamoylgonyautoxin-2/3 (C1/C2,

1 October 2024

cat. no. C1/C2-c), decarbamoylneosaxitoxin (dcneoSTX, cat. no. dcneoSTX-d), gonyautoxin-6 (GTX6, cat. no. GTX6-b), and gonyautoxin-1/4 (GTX1/4, cat. no GTX1/4-e) were purchased from the National Research Council Canada (NRC). Neosaxitoxin (neoSTX, cat. no. 41619) was purchased from Sigma-Aldrich. All the purchased toxins were lyophilized prior to making 1-5 mM stocks using MilliQ water.

Thermofluor (TF) assay

Thermofluor assays⁴³ for the STX congeners (dcSTX, GTX2/3, dcGTX2/3, GTX5, C1/C2, neoSTX, dcneoSTX, GTX6 and GTX1/4) were performed as previously described for STX^{10,23}. Briefly, two-fold serial dilutions of each toxin were prepared in 150 mM NaCl, 10 mM HEPES, pH 7.4. 20 µL samples containing 1.1 µM *Rc*Sxph or *Np*Sxph, 5x SYPRO Orange dye (Sigma-Aldrich, cat. no. S5692, stock concentration 5,000x), 0-20 µM toxin, 150 mM NaCl, 10 mM HEPES, pH 7.4 were set up in 96-well PCR plates (Bio-Rad, cat. no. MLL9601), sealed with a microseal B adhesive sealing film (Bio-Rad, cat. no. MSB1001) and centrifuged (1 min, 230xg) prior to thermal denaturation using a CFX Connect Thermal Cycler (Bio-Rad, cat. no. 1855201). Fluorescence was measured using the HEX channel (excitation λ =515-535 nm, emission λ =560-580 nm). Samples were incubated at 25°C for 2 min followed by a temperature gradient from 25°C to 95°C at 0.2°C min⁻¹, and final incubation at 95°C for 1 min.

For the toxin dose-response curves, Sxph melting temperatures (Tms) in the presence of varied toxin concentrations were calculated by fitting the denaturation curves using a Boltzmann function in GraphPad Prism (GraphPad Software) using the equation $F=F_{min}+(F_{max}-F_{min})/(1+exp((Tm-T)/C))$, where F is the fluorescence intensity at temperature T, F_{min} and F_{max} are the fluorescence intensities before and after the denaturation transition, respectively, Tm is the midpoint temperature of the transition, and C is the slope at Tm. Δ Tms for Sxph in the absence (Tm_{Sxph}) and presence (Tm_{Sxph+toxin}) of different toxin concentrations were calculated using the following equation: Δ Tm=Tm_{Sxph+toxin}-Tm_{Sxph}.

Fluorescence polarization competition (FPc) assay

Fluorescence polarization competition (FPc) assays²⁸ were performed using 100 µL total reaction volume per well and final concentrations of the fluorescent ligand and Sxph of: for *Rc*Sxph, 1 nM F-STX and 12 nM *Rc*Sxph; for *Np*Sxph, 0.5 nM F-STX and 1.6 nM *Np*Sxph. For the displacement experiments, two-fold serial dilutions of unlabeled toxins (STX, dcSTX, GTX2/3, dcGTX2/3, GTX5, C1/C2, neoSTX, dcneoSTX, GTX6 and GTX1/4) were prepared using in a buffer of 150 mM NaCl, 10 mM HEPES, pH 7.4 at the following concentration ranges: STX and GTX5,

1 October 2024

0-875 nM: dcSTX and GTX2/3. dcGTX2/3. C1/C2. 0-7 µM: GTX6. 0-14 µM: and neoSTX. dcneoSTX, and GTX1/4, 0-28 µM. Samples containing 2 nM F-STX, 24 nM RcSxph, and 150 mM NaCl, 10 mM HEPES, pH 7.4 or 1 nM F-STX, 3.2 nM NpSxph, and 150 mM NaCl, 10 mM HEPES, pH 7.4 were incubated for 1 h at room temperature $(23 \pm 2^{\circ}C)$ protected from light. Subsequently, 50 µL the equilibrated Sxph:F-STX solution was mixed with 50 µL of each toxin dilution series in 96-well black flat-bottomed polystyrene microplates (Greiner Bio-One, cat. no. 655900). The plate was sealed with an aluminum foil sealing film AlumaSeal II (Excel Scientific, cat. no. AF-100) and incubated at room temperature (23 ± 2°C) for 2 h to attain equilibrium. Measurements were performed at 25°C on a Synergy H1 microplate reader (BioTek) using the polarization filter setting (excitation λ =485 nm, emission λ =528 nm). Data were normalized using the following equation ²⁸: P = ($P_{Sxph:toxin} - P_{Toxin}$)/($P_{Ctrl} - P_{Toxin}$), where P is the polarization measured at a given toxin concentration, P_{Sxph:toxin} is the polarization of Sxph:toxin mixture, P_{Toxin} is the polarization of toxin in the absence of Sxph, and P_{Ctrl} is the maximum polarization of Sxph in the absence of unlabeled toxin. The concentrations of the competing toxin causing displacement of 50% of bound F-STX (IC₅₀) were calculated by fitting normalized fluorescence polarization as a function of toxin concentration using the nonlinear regression analysis in GraphPad Prism (GraphPad Software).

Radioligand receptor binding assay (RBA)

Radioligand receptor binding assay (RBA) analysis for PSTs followed established protocols^{12,13} with modifications. Tritiated saxitoxin ([³H]STX) was provided by American Radiolabeled Chemicals (ARC, St. Louis, MO; cat no. ARK0101). Saxitoxin dihydrochloride (STX-diHCl, cat. no. NIST-8642a) was purchased from the National Institute of Standards and Technology. All other unlabeled toxins were Certified Reference Material (CRM) purchased from the NRC Canada (cat. no. dcSTX-c, GTX2/3-d, dcGTX2/3-c, GTX5-d, C1/C2-c, dcneoSTX-d, GTX6-b, GTX1/4-d, neoSTX-d).

Assays were performed in a Multiscreen 96-well GF/B microtiter filtration plate (Millipore, cat. no. MSFBN6B) with 210 µL total per well of the following: 35 µL assay buffer (20 mM HEPES, 100 mM NaCl, 1 mM EDTA, pH 7.4), 35 µL unlabeled toxin, 35 µL [³H]STX, and 105 µL receptor solution (rat brain homogenate, *Np*Sxph, *Rc*Sxph, or *Rc*Sxph E540A). Serial dilutions of unlabeled toxins were prepared in either 0.003 N hydrochloric acid (Fisher Chemical) or 20 µM acetic acid (J.T. Baker) based on the stock storage solution with the following in-well concentration ranges: STXdiHCl, dcSTX, GTX2/3, GTX5, GTX6, 0-1 µM/well; dcneoSTX, 0-5 µM; and GTX1/4, neoSTX, dcGTX2/3, C1/C2, 0-10 µM. The [³H]STX solution was prepared in cold assay buffer at a concentration of ~1 nM. Rat brain (Hilltop Lab Animals, Scottdale, PA) homogenate was prepared

1 October 2024

according to previously published protocols^{12, 13}, then was diluted in cold assay buffer to a concentration of ~30 µg/well in assay. Saxiphilin proteins (*Np*Sxph, *Rc*Sxph, and *Rc*Sxph E540A) were prepared in cold assay buffer at a concentration of 0.14 µg/well (~7.2 x 10⁻³ µM/well). Following a 1 h incubation at 4°C, wells were vacuum filtered and washed twice with 100 µL of cold assay buffer. Scintillation cocktail (Sigma Ultima Gold cat no. L8286; 50 µL/well) was then added and incubated at room temperature for 30 minutes. Radioactivity was measured (1 min/well) using a Perkin Elmer 2450 Microbeta microplate scintillation counter. Data were normalized using B/B_{max}, where B represents the bound [³H]STX in the sample and B_{max} represents the maximum binding of [³H]STX in the absence of competing unlabeled STX. The half maximal effective concentrations (EC₅₀) were calculated by fitting normalized counts per minute (CPM) as a function of toxin concentration using the nonlinear regression sigmoidal doseresponse (variable slope) curve fit with weight by 1/Y^2 and bottoms constrained to a shared value for all data sets in GraphPad Prism.

Isothermal titration calorimetry (ITC)

ITC measurements were performed at 25°C using a MicroCal PEAQ-ITC calorimeter (Malvern Panalytical) as described previously^{10,23}. *Np*Sxph was purified using a final SEC step in 150 mM NaCl, 10 mM HEPES, pH 7.4. 1 mM or 2 mM toxin stock in MilliQ water was diluted with the SEC buffer to prepare 100 μ M or 200 μ M toxin solutions having a final buffer composition of 135 mM NaCl, 9 mM HEPES, pH 7.4. To match buffers between the Sxph and toxin solutions, the purified protein samples were diluted to 10 μ M or 20 μ M with MilliQ water to reach a buffer concentration of 135 mM NaCl, 9 mM HEPES, pH 7.4. To match buffers between the Sxph and toxin solutions, the purified protein samples were diluted to 10 μ M or 20 μ M with MilliQ water to reach a buffer concentration of 135 mM NaCl, 9 mM HEPES, pH 7.4. Protein samples were filtered through a 0.22 μ m spin filter (Millipore, cat. no. UFC30GV00) before loading into the sample cell and titrated with toxin (100 μ M for GTX5, dcSTX, GTX2/3 and C1/C2, and 200 μ M for dcGTX2/3) using a schedule of 0.4 μ L titrant injection followed by 35 injections of 1 μ L for GTX5 and dcSTX, 18 injections of 2 μ L for dcGTX2/3, and 24 injections of 1.5 μ L for GTX2/3. The calorimetric experiment settings were: reference power, 5 μ cal/s; spacing between injections, 150 s; stir speed 750 rpm; and feedback mode, high. Data analysis was performed using MicroCal PEAQ-ITC Analysis Software (Malvern Panalytical) using a fitted offset for correcting for the heat of dilution and the single binding site model.

Crystallization and structure determination

*Np*Sxph was purified using a final SEC step in 30 mM NaCl, 10 mM HEPES, pH 7.4 as previously described¹⁰, and concentrated to 30-40 mg mL⁻¹ using a 50-kDa cutoff Amicon Ultra centrifugal filter unit (Millipore, cat. no. UFC505096). Toxin stocks of dcSTX, GTX2/3, dcGTX2/3, GTX5 and

1 October 2024

C1/C2 for co-crystallization experiments were prepared at 5 mM. For co-crystallization of *Np*Sxph with STX congeners, *Np*Sxph was mixed with each toxin in a molar ratio of 1:1.2 *Np*Sxph:toxin. Samples were incubated on ice for 1 hour before setting up crystallization trays. *Np*Sxph:dcSTX, *Np*Sxph:GTX2/3, *Np*Sxph:dcGTX2/3, *Np*Sxph:GTX5, *Np*Sxph:C1/C2 crystals were obtained by hanging drop vapor diffusion at 4°C from 400 nl drops (1:1 (v/v) ratio of protein and precipitant) set with Mosquito crystal (SPT Labtech) using 20-25% (v/v) PEG 400, 4-5% (w/v) PGA-LM, 100-200 mM sodium acetate, pH 5.0. All *Np*Sxph:toxin crystals were harvested and flash-frozen in liquid nitrogen without additional cryoprotectant.

X-ray datasets for all *Np*Sxph:STX congener complexes were collected at 100K at the Advanced Light Source (ALS) 8.3.1 beamline (Berkeley, CA), processed with XDS⁴⁴ and scaled and merged with Aimless⁴⁵. The *Np*Sxph:toxin structures were solved by molecular replacement using the apo-*Np*Sxph structure (PDB:8D6G) as a search model in Phaser from PHENIX⁴⁶. The electron density map and the model were manually checked in COOT⁴⁷ and iterative refinement was performed using phenix.refine⁴⁶. The quality of all models was assessed using MolProbity⁴⁸ and refinement statistics.

Molecular dynamics simulations

Atomic coordinates of NpSxph structures with STX (PDB:8D6M)¹⁰, dcSTX (PDB:8V68), GTX2 (PDB:8V69), dcGTX2 (PDB:8V65), GTX5 (PDB:8V66), and C1 (PDB:8V67), were used alongside those of crystallographic waters to prepare the corresponding topology and coordinate files using CHARMM-GUI⁴⁹. Missing residues in the protein structures (Lys174, Arg175, Lys645, Ala646) were modeled using Prime (Schrödinger Suite 2024.1)⁵⁰. Residues from the C-terminal 3C protease cleavage sites (Ser827, Asn828, Ser829) were removed, leaving the C-termini negatively charged, while the N-termini were capped with an acetyl group. All residues were maintained in their dominant protonation states at pH 7.0, except for His71, His125, Glu211, Glu214. Asp267. His337. Asp392. and His425. which were protonated. Disulfide bonds were introduced between the following cysteine pairs: Cys10-Cys45, Cys20-Cys36, Cys27-Cys418, Cys91-Cys113, Cys124-Cys131, Cys133-Cys155. Cys163-Cys185, Cys196-Cys203, Cys205-Cys227, Cys235-Cys826, Cys259-Cys342, Cys304-Cys317, Cys314-Cys325, Cys477-Cys509, Cys487-Cys500, Cys534-Cys821, Cys370-Cys384, Cys552-Cys781, Cys589-Cys667, Cys623-Cys637, Cys634-Cys650, and Cys707-Cys721.

A water box, extending 10 Å from the protein surface, was built around the protein. Random water molecules were replaced with sodium and chlorine ions to neutralize the systems and achieve a physiological concentration of 0.121 M NaCl. The AMBER force field ff19SB⁵¹ was employed for

1 October 2024

protein residue parameterization, and the OPC water mode I^{52} was applied. Ligands were parameterized based on crystal structure coordinates using Sage (OpenFF 2.1.0)⁵³. The topology and coordinate files for the system and ligands were manually combined, and overlapping water molecules were removed to generate the final systems. The final systems had dimensions of 122 × 122 × 122 Å³ and contained approximately 223,600 atoms, including 52,600 water molecules, 119 sodium ions, and 125 chloride ions.

Simulations were carried out using GROMACS 2024.2⁵⁴. Each simulation system underwent energy minimization, followed by equilibration in the constant temperature, constant volume (NVT) ensemble for 1 ns, with harmonic restraints of 1.0 kcal·mol⁻¹·Å² on the ligand atoms and protein heavy atoms, and 0.1 kcal·mol⁻¹·Å⁻² on protein hydrogen atoms. Five independent 1-µs production runs were conducted for each toxin-bound *Np*Sxph complex in the constant temperature and constant pressure (NPT) ensemble. Temperature control was implemented with a velocity-rescaling thermostat⁵⁵, using a 1.0 ps time constant and a reference temperature of 303.15 K, while pressure control used a stochastic cell rescaling isotropic barostat⁵⁶ with a 5.0 ps time constant, a reference pressure of 1 bar, and a compressibility of 4.5 × 10⁻⁵ bar⁻¹. The LINCS algorithm⁵⁷ was used to constrain the bond lengths of hydrogen atoms. Periodic boundary conditions were applied, with a 10.0 Å cutoff for Lennard-Jones and short-range electrostatic interactions. Long-range electrostatics were computed using the particle mesh Ewald (PME)⁵⁸ method with a fourth-order interpolation scheme and fast Fourier transform (FFT) grid spacing of 1.25 Å. A continuum model correction was applied to account for long-range van der Waals interactions. The equations of motion were integrated with a 2 fs time step.

The root-mean-square deviation (RMSD) of the binding pose for each toxin over time was calculated across the five replicas for each system using MDTraj $1.9.4^{59}$. The initial 10 ns of simulation time from each trajectory were excluded from the analysis. For this analysis, snapshots in the trajectories were aligned based on the C α atoms of the protein that were within 15 Å of the ligand's center of mass, considering all trajectories. The stability of the ligand core, assessed by the RMSD of atoms shared by all ligands relative to their initial structures, was compared across the different ligands. The conformational plasticity of the α 6C1- β 6C1 loop (residues Glu784-Asp791) was assessed by calculating the root mean square fluctuation (RMSF) of C α atoms with MDTraj 1.9.4⁵⁹, with error bars calculated as the standard deviation of the RMSF over the 5 trajectories for each toxin-bound *Np*Sxph complex.

Structural interaction fingerprints (SIFts) for each toxin-bound *Np*Sxph complex were calculated using PLIP v2.3.0⁶⁰ with default settings. The initial analysis focused on identifying salt bridges,

1 October 2024

hydrogen bonds, water-bridged hydrogen bonds, π -cation interactions, hydrophobic interactions, and π -stacking interactions. However, due to the low occurrence of π -cation, hydrophobic, and π -stacking interactions (present in less than 0.1% of all frames for any ligand-bound *Np*Sxph complex), these were excluded from further analysis. To simplify the representation of the remaining interactions, ligand moieties were categorized by their functional groups: carbamate (specific to GTX5, STX, GTX2/3, and C1/2), C12 hydrate (common to all toxins), C11 sulfate (specific to GTX2/3, C1/C2, and dcGTX2/3) *N*-sulfate (specific to GTX5 and C1/C2) and the STX core (encompassing the remaining atoms in all ligands). The protein residues involved in interactions were grouped, and their contributions to the fraction of the frames where interactions were present were summed.

The occupancy of Wat541, Wat795, and Wat786 near residues Glu541, Asp795, and Asp786 respectively, was determined by calculating the minimum distance between the water molecules and the C δ or C γ atoms of Glu541 or Asp795, respectively. Due to the flexibility of the α 6C1- β 6C1 loop, the occupancy of Wat786 near Asp786 was only detected when the toxin N7 was within ~4.2 Å of the C γ of Asp786. Therefore, it was calculated as a function of the minimum distance between the toxin N7 and the C γ of Asp786. Subsequently, the water occupancy probability for each trajectory was assessed as the fraction of frames where the calculated minimum distances were less than 4.5 Å. The results are presented as mean values along with the 25% and 75% quantiles.

GTX2/3 synthesis

The preparation of gonyautoxin 2/3 (GTX 2/3) follows previously published work^{33,61}. Synthesis of this toxin relies on a key oxidative dearomatization of a pyrrole ring to form the bis-guanidine tricyclic core that is common to all paralytic shellfish poisons. Subsequent functional group interconversion steps enable installation of the C11-alcohol and C12-ketone groups. A final step sequence involving reductive deprotection to liberate the two protected guanidine groups and C11-alcohol sulfation delivers the natural product as a single isomer, gonyautoxin 3. However, upon standing in buffered solution, epimizeration of the C11-sulfate occurs to give a 3:1 equilibrium mixture of GTX 2 and 3, consistent with the original isolation report. The structure of the synthetic material was validated by ¹H NMR and mass spectrometry, and its potency was confirmed through whole-cell electrophysiology recordings against Na_V1.4 expressed in CHO cells.

Automated patch-clamp electrophysiology

1 October 2024

Chinese hamster ovary (CHO) cells stably expressing the α -subunit of human Na_V1.4 sodium channel (*Hs*Na_V1.4) (B'SYS GmbH, cat. no. BSYS-NaV1.4-CHO-C) were cultured at 37°C in a humidified incubator with 5% CO₂ in complete culture medium containing: Ham's F-12 medium with GlutaMAX (Gibco, cat. no. 31765035) supplemented with 9% (v/v) heat inactivated fetal bovine serum (HI FBS) (Gibco, cat. no. 16140071), and antibiotics (0.9% (v/v) penicillin/streptomycin solution (Gibco, cat. no. 15-140-122) and 100 µg/mL Hygromycin B (Sigma-Aldrich, cat. no. 10843555001)).

Whole-cell patch-clamp experiments were performed using a QPatch Compact (Sophion Bioscience). The extracellular solution (ECS, saline) contained 2 mM CaCl₂, 1 mM MgCl₂, 10 mM HEPES, 4 mM KCl, 145 mM NaCl, 10 mM glucose (pH 7.4 with NaOH), and osmolarity adjusted to 305 mOsm/L with sucrose. The intracellular solution (ICS) contained 140 mM CsF, 1 mM/5 mM ethylene glycol-bis(β -aminoethyl ether)-N,N,N',N'-tetraacetic acid (EGTA)/CsOH, 10 mM HEPES, 10 mM NaCl (pH 7.3 with 3M CsOH), and osmolarity adjusted to 320 mOsm/L with sucrose. All solutions were kept at room temperature (23 ± 2°C) before application to the CHO-*Hs*Na_V1.4 cells. Cells were dissociated using Detachin (AMSBIO, cat. no. T100100) and kept in serum-free medium (Sigma-Aldrich, cat. no. C5467) supplemented with 25 mM HEPES and 0.04 mg/mL soybean trypsin inhibitor until recording. Cells were washed and resuspended in the ECS to reach a cell density of 4-6 x 10⁶ cells/mL shortly prior application to the QPatch Compact.

Sodium currents were elicited by a 60 ms depolarization step from -120 mV to 0 mV, with a holding potential of -120 mV and a sweep-to-sweep interval duration of 1.64 s. All experiments were conducted at room temperature (23 ± 2°C) using single-hole QPlates (Sophion Bioscience, cat. no. SB0201).

To determine toxin dose-response curves, toxin solutions were prepared in a 3-fold serial dilution series in ECS. Cumulative concentration-response experiments for STX, dcSTX, GTX2/3, dcGTX2/3, GTX5, and C1/C2 were performed by applying increasing toxin concentrations to each CHO-*Hs*Na_V1.4 cell. Peak currents were sampled at 10 kHz and filtered at 333 Hz using an 8th order Bessel filter with leak subtraction applied. IC₅₀ values for each toxin were calculated by fitting the dose-response curves (normalized peak current (I_x/I₀) as a function of toxin concentration) using the following equation: $I_x/I_0=(I_{max}-I_{min})/(1+x/IC_{50})$, where I_x is the current amplitude at the toxin concentration *x*, I₀ is the current amplitude in the latest saline period before toxin application, and I_{max} and I_{min} are the maximum and minimum peak current amplitudes, respectively, and IC₅₀ is the half-maximal inhibitory concentration.

1 October 2024

For the toxin neutralization experiment, NpSxph protein stock solution a buffer of 150 mM NaCl and 10 mM HEPES, pH 7.4 was diluted in ECS to 25 µM for the experiments with C1/C2 and GTX5, and to 5 µM for experiments with STX, dcSTX, GTX2/3, and dcGTX2/3. NpSxph was incubated with each toxin using 5:1 [NpSxph]:[toxin] molar ratio for at least 30 min prior application to the CHO-HsNav1.4 cells. Toxin solutions and NpSxph:toxin mixture were prepared in ECS. Baseline sodium currents were first recorded in saline, and then using toxin concentrations sufficient to block ~90% peak current as established from the dose-response experiments: 50 nM for STX, 200 nM for dcSTX, 60 nM for GTX2/3, 300 nM for dcGTX2/3, 1 µM for C1/C2, and 2.7 µM for GTX5. Following toxin wash-out, NpSxph:toxin mixture was applied to assess the effect of NpSxph on the toxin response. Following the NpSxph:toxin wash-out, toxin at the desired concentration (50 nM for STX, 200 nM for dcSTX, 900 nM for GTX2/3, 300 nM for dcGTX2/3, 1 µM for C1/C2, and 2.7 µM for GTX5) was applied to assess channel block. Peak currents were sampled at 25 kHz and filtered at 8333 Hz using an 8th order Bessel filter with leak subtraction applied. Normalized current was determined by using the following equation: $I = (I_{\text{Sxph:toxin}} -$ I_{Toxin})/(I_{ctrl} - I_{Toxin}), where I_{Sxph:toxin} is the current after application of NpSxph:toxin mixture, I_{Toxin} is the current after toxin application, and I_{Ctrl} is the basal current recorded in saline.

All data analyses were performed using the Sophion Analyzer Software (Sophion Bioscience) and GraphPad Prism (GraphPad Software).

Two-electrode voltage clamp

Two-electrode voltage clamp (TEVC) recordings were performed as described previously¹⁹. In brief, a pcDNA3.1+ vector containing *Phyllobates terribilis* SCN4A (*Pt*Na_V1.4) DNA (GenBank: MZ545381.1)¹⁹ was linearized with *Xba*I, and cRNA subsequently synthesized using mMESSAGE mMACHINE T7 Transcription Kit (Invitrogen, CA, USA). *Xenopus laevis* oocytes (harvested under UCSF IACUC protocol AN193390-011) were injected with 4–8 ng *Pt*Na_V1.4 and recordings were performed 1–2 days post injection at room temperature ($23 \pm 2^{\circ}$ C). Oocytes were impaled with borosilicate recording microelectrodes (0.3–2.0 M Ω resistance), backfilled with 3M KCI. Sodium currents were recorded using ND96 solution containing the following: 96 mM NaCl; 1 mM CaCl₂; 1 mM MgCl₂; 2 mM KCl; 5 mM HEPES (pH 7.5 with NaOH). Data were acquired using an Axoclamp 900B amplifier (Molecular Devices, CA, USA) controlled by pClamp software (v10.9, Molecular Devices), digitized at 10 kHz using an Axoclamp 1550B digitizer (Molecular Devices).

The effects of toxins and Sxph-toxin combinations on $PtNa_V1.4$ were assessed using a 60-ms depolarization step from -120 to 0 mV, with holding potential of -120 mV and sweep-to-sweep duration of 10 s. Leak currents were subtracted using a P/4 protocol during data acquisition. The

1 October 2024

half maximal inhibitory concentration (IC_{50}) values for STX congeners against $PtNa_V1.4$ were determined by perfusing increasing concentrations of toxin in series and calculated using nonlinear regression analysis in Prism (v10.1, GraphPad, MA, USA). Toxin concentrations required to inhibit 90% of the current (IC_{90}) were calculated from the determined IC_{50} and Hill

slope (*H*) according to the following equation: $IC_x = \left(\frac{x}{100-x}\right)^{1/H} \times IC_{50}$.

Sxph rescue experiments were conducted by first recording baseline currents in ND96, then perfusing sufficient toxin to give ~90% block (100 nM STX, 200 nM GTX2/3, or 800 nM dcSTX). Sxph was then applied directly into the 1-mL recording chamber. For all saxiphilin:toxin ratios, the concentration of Sxph stock solution was adjusted such that the added Sxph solution was less than 1% of the total recording chamber volume. Rescue responses were normalized according to the following equation: $I_{norm} = \frac{(I_x - I_{min})}{(I_{max} - I_{min})}$, where I_x is the current following addition of Sxph at ratio *x*, I_{min} is the current following toxin block, and I_{max} was the maximal current observed in recording solution alone. Toxin rescue kinetics were modelled using non-linear regression of normalized currents over time, and quality of model fit was assessed using Akaike's Information Criterion, corrected for small sample size (AICc). Data analysis was performed using Clampfit (v11.0, Molecular Devices) and Prism (v10.1, GraphPad).

1 October 2024

Supplementary material

Supplementary material contains Figures S1-S8, Tables S1-S5, Movie S1, and Supplementary References.

1 October 2024

References

- 1. Wiese, M., D'Agostino, P.M., Mihali, T.K., Moffitt, M.C. & Neilan, B.A. Neurotoxic alkaloids: saxitoxin and its analogs. *Mar Drugs* **8**, 2185-211 (2010).
- 2. Munday, R., Thomas, K., Gibbs, R., Murphy, C. & Quilliam, M.A. Acute toxicities of saxitoxin, neosaxitoxin, decarbamoyl saxitoxin and gonyautoxins 1&4 and 2&3 to mice by various routes of administration. *Toxicon* **76**, 77-83 (2013).
- 3. Duran-Riveroll, L.M. & Cembella, A.D. Guanidinium Toxins and Their Interactions with Voltage-Gated Sodium Ion Channels. *Mar Drugs* **15**(2017).
- 4. Alonso, E., Alfonso, A., Vieytes, M.R. & Botana, L.M. Evaluation of toxicity equivalent factors of paralytic shellfish poisoning toxins in seven human sodium channels types by an automated high throughput electrophysiology system. *Arch Toxicol* **90**, 479-88 (2016).
- 5. Thottumkara, A.P., Parsons, W.H. & Du Bois, J. Saxitoxin. *Angew Chem Int Ed Engl* **53**, 5760-84 (2014).
- 6. Anderson, D.M. et al. Evidence for massive and recurrent toxic blooms of Alexandrium catenella in the Alaskan Arctic. *Proc Natl Acad Sci U S A* **118**(2021).
- 7. Dai, Y. et al. Coastal phytoplankton blooms expand and intensify in the 21st century. *Nature* **615**, 280-284 (2023).
- 8. Shen, H., Liu, D., Wu, K., Lei, J. & Yan, N. Structures of human Nav1.7 channel in complex with auxiliary subunits and animal toxins. *Science* **363**, 1303-1308 (2019).
- 9. Shen, H. et al. Structural basis for the modulation of voltage-gated sodium channels by animal toxins. *Science* **362**(2018).
- 10. Chen, Z. et al. Definition of a saxitoxin (STX) binding code enables discovery and characterization of the anuran saxiphilin family. *Proc Natl Acad Sci U S A* **119**, e2210114119 (2022).
- 11. Yen, T.-J., Lolicato, M., Thomas-Tran, R., Du Bois, J. & Minor, D.L., Jr. Structure of the Saxiphilin:saxitoxin (STX) complex reveals a convergent molecular recognition strategy for paralytic toxins. *Sci Adv* **5**(2019).
- 12. Van Dolah, F.M., Fire, S.E., Leighfield, T.A., Mikulski, C.M. & Doucette, G.J. Determination of paralytic shellfish toxins in shellfish by receptor binding assay: collaborative study. *J* AOAC Int **95**, 795-812 (2012).
- 13. International, A. Official Method 2011.27. in *Official methods of analysis of AOAC International* (ed. Latimer, G.W.) (AOAC International, Gaithersburg, MD, 2012).
- 14. Llewellyn, L.E. Saxitoxin, a toxic marine natural product that targets a multitude of receptors. *Nat Prod Rep* **23**, 200-22 (2006).
- 15. Dillon, M. et al. Current Trends and Challenges for Rapid SMART Diagnostics at Point-of-Site Testing for Marine Toxins. *Sensors (Basel)* **21**(2021).
- 16. Gawankar, S., Masten, S.J. & Lahr, R.H. Review of the occurrence, treatment technologies, and detection methods for saxitoxins in freshwaters. *J Water Health* **22**, 1472-1490 (2024).
- 17. Mahar, J., Lukacs, G.L., Li, Y., Hall, S. & Moczydlowski, E. Pharmacological and biochemical properties of saxiphilin, a soluble saxitoxin-binding protein from the bullfrog (Rana catesbeiana). *Toxicon* **29**, 53-71 (1991).

- 18. Penzotti, J.L., Lipkind, G., Fozzard, H.A. & Dudley, S.C., Jr. Specific neosaxitoxin interactions with the Na+ channel outer vestibule determined by mutant cycle analysis. *Biophys J* **80**, 698-706 (2001).
- 19. Abderemane-Ali, F. et al. Evidence that toxin resistance in poison birds and frogs is not rooted in sodium channel mutations and may rely on "toxin sponge" proteins. *J Gen Physiol* **153**(2021).
- 20. Prinzmetal, M., Sommer, H. & Leake, C.D. The pharmacological action of "mussel poison". *J Pharmacol Exp Ther* **46**, 63-73 (1932).
- 21. Kao, C.Y. & Fuhrman, F.A. Differentiation of Actions of Tetrodotoxin and Saxitoxin. *Toxicon* **5**, 25-34 (1967).
- 22. Llewellyn, L.E., Bell, P.M. & Moczydlowski, E.G. Phylogenetic survey of soluble saxitoxinbinding activity in pursuit of the function and molecular evolution of saxiphilin, a relative of transferrin. *Proc Biol Sci* **264**, 891-902 (1997).
- 23. Chen, Z., Zakrzewska, S., Hajare, H.S., Du Bois, J. & Minor, D.L., Jr. Expression, purification, and characterization of anuran saxiphilins using thermofluor, fluorescence polarization, and isothermal titration calorimetry. *STAR Protoc* **5**, 102792 (2023).
- 24. Caty, S.N. et al. Molecular physiology of chemical defenses in a poison frog. *J Exp Biol* **222**(2019).
- 25. O'Connell, L.A. et al. Rapid toxin sequestration modifies poison frog physiology. *J Exp Biol* **224**(2021).
- 26. Alvarez-Buylla, A. et al. Binding and sequestration of poison frog alkaloids by a plasma globulin. *Elife* **12**(2023).
- 27. Morabito, M.A. & Moczydlowski, E. Molecular cloning of bullfrog saxiphilin: a unique relative of the transferrin family that binds saxitoxin. *Proc Natl Acad Sci U S A* **91**, 2478-82 (1994).
- 28. Rossi, A.M. & Taylor, C.W. Analysis of protein-ligand interactions by fluorescence polarization. *Nat Protoc* **6**, 365-87 (2011).
- 29. Ruberu, S.R. et al. Receptor binding assay for the detection of paralytic shellfish poisoning toxins: comparison to the mouse bioassay and applicability under regulatory use. *Food Addit Contam Part A Chem Anal Control Expo Risk Assess* **35**, 144-158 (2018).
- Llewellyn, L.E. & Moczydlowski, E.G. Characterization of saxitoxin binding to saxiphilin, a relative of the transferrin family that displays pH-dependent ligand binding. *Biochemistry* 33, 12312-22 (1994).
- 31. Fox, J.M., Zhao, M., Fink, M.J., Kang, K. & Whitesides, G.M. The Molecular Origin of Enthalpy/Entropy Compensation in Biomolecular Recognition. *Annu Rev Biophys* **47**, 223-250 (2018).
- 32. Chodera, J.D. & Mobley, D.L. Entropy-enthalpy compensation: role and ramifications in biomolecular ligand recognition and design. *Annu Rev Biophys* **42**, 121-42 (2013).
- 33. Mulcahy, J.V. & Du Bois, J. A stereoselective synthesis of (+)-gonyautoxin 3. *Journal of the American Chemical Society* **130**, 12630-+ (2008).
- 34. Moczydlowski, E., Hall, S., Garber, S.S., Strichartz, G.S. & Miller, C. Voltage-dependent blockade of muscle Na+ channels by guanidinium toxins. *J Gen Physiol* **84**, 687-704 (1984).

- 35. Choudhary, G., Shang, L., Li, X. & Dudley, S.C., Jr. Energetic localization of saxitoxin in its channel binding site. *Biophys J* **83**, 912-9 (2002).
- 36. Thomas-Tran, R. & Du Bois, J. Mutant cycle analysis with modified saxitoxins reveals specific interactions critical to attaining high-affinity inhibition of hNaV1.7. *Proc Natl Acad Sci U S A* **113**, 5856-61 (2016).
- 37. Walker, J.R. et al. Marked difference in saxitoxin and tetrodotoxin affinity for the human nociceptive voltage-gated sodium channel (Nav1.7) [corrected]. *Proc Natl Acad Sci U S A* **109**, 18102-7 (2012).
- 38. Penzotti, J.L., Fozzard, H.A., Lipkind, G.M. & Dudley, S.C., Jr. Differences in saxitoxin and tetrodotoxin binding revealed by mutagenesis of the Na+ channel outer vestibule. *Biophys J* **75**, 2647-57 (1998).
- 39. AOAC. Official Method 959.08 paralytic shellfish poison, biological method. in *Official Methods of Analysis*, Vol. 1 (Arlington, VA, 1999).
- 40. Sommer, H. & Meyer, K.F. Paralytic shell-fish poisoning. *Arch. Pathol.* 24, 560-598 (1937).
- 41. Andresen, B.M. & Du Bois, J. De novo synthesis of modified saxitoxins for sodium ion channel study. *J Am Chem Soc* **131**, 12524-5 (2009).
- 42. Ondrus, A.E. et al. Fluorescent saxitoxins for live cell imaging of single voltage-gated sodium ion channels beyond the optical diffraction limit. *Chem Biol* **19**, 902-12 (2012).
- 43. Huynh, K. & Partch, C.L. Analysis of protein stability and ligand interactions by thermal shift assay. *Curr Protoc Protein Sci* **79**, 28 9 1-14 (2015).
- 44. Kabsch, W. Xds. Acta Crystallogr D Biol Crystallogr 66, 125-32 (2010).
- 45. Evans, P.R. & Murshudov, G.N. How good are my data and what is the resolution? *Acta Crystallogr D Biol Crystallogr* **69**, 1204-14 (2013).
- 46. Adams, P.D. et al. PHENIX: a comprehensive Python-based system for macromolecular structure solution. *Acta Crystallogr D Biol Crystallogr* **66**, 213-21 (2010).
- 47. Emsley, P. & Cowtan, K. Coot: model-building tools for molecular graphics. *Acta Crystallogr D Biol Crystallogr* **60**, 2126-32 (2004).
- 48. Williams, C.J. et al. MolProbity: More and better reference data for improved all-atom structure validation. *Protein Sci* **27**, 293-315 (2018).
- 49. Jo, S., Kim, T., Iyer, V.G. & Im, W. CHARMM-GUI: a web-based graphical user interface for CHARMM. *J Comput Chem* **29**, 1859-65 (2008).
- 50. Jacobson, M.P. et al. A hierarchical approach to all-atom protein loop prediction. *Proteins* **55**, 351-67 (2004).
- 51. Tian, C. et al. ff19SB: Amino-Acid-Specific Protein Backbone Parameters Trained against Quantum Mechanics Energy Surfaces in Solution. *J Chem Theory Comput* **16**, 528-552 (2020).
- 52. Izadi, S., Anandakrishnan, R. & Onufriev, A.V. Building Water Models: A Different Approach. *J Phys Chem Lett* **5**, 3863-3871 (2014).
- 53. Boothroyd, S. et al. Development and Benchmarking of Open Force Field 2.0.0: The Sage Small Molecule Force Field. *J Chem Theory Comput* **19**, 3251-3275 (2023).
- 54. Abraham, M.J. et al. GROMACS: High performance molecular simulations through multilevel parallelism from laptops to supercomputers. *SoftwareX* **1-2**, 19-25 (2015).

- 55. Bussi, G., Donadio, D. & Parrinello, M. Canonical sampling through velocity rescaling. *J Chem Phys* **126**, 014101 (2007).
- 56. Bernetti, M. & Bussi, G. Pressure control using stochastic cell rescaling. *J Chem Phys* **153**, 114107 (2020).
- 57. Hess, B., Bekker, H., Berendsen, H.J.C. & Fraaije, J.G.E.M. LINCS: A linear constraint solver for molecular simulations. *Journal of Computational Chemistry* **18**, 1463-1472 (1997).
- 58. Darden, T., York, D. & Pedersen, L. Particle Mesh Ewald an N.Log(N) Method for Ewald Sums in Large Systems. *Journal of Chemical Physics* **98**, 10089-10092 (1993).
- 59. McGibbon, R.T. et al. MDTraj: A Modern Open Library for the Analysis of Molecular Dynamics Trajectories. *Biophys J* **109**, 1528-32 (2015).
- 60. Salentin, S., Schreiber, S., Haupt, V.J., Adasme, M.F. & Schroeder, M. PLIP: fully automated protein-ligand interaction profiler. *Nucleic Acids Res* **43**, W443-7 (2015).
- 61. Mulcahy, J.V., Walker, J.R., Merit, J.E., Whitehead, A. & Du Bois, J. Synthesis of the Paralytic Shellfish Poisons (+)-Gonyautoxin 2, (+)-Gonyautoxin 3, and (+)-11,11-Dihydroxysaxitoxin. *J Am Chem Soc* **138**, 5994-6001 (2016).

1 October 2024

Table 1 NpSxph:Toxin thermodynamic binding parameters								
		N (sites)	Kd (nM)	∆H (kcal mol⁻¹)	ΔS (cal mol ⁻¹ K ⁻¹)	∆G (kcal mol-1)	∆∆G (kcal mol-1)	n
NpSxph	STX*	0.92 ± 0.02	2.5 ± 0.1	-18.7 ± 0.2	-23.2 ± 0.8	-11.8 ± 0.1	-	2
	dcSTX	0.81 ± 0.01	11.3 ± 4.0	-17.8 ± 0.4	-23.3 ± 2.0	-10.9 ± 0.2	0.9	2
	GTX2/3	1.11 ± 0.15	78.2 ± 18.1	-11.9 ± 1.3	-7.3 ± 4.1	-9.7 ± 0.1	2.1	2
	dcGTX2/3	1.00 ± 0.13	295 ± 74	-12.4 ± 0.0	-11.7 ± 0.4	-8.9 ± 0.1	2.9	2
	GTX5	0.90 ± 0.06	1.8 ± 0.1	-18.0 ± 0.8	-20.2 ± 3.5	-12.0 ± 0.3	-0.2	2
	C1/C2	0.86 ± 0.10	274 ± 71	-14.3 ± 2.2	-17.7 ± 0.8	-9.0 ± 0.2	2.8	2

'N' is the number of binding sites.

Kd denotes the dissociation constant.

 $\Delta \Delta G = \Delta G_{\text{Toxin}} - \Delta G_{\text{STX}}$

'n' is the number of observations.

* data from ¹⁰

Errors are S.D.

1 October 2024

Figure 1

Zakrzewska et al.



Figure 1 TF assays reveal Sxph gonyautoxin binding preferences. **A**, STX core showing modification sites R1 (magenta), R2 (yellow), and R3 (green). **B**, and **C**, Exemplar TF assay results for **B**, *Rc*Sxph and **C**, *Np*Sxph in the presence of STX, GTX5, C1/C2, dcGTC2/3, GTX2/3, or neoSTX at 0 nM (black) 19.5 nM (blue), 625 nM (cyan), 5,000 nM (orange), and 20,000 nM (red) of each toxin. **D**, and **E**, Concentration dependence of Δ Tm for **D**, *Rc*Sxph and **E**, *Np*Sxph for STX (black circles), dcSTX (blue circles), GTX5 (green triangles), GTX2/3 (orange inverted triangles), dcGTX2/3 (inverted magenta triangles), C1/C2 (purple diamonds), neoSTX (dark red squares). Error bars are S.E.M.



Figure 2 FP and RBA studies reveal Sxph gonyautoxin binding affinities. A, and B, Exemplar competition FP assays for A, *Rc*Sxph and B, *Np*Sxph for STX (black circles), dcSTX (blue circles), GTX5 (green triangles), GTX2/3 (orange inverted triangles), dcGTX2/3 (inverted magenta triangles), C1/C2 (purple diamonds), and neoSTX (dark red squares). C, and D, Comparisons of Δ Tm and $\Delta\Delta$ G values for C, *Rc*Sxph and D, *Np*Sxph. (line y = -0.7564x + 3.6 R² = 0.9668 and y =-0.7679x + 3.193 R² = 0.9485, respectively). E, Comparison of $\Delta\Delta$ G relative to STX for binding

1 October 2024

of the indicated toxins to *Np*Sxph and *Rc*Sxph (line, $y = 1.356x + 0.08794 R^2 = 0.9417$). Grey dashed line shows x=y. **F**, and **G**, Exemplar RBA curves for **F**, *Rc*Sxph and **B**, *Np*Sxph for STX (black circles), dcSTX (blue circles), GTX5 (green triangles), GTX2/3 (orange inverted triangles), dcGTX2/3 (inverted magenta triangles), C1/C2 (purple diamonds), and neoSTX (dark red squares) where B/B_{max} represents the bound [³H]STX in the sample/maximum binding of [³H]STX in the absence of competing unlabeled STX. **H**, and **I**, Comparison of binding free energies measured by FP and RBA assays for **H**, *Rc*Sxph and **I**, *Np*Sxph. (line y= 0.7331x - 2.878 R² = 0.9755 and y = 0.89314x - 0.8046 R² = 0.9604, respectively). Grey dashed line shows x=y. Error bars are S.E.M.



```
Zakrzewska et al.
```



Figure 3 ITC studies reveal *Np*Sxph STX congener binding thermodynamic properties. **A-E**, Exemplar ITC isotherms for the titration of **A**, 100 μ M dcSTX into 11.4 μ M *Np*Sxph **B**, 100 μ M GTX2/3 into 11.3 μ M *Np*Sxph **C**, 200 μ M dcGTX2/3 into 22.4 μ M *Np*Sxph **D**, 100 μ M GTX5 into 12.4 μ M *Np*Sxph and **E**, 100 μ M C1/C2 into 11.9 μ M *Np*Sxph. Kd and Δ H values are indicated. Toxin structures are shown. **F**, Comparison of Δ G_{ITC} and Δ G_{FP} for *Np*Sxph binding to each of the indicated toxins (line y =1.287x + 2.263, R² = 0.9287). Grey dashed line shows x=y.





1 October 2024

congener structures. Structures are grouped by STX modification type: Carbamoyl modification, dcSTX (yellow) and GTX5(B1) (magenta); Carbamoyl and C11 modification, dcGTX2 (cyan) and C1 (marine); and C11 modification, GTX2 (light pink). **H**, Comparison of the STX congeners from 'F' with STX from the *Np*Sxph:STX complex (PDB:8D6M)¹⁰. STX congener colors are as in 'F'. Modification sites R1 (carbamoyl) (magenta) and R3 (C11) (green) are highlighted. **I**, Electrostatic surface potentials for *Np*Sxph:STX complex (PDB:8D6M)¹⁰ and the indicated *Np*Sxph:STX congener complexes. STX and STX congeners are shown in space filling.



Figure 5 *Np***Sxph rescues** *Pt***N**_a**v1.4 block by diverse STX congeners. A** and **B**, Exemplar TEVC recordings of *Pt***N**_a**v1.4** before (Ctrl, black) and after (red) application of 100 nM STX, and after application of Sxph in the presence of 100 nM STX at the indicated [Sxph]:[STX] molar ratios for **A**, *Np*Sxph (teal) and **B**, *Rc*Sxph (orange). **C**, [*Np*Sxph]:[STX] (teal) and [*Rc*Sxph]:[STX] (orange) dose-response curves in the presence of 100 nM STX. **D**, Exemplar TEVC recordings of *Pt*Nav1.4 expressed in *X. laevis* oocytes in the presence of the indicated STX congener (dcSTX, GTX5, and GTX2/3) concentrations. Inset shows the voltage pulse protocol. **E**, Toxin response curves for STX (black circles, from Abderemane-Ali et al., 2021), GTX2/3 (purple triangles), dcSTX (maroon squares), and GTX5 (magenta inverted triangles). **F** and **G**, Exemplar *Pt*Nav1.4 responses to application of **F**, 200 nM GTX2/3 (red) (30 minutes post application), **G**, 800 nM dcSTX (red) (one minute post application), and 3:1 [*Np*Sxph]:[toxin] application (teal). For C and E, data represents mean with S.E.M. from 6–8 oocytes.



Figure 6 *Np***Sxph rescues** *Hs***Na**_V**1.4** from block by diverse STX congeners. **A**, Exemplar whole-cell patch clamp recordings of *Hs*Na_V1.4 expressed in CHO cells in the presence of the indicated STX congener (STX, dcSTX, GTX2/3, dcGTX2/3, GTX5, and C1/C2) concentrations. **B**, Toxin dose-response curves for STX (black circles), dcSTX (blue circles), GTX5 (green triangles), GTX2/3 (orange inverted triangles), dcGTX2/3 (magenta inverted triangles), and C1/C2 (purple diamonds). **C**, Exemplar whole-cell patch clamp *Hs*Na_V1.4 responses in the absence (Ctrl, black) and presence (red) of 50 nM STX, 200 nM dcSTX, 60 nM GTX2/3, 300 nM dcGTX2/3, 1

1 October 2024

 μ M C1/C2, and 2.7 μ M GTX5, and then after 5:1 [*Np*Sxph]:[Toxin] application (teal). **D**, Exemplar whole-cell patch clamp time course showing *Hs*Na_V1.4 peak currents from 'C' after application of STX congener (50 nM STX (grey), 200 nM dcSTX (blue), 60 nM GTX2/3 (orange), 300 nM dcGTX2/3 (pink), 1 μ M C1/C2 (purple), and 2.7 μ M GTX5 (green)), followed by toxin wash-out, then application of 5:1 [*Np*Sxph]:[Toxin] mixture (teal). Following the [*Np*Sxph]:[Toxin] wash-out, toxin at the desired concentration (50 nM STX (grey), 200 nM dcSTX (blue), 60 nM GTX2/3 (orange), 300 nM dcGTX2/3 (pink), 1 μ M C1/C2 (purple), and 2.7 μ M GTX5 (green)) was applied. **E**, Fraction recovery for 5:1 *Np*Sxph:toxin. Colors as in 'D'. For B and E, error bars are S.E.M.



Figure 7 Structural comparison of STX binding site in Nav1.7 and *Np***Sxph. A**, Top view of the STX binding site in Nav1.7 structure in complex with STX (PDB:6J8G)⁸. **Left**, Comparison of STX (orange) with C1 (blue) shown in stick models. Modification sites for C1 are highlighted, sulfate on the carbamate (Cb, red) and sulfate on the C11 position (C11, blue). **Right**, C1 shown in space filling. **B**, Schematic of C1 model from 'A'. Clash site in DIII is shaded. **C**, Toxin binding site in *Np*Sxph:C1 complex (PDB:8V67). C1 shown in space filling. **D**, Schematic of *Np*Sxph:C1 complex from 'C'.