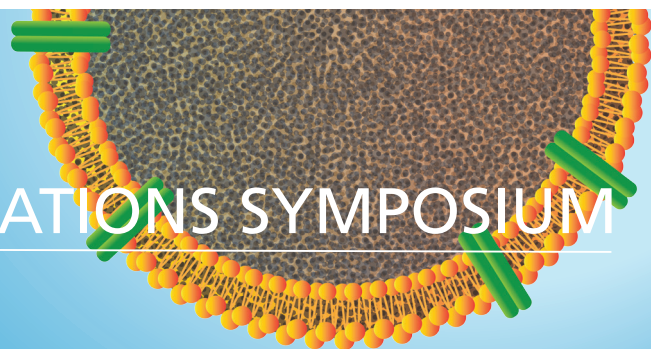


# ION CHANNEL MODULATIONS SYMPOSIUM

5 & 6 November 2022

Venue: New York University (the Schwartz Lecture Hall and Lobby),  
Entrance 435 E. 30<sup>th</sup> Science Building, USA



## SPONSORSHIPS

The Ion Channel Modulation Symposium (USA) is a joint event hosted by **Sophion Bioscience** and **New York University**. The event will be held at New York University (the Schwartz Lecture Hall and Lobby), Entrance 435 E. 30<sup>th</sup> Science Building, USA on the **5 & 6 of November 2022**.

We have designed three sponsorship packages:

### Platinum Sponsorship (Cost \$5,500):

- Company name and logo on the ICMS2022 USA website and circulated materials
- A 'table top' exhibition space
- 'Zap Stands' / roll-ups (one per stand) can be used behind the tables (not 'nomadic' style)
- Full symposium attendance for two persons
- Invitation to the Symposium reception

### Gold Sponsorship (Cost \$3,200):

- Company name and logo on the ICMS2022 USA website and circulated materials
- A 'table top' exhibition space
- 'Zap Stands' / roll-ups (one per stand) can be used behind the tables (not 'nomadic' style)
- Full symposium attendance for one person
- Invitation to the Symposium reception

### Silver Sponsorship (Cost \$550):

- Company name and logo on the ICMS2022 USA website
- Option to have literature available to the conference attendees on a shared table top in the exhibition area

The number of sponsorship opportunities for the ICMS2022 USA is limited to so please apply early to avoid disappointment.

Please fill in the below form and send it by mail to [duncan.jarman@sophion.com](mailto:duncan.jarman@sophion.com). Once the application has been approved we will inform you by mail. We will need a purchase order from you in order for us to issue an invoice. Your sponsorship is not secured until we have received the your payment. We reserve the right to reject any sponsorship applicant.

Company	
Address 1	
City and Post code	
Contact person	
Mail address contact person	
Phone no.	
Type of sponsorship	Platinum: <input type="checkbox"/> Gold: <input type="checkbox"/> Silver: <input type="checkbox"/>
PO number:	
Signature & date:	