

Application Report

Pharmacological evaluation of GABA_A receptor subtypes on Qube 384

High-throughput screening and cumulative concentration-response relationship of a GABA_A($\alpha_5\beta_3\gamma_2$) cell line. Characterization of GABA_A receptor subtypes

Summary

Studies of GABA_A ion channels using the automated patch clamp platform Qube 384 with focus on:

- Short ligand exposure with repetitive stimulations with EC₅₀ concentrations of GABA
- Effects of agonists, antagonists and modulators
- Cumulative and non-cumulative concentration-response relationships
- Characterizing the pharmacological properties of four cell lines expressing different GABA_A subtypes

Introduction

γ -aminobutyric acid (GABA) is the major inhibitory neurotransmitter of the central nervous system and pathophysiological changes in GABA signalling is the leading cause in a large group of neurological and psychiatric disorders including epilepsy, schizophrenia and depression. Consequently, manipulation of the GABA signalling holds a great therapeutic potential¹⁻⁴.

Ionotropic GABA receptors consist of 5 membrane-spanning subunits^{5,6}, of which 19 different have been identified in humans (α_{1-6} , β_{1-3} , γ_{1-3} , δ , ϵ , θ , π , ρ_{1-3}). Receptors in different cell types have different subunit composition and differ both in pharmacology and subcellular location⁷. Due to the heterogeneity of GABA_A receptors, they can perform different inhibitory tasks: Positive modulation of GABA_A receptors can be anticonvulsant, hypnotic, anaesthetic and anxiolytic, whereas negative modulation can enhance cognition but also be anxiogenic and proconvulsant¹⁻⁴.

Here we use four GABA_A/HEK293 cell lines and a range of tool compounds to demonstrate two different approaches to GABA_A

receptor evaluation on Qube. We show 1) a cumulative dose-response experiment with a GABA_A($\alpha_5\beta_3\gamma_2$) cell line and 2) a GABA_A receptor subtype screen, using a single compound plate layout fitting various receptor subtypes. This compound plate enables evaluation of both non-cumulative concentration-response relationships of GABA and bicuculline (competitive antagonist) and the potentiation with diazepam (positive allosteric modulator), simultaneously.

Results and discussion

Short agonist exposure, GABA_A($\alpha_5\beta_3\gamma_2$)

Desensitization due to prolonged or repeated agonist exposure poses a challenge when studying GABA_A receptors in an in vitro setting. Therefore, the effect of GABA on the GABA_A($\alpha_5\beta_3\gamma_2$) receptor was evaluated on Qube, employing the stacked delivery feature, where both the GABA-containing solution and the wash-out solution are stacked in the pipette. In this way, the exposure time is reduced to less than one second.

With the stacked delivery feature, consecutive applications of GABA (12 μ M) could be made without significant rundown (Figure 1A). 12 μ M GABA, which is close to the EC₅₀ value, elicited on average 19.6 nA (\pm 5.7 nA, SD) current pr. site using a multihole (x10 patch holes) QChip (see Figure 1B for a QChip view).

The success rate was up to 88% per plate with the following criteria:

Resistance > 100 M Ω per cell

Capacitance > 5 pF per cell

Current amplitude at 12 μ M GABA > 300 pA per cell

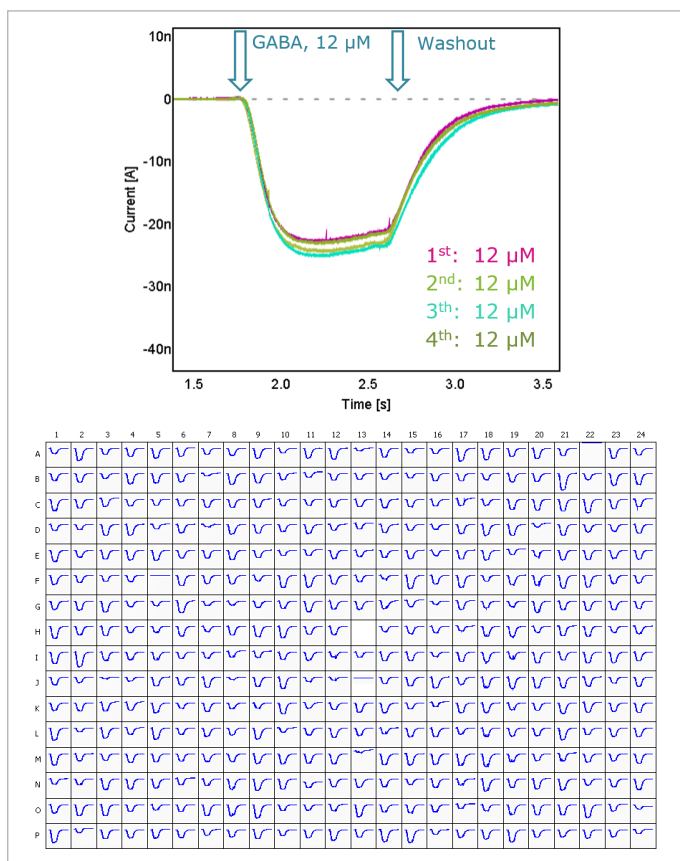


Fig. 1: Top: 4 consecutive applications of 12 μM GABA using the stacked delivery feature. Bottom: Plate view of a QChip, showing the 384 individual responses to 12 μM GABA.

Cumulative GABA concentration response

The concentration-response relationship of GABA on the $\text{GABA}_A(\alpha_5\beta_3\gamma_2)$ receptor was evaluated on Qube. GABA was applied in increasing concentrations (3-fold dilution from 400 μM , Figure 2). The EC_{50} value for the cumulative concentration response was found to be 10.9 μM (CI95%: 10.0 to 12.4 μM) and the Hill slope was 1.5 (± 0.1 , SD).

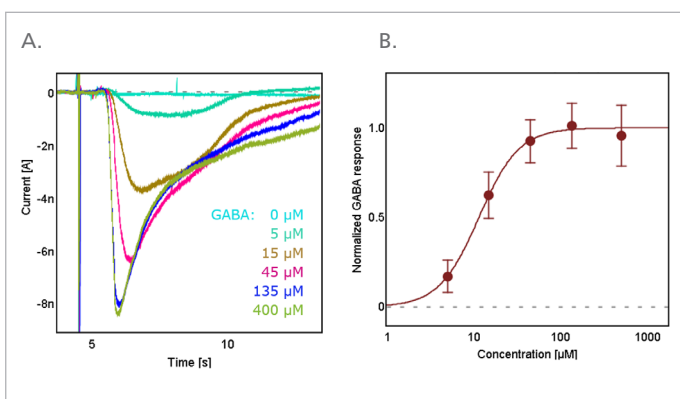


Fig. 2: Cumulative concentration-response relationship of GABA binding to the $\text{GABA}_A(\alpha_5\beta_3\gamma_2)$ receptor. A) Typically recorded currents in response to increasing concentrations of GABA. B) Peak current (normalized to highest value in experiment, average \pm SD, $n=192$) as a function of GABA concentration for the whole QChip.

GABA_A receptor subtype screen

The GABA_A receptor subtypes differ both in pharmacology and physiological function. We designed a compound plate layout (Figure 3) for the screen, using one compound plate, four QChips and four cell lines to evaluate the pharmacological properties of four GABA_A receptor types ($\alpha_5\beta_3\gamma_2$, $\alpha_1\beta_2\gamma_2$, $\alpha_2\beta_3\gamma_2$ and $\alpha_4\beta_3\delta$). The experiment results included non-cumulative GABA and bicuculline (competitive antagonist) concentration-response experiments. In addition, the potentiation of diazepam (positive allosteric modulator) was evaluated at different concentrations.

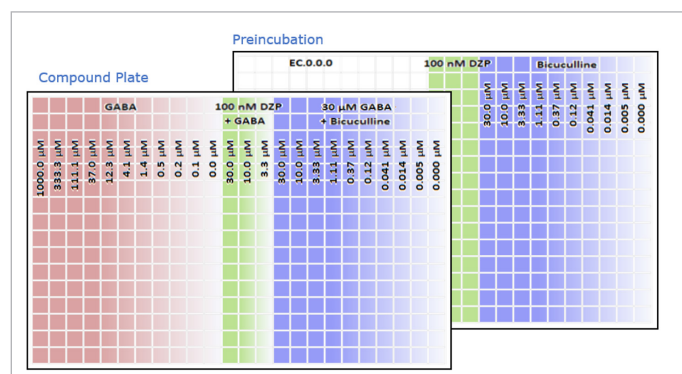


Fig. 3: Compound plate layout: 100 nM diazepam and varying concentrations of bicuculline were washed in (pre-incubation) prior to GABA application (compound plate).

The results of the screen are displayed in Figure 4-7 for the four different receptor types ($\alpha_5\beta_3\gamma_2$, $\alpha_1\beta_2\gamma_2$, $\alpha_2\beta_3\gamma_2$ and $\alpha_4\beta_3\delta$).

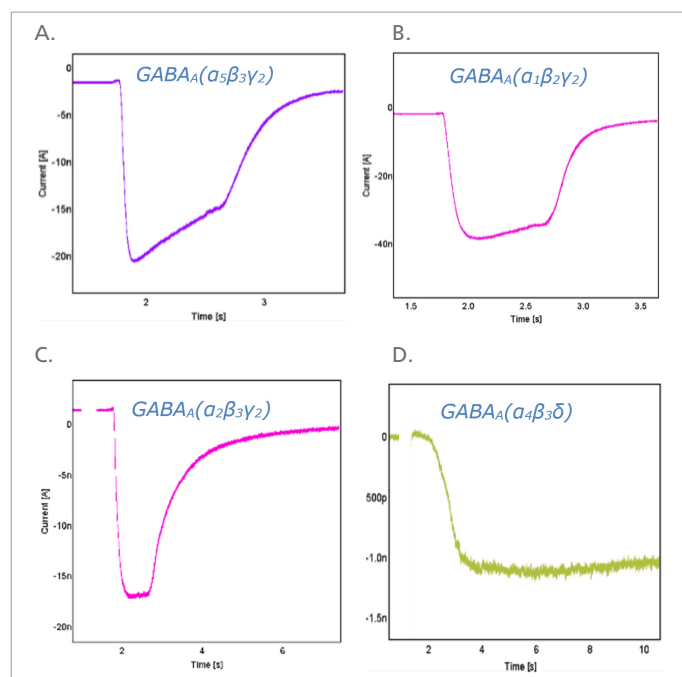


Fig. 4: GABA_A receptor kinetics. Typical recordings from one multihole site (10 cells/site). The effect of GABA was evaluated employing the stacked delivery feature, with an exposure time of less than one second (with exception from $\text{GABA}_A(\alpha_4\beta_3\delta)$ where a regular and not stacked delivery feature was employed due to the slow kinetics of δ -subunit containing GABA_A receptors). A) $\text{GABA}_A(\alpha_5\beta_3\gamma_2)$: On average, 111 μM GABA elicited a 26.3 nA (± 8.5 nA) response (peak current). B) $\text{GABA}_A(\alpha_1\beta_2\gamma_2)$: On average, 111 μM GABA elicited a 31.6 nA (± 8.3) response (peak current). C) $\text{GABA}_A(\alpha_2\beta_3\gamma_2)$: On average, 1 mM GABA elicited a 9.6 nA (± 4.1) response (peak current). D) $\text{GABA}_A(\alpha_4\beta_3\delta)$: On average, 111 μM GABA elicited a 0.54 nA (± 0.17) response (peak current). Deviations are \pm SD.

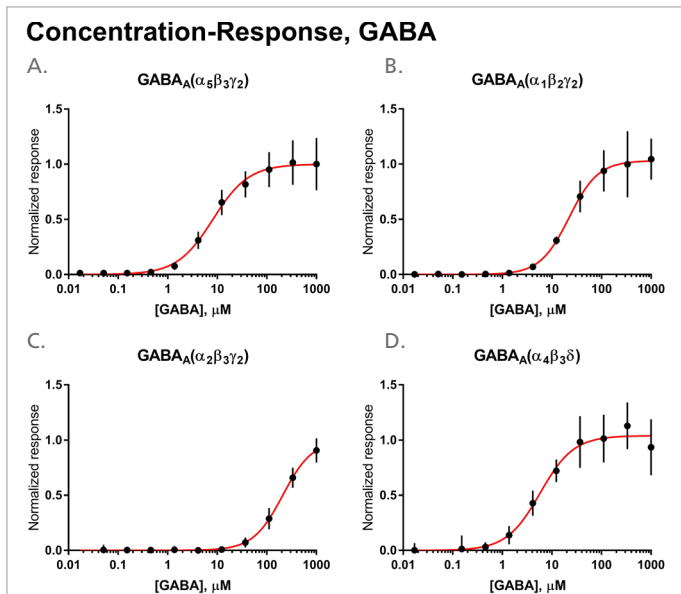


Fig. 5: GABA concentration-response of the four GABA_A receptor subtypes. A) For GABA_A($\alpha_5\beta_3\gamma_2$) the EC₅₀ value was 10.4 μ M (CI95%: 8.5 to 13.7 μ M) and the Hill slope was 1.2 (± 0.07). B) For GABA_A($\alpha_1\beta_2\gamma_2$) the EC₅₀ value was 22.1 μ M (CI95%: 19.6 to 25.0 μ M) and the Hill slope was 1.5 (± 0.1). C) For GABA_A($\alpha_2\beta_3\gamma_2$) the EC₅₀ value was 0.21 mM (CI95%: 0.19 to 0.23 μ M) and the Hill slope 1.45 (± 0.08). D) For GABA_A($\alpha_4\beta_3\delta$) the EC₅₀ value was 5.7 μ M (CI95%: 4.5 to 7.3 μ M) and the Hill slope was 1.2 (± 0.2). Error bars: \pm SD.

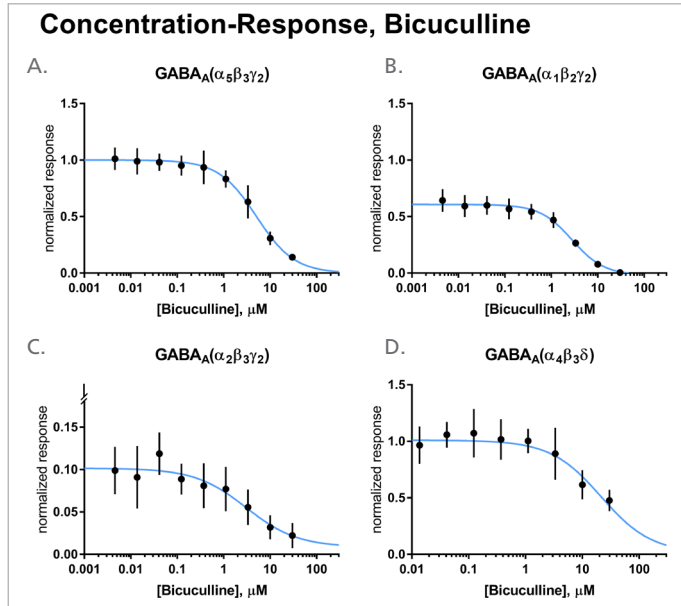


Fig. 6: Bicuculline concentration-response relationship of four GABA_A receptor subtypes. Prior to the exposure to 30 μ M GABA, the cells were preincubated with the test concentration of bicuculline. The response is normalized to the max GABA current. A) For GABA_A($\alpha_5\beta_3\gamma_2$) the IC₅₀ value at 30 μ M GABA was 5.1 μ M (CI95%: 4.4 to 5.8 μ M) and the Hill slope was -1.1 (± 0.06). B) For GABA_A($\alpha_1\beta_2\gamma_2$) the IC₅₀ value at 30 μ M GABA was 2.7 μ M (CI95%: 2.5 to 3.1 μ M) and the Hill slope was -1.3 (± 0.08). C) For GABA_A($\alpha_2\beta_3\gamma_2$) the IC₅₀ value at 30 μ M GABA was 3.8 μ M (CI95%: 2.4 to 5.7 μ M) and the Hill slope was -0.63 (± 0.08). D) For GABA_A($\alpha_4\beta_3\delta$) the IC₅₀ value at 30 μ M GABA was 16 μ M (CI95%: 12 to 21 μ M) and the Hill slope was -1.1 (± 0.1). Error bars: \pm SD.

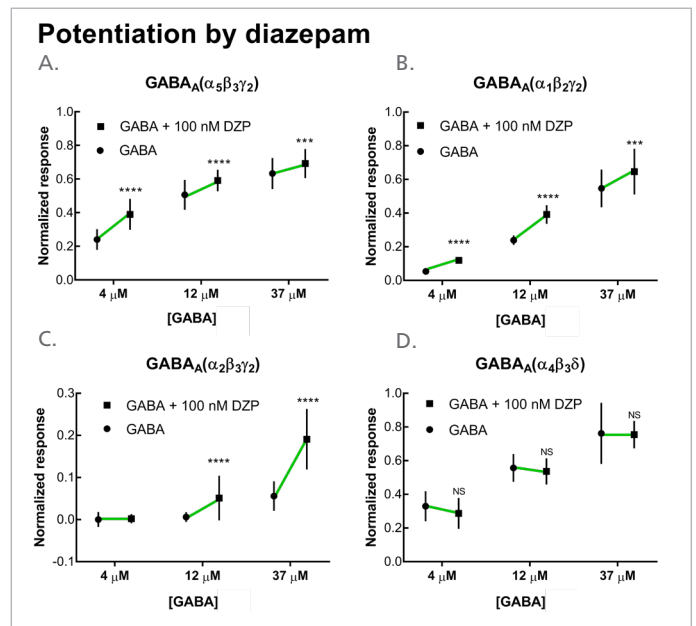


Fig. 7: Potentiation by diazepam. Diazepam is a benzodiazepine, binding to the benzodiazepine binding site which is located between the α and the γ subunit of the GABA_A receptors. A diazepam effect can hence confirm a probe γ subunit expression. A) Diazepam (100 nM) potentiated the GABA_A($\alpha_5\beta_3\gamma_2$) response: 4 μ M: 162% \pm 7%, $p < 0.0001$, 12 μ M: 117% \pm 3%, $p < 0.0001$ and 37 μ M: 109% \pm 3%, $p < 0.05$. B) Diazepam (100 nM) potentiated the GABA_A($\alpha_1\beta_2\gamma_2$) response: 4 μ M: 222% \pm 7%, $p < 0.0001$, 12 μ M: 164% \pm 5%, $p < 0.0001$ and 37 μ M: 116% \pm 6%, $p < 0.01$. C) Diazepam (100 nM) potentiated the GABA_A($\alpha_2\beta_3\gamma_2$) response. 4 μ M GABA did however not induce a measurable response and the response to 12 μ M GABA was only measurable in the presence of diazepam. The response to 37 μ M GABA was increased by 100nM Diazepam: 341% \pm 27%, $p < 0.0001$. D) Diazepam (100 nM) did not potentiate the GABA response. A δ -subunit-containing GABA_A receptor does not contain a benzodiazepine binding site, hence the lack of diazepam effect. Error bars: \pm SD.

Methods

The four cell lines expressing human GABA_A receptors were cultured according to the suppliers' description. ($\alpha_5\beta_3\gamma_2$)/HEK 293 was kindly supplied by Charles River Laboratories, Cleveland, OH, ($\alpha_5\beta_3\gamma_2$)/HEK 293 and ($\alpha_5\beta_3\gamma_2$)/HEK 293 was kindly supplied by SB Drug Discovery, Glasgow, UK and ($\alpha_5\beta_3\gamma_2$)/CHO was kindly supplied by B'SYS, Witterswil, CH.

All experiments were carried out at ambient temperature using Qube 384 multi-hole consumables and patched using a standard whole cell protocol.

Data analysis was performed using the Sophion Assay Software and GraphPad Prism 7.03 (GraphPad Software Inc.).

References:

1. Foster, A. & Kemp, J. Glutamate- and GABA-based CNS therapeutics. *Curr. Opin. Pharmacol.* 6, 7–17 (2006).
2. Johnston, G. A. R. GABA(A) receptor channel pharmacology. *Curr. Pharm. Des.* 11, 1867–85 (2005).
3. Watanabe, M., Maemura, K., Kanbara, K., Tamayama, T. & Hayasaki, H. GABA and GABA receptors in the central nervous system and other organs. *Int. Rev. Cytol.* 213, 1–47 (2002).
4. Chapouthier, G. & Venault, P. A pharmacological link between epilepsy and anxiety? *Trends Pharmacol. Sci.* 22, 491–3 (2001).
5. Kaila, K., Voipio, J., Paalasmaa, P., Pasternack, M. & Deisz, R. A. The role of bicarbonate in GABA_A receptor-mediated IPSPs of rat neocortical neurones. *J. Physiol.* 464, 273–89 (1993).
6. Farrant, M. & Kaila, K. The cellular, molecular and ionic basis of GABA(A) receptor signalling. *Prog. Brain Res.* 160, 59–87 (2007).
7. Barnard, E. A. et al. International Union of Pharmacology. XV. Subtypes of gamma-aminobutyric acidA receptors: classification on the basis of subunit structure and receptor function. *Pharmacol. Rev.* 50, 291–313 (1998).

Author:

Kim Boddum , Application scientist

Sophion Bioscience A/S, Baltorpevej 154, 2750 Ballerup, Denmark
Phone: +45 4460 8800, E-mail: info@sophion.com

sophion.com