

## Application Report

# Pharmacological evaluation of GABA<sub>A</sub> receptor subtypes on Qube 384

High-throughput screening and cumulative concentration-response relationship of a GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) cell line. Characterization of GABA<sub>A</sub> receptor subtypes

### Summary

Studies of GABA<sub>A</sub> ion channels using the automated patch clamp platform Qube 384 with focus on:

- Short ligand exposure with repetitive stimulations with EC<sub>50</sub> 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<sub>A</sub> subtypes

### Introduction

$\gamma$ -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 potential<sup>1-4</sup>.

Ionotropic GABA receptors consist of 5 membrane-spanning subunits<sup>5,6</sup>, of which 19 different have been identified in humans ( $\alpha_{1-6}$ ,  $\beta_{1-3}$ ,  $\gamma_{1-3}$ ,  $\delta$ ,  $\epsilon$ ,  $\theta$ ,  $\pi$ ,  $\rho_{1-3}$ ). Receptors in different cell types have different subunit composition and differ both in pharmacology and subcellular location<sup>7</sup>. Due to the heterogeneity of GABA<sub>A</sub> receptors, they can perform different inhibitory tasks: Positive modulation of GABA<sub>A</sub> receptors can be anticonvulsant, hypnotic, anaesthetic and anxiolytic, whereas negative modulation can enhance cognition but also be anxiogenic and proconvulsant<sup>1-4</sup>.

Here we use four GABA<sub>A</sub>/HEK293 cell lines and a range of tool compounds to demonstrate two different approaches to GABA<sub>A</sub>

receptor evaluation on Qube. We show 1) a cumulative dose-response experiment with a GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) cell line and 2) a GABA<sub>A</sub> 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<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ )

Desensitization due to prolonged or repeated agonist exposure poses a challenge when studying GABA<sub>A</sub> receptors in an in vitro setting. Therefore, the effect of GABA on the GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) 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  $\mu$ M) could be made without significant rundown (Figure 1A). 12  $\mu$ M GABA, which is close to the EC<sub>50</sub> 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 $\Omega$  per cell

Capacitance > 5 pF per cell

Current amplitude at 12  $\mu$ M GABA > 300 pA per cell

## GABA<sub>A</sub> receptor subtype screen

The GABA<sub>A</sub> 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<sub>A</sub> 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.

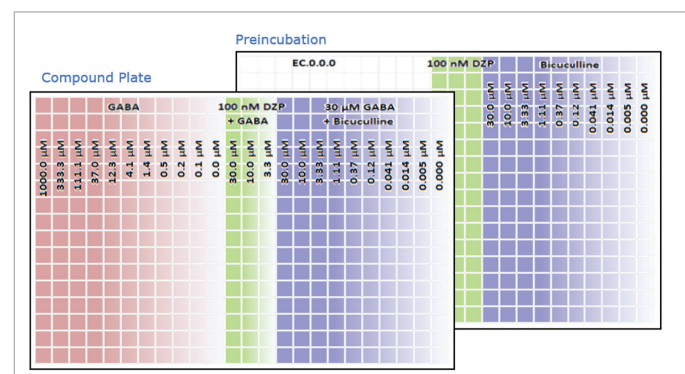


Fig. 3: Compound plate layout: 100 nM diazepam and varying concentrations of bicuculline were washed in (pre-incubation) prior to GABA application (compound plate).

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  $\alpha_4\beta_3\delta$ ).

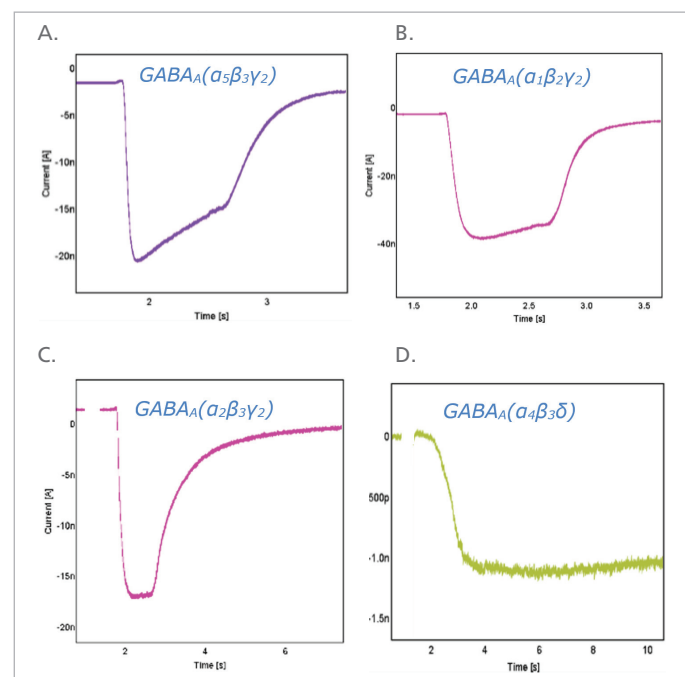


Fig. 4: GABA<sub>A</sub> receptor kinetics. Typical recordings from one multihole site (10 cells/site). The effect of GABA was evaluated employing the stacked delivery feature, with an exposure time of less than one second (with exception from GABA<sub>A</sub> ( $\alpha_4\beta_3\delta$ ) where a regular and not stacked delivery feature was employed due to the slow kinetics of  $\delta$ -subunit containing GABA<sub>A</sub> receptors).

A) GABA<sub>A</sub> ( $\alpha_5\beta_3\gamma_2$ ): On average, 111  $\mu$ M GABA elicited a 26.3 nA ( $\pm$  8.5 nA) response (peak current). B) GABA<sub>A</sub> ( $\alpha_1\beta_2\gamma_2$ ): On average, 111  $\mu$ M GABA elicited a 31.6 nA ( $\pm$  8.3) response (peak current). C) GABA<sub>A</sub> ( $\alpha_2\beta_3\gamma_2$ ): On average, 1 mM GABA elicited a 9.6 nA ( $\pm$  4.1) response (peak current). D) GABA<sub>A</sub> ( $\alpha_4\beta_3\delta$ ): On average, 111  $\mu$ M GABA elicited a 0.54 nA ( $\pm$  0.17) response (peak current). Deviations are  $\pm$  SD.

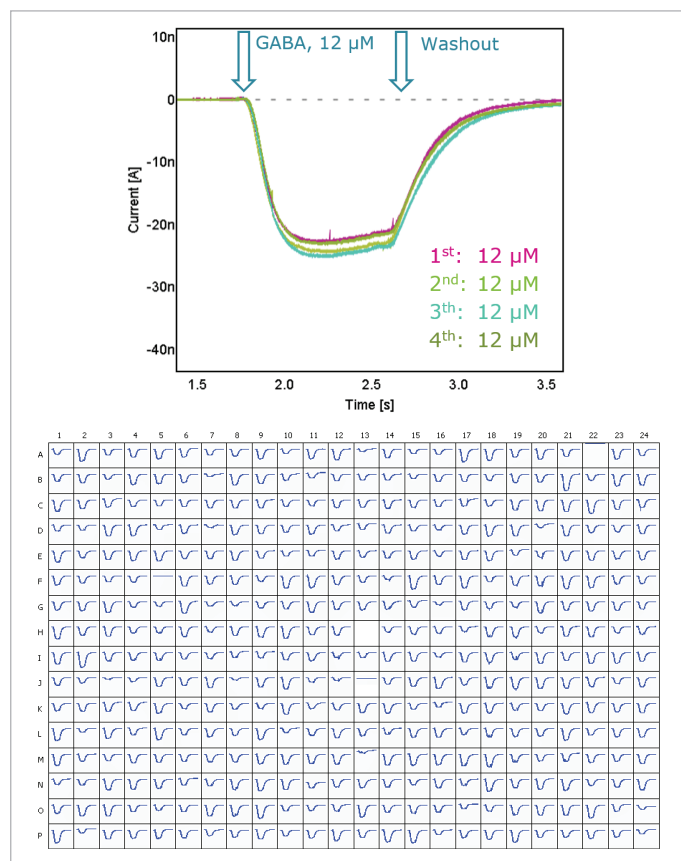


Fig. 1: Top: 4 consecutive applications of 12  $\mu$ M GABA using the stacked delivery feature. Bottom: Plate view of a QChip, showing the 384 individual responses to 12  $\mu$ M GABA.

## Cumulative GABA concentration response

The concentration-response relationship of GABA on the GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) receptor was evaluated on Qube. GABA was applied in increasing concentrations (3-fold dilution from 400  $\mu$ M, Figure 2). The EC<sub>50</sub> value for the cumulative concentration response was found to be 10.9  $\mu$ M (CI95%: 10.0 to 12.4  $\mu$ M) and the Hill slope was 1.5 ( $\pm$ 0.1, SD).

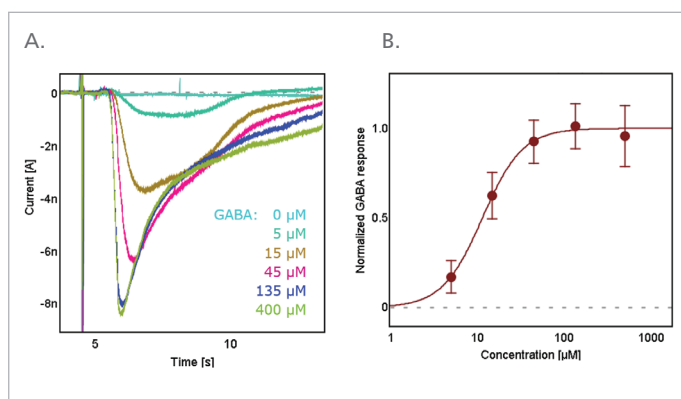


Fig. 2: Cumulative concentration-response relationship of GABA binding to the GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) receptor. A) Typically recorded currents in response to increasing concentrations of GABA. B) Peak current (normalized to highest value in experiment, average  $\pm$  SD, n=192)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

## Concentration-Response, GABA

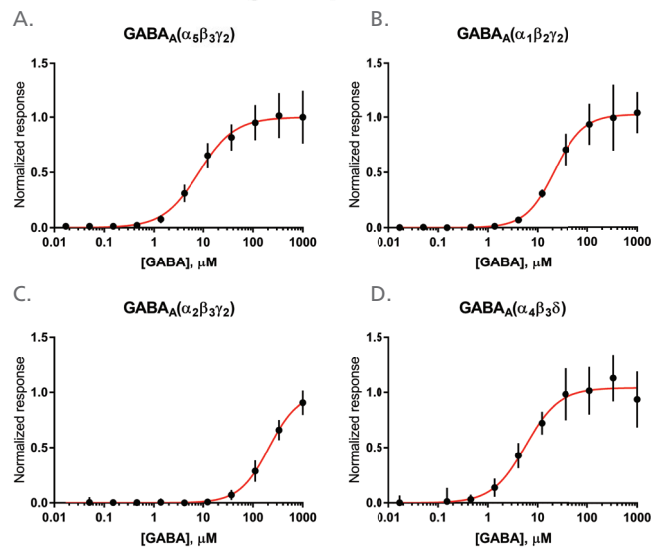


Fig. 5: GABA concentration-response of the four GABA<sub>A</sub> receptor subtypes. A) For GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) the EC<sub>50</sub> value was 10.4  $\mu$ M (CI95%: 8.5 to 13.7  $\mu$ M) and the Hill slope was 1.2 ( $\pm 0.07$ ). B) For GABA<sub>A</sub>( $\alpha_1\beta_2\gamma_2$ ) the EC<sub>50</sub> value was 22.1  $\mu$ M (CI95%: 19.6 to 25.0  $\mu$ M) and the Hill slope was 1.5 ( $\pm 0.1$ ). C) For GABA<sub>A</sub>( $\alpha_2\beta_3\gamma_2$ ) the EC<sub>50</sub> value was 0.21 mM (CI95%: 0.19 to 0.23  $\mu$ M) and the Hill slope 1.45 ( $\pm 0.08$ ). D) For GABA<sub>A</sub>( $\alpha_4\beta_3\delta$ ) the EC<sub>50</sub> value was 5.7  $\mu$ M (CI95%: 4.5 to 7.3  $\mu$ M) and the Hill slope was 1.2 ( $\pm 0.2$ ). Error bars:  $\pm$  SD.

## Concentration-Response, Bicuculline

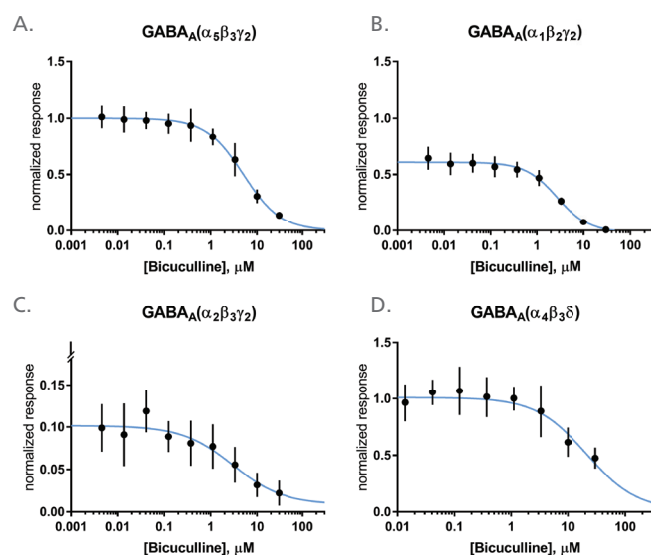


Fig. 6: Bicuculline concentration-response relationship of four GABA<sub>A</sub> receptor subtypes. Prior to the exposure to 30  $\mu$ M GABA, the cells were preincubated with the test concentration of bicuculline. The response is normalized to the max GABA current. A) For GABA<sub>A</sub>( $\alpha_5\beta_3\gamma_2$ ) the IC<sub>50</sub> value at 30  $\mu$ M GABA was 5.1  $\mu$ M (CI95%: 4.4 to 5.8  $\mu$ M) and the Hill slope was -1.1 ( $\pm 0.06$ ). B) For GABA<sub>A</sub>( $\alpha_1\beta_2\gamma_2$ ) the IC<sub>50</sub> value at 30  $\mu$ M GABA was 2.7  $\mu$ M (CI95%: 2.5 to 3.1  $\mu$ M) and the Hill slope was -1.3 ( $\pm 0.08$ ). C) For GABA<sub>A</sub>( $\alpha_2\beta_3\gamma_2$ ) the IC<sub>50</sub> value at 30  $\mu$ M GABA was 3.8  $\mu$ M (CI95%: 2.4 to 5.7  $\mu$ M) and the Hill slope was -0.63 ( $\pm 0.08$ ). D) For GABA<sub>A</sub>( $\alpha_4\beta_3\delta$ ) the IC<sub>50</sub> value at 30  $\mu$ M GABA was 16  $\mu$ M (CI95%: 12 to 21  $\mu$ M) and the Hill slope was -1.1 ( $\pm 0.1$ ). Error bars:  $\pm$  SD.

## Potentialiation by diazepam

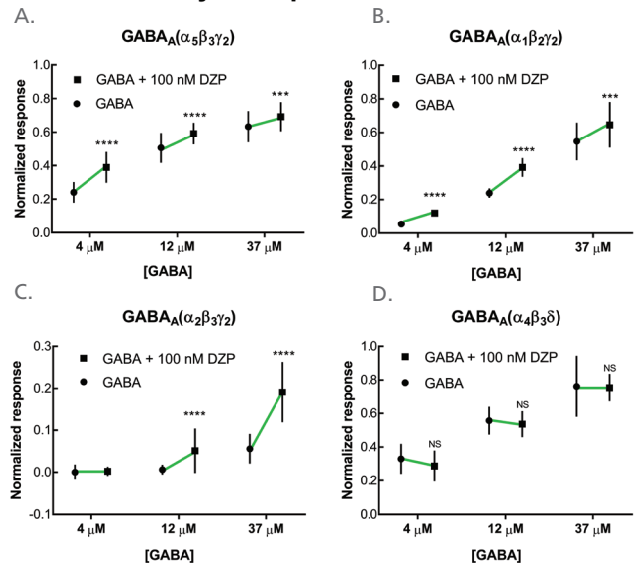


Fig. 7: Potentiation by diazepam. Diazepam is a benzodiazepine, binding to the benzodiazepine binding site which is located between the  $\alpha$  and the  $\gamma$  subunit of the GABA<sub>A</sub> receptors. A diazepam effect can hence confirm a proper  $\gamma$  subunit expression. A) Diazepam (100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

## Methods

The four HEK cell lines expressing human GABA<sub>A</sub> receptors were cultured according to the suppliers' description. (α<sub>5</sub>β<sub>3</sub>γ<sub>2</sub>)/HEK 293 was kindly supplied by Charles River Laboratories, Cleveland, OH, (α<sub>5</sub>β<sub>3</sub>γ<sub>2</sub>)/HEK 293 and (α<sub>5</sub>β<sub>3</sub>γ<sub>2</sub>)/HEK 293 was kindly supplied by SB Drug Discovery, Glasgow, UK and (α<sub>5</sub>β<sub>3</sub>γ<sub>2</sub>)/HEK 293 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<sub>A</sub> 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