

Activation and inhibition of GABA_A $\alpha 5\beta 3\gamma 2$ receptors on the Port-a-Patch

The electrophysiology team at Nanion Technologies GmbH, Munich. Cells were kindly supplied by Merck Millipore.

Summary

Gamma aminobutyric acid type A (GABA_A) receptors are the most important inhibitory neurotransmitter receptors in the mammalian central nervous system (CNS). They are opened by GABA allowing the passage of chloride ions across the membrane. GABA_A channels are modulated by a variety of different drugs including benzodiazepines, barbiturates, neuroactive steroids, anesthetics, and convulsants¹. The receptors are heteropentameric and depending on the subunit combination, they exhibit different electrophysiological and pharmacological properties². Six α -, three β -, three γ -, one δ -, one ϵ -, one π -, one θ - and three ρ -subunits have been cloned, including splice variants of some of these subunits^{3,4}. Functional GABA_A receptors typically assemble with two α , two β , and one γ subunit, with alternating α and β subunits connected by a γ subunit¹. GABA_A receptors play a critical role in regulating excitability of the brain, anxiety, vigilance, as well as learning and memory². The $\alpha 5$ subunit is highly expressed in the hippocampus and olfactory bulb and expressed in low levels in other brain regions including the cortex, subiculum, hypothalamus, sympathetic preganglionic neurons, and amygdala⁵. GABA_A receptors containing the $\alpha 5$ subunit cluster at both extrasynaptic sites as well as synaptic sites thus contributing to tonic currents and synaptic GABA-ergic neurotransmission⁵. $\alpha 5$ -containing receptors exhibit unique physiology and pharmacology and they are potential pharmacological targets for the treatment of neurodevelopmental disorders, depression, schizophrenia, and mild cognitive impairment.

The Port-a-Patch with External Perfusion System was used to record $\alpha 5\beta 3\gamma 2$ receptors expressed in HEK293 cells.

Results

GABA_A receptors with the subunit combination $\alpha 5\beta 3\gamma 2$ were activated by increasing concentrations of GABA (Figure 1). Currents started to activate at concentrations above 3 μ M and reached maximum at a concentration of 30 μ M. A concentration response curve was constructed for an average of 8 cells (Figure 1B). The EC₅₀ was calculated to be $10.8 \pm 0.2 \mu$ M (n = 8), in good agreement with the literature for cloned GABA_A receptors expressing the $\alpha 5$ -subunit⁶.

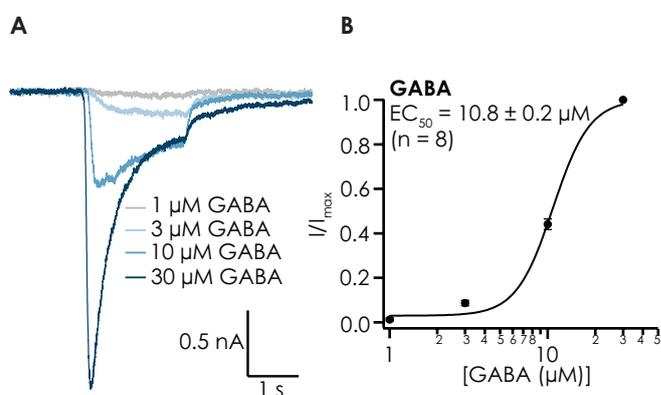


Figure 1: **A** Whole cell current responses from HEK293 cells expressing $\alpha 5\beta 3\gamma 2$ by increasing concentrations of GABA. Shown are current responses from an example cell with perfusion of external GABA at the concentrations indicated for 2 s. Following this, GABA was washed away by external recording solution. Holding potential was -80 mV. **B** Concentration response curve for an average of 8 cells was constructed and is shown. EC₅₀ = $10.8 \pm 0.2 \mu$ M (n = 8).

Application Note

GABA_A currents mediated by $\alpha 5\beta 3\gamma 2$ could be repetitively activated by 1 mM GABA (Figure 2). A small decrease in peak amplitude was observed between the first and second application of GABA (14% reduction from -525 ± 91 pA to -455 ± 65 pA). Following this, repeated application of GABA resulted in identical peak current amplitudes (< 3% difference compared with the previous amplitude).

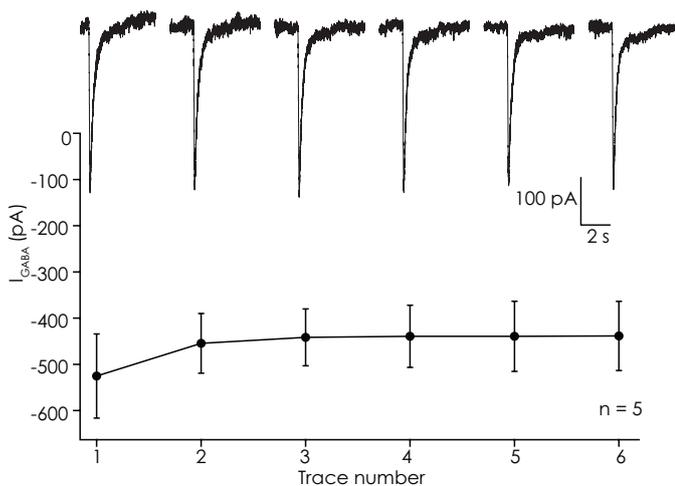


Figure 2: Traces from an example cell showing repetitive activation by 1 mM GABA (top) and the corresponding peak amplitudes for an average of 5 cells (bottom).

GABA_A receptors with the subunit combination $\alpha 5\beta 3\gamma 2$ were activated by 30 μ M GABA and blocked by the competitive antagonist, bicuculline, when co-applied with GABA (Figure 3).

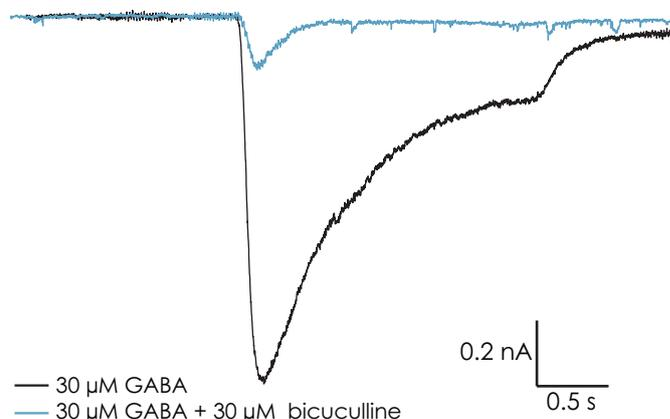


Figure 3: $\alpha 5\beta 3\gamma 2$ GABA_A receptors activated by 30 μ M GABA were blocked by 30 μ M bicuculline when co-applied.

In conclusion, GABA_A receptors with the subunit combination $\alpha 5\beta 3\gamma 2$ were reliably recorded on the Port- α -Patch with External Perfusion System. Perfusion of GABA was triggered automatically by the amplifier and repetitive activation of GABA_A receptors was possible using a high concentration of GABA. In this way, both inhibitors, such as bicuculline, and other modulators (not shown) could be co-applied and characterized using the Port- α -Patch.

References

1. Sieghart, W. 2006. *Advances in Pharmacology*. 54(5): 231-263
2. Sarto-Jackson, I. & Sieghart, W. 2008. *Mol. Memb. Biol* 25(4): 302-310
3. Simon, J., et al., 2004. *JBC*. 279(40): 41422-41435
4. Olsen, R.W. & Sieghart, W. 2009. *Neuropharm*. 56(1): 141-148
5. Jacob, T.C. 2019. *Front. Mol. Neurosci*. 12:179. doi: 10.3389/fnmol.2019.00179
6. Petroski, R.E., et al. 2006. *JPET*. 317(1): 369-377

Methods

Cells

HEK293 cells stably expressing $\alpha 5\beta 3\gamma 2$ GABA_A receptors were kindly provided by Merck Millipore.

Cell culture

HEK293 cells stably expressing $\alpha 5\beta 3\gamma 2$ GABA_A receptors were cultured using standard culture conditions for the Port- α -Patch.

Electrophysiology

Whole cell patch clamp recordings were conducted according to Nanion's standard procedure for the Port- α -Patch. Currents were elicited by 2 s perfusion of GABA at different concentrations. A continuous holding potential of -80 mV was used. Bicuculline was pre-incubated for 2 mins prior to co-application with GABA.