

Orphan receptor GPR158 serves as a metabotropic glycine receptor: mGlyR

Thibaut Laboute^{1,4}, Stefano Zucca¹, Matthew Holcomb², Dipak N. Patil^{1†}, Chris Garza², Brittany A. Wheatley³, Raktim N. Roy³, Stefano Forli², Kirill A. Martemyanov^{1*}

¹Department of Neuroscience, UF Scripps Biomedical Research, Jupiter, FL 33458, USA.

²Department of Integrative Structural and Computational Biology, The Scripps Research Institute, La Jolla, CA 92037 USA.

³Department of Integrative Structural and Computational Biology, UF Scripps Biomedical Research, Jupiter, FL 33458, USA.

⁴UMR1253, iBrain, Université de Tours, INSERM, Tours, France

*Corresponding author. Email: kmartemyanov@ufl.edu

†Present address: Lilly Biotechnology Center, Eli Lilly and Company, 10290 Campus Point Dr., San Diego, CA 92121, USA.

Glycine is a major neurotransmitter involved in several fundamental neuronal processes. The identity of the metabotropic receptor mediating slow neuromodulatory effects of glycine is unknown. We identified an orphan G protein-coupled receptor, GPR158, as a metabotropic glycine receptor (mGlyR) and validated its pharmacology using Promega Bright NanoLuc® BRET technology. We used a live-cell assay to monitor GAP activity by following kinetics of G protein deactivation. In this assay, activation of G proteins by GPCR stimulation generates the BRET signal upon interaction of liberated Venus-Gbg subunits with the masGRK3CT-NanoLuc® reporter illuminated with Nano-Glo®. This signal is quenched when Gα deactivation is triggered by GPCR antagonism and recombines with Venus-Gbg to form inactive heterotrimer. Glycine and a related modulator, taurine, directly bind to a Cache domain of GPR158, and this event inhibits the activity of the intracellular signaling complex regulator of G protein signaling 7/G protein b5 (RGS7-Gb5), which is associated with the receptor. Glycine signals through mGlyR to inhibit production of the second messenger adenosine 3',5'-monophosphate. We further show that glycine, but not taurine, acts through mGlyR to regulate neuronal excitability in cortical neurons. These results identify a major neuromodulatory system involved in mediating metabotropic effects of glycine, with implications for understanding cognition and affective states.

This discovery is now published in Science,

<https://www.science.org/doi/10.1126/science.add7150>