

Anti-phospho-GABABR2 (Ser783)

Rabbit Polyclonal Antibody

Catalog # **G803-465BR**

Lot # Z2014-8

Cited Applications

WB, IF

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the GABABR2 protein phosphorylated at serine 783

Cross Reactivity

Rat, Mouse, Human, Non-Human Primate, Chicken, Bovine and Xenopus

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser 783 of GABABR2

Formulation

10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

Stability

1yr at -20°C from date of shipment

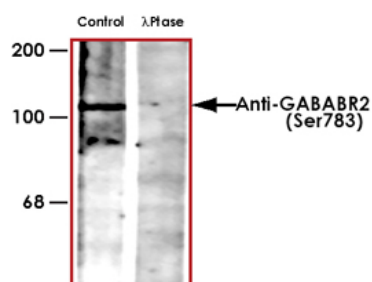
Scientific Background

Gamma-aminobutyric acid (GABA) is the primary inhibitory neurotransmitter in the central nervous system. There are two major classes of GABA receptors: the GABAA and the GABAB subtype of receptors. GABAB receptors are heterodimeric G protein-coupled receptors that mediate slow synaptic inhibition in the central nervous system. It has recently been demonstrated that AMPK binds directly to GABAB receptors and phosphorylates S783 in the cytoplasmic tail of the R2 subunit and that S783 plays a critical role in enhancing neuronal survival after ischemia as phosphorylation of S783 is evident in many brain regions and is increased dramatically after ischemic injury to the brain.

References

1. Kuramoto N, Wilkins ME, Fairfax BP, et al, Phospho-dependent functional modulation of GABA(B) receptors by the metabolic sensor AMP-dependent protein kinase. *Neuron* 53:233-247. 2007.

Sample Data



Western blot of rat synaptic membrane showing specific immunolabeling of the ~102 k GABABR2 protein phosphorylated at Ser783 (control). The phosphospecificity of this labeling is shown in the second lane (lambda-phosphatase: λ-Ptase). The blot is identical to the control except that it was incubated in λ-Ptase (1200 units for 30 min) before being exposed to the Anti-phospho-GABABR2 (Ser783). The immunolabeling is completely eliminated by treatment with λ-Ptase.

Anti-phospho-GABABR2 (Ser783)

Rabbit Polyclonal Antibody

Catalog #	G803-465BR
Lot #	Z2014-8
Purification	Affinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com

www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.