

PRODUCT INFORMATION

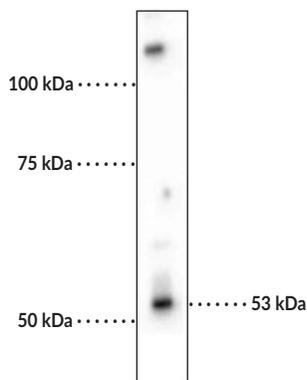


GABA_A Receptor β_3 Subunit Polyclonal Antibody Item No. 29275

Overview and Properties

Contents:	This vial contains 100 μ l of affinity-purified rabbit polyclonal antibody.
Synonyms:	GABRB3, Gamma-aminobutyric Acid Receptor Subunit β_3
Immunogen:	Fusion protein from the cytoplasmic loop of the β_3 subunit of the rat GABA _A receptor
Molecular Weight:	~53 kDa
Species Reactivity:	(+) Mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 μ g/ml BSA, and 50% glycerol
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:300 and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



WB of rat brain lysate showing specific immunolabeling of the ~53 kDa β_3 -subunit of the GABA_A receptor.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/23/2020

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

GABA_A receptors are ligand-gated chloride channels that mediate the effects of the inhibitory neurotransmitter GABA in the CNS.^{1,2} They are postsynaptic heteropentameric receptors that contain protein subunits from the following isoforms: α_{1-6} , β_{1-4} , γ_{1-3} , δ , ϵ , π , θ , and ρ_{1-3} , arranged around a central pore. Phasic inhibitory synaptic transmission is regulated by $\alpha_1\beta_2\gamma_2$ subunit-containing GABA_A receptors, the major isoform found in the brain.^{2,3} The β subunit of GABA_A receptors interfaces with an α subunit to form the GABA binding site that initiates GABA-induced action potentials and forms the benzodiazepine binding site with the γ subunit. β_3 subunit-containing GABA_A receptors are widely expressed in the cerebral cortex, cerebellum, olfactory bulb, and hippocampus.⁴ Phosphorylation of the β_3 subunit by PKA or PKC inhibits binding of the β_3 subunit with the clathrin adaptor protein AP2 and reduces GABA_A receptor endocytosis.⁵ Mutations in *GABRB3*, which encodes the β_3 subunit isoform, have been found in patients with childhood absence epilepsy (CAE), infantile spasms (IS), and Lennox-Gastaut syndrome (LGS).² Cayman's GABA_A Receptor β_3 Subunit Polyclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes the GABA_A receptor β_3 subunit at approximately 53 kDa from rat and mouse samples.

References

1. Crestani, F. and Rudolph, U. Behavioral functions of GABA_A receptor subtypes - the Zurich experience. *Adv. Pharmacol.* **72**, 37-51 (2015).
2. Hirose, S. Mutant GABA_A receptor subunits in genetic (idiopathic) epilepsy. *Prog. Brain Res.* **213**, 55-85 (2014).
3. Wongsamitkul, N., Maldifassi, M.C., Simeone, X., et al. α subunits in GABA_A receptors are dispensable for GABA and diazepam action. *Sci. Rep.* **7(1)**, 15498 (2017).
4. Miralles, C.P., Li, M., Mehta, A.K., et al. Immunocytochemical localization of the β_3 subunit of the γ -aminobutyric acid_A receptor in the rat brain. *J. Comp. Neurol.* **413(4)**, 535-548 (1999).
5. Vithlani, M., Moss, S.J., and Terunuma, M. The dynamic modulation of GABA_A receptor trafficking and its role in the formation of inhibitory synapses. *Physiol. Rev.* **91(3)**, 1009-1022 (2011).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM