

PRODUCT INFORMATION



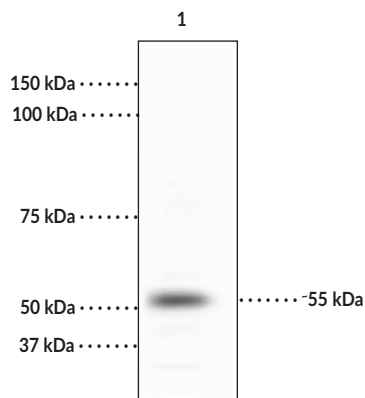
GABA_A Receptor β_1 Subunit Polyclonal Antibody

Item No. 29273

Overview and Properties

Contents:	This vial contains 100 μ l of affinity-purified polyclonal antibody from pooled serum.
Synonyms:	GABA _A Receptor Subunit β_1 , GABRB2, Gamma-aminobutyric Acid (GABA) A Receptor, β_1 , Gamma-aminobutyric Acid Receptor Subunit β_1
Immunogen:	Fusion protein from the cytoplasmic loop of the β_1 subunit of the rat GABA _A receptor
Molecular Weight:	~55 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:	Western blot (WB); the recommended starting dilution is 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Mouse whole brain lysates

WB of mouse whole brain lysates showing specific immunolabeling of the ~55 kDa β_1 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, 12/10/2020

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. Phosphorylation of the β_1 subunit by PKA or PKC inhibits binding of the β_1 subunit with the clathrin adaptor protein AP2 and reduces GABA_A receptor endocytosis.⁴ Mice expressing mutations in *Gabrb1*, which encodes the β_1 subunit isoform, have increased GABA_A receptor-mediated tonic inhibition in the nucleus accumbens and increased alcohol consumption compared to wild-type mice.⁵ SNPs in *GABRB1* are associated with increased impulsivity and reward sensitivity in human adolescents.⁶ Cayman's GABA_A Receptor β_1 Polyclonal Antibody can be used for Western blot (WB) applications. The antibody recognizes the GABA_A receptor β_1 subunit at approximately 55 kDa from mouse and rat samples.

References

1. Crestani, F. and Rudolph, U. Behavioral functions of GABA_A receptor subtypes - the Zurich experience. *Diversity and Functions of GABA receptors: A tribute to Hanns Möhler, part A*. Rudolph, U., editor, 1st edition, Academic Press (2015).
2. Hirose, S. Mutant GABA_A receptor subunits in genetic (idiopathic) epilepsy. *Genetics of epilepsy*. Steinlein, O., editor, 1st edition, Elsevier (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. 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).
5. Anstee, Q.M., Knapp, S., Maguire, E.P., et al. Mutations in the *GABRB1* gene promote alcohol consumption through increased tonic inhibition. *Nat. Commun.* **4**, 2816 (2013).
6. Duka, T., Nikolaou, K., King, S.L., et al. *GABRB1* single nucleotide polymorphism associated with altered brain responses (but not performance) during measures of impulsivity and reward sensitivity in human adolescents. *Front. Behav. Neurosci.* **11**(24), (2017).

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