

PRODUCT INFORMATION



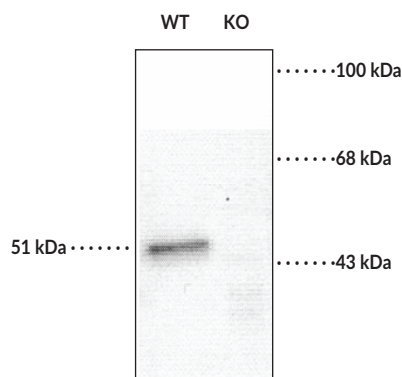
GABA_A Receptor α_1 Subunit (N-Term) Polyclonal Antibody

Item No. 29267

Overview and Properties

Contents:	This vial contains 100 μ l of affinity-purified rabbit polyclonal antibody.
Synonyms:	γ -Aminobutyric Acid Receptor Subunit α_1 , GABA _A Receptor Subunit α_1 , GABRA1
Immunogen:	Peptide from the N-terminal region of the α_1 subunit of the rat GABA _A receptor conjugated to keyhole limpet hemocyanin (KLH)
Molecular Weight:	~51 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 1:100 and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



WB of mouse forebrain lysates from wild type (WT) and α_1 -knockout animals showing specific immunolabeling of the ~51 kDa α_1 -subunit of the GABA_A receptor. The labeling was absent from a lysate prepared from α_1 -knockout animals.

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, 06/22/2020

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 α subunit of GABA_A receptors interfaces with a β subunit to form the GABA binding site that initiates GABA-induced action potentials and forms the benzodiazepine binding site with the γ subunit. Mutations in *GABRA1*, which encodes the α_1 subunit isoform, are found in patients with autosomal dominant juvenile myoclonic epilepsy (JME), infantile spasms, and childhood absence epilepsy (CAE).² *GABRA1* expression is upregulated in ipsilateral, but not contralateral, perilesional tissue in a rat model of cortical ischemia-reperfusion injury.⁴ Cayman's GABA_A Receptor α_1 Subunit (N-Term) Polyclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes the GABA_A receptor α_1 subunit at approximately 51 kDa from mouse and rat 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. Neumann-Haefelin, T., Bosse, F., Redecker, C., *et al.* Upregulation of GABA_A-receptor α_1 - and α_2 -subunit mRNAs following ischemic cortical lesions in rats. *Brain Res.* **816(1)**, 234-237 (1999).

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