

PRODUCT INFORMATION

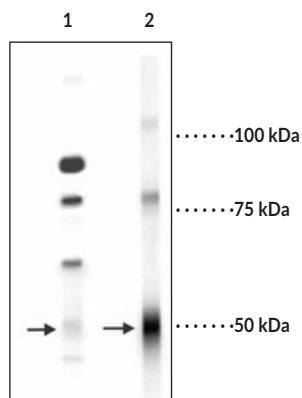


GABA_A Receptor δ subunit (N-Term) Polyclonal Antibody Item No. 10600

Overview and Properties

Contents:	This vial contains 100 μ l of affinity-purified antibody.
Synonym:	γ -Aminobutyric Acid A Receptor
Immunogen:	Fusion protein from the N-terminal region of the δ subunit of rat GABA _A receptor
Molecular Weight:	50 kDa
Species Reactivity:	(+) Mouse and rat; other species not tested
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:	Immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:500, 1:200, and 1:1,000, respectively. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Mouse whole brain lysate
Lane 2: Mouse synaptic plasma membrane lysate

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, 04/10/2020

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

γ -Aminobutyric acid (GABA) is the primary inhibitory neurotransmitter in the central nervous system, causing a hyperpolarization of the membrane through the opening of a Cl^- channel associated with the GABA_A receptor (GABA_A -R) subtype. GABA_A -Rs are important therapeutic targets for a range of sedative, anxiolytic, and hypnotic agents and are implicated in several diseases including epilepsy, anxiety, depression, and substance abuse. The GABA_A -R is a multimeric subunit complex. To date six α 's, four β 's, and four γ 's, plus alternative splicing variants of some of these subunits, have been identified. Injection in oocytes or mammalian cell lines of cRNA coding for α and β subunits results in the expression of functional GABA_A -Rs sensitive to GABA. However, co-expression of a γ subunit is required for benzodiazepine modulation. The various effects of the benzodiazepines in brain may also be mediated *via* different α subunits of the receptor. More recently there have been a number of studies demonstrating that the δ subunit of the receptor may affect subunit assembly and may also confer differential sensitivity to neurosteroids and to ethanol.

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