

PRODUCT DATA SHEET



CATALOG NO.: SA-516

LOT NO.: temp

PRODUCT: Anti-GABA_A Receptor, α_5 -Subunit, C-terminus

IMMUNOGEN: Recombinant protein from the cytosolic loop of the α_5 -Subunit of rat GABA_A Receptor.

SPECIFICITY: Recognizes the 55 kDa α_5 -Subunit of the GABA_A Receptor protein. Recognizes human, mouse, rat, canine, zebrafish and bovine subunit α_5 of the GABA_A receptor.

APPLICATIONS: WB, 1:1000

FORM / PURIFICATION: Affinity purified rabbit polyclonal antibody in 10 mM HEPES (pH 7.5), 150 mM NaCl, 10% BSA and 50% glycerol.

STORAGE: -20°C.

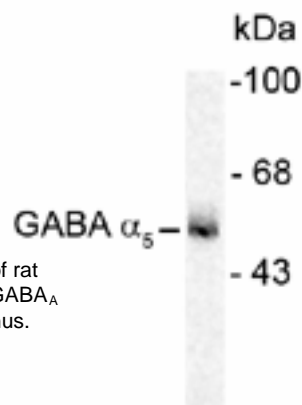


Figure: Western Blot analysis of rat hippocampus lysate with Anti-GABA_A Receptor, α_5 -Subunit, C-terminus.

REFERENCES:

1. NJ Brandon *et al.* *Mol Cell Neurosci.* 2003 **22** 87
2. RM McKernan *et al.* *Nat Neurosci* 2000 **3** 587
3. W Ogris *et al.* *Biochem Pharmacol.* 2004 **68** 1621
4. A Pörtl *et al.* *J Neurochem* 2003 **87** 1444

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.