

GluR8 Antibody [#C0212]

Catalog Number: C0212-1, C0212-2, C0212

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No.: O00222

All Names: GPRC1H, MGLUR8, mGlu8, GLUTAMATE RECEPTOR, METABOTROPIC, 8, glutamate receptor, metabotropic 8 precursor

Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Storage/Stability: Store at -20°C/1 year

Immunogen: The antiserum was produced against synthesized peptide derived from human GluR8.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity GluR8 antibody detects endogenous levels of total GluR8 protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:300~1:1000 IHC: 1:50~1:100 ELISA: 1:10000

References: Patrick. M. Downey, J Biomol Screen, Dec 2005; 10: 841 - 848.
RJ Wenthold, J. Biol. Chem., Jan 1994; 269: 1332 - 1339.
Toshiaki Watanabe, Cancer Res., Oct 2006; 66: 9804 - 9808.
V. Bruno, J. Neurosci., Sep 2000; 20: 6413.