

## Glutamate receptor 2 (Ab-880) Antibody [#2B7096]

- Catalog Number:** 2B7096-1, 2B7096-2, 2B7096
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P42262
- All Names:** GRIA2, GluR-2, GluR-B, GluR-K2, Glutamate receptor 2 precursor, Glutamate receptor ionotropic, AMPA 2
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), 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 non-phosphopeptide derived from human Glutamate receptor 2 around the phosphorylation site of serine 880 (I-E-S-V-K).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Glutamate receptor 2 (Ab-880) antibody detects endogenous levels of total Glutamate receptor 2 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000    IHC: 1:50~1:100  
ELISA: 1:20000
- References:** Kenneth J. Seidenman, J. Neurosci., Oct 2003; 23: 9220.  
Baosong Liu, J. Neurosci., May 2006; 26: 5309 - 5319.  
Jie Fu, J. Neurosci., Aug 2003; 23: 7592.  
Jose L. Perez, J. Neurosci., Aug 2001; 21: 5417.