

Ephrin-B3 Antibody [#C0182]

- Catalog Number:** C0182-1, C0182-2, C0182
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : Q15768
All Names: EFL6, EPLG8, LERK8, Ephrin B3, LIGAND OF EPH-RELATED KINASE 8, eph-related receptor tyrosine kinase ligand 8, ephrin-B3 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 Ephrin-B3.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Ephrin-B3 antibody detects endogenous levels of total Ephrin-B3 protein.
Reactivity: Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Mitsutoshi Nakada, Cancer Res., Sep 2006; 66: 8492 - 8500.
Klas Kullander, Genes & Dev., Apr 2001; 15: 877.
M. Douglas Benson, PNAS, Jul 2005; 102: 10694 - 10699.
Stephanie R. Kadison, J. Neurosci., Aug 2006; 26: 8909 - 8914.