

Ephrin-A4 Antibody [#C0181]

- Catalog Number:** C0181-1, C0181-2, C0181
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P52798
- All Names:** EFL4, EPLG4, LERK4, eph-related receptor tyrosine kinase ligand 4, ephrin A4 isoform a, ephrin A4 isoform b, ephrin A4 isoform c, ligand of eph-related kinase 4
- 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-A4.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Ephrin-A4 antibody detects endogenous levels of total Ephrin-A4 protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000 ELISA: 1:5000
- References:** Hans-Christian Aasheim, Blood, Jan 2000; 95: 221 - 230
L. Scott Janis, J. Neurosci., Jun 1999; 19: 4962.
Amy E. Merrill, Hum. Mol. Genet., Apr 2006; 15: 1319 - 1328.
Katrin Mackarehstschian, Cereb Cortex, Sep 1999; 9: 601 - 610.