

PDGFB Antibody [#C0299]

- Catalog Number:** C0299-1, C0299-2, C0299
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P01127
- All Names:** PDGF2, ONCOGENE SIS, PDGF B CHAIN, PLATELET-DERIVED GROWTH FACTOR B CHAIN, SIMIAN SARCOMA VIRAL ONCOGENE HOMOLOG, SSV PDGFB/COL1A1 FUSION GENE, V-SIS PLATELET-DERIVED GROWTH FACTOR BETA POLYPEPTIDE
- 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 PDGFB.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** PDGFB antibody detects endogenous levels of total PDGFB protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 ELISA: 1:10000
- References:** Li Fang, Arterioscler. Thromb. Vasc. Biol., Apr 2004; 24: 787 - 792.
EJ Battegay, J. Cell Biol., May 1994; 125: 917.
RA Seifert, J. Biol. Chem., Feb 1993; 268: 4473 - 4480.
Jason Liu, Am J Physiol Lung Cell Mol Physiol, May 1998; 274: 702.