

## SHP-2 (Ab-580) Antibody [#B0028]

- Catalog Number:** B0028-1, B0028-2, B0028
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q06124
- All Names:** EC 3.1.3.48, PTN11, PTP-1D, PTP-2C, PTP2C, PTPN11, Protein-tyrosine phosphatase 2C, Protein-tyrosine phosphatase, non-receptor type 11, SH-PTP2, SH-PTP3, SHPTP2
- 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 SHP-2 around the phosphorylation site of tyrosine 580 (R-V-Y<sup>P</sup>-E-N).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** SHP-2 (Ab-580) antibody detects endogenous levels of total SHP-2 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1000 IHC: 1:50~100  
ELISA: 1:20000
- References:** Minako Yamaoka-Tojo, *Arterioscler. Thromb. Vasc. Biol.*, Sep 2006; 26: 1991 - 1997.  
Liang Liang, *Endocrinology*, Sep 2000; 141: 3328.  
Tony Avril, *J. Immunol.*, Dec 2004; 173: 6841 - 6849.  
Laura A. Maile, *Mol. Biol. Cell*, Sep 2003; 14: 3519.