

PTEN (Ab-380/382/383) Antibody [#B7203]

- Catalog Number:** B7203-1, B7203-2, B7203
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), 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 PTEN around the phosphorylation site of Serine 380 and Threonine 382/383.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** PTEN (Ab-380/382/383) Antibody detects endogenous levels of total PTEN protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC:1:50~100 ELISA: 1:20000
- Swiss-Prot No. :** P60484
- References:** Al-Khouri AM, et al. (2005) J Biol Chem. 280(42):35195-35202.
Torres J, et al. (2001) J Biol Chem. 276(2): 993-998.
Vazquez F, et al. (2000) Mol Cell Biol. 20(14): 5010-5018.