

p56Dok-2 (Ab-299) Antibody [#B7059]

- Catalog Number:** B7059-1, B7059-2, B7059
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** O60496
- All Names:** DOK2, Docking protein 2
- 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 non-phosphopeptide derived from human p56Dok-2 around the phosphorylation site of tyrosine 299 (G-E-Y^P-A-V).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** p56Dok-2 (Ab-299) antibody detects endogenous levels of total p56Dok-2 protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
- References:** Feng Cong, et.al. (1999) Mol. Cell. Biol; 19: 8314-8325.
Serge Lemay, et.al. (2000) Mol. Cell. Biol; 20: 2743-2754.
Ute Schaeper, et.al.(2000) J. Cell Biol; 149: 1419.
Miyuki Honma, et.al. (2006) Genes Cells; 11: 143-151.