

VASP (Ab-238) Antibody [#B7250]

- Catalog Number:** B7250-1, B7250-2, B7250
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- 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 VASP around the phosphorylation site of serine 238.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** VASP (Ab-238) antibody detects endogenous levels of total VASP protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 ELISA: 1:20000
- Swiss-Prot No. :** P50552
- References:** Wang HG, et al. (2005) euron; 45(3): 389-403.
Melichar VO, et al. (2004) Proc Natl Acad Sci U S A; 101(47): 16671-6.
Jensen BO, et al. (2004) Blood; 104(9): 2775-82.
Zhuang S, et al. (2004) J Biol Chem; 279(11): 10397-407.
Smolenski A, et al. (1998) J Biol Chem; 273(32): 20029-35.
Butt E, et al. (1994) J Biol Chem; 269(20): 14509-17.