

## Ezrin (Ab-353) Antibody [#B7080]

- Catalog Number:** B7080-1, B7080-2, B7080
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P15311
- All Names:** Cytovillin, EZRI, VIL2, Villin 2, p81
- 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 Ezrin around the phosphorylation site of tyrosine 353 (Q-D-Y<sup>P</sup>-E-E).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Ezrin (Ab-353) antibody detects endogenous levels of total Ezrin protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:500~1:1000      ELISA: 1:20000
- References:** Zhao H, et al. (2004) Proc Natl Acad Sci U S A 101 (25): 9485-90.  
Wang Q, et al. (2003) J Biol Chem 278(48): 47731-43.  
Gautreau A, et al. (1999) Proc Natl Acad Sci U S A: 96(13): 7300-5.  
Crepaldi T, et al. (1997) J Cell Biol: 28; 138(2): 423-34.