

Ezrin (Ab-146) Antibody [#B8030]

Catalog Number:	B8030
Concentration:	1mg/ml
Swiss-Prot No.:	P15311
Other Names:	Cytovillin; EZRI; p81; VIL2; Villin 2
Storage/Stability:	Store at -20°C/1 year
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.
Immunogen:	The antiserum was produced against synthesized peptide derived from Internal of human Ezrin.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Ezrin (Ab-146) antibody detects endogenous levels of total Ezrin protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Gould K.L., EMBO J. 8:4133-4142(1989). Bechtel S., BMC Genomics 8:399-399(2007). Mungall A.J., Nature 425:805-811(2003).