

VAV2 (Ab-142) Antibody [#B1241]

Catalog Number:	B1241
Concentration:	1mg/ml
Swiss-Prot No.:	P52735
Other Names:	Protein vav-2, VAV2
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 non-phosphopeptide derived from human VAV2 around the phosphorylation site of tyrosine 142 (D-V-Y ^P -R-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	VAV2 (Ab-142) antibody detects endogenous levels of total VAV2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Henske E.P., Ann. Hum. Genet. 59:25-37(1995). Tamas P., J. Biol. Chem. 278:5163-5171(2003). RIKEN structural genomics initiative (RSGI); Submitted (APR-2007) to the PDB data bank.