

VAV1 Antibody [#B0593]

Catalog Number:	B0593
Concentration:	1mg/ml
Swiss-Prot No.:	P15498
Other Names:	VAV, Vav proto-oncogene
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 VAV1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	VAV1 antibody detects endogenous levels of total VAV1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Coppola J., Cell Growth Differ. 2:95-105(1991). Katzav S., Mol. Cell. Biol. 11:1912-1920(1991). Ramos-Morales F., Oncogene 11:1665-1669(1995).