

WAVE1 (Ab-125) Antibody [#B0598]

Catalog Number:	B0598
Concentration:	1mg/ml
Swiss-Prot No.:	Q92558
Other Names:	KIAA0269, SCAR1, Verprolin homology domain-containing protein 1, WAS1, WASF1, WASP-family protein member 1, Wiskott-Aldrich syndrome protein family member 1
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 WAVE1 around the phosphorylation site of tyrosine 125 (E-T-Y ^P -D-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	WAVE1 (Ab-125) antibody detects endogenous levels of total WAVE1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	John P. Dahl, J. Neurosci., Apr 2003; 23: 3343. Aiwu Cheng, J. Neurosci., Feb 2007; 27: 1519 - 1528. Koji Eto, Blood, Jul 2007; 10.1182/blood-2007-04-085860. Atsushi Oda, Blood, Apr 2005; 105: 3141 - 3148.