

VEGFR2 (Ab-1175) Antibody [#B7252]

Catalog Number:	B7252
Concentration:	1mg/ml
Swiss-Prot No.:	P35968
Other Names:	EC 2.7.10.1; FLK1; KDR; kinase insert domain receptor; protein-tyrosine kinase receptor Flk-1; vascular endothelial growth factor receptor 2; VGFR2; VGR2
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 VEGFR2 around the phosphorylation site of tyrosine 1175 (K-D-Y ^P -I-V).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	VEGFR2 (Ab-1175) antibody detects endogenous levels of total VEGFR2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Yin L.Y., Submitted (DEC-1997) to the EMBL/GenBank/DDBJ databases. Yu Y., Submitted (MAY-1998) to the EMBL/GenBank/DDBJ databases. Terman B.I., Oncogene 6:1677-1683(1991).