

## VEGFR1 Antibody [#B0594]

<b>Catalog Number:</b>	B0594
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P17948
<b>Other Names:</b>	EC 2.7.1.112, EC 2.7.10.1, FLT, FLT1, FRT, Flt-1, Fms-like tyrosine kinase 1, Tyrosine-protein kinase FRT, Tyrosine-protein kinase receptor FLT, VGFR1, Vascular endothelial growth factor receptor 1 precursor, Vascular permeability factor receptor
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human VEGFR1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	VEGFR1 antibody detects endogenous levels of total VEGFR1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 90%, Positives = 100%); Rat (Identities = 90%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	G. S. Robinson, FASEB J, Mar 2001; 10.1096/fj.00-0598fje. Donald P. Lesslie, AACR Meeting Abstracts, Apr 2005; 2005: 1305. Jack L. Arbiser, Am. J. Pathol., Apr 2000; 156: 1469. Feilim Mac Gabhann, Am J Physiol Heart Circ Physiol, Jan 2007; 292: H459 - H474.