

## EDD Antibody [#C0354]

<b>Catalog Number:</b>	C0354
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O95071
<b>Other Names:</b>	CAK, CD167a antigen, Cell adhesion kinase, Discoidin receptor tyrosine kinase, EC 2.7.1.112, EC 2.7.10.1, EDDR1, Epithelial discoidin domain receptor 1 precursor, HGK2, Protein-tyrosine kinase RTK 6, RTK6, TRK E, TRKE, Tyrosine kinase DDR, Tyrosine- protein kinase CAK, DDR1
<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 EDD.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	EDD antibody detects endogenous levels of total EDD protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	Michelle J. Henderson, J. Biol. Chem., Dec 2006; 281: 39990 - 40000. Lisa N. Kinch, Protein Sci., Feb 2005; 14: 360 - 367. Michelle J. Henderson, J. Biol. Chem., Jul 2002; 277: 26468 - 26478. Darren N. Saunders, Mol. Cell. Biol., Aug 2004; 24: 7225 - 7234.