

GRAK Antibody [#C13060]

Catalog Number:	C13060
Concentration:	1mg/ml
Swiss-Prot No.:	P49863
Other Names:	Granzyme K [Precursor], EC 3.4.21.-, Granzyme-3, NK-tryptase-2, NK-TRYP-2, Fragmentin-3, GZMK, TRYP2
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 GRAK.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRAK antibody detects endogenous levels of total GRAK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Przetak M.M., FEBS Lett. 364:268-271(1995). Sayers T.J., J. Leukoc. Biol. 59:763-768(1996). Hameed A., J. Immunol. 141:3142-3147(1988).