

GRAA Antibody [#C13058]

Catalog Number:	C13058
Concentration:	1mg/ml
Swiss-Prot No.:	P12544
Other Names:	Granzyme A , Granzyme A , Cytotoxic T-lymphocyte proteinase 1, Hanukkah factor, H factor, HF, Granzyme-1, CTL tryptase, Fragmentin-1, GZMA, CTLA3, HFSP
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 GRAA.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRAA antibody detects endogenous levels of total GRAA protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	The MGC Project Team, e Res. 14:2121-2127(2004). Bell J.K., Nat. Struct. Biol. 10:527-534(2003). Hink-Schauer C., Nat. Struct. Biol. 10:535-540(2003).