

CST8 Antibody [#C15271]

Catalog Number:	C15271
Concentration:	1mg/ml
Swiss-Prot No.:	O60676
Other Names:	Cystatin-8; Cystatin-related epididymal spermatogenic protein
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 C-terminal of human CST8.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CST8 antibody detects endogenous levels of total CST8 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Deloukas P., Nature 414:865-871(2001). Cornwall G.A., Mol. Reprod. Dev. 41:37-46(1995). Ramachandran P., J. Proteome Res. 5:1493-1503(2006).