

CLUS Antibody [#C13035]

Catalog Number:	C13035
Concentration:	1mg/ml
Swiss-Prot No.:	P10909
Other Names:	Clusterin [Precursor]; Complement-associated protein SP-40,40; Complement cytolysis inhibitor; CLI; NA1/NA2; Apolipoprotein J; Apo-J; Testosterone-repressed prostate message 2; TRPM-2; Clusterin beta chain (ApoJalpha) (Complement cytolysis inhibitor a chain); Clusterin alpha chain (ApoJbeta) (Complement cytolysis inhibitor b chain); CLU; APOJ, CLI
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 CLUS.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CLUS antibody detects endogenous levels of total CLUS 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:	Kirschbaum L., EMBO J. 8:711-718(1989). de Silva H.V., Biochemistry 29:5380-5389(1990). Wong P., Eur. J. Biochem. 221:917-925(1994).