

Collagen XXV Antibody [#C12227]

Catalog Number:	C12227
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BXS0
Other Names:	Collagen alpha-1(XXV) chain, CLAC-P, Alzheimer disease amyloid-associated protein, AMY, Collagen-like Alzheimer amyloid plaque component, COPA1
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 Collagen XXV.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen XXV antibody detects endogenous levels of total Collagen XXV protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Hashimoto T., EMBO J. 21:1524-1534(2002). Soederberg L., J. Neuropathol. Exp. Neurol. 62:1108-1117(2003). Soederberg L., J. Biol. Chem. 280:1007-1015(2005).