

SEMA4A Antibody [#C11691]

Catalog Number:	C11691
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H3S1
Other Names:	Semaphorin-4A [Precursor], Semaphorin-B, Sema B, SEMA4A, SEMAB, SEMB, UNQ783/PRO1317
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 SEMA4A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SEMA4A antibody detects endogenous levels of total SEMA4A protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	Clark H.F., Genome Res. 13:2265-2270(2003). The MGC Project Team, Genome Res. 14:2121-2127(2004). Abid A., J. Med. Genet. 43:378-381(2006).