

STMN4 Antibody [#C18953]

Catalog Number:	C18953
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H169
Other Names:	RB3; stathmin-4; stathmin-like 4; stathmin-like B3; stathmin-like-protein RB3
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 STMN4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	STMN4 antibody detects endogenous levels of total STMN4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Coadwell W.J., Submitted (JAN-2001) to the EMBL/GenBank/DDBJ databases. Wiemann S., Genome Res. 11:422-435(2001). The MGC Project Team, Genome Res. 14:2121-2127(2004).