

SFRS7 Antibody [#C18943]

Catalog Number:	C18943
Concentration:	1mg/ml
Swiss-Prot No.:	Q16629
Other Names:	SFR7; Splicing factor 9G8; Splicing factor, arginine/serine-rich 7
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 SFRS7.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SFRS7 antibody detects endogenous levels of total SFRS7 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Cavaloc Y., EMBO J. 13:2639-2649(1994). Popielarz M., J. Biol. Chem. 270:17830-17835(1995). The MGC Project Team, Genome Res. 14:2121-2127(2004).