

hnRNP L Antibody [#C10375]

Catalog Number:	C10375
Concentration:	1mg/ml
Swiss-Prot No.:	P14866
Other Names:	HNRPL, HNRPL;, heterogeneous nuclear ribonucleoprotein L, hnRNP-L
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 N-terminal of human hnRNP L.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	hnRNP L antibody detects endogenous levels of total hnRNP L protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Pinol-Roma S., J. Cell Biol. 109:2575-2587(1989). Ito M., Cancer Res. 61:2038-2046(2001). Rasmussen H.H., Electrophoresis 13:960-969(1992).