

ELAV2/4 Antibody [#C10486]

Catalog Number:	C10486
Concentration:	1mg/ml
Swiss-Prot No.:	Q12926/P26378
Other Names:	ELAV-like protein 4, Paraneoplastic encephalomyelitis antigen HuD, Hu-antigen D, ELAVL4, HUD, PNEM
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 ELAV2/4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ELAV2/4 antibody detects endogenous levels of total ELAV2/4 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
References:	Szabo A., Cell 67:325-333(1991). Behrends U., Int. J. Cancer 100:669-677(2002). Gregory S.G., Nature 441:315-321(2006).