

YTHDF1 Antibody [#C11643]

Catalog Number:	C11643
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BYJ9
Other Names:	YTH domain family protein 1, Dermatomyositis associated with cancer putative autoantigen 1, DACA-1, YTHDF1, C20orf21
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 YTHDF1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	YTHDF1 antibody detects endogenous levels of total YTHDF1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Ota T., Nat. Genet. 36:40-45(2004). Deloukas P., Nature 414:865-871(2001). The MGC Project Team, Genome Res. 14:2121-2127(2004).