

DCP1A Antibody [#C11740]

Catalog Number:	C11740
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NPI6
Other Names:	mRNA-decapping enzyme 1A, EC 3.-.-., Transcription factor SMIF, Smad4-interacting transcriptional co-activator, DCP1A, SMIF
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 DCP1A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	DCP1A antibody detects endogenous levels of total DCP1A protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Lykke-Andersen J., Mol. Cell. Biol. 22:8114-8121(2002). Bai R.-Y., Nat. Cell Biol. 4:181-190(2002). Ota T., Nat. Genet. 36:40-45(2004).