

ELOA2 Antibody [#C13048]

Catalog Number:	C13048
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IYF1/Q8NG57
Other Names:	RNA polymerase II transcription factor SIII subunit A2, Elongin-A2, Elongin-A2; EloA2, Transcription elongation factor B polypeptide 3B, TCEB3B, TCEB3L
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 ELOA2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ELOA2 antibody detects endogenous levels of total ELOA2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:5000
References:	Aso T., J. Biol. Chem. 275:6546-6552(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004). Sjoeblohm T., Science 314:268-274(2006).