

## EIF3K Antibody [#C15707]

<b>Catalog Number:</b>	C15707
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9UBQ5
<b>Other Names:</b>	Eukaryotic translation initiation factor 3 subunit K, Eukaryotic translation initiation factor 3 subunit 12; eIF-3 p25; eIF-3 p28; eIF3k; Muscle-specific gene M9 protein; PLAC-24
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human EIF3K.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	EIF3K antibody detects endogenous levels of total EIF3K protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Mayeur G.L., Eur. J. Biochem. 270:4133-4139(2003). Nawa G., Submitted (OCT-1998) to the EMBL/GenBank/DDBJ databases. Ye M., Submitted (JUL-1998) to the EMBL/GenBank/DDBJ databases.