

EIF2AK1 Antibody [#C15698]

Catalog Number:	C15698
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BQI3
Other Names:	E2AK1; EC 2.7.11.1; EIF2AK1; eukaryotic translation initiation factor 2 alpha kinase 1; HCR; heme-controlled repressor; heme-regulated eukaryotic initiation factor eIF-2-alpha kinase; heme-regulated inhibitor; hemin-sensitive initiation factor-2 alpha kinase; kinase HRI
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 EIF2AK1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	EIF2AK1 antibody detects endogenous levels of total EIF2AK1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 94%, Positives = 94%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Hwang S.-Y., Mol. Cells 10:584-591(2000). Cannon G., Submitted (AUG-1999) to the EMBL/GenBank/DDBJ databases. Nagase T., DNA Res. 7:65-73(2000).