

## GUF1 Antibody [#C16021]

<b>Catalog Number:</b>	C16021
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8N442
<b>Other Names:</b>	GTP-binding protein GUF1 homolog
<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 GUF1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GUF1 antibody detects endogenous levels of total GUF1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 84%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Ota T., Nat. Genet. 36:40-45(2004). The MGC Project Team, Genome Res. 14:2121-2127(2004). Vasilescu J., J. Proteome Res. 6:298-305(2007).