

## IFM3 Antibody [#C12462]

<b>Catalog Number:</b>	C12462
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q01628
<b>Other Names:</b>	IFM3; INIB; interferon-induced transmembrane protein 3; interferon-inducible protein 1-8U
<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 IFM3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	IFM3 antibody detects endogenous levels of total IFM3 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Lewin A.R., Eur. J. Biochem. 199:417-423(1991). The MGC Project Team; Genome Res. 14:2121-2127(2004).