

## ELOVL3 Antibody [#C15619]

<b>Catalog Number:</b>	C15619
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9HB03
<b>Other Names:</b>	Elongation of very long chain fatty acids protein 3, Cold-inducible glycoprotein of 30 kDa
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 ELOVL3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ELOVL3 antibody detects endogenous levels of total ELOVL3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100   ELISA: 1:20000
<b>References:</b>	The MGC Project Team; Genome Res. 14:2121-2127(2004). Semina E.V., Submitted (AUG-2000) to the EMBL/GenBank/DDBJ databases.