

## PMS1 Antibody [#C13103]

<b>Catalog Number:</b>	C13103
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P54277
<b>Other Names:</b>	DNA mismatch repair protein PMS1, PMSL1
<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 PMS1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PMS1 antibody detects endogenous levels of total PMS1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100   ELISA: 1:10000
<b>References:</b>	Nicolaides N.C., Nature 371:75-80(1994). Rieder M.J., Submitted (APR-2003) to the EMBL/GenBank/DDBJ databases. Wang Q., Hum. Genet. 105:79-85(1999).