

## MMP-10 Antibody [#C0263]

<b>Catalog Number:</b>	C0263
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P09238
<b>Other Names:</b>	STMY2, STROMELYSIN II, matrix metalloproteinase 10 (stromelysin 2)
<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 MMP-10.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	MMP-10 antibody detects endogenous levels of total MMP-10 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 81%, Positives = 100%); Rat (Identities = 81%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Muller D., Biochem. J. 253:187-192(1988). Kalnine N., Submitted (OCT-2004) to the EMBL/GenBank/DDBJ databases. The MGC Project Team; Genome Res. 14:2121-2127(2004). Livingston R.J., Submitted (SEP-2004) to the EMBL/GenBank/DDBJ databases.