

## GSTT1/4 Antibody [#C15930]

<b>Catalog Number:</b>	C15930
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P30711/A8MPT4
<b>Other Names:</b>	EC 2.5.1.18; glutathione S-transferase theta 1; glutathione transferase T1-1; GST class-theta-1
<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 GSTT1/4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	GSTT1/4 antibody detects endogenous levels of total GSTT1/4 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Pemble S., Biochem. J. 300:271-276(1994). Jemth P., Arch. Biochem. Biophys. 348:247-254(1997). Sprenger R., Pharmacogenetics 10:557-565(2000).