

## PXMP4 Antibody [#C17633]

<b>Catalog Number:</b>	C17633
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9Y6I8
<b>Other Names:</b>	24 kDa peroxisomal intrinsic membrane protein; PMP24
<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 N-terminal of human PXMP4.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PXMP4 antibody detects endogenous levels of total PXMP4 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Reguenga C., Biochim. Biophys. Acta 1445:337-341(1999). Wiemann S., Genome Res. 11:422-435(2001). Deloukas P., Nature 414:865-871(2001).