

POM121 Antibody [#C17074]

Catalog Number:	C17074
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y2N3/Q96HA1/A6NF01/A8CG34
Other Names:	nuclear envelope pore membrane protein POM 121; NUP121; P145; PO121; POM121 membrane glycoprotein
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human POM121.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	POM121 antibody detects endogenous levels of total POM121 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ishikawa K., DNA Res. 5:169-176(1998). Ohara O., Submitted (JAN-2004) to the EMBL/GenBank/DDBJ databases. Ota T., Nat. Genet. 36:40-45(2004).