

PITPNB Antibody [#C17679]

Catalog Number:	C17679
Concentration:	1mg/ml
Swiss-Prot No.:	P48739
Other Names:	Phosphatidylinositol transfer protein beta isoform; PI-TP-beta; PI-TPbeta; PIPNB; PPI2; PtdIns transfer protein beta; PtdInsTP
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 PITPNB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PITPNB antibody detects endogenous levels of total PITPNB protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%); Rat (Identities = 92%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Tanaka S., Biochim. Biophys. Acta 1259:199-202(1995). Collins J.E., Genome Biol. 5:RESEARCH84.1-RESEARCH84.11(2004). Dunham I., Nature 402:489-495(1999).