

PEX11B Antibody [#C17630]

Catalog Number:	C17630
Concentration:	1mg/ml
Swiss-Prot No.:	O96011
Other Names:	peroxin-11B; peroxisomal biogenesis factor 11B; PEX11-BETA; PEX11beta; PX11B
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 PEX11B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PEX11B antibody detects endogenous levels of total PEX11B protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Abe I., Biochem. Biophys. Res. Commun. 252:529-533(1998). Schrader M., J. Biol. Chem. 273:29607-29614(1998). Zuo D., Submitted (JUN-2004) to the EMBL/GenBank/DDBJ databases.