

Rabphilin 3A (Ab-237) Antibody [#B0563]

Catalog Number:	B0563
Concentration:	1mg/ml
Swiss-Prot No.:	P47708
Other Names:	RP3A, RPH3A, exophilin 1
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 non-phosphopeptide derived from human rabphilin 3A around the phosphorylation site of serine 234 (R-A-S ^P -E-A).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Rabphilin 3A (Ab-237) antibody detects endogenous levels of total rabphilin 3A protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	H Shirataki, Mol. Cell. Biol., Apr 1993; 13: 2061 - 2068. N Masumoto, J. Cell Biol., Dec 1996; 135: 1741. M.E. Burns, J. Gen. Physiol., Feb 1998; 111: 243.