

RGS16 Antibody [#B1232]

Catalog Number:	B1232
Concentration:	1mg/ml
Swiss-Prot No.:	O15492
Other Names:	A28-RGS14P, RGS-R, RGSG, RGSR, Regulator of G-protein signaling 16, Retinally abundant regulator of G-protein signaling
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 RGS16.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RGS16 antibody detects endogenous levels of total RGS16 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Buckbinder L., Proc. Natl. Acad. Sci. U.S.A. 94:7868-7872(1997). Snow B.E., Gene 206:247-253(1998). Puhl H.L. III, Submitted (MAR-2002) to the EMBL/GenBank/DDBJ databases.