

Rho/Rac Guanine Nucleotide Exchange Factor 2 (Ab-885) Antibody [#B1233]

Catalog Number:	B1233
Concentration:	1mg/ml
Swiss-Prot No.:	Q8TDA3
Other Names:	GEF-H1, GEFH1, Guanine nucleotide exchange factor GEF-H1, proliferating cell nucleolar antigen p40, rho/rac guanine nucleotide exchange factor 2
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 Rho/Rac Guanine Nucleotide Exchange Factor 2 around the phosphorylation site of serine 885 (R-R-S ^P -L-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Rho/Rac Guanine Nucleotide Exchange Factor 2 (Ab-885) antibody detects endogenous levels of total Rho/Rac Guanine Nucleotide Exchange Factor 2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Krendel M., Nat. Cell Biol. 4:294-301(2002). Mural R.J., Submitted (SEP-2005) to the EMBL/GenBank/DDBJ databases. Hall R., Submitted (FEB-2008) to the EMBL/GenBank/DDBJ databases.