

## ARHGDI Antibody [#B1189]

<b>Catalog Number:</b>	B1189
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P52565
<b>Other Names:</b>	GDIR, Rho GDI 1, Rho GDP-dissociation inhibitor 1, Rho-GDI alpha
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human ARHGDI.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ARHGDI antibody detects endogenous levels of total ARHGDI protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 90%, Positives = 90%); Rat (Identities = 90%, Positives = 90%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Leffers H., Exp. Cell Res. 209:165-174(1993). Maeda A., Submitted (APR-1993) to the EMBL/GenBank/DDBJ databases. Chuang T.H., Submitted (JUL-1992) to the EMBL/GenBank/DDBJ databases.