

BAI2 Antibody [#G214]

Catalog Number:	G214
Concentration:	1mg/ml
Swiss-Prot No.:	O60241
Other Names:	Brain-specific angiogenesis inhibitor 2;BAI2
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 BAI2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BAI2 antibody detects endogenous levels of total BAI2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 86%, Positives = 93%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Shiratsuchi T., Cytogenet. Cell Genet. 79:103-108(1997). Suwa M., Submitted (JUL-2001) to the EMBL/GenBank/DDBJ databases. Gregory S.G., Nature 441:315-321(2006).