

BRI3B Antibody [#C11418]

Catalog Number:	C11418
Concentration:	1mg/ml
Swiss-Prot No.:	Q8WY22
Other Names:	BRI3-binding protein, I3-binding protein, Cervical cancer 1 proto-oncogene-binding protein KG19, HCCR-1, BRI3BP, KG19
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 BRI3B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BRI3B antibody detects endogenous levels of total BRI3B protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Lin L., Biochem. Genet. 39:369-377(2001). The MGC Project Team, Genome Res. 14:2121-2127(2004). Kim J.W., Submitted (FEB-2002) to the EMBL/GenBank/DDBJ databases.