

BRS3 Antibody [#G217]

Catalog Number:	G217
Concentration:	1mg/ml
Swiss-Prot No.:	P32247
Other Names:	Bombesin receptor subtype-3; BRS-3
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 BRS3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BRS3 antibody detects endogenous levels of total BRS3 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Fathi Z., J. Biol. Chem. 268:5979-5984(1993). Gorbulev V., FEBS Lett. 340:260-264(1994). Ross M.T., Nature 434:325-337(2005).