

BRMS1 Antibody [#C11729]

Catalog Number:	C11729
Concentration:	1mg/ml
Swiss-Prot No.:	Q9HCU9
Other Names:	Breast cancer metastasis-suppressor 1, BRMS1
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 N-terminal of human BRMS1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	BRMS1 antibody detects endogenous levels of total BRMS1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Seraj M.J., Cancer Res. 60:2764-2769(2000). The MGC Project Team, Genome Res. 14:2121-2127(2004). Janneau J.-L., . Clin. Endocrinol. Metab. 87:5336-5339(2002).