

CRFR2 Antibody [#G228]

Catalog Number:	G228
Concentration:	1mg/ml
Swiss-Prot No.:	Q13324
Other Names:	Corticotropin-releasing factor receptor 2 [Precursor]; CRF-R 2; CRF2; Corticotropin-releasing hormone receptor 2; CRH-R 2; CRHR2; CRF2R, CRH2R
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 CRFR2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CRFR2 antibody detects endogenous levels of total CRFR2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Liaw C.W., Endocrinology 137:72-77(1996). Kostich W.A., Mol. Endocrinol. 12:1077-1085(1998). Hillier L.W., Nature 424:157-164(2003).