

CNR1 Antibody [#G226]

Catalog Number:	G226
Concentration:	1mg/ml
Swiss-Prot No.:	P21554
Other Names:	CANN6; cannabinoid receptor 1; CB-R; CB1; CB1R; CNR
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 CNR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CNR1 antibody detects endogenous levels of total CNR1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Gerard C., Nucleic Acids Res. 18:7142-7142(1990). Gerard C., Biochem. J. 279:129-134(1991). Shire D., J. Biol. Chem. 270:3726-3731(1995).