

GRIN2D Antibody [#C15922]

Catalog Number:	C15922
Concentration:	1mg/ml
Swiss-Prot No.:	O15399
Other Names:	EB11; estrogen receptor binding CpG island; glutamate [NMDA] receptor epsilon 4; glutamate receptor, ionotropic, N-methyl D-aspartate 2D; GRIN2D; N-methyl D-aspartate receptor 2D; NMDE4; NR2D
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 GRIN2D.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRIN2D antibody detects endogenous levels of total GRIN2D protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Hess S.D., J. Neurochem. 70:1269-1279(1998). Grimwood J., Nature 428:529-535(2004). Sjoebloom T., Science 314:268-274(2006).