

JIP3 Antibody [#C11903]

Catalog Number:	C11903
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UPT6
Other Names:	C-jun-amino-terminal kinase interacting protein 3, JIP-3, JIP3, JNK MAP kinase scaffold protein 3, JNK-interacting protein 3, JNK/SAPK-associated protein 1, JSAP1, KIAA1066, MAPK8IP3, Mitogen- activated protein kinase 8-interacting protein 3, SYD2, Sunday driver 2
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 JIP3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	JIP3 antibody detects endogenous levels of total JIP3 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Hattori A., DNA Res. 7:357-366(2000). Martin J., Nature 432:988-994(2004). Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004).