

MAFF Antibody [#C11887]

Catalog Number:	C11887
Concentration:	1mg/ml
Swiss-Prot No.:	Q9ULX9
Other Names:	Transcription factor MafF, V-maf musculoaponeurotic fibrosarcoma oncogene homolog F, U-Maf, MAFF
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 MAFF.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MAFF antibody detects endogenous levels of total MAFF protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Moran J.A., Biochem. J. 361:371-377(2002). Dunham I., Nature 402:489-495(1999).