

MRRF Antibody [#C18437]

Catalog Number:	C18437
Concentration:	1mg/ml
Swiss-Prot No.:	Q96E11
Other Names:	Ribosome-recycling factor, mitochondrial; RRF; Ribosome-releasing factor, mitochondrial
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 MRRF.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRRF antibody detects endogenous levels of total MRRF protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 86%, Positives = 100%); Rat (Identities = 86%, Positives = 100%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Ota T., Nat. Genet. 36:40-45(2004). Humphray S.J., Nature 429:369-374(2004). Zhang Y., Biochim. Biophys. Acta 1443:245-250(1998).