

RNUXA Antibody [#C18456]

Catalog Number:	C18456
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H814
Other Names:	PHAX; phosphorylated adapter RNA export protein; RNA U small nuclear RNA export adapter
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 RNUXA.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RNUXA antibody detects endogenous levels of total RNUXA protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Ebert L., Submitted (JUN-2004) to the EMBL/GenBank/DDBJ databases. Boulon S., Mol. Cell 16:777-787(2004).