

Ubinuclein Antibody [#C11739]

Catalog Number:	C11739
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NPG3
Other Names:	Ubiquitously expressed nuclear protein, VT4, Ubinuclein, UBN1
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 Ubinuclein.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Ubinuclein antibody detects endogenous levels of total Ubinuclein protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Aho S., J. Cell Biol. 148:1165-1176(2000). Tagami H., Cell 116:51-61(2004). Beausoleil S.A., Proc. Natl. Acad. Sci. U.S.A. 101:12130-12135(2004).