

SUPT16H Antibody [#C11966]

Catalog Number:	C11966
Concentration:	1mg/ml
Swiss-Prot No.:	Q9Y5B9
Other Names:	CDC68, Chromatin-specific transcription elongation factor 140 kDa subunit, FACT140
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 SUPT16H.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SUPT16H antibody detects endogenous levels of total SUPT16H protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Kang S.-W., Genes Cells 5:251-263(2000). Wada T., Mol. Cell 5:1067-1072(2000). Keller D.M., Mol. Cell 7:283-292(2001).