

H-NUC Antibody [#C12168]

Catalog Number:	C12168
Concentration:	1mg/ml
Swiss-Prot No.:	P30260
Other Names:	CDC27Hs, Cell division cycle protein 27 homolog, H-NUC
Storage/Stability:	Store at -20°C/1 yearH-NUC
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 H-NUC.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	H-NUC antibody detects endogenous levels of total H-NUC protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Tugendreich S., Proc. Natl. Acad. Sci. U.S.A. 90:10031-10035(1993). Chen P.L., Cell Growth Differ. 6:199-210(1995). Kraft C., EMBO J. 22:6598-6609(2003).