

ASF1B Antibody [#C11765]

Catalog Number:	C11765
Concentration:	1mg/ml
Swiss-Prot No.:	Q9NVP2
Other Names:	Histone chaperone ASF1B, Anti-silencing function protein 1 homolog B, hAsf1, hAsf1b, CCG1-interacting factor A-II, CIA-II, hCIA-II, ASF1B
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 ASF1B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	ASF1B antibody detects endogenous levels of total ASF1B protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Sillje H.H.W., Curr. Biol. 11:1068-1073(2001). ehara T., J. Biol. Chem. 278:35660-35667(2003). Ota T., Nat. Genet. 36:40-45(2004).