

## SKP1A/p19 Antibody [#C0323]

<b>Catalog Number:</b>	C0323
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P63208
<b>Other Names:</b>	EMC19, OCP-2, OCP-II, OCP2, RNA polymerase II elongation factor-like, RNA polymerase II elongation factor-like OCP2, S-phase kinase-associated 1A, S-phase kinase-associated 1A (p19A), SIII, SKP1, TCEB1L, cyclin A/CDK2-associated p19, organ of Corti 2, p19A, p19skp1, transcription elongation factor B, transcription elongation factor B (SIII), polypeptide 1-like
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from Internal of human SKP1A/p19.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	SKP1A/p19 antibody detects endogenous levels of total SKP1A/p19 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
<b>References:</b>	SILVIA MANDEL, Ann. N.Y. Acad. Sci., Aug 2005; 1053: 356 - 375. Deborah French, Blood (ASH Annual Meeting Abstracts), Nov 2004; 104: 144. Patana Teng-umnuay, J. Biol. Chem., Dec 1999; 274: 36392. Slim Sassi, Glycobiology, Apr 2001; 11: 283 - 295.