

## Ki67 Antibody [#C0250]

<b>Catalog Number:</b>	C0250
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P46013
<b>Other Names:</b>	Antigen KI-67, MKI67
<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 C-terminal of human ki67.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Ki67 antibody detects endogenous levels of total ki67protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Mitch Dowsett, J Natl Cancer Inst, Jan 2007; 99: 167 - 170. Phyllis A. Gimotty, J. Clin. Oncol, Nov 2005; 23: 8048 - 8056. Tim J. Dudderidge, Clin. Cancer Res., Apr 2005; 11: 2510 - 2517. Scott F. Sieg, J. Immunol., Nov 2003; 171: 5208.