

## p73 Antibody [#C10080]

<b>Catalog Number:</b>	C10080
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O15350
<b>Other Names:</b>	p53-like transcription factor; p53-related protein; P73 alpha protein; p73-alpha; TP73; TRP73; Tumor protein p73
<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 p73.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p73 antibody detects endogenous levels of total p73 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Kaghad M., Cell 90:809-819(1997). Mai M., Genomics 51:359-363(1998). De Laurenzi V., J. Exp. Med. 188:1763-1768(1998).