

## p97 MAPK (Ab-189) Antibody [#B0919]

<b>Catalog Number:</b>	B0919
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q16659
<b>Other Names:</b>	EC 2.7.11.24, ERK-3, ERK3, Extracellular signal- regulated kinase 3, MAP kinase isoform p97, MAPK6, MK06, Mitogen-activated protein kinase 6, PRKM4, PRKM6, kinase ERK3, p55-MAPK, p97-MAPK
<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 non-phosphopeptide derived from human p97 MAPK around the phosphorylation site of serine 189 (H-L-S <sup>P</sup> -E-G).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	p97 MAPK (Ab-189) antibody detects endogenous levels of total p97 MAPK protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Zhu A.X., Mol. Cell. Biol. 14:8202-8211(1994). Meloche S., Oncogene 13:1575-1579(1996). Saranath D., Submitted (SEP-2001) to the EMBL/GenBank/DDBJ databases.