

## Desmin (Ab-60) Antibody [#B1196]

<b>Catalog Number:</b>	B1196
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P17661
<b>Other Names:</b>	DES; DESM
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 Desmin around the phosphorylation site of serine 60 (R-T-S <sup>P</sup> -G-G).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Desmin (Ab-60) antibody detects endogenous levels of total Desmin protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Li Z., Gene 78:243-254(1989). Vicart P., Hum. Genet. 98:422-429(1996). Goldfarb L.G., Nat. Genet. 19:402-403(1998).