

## Caldesmon (Ab-789) Antibody [#B0462]

<b>Catalog Number:</b>	B0462
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q05682
<b>Other Names:</b>	CAD, CALD1, CDM, L-caldesmon, Non-muscle caldesmon
<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 Caldesmon around the phosphorylation site of serine 789 (V-T-S <sup>P</sup> -P-T).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Caldesmon (Ab-789) antibody detects endogenous levels of total Caldesmon protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Robert Eves, J. Cell Sci., May 2006; 119: 1691 - 1702. Shigeko Yamashiro, Mol. Biol. Cell, Jan 2001; 12: 239. KS Warren, J. Cell Biol., Apr 1994; 125: 359.