

## ATF-2 (Ab-472) Antibody [#B8294]

<b>Catalog Number:</b>	B8294
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P15336
<b>Other Names:</b>	Activating 2; ATF2; cAMP response element binding protein CRE- BP1; CREB2; CREBP1; Cyclic-AMP-dependent ATF-2; HB16; MXBP protein
<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 ATF-2 around the phosphorylation site of serine 472 (A-L-S <sup>P</sup> -Q-I).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ATF-2 (Ab-472) antibody detects endogenous levels of total ATF-2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Maekawa T., EMBO J. 8:2023-2028(1989). Yang L., J. Immunol. 158:2522-2525(1997). Hillier L.W., Nature 434:724-731(2005).