

CREB (Ab-121) Antibody [#B0471]

Catalog Number:	B0471
Concentration:	1mg/ml
Swiss-Prot No.:	P16220
Other Names:	CREB-1, CREB1, cAMP-response element binding protein
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human CREB around the phosphorylation site of serine 121 (T-D-S ^P -Q-K).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CREB (Ab-121) antibody detects endogenous levels of total CREB protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Lionel Groussin, J. Clin. Endocrinol. Metab., Jan 2000; 85: 345 - 354. Deepa B. Shankar, Blood (ASH Annual Meeting Abstracts), Nov 2005; 106: 367. LA Berkowitz and MZ Gilman, PNAS, Jul 1990; 87: 5258.