

Anti-CREB

Rabbit Polyclonal Antibody

Catalog # C49-63R

Lot # H200-3

Cited Applications

- Immunoprecrecipitation (2 µg)
- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the CREB protein

Cross Reactivity

- Western blot and immunoprecipitation of human and mouse CREB

CREB from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 127-131 of human CREB

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

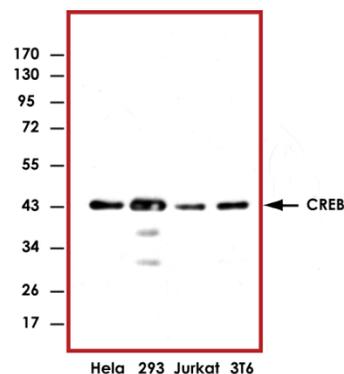
Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

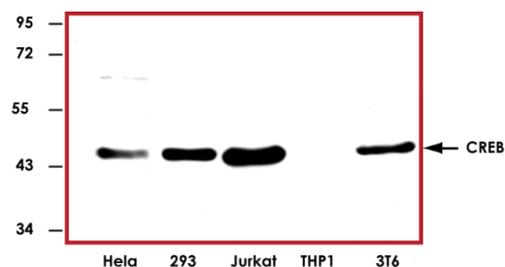
Scientific Background

CREB1 is a member of the cAMP responsive element binding protein family of transcription factors that are critical mediators of gene expression in response to extracellular signals and are essential regulators of adaptive behavior and long-term memory formation (1). Activation of CREB1 by protein kinase A-mediated phosphorylation has been implicated in the survival of mammalian cells. Mice lacking CREB in the central nervous system during development show extensive apoptosis of postmitotic neurons (2). CREB binds as homo- and heterodimers to promoters containing CRE and activator protein 1 (AP-1) sites to alter target-gene expression).

Sample Data



Representative immunoprecipitation with 2 µg of Anti-CREB using 500 µg of HeLa, HEK-293T, Jurkat, and 3T6 cell lysate.



Representative western blot with Anti-CREB (1:1000) using 30 µg of HeLa, HEK-293T, Jurkat, THP1, and 3T6 cell lysate.

References

1. Valverde, O. et al: Modulation of anxiety-like behavior and morphine dependence in CREB-deficient mice. *Neuropsychopharmacology*. 2004 Jun; 29(6): 1122-33.
2. Mantamadiotis, T. et al: Disruption of CREB function in brain leads to neurodegeneration. *Nat Genet*. 2002 May; 31(1): 47-54.

Anti-CREB

Rabbit Polyclonal Antibody

Catalog Number

C49-63R

Specific Lot Number

H200-3

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com

www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.