

PRODUCT INFORMATION



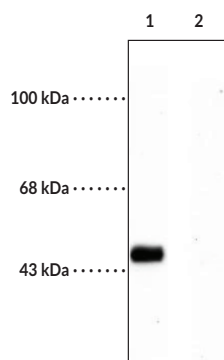
CREB (Phospho-Ser¹³³) Polyclonal Antibody

Item No. 10009181

Overview and Properties

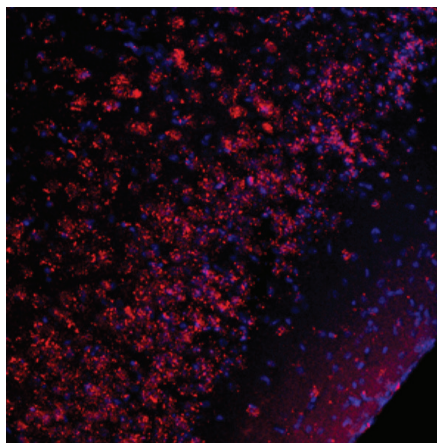
Contents:	This vial contains 100 µl of affinity-purified polyclonal antibody.
Immunogen:	Synthetic peptide from amino acid residues surrounding the phospho-Ser ¹³³ of rat CREB
Molecular Weight:	45 kDa
Cross Reactivity:	(+) Rat CREB
Species Reactivity:	(+) Mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:1,000 for IHC and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: - λ Ptase
Lane 2: + λ Ptase

WB of rat hippocampal lysate stimulated with forskolin showing specific immunolabeling of the ~45 kDa CREB phosphorylated at Ser¹³³ in the first lane (-). Phosphospecificity is shown in the second lane (+) where immunolabeling is completely eliminated by lysate treatment with λ-Ptase (800 units/1 mg protein for 30 min).



Immunofluorescent labeling of mouse piriform cortex labeled with CREB (Phospho-Ser¹³³) Polyclonal Antibody (red). Cell nuclei have been labeled with DAPI DNA stain (blue).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/29/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

It is well known that the control of gene expression involves activation of protein kinase cascades that regulate transcription factors within the nucleus.¹ The cyclic AMP response element binding protein (CREB) is one of the best characterized stimulus-induced transcription factors.² This transcription factor is a component of intracellular signaling events that regulate a wide range of biological functions, from spermatogenesis to circadian rhythms and memory.^{3,4} A variety of protein kinases including protein kinase A (PKA), mitogen-activated protein kinases (MAPKs), and Ca²⁺/calmodulin-dependent protein kinases (CaMKs) phosphorylate CREB at serine 133 (Ser¹³³), and phosphorylation of Ser¹³³ is required for CREB-mediated transcription.^{5,6}

References

1. Karin, M. and Hunter, T. Transcriptional control by protein phosphorylation: Signal transmission from the cell surface to the nucleus. *Curr. Biol.* **5**(7), 747-757 (1995).
2. Montminy, M. Transcriptional regulation by cyclic AMP. *Annu. Rev. Biochem.* **66**, 807-822 (1997).
3. Shaywitz, A.J. and Greenberg, M.E. CREB: A stimulus-induced transcription factor activated by a diverse array of extracellular signals. *Annu. Rev. Biochem.* **68**, 821-861 (1999).
4. Silva, A.J., Kogan, J.H., Frankland, P.W., et al. CREB and memory. *Annu. Rev. Neurosci.* **21**, 127-148 (1998).
5. Johannessen, M., Delghandi, M.P., and Moens, U. What turns CREB on? *Cell. Signal.* **16**(11), 1211-1227 (2004).
6. Kornhauser, J.M., Cowan, C.W., Shaywitz, A.J., et al. CREB transcriptional activity in neurons is regulated by multiple, calcium-specific phosphorylation events. *Neuron* **34**(2), 221-233 (2002).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM