



## CREB (Phospho-Ser133) Transcription Factor Assay Kit

---

Item No. 10009846

[www.caymanchem.com](http://www.caymanchem.com)

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

## TABLE OF CONTENTS

<b>GENERAL INFORMATION</b>	3	Materials Supplied
	4	Safety Data
	4	Precautions
	5	If You Have Problems
	5	Storage and Stability
	5	Materials Needed but Not Supplied
<b>INTRODUCTION</b>	6	Background
	7	About This Assay
<b>PRE-ASSAY PREPARATION</b>	9	Sample Buffer Preparation
	14	Purification of Cellular Nuclear Extracts
	15	Reagent Preparation
<b>ASSAY PROTOCOL</b>	18	Plate Set Up
	20	Performing the Assay
	24	Assay Procedure Summary
<b>ANALYSIS</b>	26	Performance Characteristics
<b>RESOURCES</b>	28	Troubleshooting
	29	References
	30	Plate Template
	31	Notes
	31	Warranty and Limitation of Remedy

## GENERAL INFORMATION

### Materials Supplied

Kit components may be stored at -20°C prior to use. For long term storage, the Positive Control should be thawed on ice, aliquoted at 25 µl/vial, and stored at -80°C. After opening kit, we recommend each kit component be stored according to the temperature listed below.

Item Number	Item	Quantity	Storage
10006880	Transcription Factor Binding Assay Buffer (4X)	1 vial/3 ml	4°C
10007472	Transcription Factor Reagent A	1 vial/120 µl	-20°C
10009849	TF CREB (Phospho-Ser <sup>133</sup> ) Positive Control	1 vial/150 µl	-80°C
10006882	Transcription Factor Antibody Binding Buffer (10X)	1 vial/3 ml	4°C
10009848	TF CREB (Phospho-Ser <sup>133</sup> ) Primary Antibody	1 vial/120 µl	-20°C
400062	Wash Buffer Concentrate (400X)	1 vial/5 ml	RT
400035	Polysorbate 20	1 vial/3 ml	RT
10009488	TF CREB (Phospho-Ser <sup>133</sup> ) Competitor dsDNA	1 vial/120 µl	-20°C
10006884	Transcription Factor Goat Anti-Rabbit HRP Conjugate	1 vial/120 µl	-20°C
10009489	TF CREB (Phospho-Ser <sup>133</sup> ) 96-Well Strip Plate	1 plate	4°C
400012	96-Well Cover Sheet	1 cover	RT
10006888	Transcription Factor Developing Solution	1 vial/12 ml	4°C
10006889	Transcription Factor Stop Solution	1 vial/12 ml	RT

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



**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.

## Precautions

**Please read these instructions carefully before beginning this assay.**

Kit components may be stored at -20°C prior to use. For long term storage, the Positive Control should be thawed on ice, aliquoted at 25 µl/vial and stored at -80°C. If the assay will be used on multiple days, we recommend each kit component be stored according to the temperatures listed in the booklet.

## If You Have Problems

### Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888  
Fax: 734-971-3641  
E-Mail: techserv@caymanchem.com  
Hours: M-F 8:00 AM to 5:30 PM EST

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

## Storage and Stability

This kit will perform as specified if stored as directed in the **Materials Supplied** section, on page 3, and used before the expiration date indicated on the outside of the box.

## Materials Needed But Not Supplied

1. A plate reader capable of measuring absorbance at 450 nm
2. Adjustable pipettes and a repeating pipettor
3. A source of UltraPure water; glass Milli-Q or HPLC-grade water is acceptable
4. 300 mM dithiothreitol (DTT)
5. Nuclear Extraction Kit available from Cayman (Item No. 10009277) or buffers for preparation of nuclear extracts (see pages 9-13)

*NOTE: The components in each kit lot have been quality assured and warranted in this specific combination only; please do not mix them with components from other lots.*

## Background

Cyclic AMP response element-binding protein (CREB) belongs to a large family of structurally related transcription factors called bZIP and includes AFT1-4, c-Fos, c-Myc, c-Jun, and C/EBP.<sup>1,2</sup> These basic region leucine zipper (bZIP) transcription factors have a DNA-binding domain and a dimerization domain that contains the leucine zipper motif. CREB proteins recognize and bind to the cAMP-responsive element promoter (CRE) site and regulate transcription of many downstream genes that play important roles in metabolic regulation, depression, and in signaling pathways that enable long term memory.<sup>3</sup>

The best characterized activator of CREB is cAMP-dependent protein kinase A (PKA)-mediated phosphorylation of CREB on Serine133 (Ser<sup>133</sup>) following G protein-coupled receptor (GPCR) stimulation by growth factors, neurotransmitters, or peptide hormones.<sup>4</sup> For example, binding of ligands to their cognate GPCR ultimately leads to an immediate increase in intracellular cAMP levels resulting in liberation of the catalytic subunits of PKA. The free catalytic subunits enter the cell nucleus where they phosphorylate CREB at Ser<sup>133</sup> activating a cascade of events that includes recruitment of CREB-binding protein (CBP) and p300 along with other cofactors in the assembly of a large transcriptional complex.<sup>6</sup>

CBP and p300 contain intrinsic histone deacetylation (HDAC) activity that assist in the activation of transcription and ultimately enable the synthesis of RNA by RNA polymerase II. Additional diverse stimuli such as hypoxia, growth factors, UV light, survival signals, and stress signals are also known activators of CREB.<sup>4</sup> These various stimuli act through a variety of complex signaling pathways that converge to phosphorylate CREB and induce transcription or inactivate the CREB transcription complex. The activated kinases able to phosphorylate CREB include (but are not limited to) PKA, AKT, MAPK-activated ribosomal S6 kinases (RSKs), and MSK1. Many of these kinases phosphorylate CREB at residues other than Ser<sup>133</sup> to initiate binding to DNA and transcription of downstream targets. Other post-translational modifications of CREB include acetylation, sumoylation, and glycosylation which add to the complexity and diversity of its transcriptional regulation in the cell.<sup>5</sup> CREB is involved in long term memory and could potentially be utilized for targeted drug therapy for memory loss in Alzheimer's disease, stroke, drug abuse, and head trauma.

## About This Assay

Cayman's CREB (Phospho-Ser<sup>133</sup>) Transcription Factor Assay is a non-radioactive, sensitive method for detecting CREB DNA binding activity in nuclear extracts and whole cell lysates. A 96-well enzyme-linked immunosorbent assay (ELISA) replaces the cumbersome radioactive electrophoretic mobility shift assay (EMSA). A specific double stranded DNA (dsDNA) consensus sequence containing the cAMP response element (CRE) is immobilized to the wells of a 96-well plate (see Figure 1, on page 8). CREB contained in a nuclear extract or whole cell lysate binds specifically to the CRE. The activated CREB transcription factor complex is detected by addition of a specific primary antibody directed against Phospho-Ser<sup>133</sup> on CREB. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm.

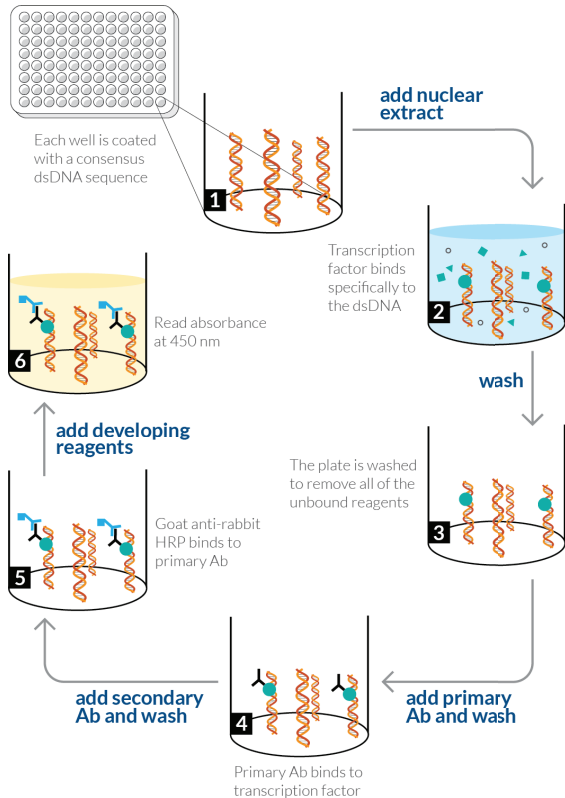


Figure 1. Schematic of the Transcription Factor Binding Assay

## PRE-ASSAY PREPARATION

### Sample Buffer Preparation

All buffers and reagents below required for preparation of Nuclear Extracts can be purchased directly from Cayman. Alternatively, Cayman's Nuclear Extraction Kit (Item No. 10009277) can be used to isolate Nuclear Proteins.

#### 1. Nuclear Extraction PBS (10X)

1.71 M NaCl, 33.53 mM KCl, 126.8 mM  $\text{Na}_2\text{HPO}_4$ , 22.04 mM  $\text{KH}_2\text{PO}_4$ , pH 7.4

#### 2. Nuclear Extraction PBS (1X)

Dilute 100 ml of 10X stock with 900 ml distilled  $\text{H}_2\text{O}$

#### 3. Nuclear Extraction Phosphatase Inhibitor Cocktail (50X)

0.5 M NaF

0.05 M  $\beta$ -glycerophosphate

0.05 M  $\text{Na}_3\text{VO}_4$

Store at  $-80^\circ\text{C}$

#### 4. Nuclear Extraction PBS/Phosphatase Inhibitor Solution (1X)

Add 200  $\mu\text{l}$  of 50X Phosphatase Inhibitor Solution to 10 ml of 1X Nuclear Extraction PBS, mix well, and keep on ice. Make fresh daily.

#### 5. Nuclear Extraction Protease Inhibitor Cocktail (100X)

10 mM AEBSF

0.5 mM Bestatin

0.2 mM Leupeptin Hemisulfate Salt

0.15 mM E-64

0.1 mM Pepstatin A

0.008 mM Aprotinin from Bovine Lung

Made in DMSO, store at -20°C

#### 6. Nuclear Extraction Hypotonic Buffer (10X)

100 mM HEPES, pH 7.5, containing 40 mM NaF, 100  $\mu$ M Na<sub>2</sub>MoO<sub>4</sub>, and 1 mM EDTA

Store at 4°C

#### 7. Complete Hypotonic Buffer (1X)

Prepare as outlined in Table 1. The phosphatase and protease inhibitors lose activity shortly after dilution; therefore any unused 1X Complete Hypotonic Buffer should be discarded.

Reagent	150 mm plate ~1.5 x 10 <sup>7</sup> cells
Hypotonic Buffer (10X)	100 $\mu$ l
Phosphatase Inhibitors (50X)	20 $\mu$ l
Protease Inhibitors (100X)	10 $\mu$ l
Distilled Water	870 $\mu$ l
Total Volume	1,000 $\mu$ l

Table 1. Preparation of Complete Hypotonic Buffer

**8. Nonidet P-40 Assay Reagent (10%)**

Nonidet P-40 or suitable substitute at a concentration of 10% (v/v) in H<sub>2</sub>O  
Store at room temperature

**9. Nuclear Extraction Buffer (2X)**

20 mM HEPES, pH 7.9, containing, 0.2 mM EDTA, 3 mM MgCl<sub>2</sub>, 840 mM NaCl, and 20% glycerol (v/v)

Store at 4°C

**10. Complete Nuclear Extraction Buffer (1X)**

Prepare as outlined in Table 2. Some of the phosphatase and protease inhibitors lose activity shortly after dilution; therefore any remaining 1X Extraction Buffer should be discarded.

Reagent	150 mm plate ~1.5 x 10 <sup>7</sup> cells
Nuclear Extraction Buffer (2X)	75 µl
Protease Inhibitors (100X)	1.5 µl
Phosphatase Inhibitors (50X)	3.0 µl
DTT (10 mM)	15 µl
Distilled Water	55.5 µl
Total Volume	150 µl

**Table 2. Preparation of Complete Nuclear Extraction Buffer**

## Purification of Cellular Nuclear Extracts

Cayman's Nuclear Extraction Kit (Item No. 10009277) can be used to isolate nuclear proteins. Alternatively, the procedure described below can be used for a 15 ml cell suspension grown in a T75 flask or adherent cells (100 mm dish 80-90% confluent) where  $10^7$  cells yields approximately 50  $\mu\text{g}$  of nuclear protein.

1. Collect  $10^7$  cells in pre-chilled 15 ml tubes.
2. Centrifuge suspended cells at  $300 \times g$  for five minutes at  $4^\circ\text{C}$ .
3. Discard the supernatant. Resuspend cell pellet in 5 ml of ice-cold 1X Nuclear Extraction PBS/Phosphatase Inhibitor Solution and centrifuge at  $300 \times g$  for five minutes at  $4^\circ\text{C}$ . Repeat one time.
4. Discard the supernatant. Add 500  $\mu\text{l}$  ice-cold 1X Complete Hypotonic Buffer. Mix gently by pipetting and transfer resuspended pellet to pre-chilled 1.5 ml microcentrifuge tube.
5. Incubate cells on ice for 15 minutes allowing cells to swell.
6. Add 100  $\mu\text{l}$  of 10% Nonidet P-40. Mix gently by pipetting.
7. Centrifuge for 30 seconds (pulse spin) at  $4^\circ\text{C}$  in a microcentrifuge. Transfer the supernatant which contains the cytosolic fraction to a new tube and store at  $-80^\circ\text{C}$ .
8. Resuspend the pellet in 100  $\mu\text{l}$  ice-cold Extraction Buffer (with protease and phosphatase inhibitors). Vortex 15 seconds at the highest setting and then gently rock the tube on ice for 15 minutes using a shaking platform. Vortex sample for 30 seconds at the highest setting and gently rock for an additional 15 minutes.
9. Centrifuge at  $14,000 \times g$  for 10 minutes at  $4^\circ\text{C}$ . The supernatant contains the nuclear fraction. Aliquot to clean chilled tubes, flash freeze, and store at  $-80^\circ\text{C}$ . Avoid freeze/thaw cycles. The extracts are ready to use in the assay.
10. Keep a small aliquot of the nuclear extract to quantitate the protein concentration.

## Reagent Preparation

### 1. Transcription Factor Antibody Binding Buffer (10X)

One vial (Item No. 10006882) contains 3 ml of a 10X stock of Transcription Factor Antibody Binding Buffer (ABB) to be used for diluting the primary and secondary antibodies. To prepare 1X ABB, dilute 1:10 by adding 27 ml of UltraPure water. Store at  $4^\circ\text{C}$  for up to two months.

### 2. Wash Buffer Concentrate (400X)

One vial (Item No. 400062) contains 5 ml of 400X Wash Buffer. Dilute the contents of the vial to a total volume of 2 liters with UltraPure water and add 1 ml of Polysorbate 20 (Item No. 400035). *NOTE: Polysorbate 20 is a viscous liquid and cannot be measured by a pipet. A positive displacement device such as a syringe should be used to deliver small quantities accurately.* A smaller volume of Wash Buffer Concentrate can be prepared by diluting the Wash Buffer Concentrate 1:400 and adding Polysorbate 20 (0.5 ml/liter of Wash Buffer). Store at  $4^\circ\text{C}$  for up to two months.



### 3. Transcription Factor Binding Assay Buffer (4X)

One vial (Item No. 10006880) contains 3 ml of a 4X stock of Transcription Factor Binding Assay Buffer (TFB). Prepare Complete Transcription Factor Binding Assay Buffer (CTFB) immediately prior to use in 1.5 ml centrifuge tubes or 15 ml conical tubes as outlined below in Table 3. This buffer is now referred to as CTFB. *NOTE: It is recommended that the CTFB be used the same day it is prepared.*

Component	Volume/ Well	Volume/ Strip	Volume/ 96-well plate
UltraPure water	73 $\mu$ l	584 $\mu$ l	7,008 $\mu$ l
Transcription Factor Binding Assay Buffer (4X)	25 $\mu$ l	200 $\mu$ l	2,400 $\mu$ l
Reagent A (Item No. 10007472)	1 $\mu$ l	8 $\mu$ l	96 $\mu$ l
300 mM DTT	1 $\mu$ l	8 $\mu$ l	96 $\mu$ l
Total Required	100 $\mu$ l	800 $\mu$ l	9,600 $\mu$ l

Table 3. Preparation of Complete Transcription Factor Binding Assay Buffer.

### 4. Transcription Factor CREB (Phospho-Ser<sup>133</sup>) Positive Control

One vial (Item No. 10009849) contains 150  $\mu$ l of unstimulated HeLa nuclear extract. This lysate is provided as a positive control for (Phospho-Ser<sup>133</sup>) CREB; it is not intended for plate to plate comparisons. The Positive Control provided is sufficient for 15 reactions and will provide a strong signal (>0.5 AU at 450 nm) when used at 10  $\mu$ l/well. When using this Positive Control, a decrease in signal may occur with repeated freeze/thaw cycles. It is recommended that the Positive Control be aliquoted at 25  $\mu$ l per vial and stored at -80°C to avoid loss in signal from repeated freeze/thaw cycles.

## Plate Set Up

There is no specific pattern for using the wells on the plate. A typical layout of CREB (Phospho-Ser<sup>133</sup>) positive control (PC), competitor dsDNA (C1), and samples of nuclear extracts (S1-S44) to be measured in duplicate is given below in Figure 2. We suggest you record the contents of each well on the template sheet provided (see page 30).

	1	2	3	4	5	6	7	8	9	10	11	12
A	S1	S1	S9	S9	S17	S17	S25	S25	S33	S33	S41	S41
B	S2	S2	S10	S10	S18	S18	S26	S26	S34	S34	S42	S42
C	S3	S3	S11	S11	S19	S19	S27	S27	S35	S35	S43	S43
D	S4	S4	S12	S12	S20	S20	S28	S28	S36	S36	S44	S44
E	S5	S5	S13	S13	S21	S21	S29	S29	S37	S37	NSB	NSB
F	S6	S6	S14	S14	S22	S22	S30	S30	S38	S38	PC	PC
G	S7	S7	S15	S15	S23	S23	S31	S31	S39	S39	Blk	Blk
H	S8	S8	S16	S16	S24	S24	S32	S32	S40	S40	C1	C1

S1-S44 - Sample Wells

NSB - Non-specific Binding Wells

PC - Positive Control Wells

Blk - Blank Wells

C1 - Competitor dsDNA Wells

Figure 2. Sample plate format

## Pipetting Hints

- Use different tips to pipette each reagent.
- Before pipetting each reagent, equilibrate the pipette tip in that reagent (*i.e.*, slowly fill the tip and gently expel the contents, repeat several times).
- Do not expose the pipette tip to the reagent(s) already in the well.

## General Information

- It is not necessary to use all the wells on the plate at one time; however a positive control should be run every time.
- For each plate or set of strips it is recommended that two Blk, two NSB, and two PC wells be included.

## Performing the Assay

### Binding of active CREB (Phospho-Ser<sup>133</sup>) to the consensus sequence:

1. Equilibrate the plate and buffers to room temperature prior to opening. Remove the plate from the foil and select the number of strips needed. The 96-well plate supplied with this kit is ready to use.

*NOTE: If you are not using all of the strips at once, place the unused strips back in the plate packet and store at 2-4°C. Be sure that the packet is sealed with the desiccant inside.*

2. Prepare the CTFB as outlined in Table 3, on page 16.
3. Add appropriate amount of reagent(s) listed below to the designated wells as follows:

**Blk** - add 100 µl of CTFB to designated wells.

**NSB** - add 100 µl of CTFB to designated wells. Do not add samples or positive control to these wells.

**C1** - Add 80 µl of CTFB prior to adding 10 µl of TF CREB (Phospho-Ser<sup>133</sup>) Competitor dsDNA (Item No. 10009488) to designated wells. Add 10 µl of control cell lysate or unknown sample.

*NOTE: Competitor dsDNA must be added prior to adding the Positive Control or nuclear extracts.*

**S1-S44** - Add 90 µl of CTFB followed by 10 µl of Nuclear Extract to designated wells. A protocol for isolation of nuclear extracts is given on page 14.

**PC** - Add 90 µl of CTFB followed by 10 µl of Positive Control to appropriate wells.

4. Use the cover provided to seal the plate. Incubate overnight at 4°C or one hour at room temperature without agitation (incubation for one hour will result in a less sensitive assay).
5. Empty the wells and wash five times with 200 µl of 1X Wash Buffer. After each wash empty the wells in the sink. After the final wash (wash #5), tap the plate on a paper towel to remove any residual Wash Buffer.

### Addition of TF CREB (Phospho-Ser<sup>133</sup>) Primary Antibody

1. Dilute the TF CREB (Phospho-Ser<sup>133</sup>) Primary Antibody (Item No. 10009848) 1:100 in 1X ABB as outlined in Table 4 below. Add 100 µl of diluted CREB (Phospho-Ser<sup>133</sup>) Primary Antibody to each well except the Blk wells.

Component	Volume/ Well	Volume/ Strip	Volume/ 96-well plate
1X ABB	99 µl	792 µl	9,504 µl
CREB (Phospho-Ser <sup>133</sup> ) Primary Antibody	1 µl	8 µl	96 µl
Total required	100 µl	800 µl	9,600 µl

**Table 4. Dilution of Primary Antibody**

2. Use the adhesive cover provided to seal the plate.
3. Incubate the plate for one hour at room temperature without agitation.
4. Empty the wells and wash each well five times with 200 µl of 1X Wash Buffer. After each wash, empty the contents of the plate into the sink. After the final wash (wash #5), tap the plate three to five times on a paper towel to remove any residual Wash Buffer.

### Addition of the Transcription Factor Goat Anti-Rabbit HRP Conjugate

1. Dilute the Transcription Factor Goat Anti-Rabbit HRP Conjugate (Item No. 10006884) 1:100 in 1X ABB as outlined in Table 5 below. Add 100  $\mu$ l of diluted Secondary Antibody to each well except the Blk wells.

Component	Volume/ Well	Volume/ Strip	Volume/ 96-well plate
1X ABB	99 $\mu$ l	792 $\mu$ l	9,504 $\mu$ l
Goat Anti-Rabbit HRP Conjugate	1 $\mu$ l	8 $\mu$ l	96 $\mu$ l
Total required	100 $\mu$ l	800 $\mu$ l	9,600 $\mu$ l

**Table 5. Dilution of Secondary Antibody**

2. Use the adhesive cover provided to seal the plate.
3. Incubate for one hour at room temperature without agitation.
4. Empty the wells and wash five times with 200  $\mu$ l of 1X Wash Buffer. After each wash, empty the contents of the plate into the sink. After the final wash (wash #5), tap the plate three to five times on a paper towel to remove any residual Wash Buffer.

### Develop and Read the Plate

1. To each well being used add 100  $\mu$ l of Transcription Factor Developing Solution (Item No. 10006888) which has been equilibrated to room temperature.
2. Incubate the plate for 15 to 45 minutes at room temperature with gentle agitation protected from light. Allow the wells to turn medium to dark blue prior to adding Transcription Factor Stop Solution (Item No. 10006889). (This reaction can be monitored by taking absorbance measurements at 655 nm prior to stopping the reactions; An OD<sub>655</sub> of 0.4-0.5 yields an OD<sub>450</sub> of approximately 1). Monitor development of sample wells to ensure adequate color development prior to stopping the reaction. *NOTE: Do not overdevelop; however PC wells may need to overdevelop to allow adequate color development in sample wells.*
3. Add 100  $\mu$ l of Stop Solution per well being used. The solution within the wells will change from blue to yellow after adding the Stop Solution.
4. Read absorbance at 450 nm within five minutes of adding the Stop Solution. Blank the plate reader according to the manufacturer's requirements using the blank wells.

## Assay Procedure Summary

*NOTE: This procedure is provided as a quick reference for experienced users. Follow the detailed procedure when initially performing the assay.*

1. Prepare CTFB as described in the Pre-Assay Preparation section, Table 3 on page 16.
2. Add 90  $\mu$ l CTFB per sample well (80  $\mu$ l if adding Competitor dsDNA), 100  $\mu$ l to Blk and NSB wells).
3. Add 10  $\mu$ l of Competitor dsDNA (optional) to appropriate wells.
4. Add 10  $\mu$ l of Positive Control to appropriate wells.
5. Add 10  $\mu$ l of sample containing activated CREB (Phospho-Ser<sup>133</sup>) to appropriate wells.
6. Incubate overnight at 4°C or one hour at room temperature without agitation.
7. Wash each well five times with 200  $\mu$ l of 1X Wash Buffer.
8. Add 100  $\mu$ l of diluted CREB (Phospho-Ser<sup>133</sup>) Primary Antibody per well (except Blk wells).
9. Incubate one hour at room temperature without agitation.
10. Wash each well five times with 200  $\mu$ l of 1X Wash Buffer.
11. Add 100  $\mu$ l of diluted Secondary Antibody (except Blk wells).
12. Incubate one hour at room temperature without agitation.
13. Wash each well five times with 200  $\mu$ l of 1X Wash Buffer.
14. Add 100  $\mu$ l of Developing Solution per well.
15. Incubate 15 to 45 minutes with gentle agitation.
16. Add 100  $\mu$ l of Stop Solution per well.
17. Measure the absorbance at 450 nm.

Steps	Reagent	Blk	NSB	PC	C1	S1-S44
1. Add reagents	CTFB	100 $\mu$ l	100 $\mu$ l	90 $\mu$ l	80 $\mu$ l	90 $\mu$ l
	Competitor dsDNA				10 $\mu$ l	
	Positive Control			10 $\mu$ l	10 $\mu$ l	
	Samples					10 $\mu$ l
2. Incubate	Cover plate and incubate overnight at 4°C or one hour at RT without agitation					
3. Wash	Wash all wells five times					
4. Add reagents	Primary Antibody		100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
5. Incubate	Cover plate and incubate one hour at RT without agitation					
6. Wash	Wash all wells five times					
7. Add reagents	Secondary Antibody		100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
8. Incubate	Cover plate and incubate one hour at RT without agitation					
9. Wash	Wash all wells five times					
10. Add reagents	Developer Solution	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
11. Incubate	Monitor development in wells					
12. Add reagents	Stop Solution	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l
13. Read	Read plate at wavelength of 450 nm					

**Table 6. Quick Protocol Guide**

## Performance Characteristics

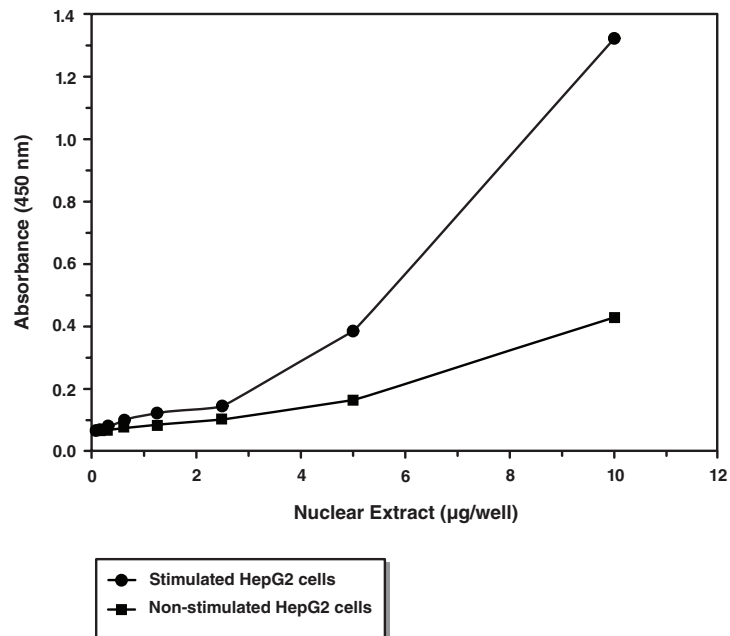


Figure 3. Assay of nuclear fractions isolated from stimulated (10  $\mu\text{M}$  Forskolin for 30 minutes) and nonstimulated HepG2 cells demonstrating activated CREB (Phospho-Ser<sup>133</sup>).

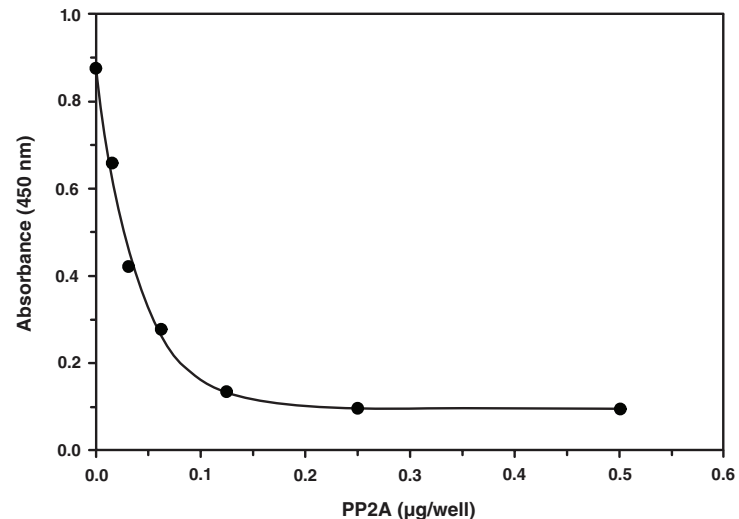


Figure 4. Assay of stimulated PC-12 nuclear extracts following treatment with PP2A. This experiment demonstrates the specificity of the CREB (Phospho-Ser<sup>133</sup>) Primary Antibody.

## RESOURCES

### Troubleshooting

Problem	Possible Causes	Recommended Solutions
No signal or weak signal in control wells	<ul style="list-style-type: none"> <li>A. Omission of key reagent</li> <li>B. Plate reader settings not correct</li> <li>C. Reagent/reagents expired</li> <li>D. Salt concentrations affected binding between DNA and protein</li> <li>E. Developing reagent used cold</li> <li>F. Developing reagent not added at correct volume</li> </ul>	<ul style="list-style-type: none"> <li>A. Check that all reagents have been added and in the correct order; perform the assay using the Positive Control</li> <li>B. Check wavelength setting on plate reader and change to 450 nm</li> <li>C. Check expiration date on reagents</li> <li>D. Reduce the amount of nuclear extract used in the assay, or reduce the amount of salt in the nuclear extracts (alternatively can perform buffer exchange)</li> <li>E. Prewarm the Developing Solution to room temperature prior to use</li> <li>F. Check pipettes to ensure correct amount of Developing Solution was added to wells</li> </ul>
High signal in all wells	<ul style="list-style-type: none"> <li>A. Incorrect dilution of antibody (too high)</li> <li>B. Improper/inadequate washing of wells</li> <li>C. Overdeveloping</li> </ul>	<ul style="list-style-type: none"> <li>A. Check antibody dilutions and use amounts outlined in instructions</li> <li>B. Follow the protocol for washing wells using the correct number of times and volumes</li> <li>C. Decrease the incubation time when using the developing reagent</li> </ul>
High background (NSB)	Incorrect dilution of antibody (too high)	Check antibody dilutions and use amounts outlined in the instructions

Problem cont.	Possible Causes cont.	Recommended Solutions cont.
Weak signal in sample wells	<ul style="list-style-type: none"> <li>A. Sample concentration is too low</li> <li>B. Incorrect dilution of antibody</li> <li>C. Salt concentrations affecting binding between DNA and protein</li> </ul>	<ul style="list-style-type: none"> <li>A. Increase the amount of nuclear extract used; loss of signal can occur with multiple freeze/thaw cycles of the sample; prepare fresh nuclear extracts and aliquot as outlined in product insert</li> <li>B. Check antibody dilutions and use amounts outlined in the instructions</li> <li>C. Reduce the amount of nuclear extract used in the assay or reduce the amount of salt in the nuclear extracts (alternatively can perform buffer exchange)</li> </ul>

### References

- Alberta, J.A., Augers, K.R., Batt, D., *et al.* Platelet-derived growth factor stimulation of monocyte chemoattractant protein-1 gene expression is mediated by transient activation of the phosphoinositide 3-kinase signal transduction pathway. *J. Biol. Chem.* **274**(43), 31,062-31,067 (1999).
- Heath, V.J., Gillespie, D.A.F., and Crouch, D.H. Inhibition of the terminal stages of adipocyte differentiation by cMyc. *Exp. Cell Res.* **254**, 91-98 (2000).
- Silva, A.J., Kogan, J.H., Frankland, P.W., *et al.* CREB and memory. *Annu. Rev. Neurosci.* **21**, 127-148 (1998).
- Johannessen, M., Delghandi, M.P., and Moens, U. What turns CREB on? *Cell. Signal.* **16**, 1211-1227 (2004).
- Pennisi, E. Opening the way to gene activity. *Science* **275**, 155-157 (1997).
- Shaywitz, A.J. and Greenberg, M.E. A stimulus-induced transcription factor activated by a diverse array of extracellular signals. *Annu. Rev. Biochem.* **68**, 821-861 (1999).

1								
2								
3								
4								
5								
6								
7								
8								
9								
10								
11								
12								
	A	B	C	D	E	F	G	H

NOTES

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.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©08/30/2016, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.



