

PRODUCT INFORMATION



TF CREB (Phospho-Ser¹³³) Primary Antibody

Item No. 10009848

Overview and Properties

Storage: -20°C (as supplied)
Stability: ≥2 years
Host: Rabbit
Applications: CREB (Phospho-Ser¹³³) Transcription Factor Assay Kit

Laboratory Procedures

This antibody has been specifically formulated to be used in the CREB (Phospho-Ser¹³³) Transcription Factor Assay Kit (Item No. 10009846). Prepare the primary antibody diluting 1:100 in 1X Antibody Binding Buffer (ABB) as outlined below. Add 100 µl of diluted CREB (Phospho-Ser¹³³) antibody to each well except the Blank (Blk) wells. Discard any unused diluted primary antibody.

| Component | Volume/Well | Volume/Strip | Volume/96-Well Plate |
|--|-------------|--------------|----------------------|
| UltraPure Water | 89 µl | 712 µl | 8,544 µl |
| Transcription Factor Antibody Binding Buffer (10X) | 10 µl | 80 µl | 960 µl |
| Transcription Factor CREB (Phospho-Ser ¹³³) Primary Antibody | 1 µl | 8 µl | 96 µl |
| Total Required | 100 µl | 800 µl | 9,600 µl |

CAUTIONS

For laboratory research use only. This reagent is not for human or veterinary use. This product has been tested and formulated to work exclusively with the CREB (Phospho-Ser¹³³) Transcription Factor Assay Kit from Cayman (Item No. 10009846). This product may not perform as described if used with other assay reagents or protocols.

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, 02/17/2022

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