

**SREBP-2 Cell-Based  
Translocation Assay Kit**

Item No. 10009239

## TABLE OF CONTENTS

<b>GENERAL INFORMATION</b>	3	Materials Supplied
	4	Precautions
	4	If You Have Problems
	4	Storage and Stability
	4	Materials Needed but Not Supplied
<b>INTRODUCTION</b>	5	Background
	5	About This Assay
<b>ASSAY PROTOCOL</b>	6	Treatment of the Cells
	7	Immunofluorescent Staining Procedure
<b>PERFORMANCE CHARACTERISTICS</b>	8	Cell Staining
<b>RESOURCES</b>	9	Troubleshooting
	10	References
	10	Related Products
	11	Warranty and Limitation of Remedy
	12	Notes

## GENERAL INFORMATION

### Materials Supplied

Kit will arrive packaged as a -20°C kit. For best results, remove components and store as stated below.

Item Number	Item	100 Tests Quantity	Storage
10009899	Cell-Based Assay Fixative	1 vial	Room Temperature
10009900	SREBP-2 Cell-Based Assay Primary Antibody	1 vial	-20°C
10009906	Cell-Based Assay Blocking Solution	1 vial	4°C
10009907	DyLight™ 549-Conjugated Goat Anti-Rabbit Secondary Antibody	1 vial	-20°C
10009869	Cell-Based Assay U-18666A	1 vial	-20°C

*NOTE: DyLight™ 549 is a product of Thermo Fisher Scientific.*

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 975-3999. We cannot accept any returns without prior authorization.



**WARNING: This product is for laboratory research use only: not for administration to humans. Not for human or veterinary diagnostic or therapeutic use.**

## Precautions

Please read these instructions carefully before beginning this assay.  
For research use only. Not for human or diagnostic use.

## 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:** techserve@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 and used before the expiration date indicated on the outside of the box.

## Materials Needed But Not Supplied

1. A 6-, 12-, 24-, or 96-well plate.
2. Raw 264.7 cells (can be obtained from ATCC) or any cells in which translocation of SREBP-2 is active.
3. Immunocytochemical staining buffer, TBS or PBS, pH 7.4.
4. Triton-X 100.
5. A fluorescence microscope equipped with a filter capable of measuring excitation and emission at 550 and 568 nm, respectively.

## INTRODUCTION

### Background

Lipid homeostasis in vertebrate cells is regulated by a family of basic helix-loop-helix (HLH) transcription factors called sterol regulatory element-binding proteins (SREBPs). SREBPs directly activate the expression of over 30 genes involved in both the synthesis and uptake of cholesterol, fatty acids, triglycerides, and phospholipids.<sup>1,2</sup> Each SREBP, including SREBP-1a, SREBP-1c, and SREBP-2, consists of three domains: an amino-terminal transcription factor domain of ~480 amino acids, a middle hydrophobic region of ~80 amino acids containing two hydrophobic transmembrane segments, and a carboxy-terminal regulatory domain of ~590 amino acids.<sup>3</sup> SREBP-2 regulates cholesterol synthesis by activating the transcription of genes for HMG-CoA reductase and other enzymes of the cholesterol synthetic pathway.<sup>4</sup> It is also involved in activating genes required to generate NADPH, which is consumed at multiple stages in cholesterol biosynthesis. SREBP-2 is ubiquitously expressed.<sup>5</sup> Under basal conditions SREBP is bound to endoplasmic reticulum membranes as a glycosylated precursor protein and SREBP cleavage-activating protein (SCAP) is bound to sterol. Upon depletion of cholesterol, SCAP becomes activated and escorts SREBP to the Golgi where it is proteolytically cleaved by site 1 protease and site 2 protease, S1P and S2P, respectively.<sup>4,6</sup> The active transcription factor consisting of the NH<sub>2</sub>-terminal domain, designated as nuclear SREBP (nSREBP), is translocated into the nucleus to stimulate transcription of genes involved in the uptake and synthesis of cholesterol. This protein activation characteristic makes it possible to study modulators of SREBP-2 through sub-cellular localization of the protein using immunocytochemical staining with a specific antibody.<sup>7</sup>

Reduction of circulating cholesterol and modulation of lipid biosynthesis have important clinical implications for many diseases including obesity, type 2 diabetes, and atherosclerosis. Thus, identification of SREBP-2 activators/inhibitors can be highly relevant in the search for cures to these diseases.

### About This Assay

Cayman's SREBP-2 Cell-Based Translocation Assay Kit provides highly specific SREBP-2 primary antibody together with a DyLight™ (trademarked by Thermo Fisher Scientific.) conjugated secondary antibody in a ready to use format. Also included as a positive control is a cholesterol trafficking inhibitor, U-18666A, which has been shown to activate SREBP-2 translocation into nuclei by scientists at Cayman.

## Treatment of the Cells

The following protocol is designed for a 96-well plate. Adjust volumes accordingly for other plate sizes.

1. Seed wells of a 96-well plate with  $3 \times 10^4$  cells/well. Grow cells overnight.
2. The next day, treat cells with or without experimental compounds for 48-72 hours, or for the period of time used in your typical experimental protocol. Cell-Based Assay U-18666A (Item No. 10009869), a cholesterol transport inhibitor, is included as a positive control.
3. After the designated incubation period, examine the effect of testing compounds on SREBP-2 activation by the following immunofluorescent staining procedure in section B.

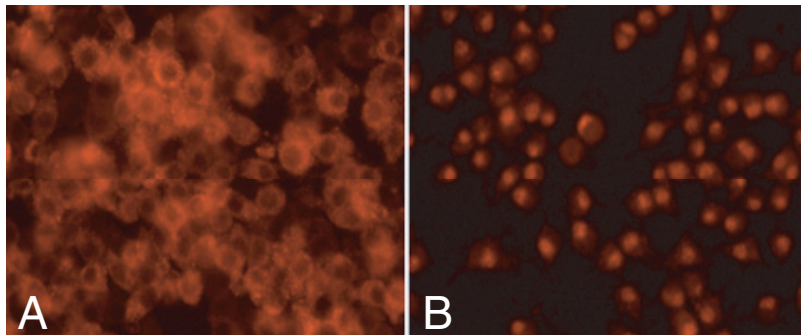
## Immunofluorescent staining procedure

**NOTE: Perform all steps at room temperature.**

The reagent volumes provided below are for use on a 96-well plate. Adjust volumes accordingly for other plate sizes.

1. Remove most of the culture medium from the wells.
2. Wash cells briefly with TBS, pH 7.4.
3. Fix the cells with Cell-Based Assay Fixative Solution (Item No. 10009899) for 10 minutes.
4. Wash the cells three times with TBS containing 0.1% Triton-X 100 (TBST) for five minutes each.
5. Incubate the cells with Cell-Based Assay Blocking Solution (Item No. 10009906) for 30 minutes.
6. Incubate the cells for one hour with 100  $\mu$ l of SREBP-2 Cell-Based Assay Primary Antibody (Item No. 10009900) diluted 1:100 in TBST.
7. Wash the cells three times with TBST for five minutes each.
8. Incubate the cells in the dark for one hour with 100  $\mu$ l of DyLight™ 549-Conjugated Goat Anti-Rabbit Secondary Antibody (Item No. 10009907) diluted 1:100 in TBST.
9. Wash the cells three times with TBST for five minutes each.
10. Examine the staining under a fluorescent microscope with a filter capable of measuring excitation and emission at 550 and 568 nm, respectively. Store the plate at 4°C in the dark for later analysis.

Cell Staining



**Figure 1: Translocation of SREBP-2 into nuclei by U-18666A.** Raw 264.7 cells were seeded in a 96-well plate at a density of  $3 \times 10^4$  cells/well and cultured overnight. The next day, cells were treated with DMSO (vehicle) or  $24 \mu\text{M}$  U-18666A for 72 hours. *Panel A:* Cells treated with DMSO alone demonstrate cytoplasmic localization of SREBP-2, indicating that most of cells have inactive protein. *Panel B:* U-18666A treatment for three days induced SREBP-2 translocation into the nuclei, indicating that blockage of cholesterol transport in these cells activates the protein.

Troubleshooting

Problem	Possible Causes	Recommended Solutions
No signal in any wells	Omission of key reagent	Check that all reagents have been added and were added in the correct order
High signal in all wells	Overgrowth of cells	Make sure to plate cells at low density before starting treatment

## References

1. Osborne, T.F. Sterol regulatory element-binding proteins (SREBP's): Key regulators of nutritional homeostasis and insulin action. *J. Biol. Chem.* **275**(42), 32379-32382 (2000).
2. Lin, J., Yang, R., Tarr, P.T., *et al.* Hyperlipidemic effects of dietary saturated fats mediated through PGC-1 $\beta$  coactivation of SREBP. *Cell* **120**, 261-273 (2005).
3. Brown, M.S. and Goldstein, J.L. The SREBP pathway: Regulation of cholesterol metabolism by proteolysis of a membrane-bound transcription factor. *Cell* **89**, 331-340 (1997).
4. Sakai, J., Nohturfft, A., Goldstein, J.L., *et al.* Cleavage of sterol regulatory element-binding proteins (SREBPs) at site-1 requires interaction with SREBP cleavage-activating protein. Evidence from *in vivo* competition studies. *J. Biol. Chem.* **273**(10), 5785-5793 (1998).
5. Iizuka, K., Bruick, R.K., Liang, G., *et al.* Deficiency of carbohydrate response element-binding protein (ChREBP) reduces lipogenesis as well as glycolysis. *Proc. Natl. Acad. Sci. USA* **101**(19), 7281-7286 (2004).
6. Matsuda, M., Korn, B.S., Hammer, R.E., *et al.* SREBP cleavage-activating protein (SCAP) is required for increased lipid synthesis in liver induced by cholesterol deprivation and insulin elevation. *Genes & Development* **15**, 1206-1216 (2001).
7. Smith, L.H., Petrie, M.S., Morrow, J.D., *et al.* The sterol response element binding protein regulates cyclooxygenase-2 gene expression in endothelial cells. *J. Lipid Res.* **46**, 862-871 (2005).

## Related Products

Adipogenesis Assay Kit - Item No. 10006908  
ChREBP Cell-Based Translocation Assay Kit - Item No. 10010060  
Glucose Assay Kit - Item No. 10009582  
LDL Receptor Polyclonal Antibody - Item No. 10007665  
p38 MAPK (Phospho-Thr<sup>180</sup>/Tyr<sup>182</sup>) Phosphorylation/Translocation  
Cell-Based Assay Kit - Item No. 10010374  
PCSK9 (human) Polyclonal Antibody - Item No. 10007185  
PCSK9 (murine) Polyclonal Antibody - Item No. 10008811  
PCSK9 Western Ready Control - Item No. 10009567  
SREBP-2 Polyclonal Antibody - Item No. 10007663  
SREBP-2 Transcription Factor Assay Kit - Item No. 10007819  
SREBP-2 Western Ready Control - Item No. 10009749

## Warranty and Limitation of Remedy

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery. Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence. This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

**For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.**

## NOTES

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.

©06/22/2011, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.