

PRODUCT INFORMATION

Transcription Factor SREBP 96-Well Strip Plate

Item No. 10008858



Overview and Properties

Contents:	A flat-bottom 96-well plate coated with dsDNA consensus sequence specific for SREBP.
Storage:	-4°C
Stability:	≥3 years
Application:	This plate is designed for use in Cayman's SREBP Transcription Factor Assay Kits (Item Nos. 10007819 and 10010854). The plate is ready to use as supplied.

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 SREBP Transcription Factor Assay Kits (Item Nos. 10007819 and 10010854). 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, 10/20/2021

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