

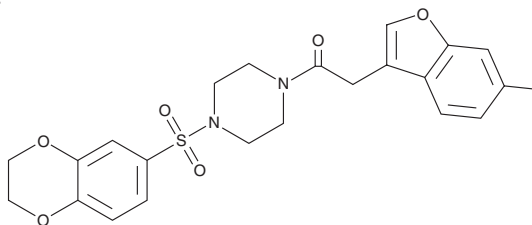
PRODUCT INFORMATION



SR-4559

Item No. 18654

CAS Registry No.: 847939-23-5
Formal Name: 1-[4-[(2,3-dihydro-1,4-benzodioxin-6-yl)sulfonyl]-1-piperazinyl]-2-(6-methyl-3-benzofuranyl)-ethanone
Synonym: Perilipin 1/5 Inhibitor 1
MF: C₂₃H₂₄N₂O₆S
FW: 456.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SR-4559 is supplied as a solid. A stock solution may be made by dissolving the SR-4559 in the solvent of choice, which should be purged with an inert gas. SR-4559 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SR-4559 is an inhibitor of the protein-protein interaction between α/β -hydrolase domain-containing protein 5 (ABHD5) and perilipin 1 (PLIN1; IC₅₀ = 510 nM).¹ It induces fatty acid release in primary mouse brown adipocytes (EC₅₀ = ~4 μ M).

Reference

1. Sanders, M.A., Madoux, F., Mladenovic, L., *et al.* Endogenous and synthetic ABHD5 ligands regulate ABHD5-perilipin interactions and lipolysis in fat and muscle. *Cell Metab.* **22(5)**, 851-860 (2015).

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, 09/17/2025

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