

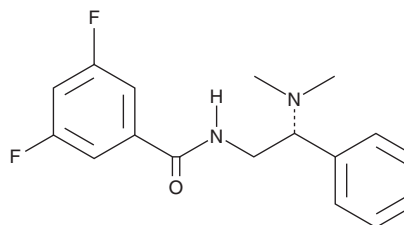
PRODUCT INFORMATION



(R)-HTS-3

Item No. 37563

CAS Registry No.: 2771208-83-2
Formal Name: N-[(2R)-2-(dimethylamino)-2-phenylethyl]-3,5-difluoro-benzamide
MF: C₁₇H₁₈F₂N₂O
FW: 304.3
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

(R)-HTS-3 is an inhibitor of lysophosphatidylcholine acyltransferase 3 (LPCAT3; IC₅₀ = 0.09 μM), an enzyme involved in the synthesis of C20:4-containing phospholipids.¹ It reduces the incorporation of arachidonic acid-d₈ (Item No. 390010) into phospholipids when used at concentrations of 1, 3, and 10 μM. (R)-HTS-3 (10 μM) reduces RSL3-induced ferroptosis in HT-1080 and 786-O cancer cells.

Reference

1. Reed, A., Ichu, T.-A., Milosevich, N., *et al.* LPCAT3 inhibitors remodel the polyunsaturated phospholipid content of human cells and protect from ferroptosis. *ACS Chem. Biol.* **17**(6), 1607-1618 (2022).

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, 10/12/2022

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