

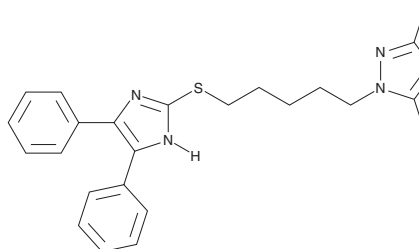
PRODUCT INFORMATION



RP 70676

Item No. 39529

CAS Registry No.: 136609-26-2
Formal Name: 1-[5-[(4,5-diphenyl-1H-imidazol-2-yl)thio]pentyl]-3,5-dimethyl-1H-pyrazole
MF: C₂₅H₂₈N₄S
FW: 416.6
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

RP 70676 is supplied as a solid. A stock solution may be made by dissolving the RP 70676 in the solvent of choice, which should be purged with an inert gas. RP 70676 is soluble in acetonitrile and DMSO.

Description

RP 70676 is an inhibitor of acyl-coenzyme A:cholesterol acyltransferase (ACAT; IC₅₀ = 44 nM for the rabbit enzyme).¹ It reduces plasma and arterial cholesterol levels in rabbits when administered at a dose of 1 mg/kg.

Reference

1. Ashton, M.J., Bridge, A.W., Bush, R.C., *et al.* RP 70676: A potent systemically available inhibitor of acyl-CoA:cholesterol O-acyl transferase (ACAT). *Bioorg. Med. Chem. Lett.* **2(5)**, 375-380 (1992).

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, 11/15/2023

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