

PRODUCT INFORMATION



GSK1940029

Item No. 36704

CAS Registry No.: 1150701-66-8
Formal Name: 1-[(3,4-dichlorophenyl)methyl]-N-[4-(hydroxymethyl)phenyl]-5-methyl-1H-1,2,3-triazole-4-carboxamide

MF: C₁₈H₁₆Cl₂N₄O₂

FW: 391.3

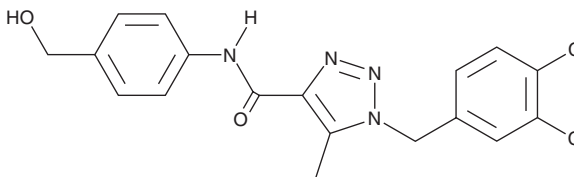
Purity: ≥98%

UV/Vis.: λ_{max}: 268 nm

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

GSK1940029 is supplied as a solid. A stock solution may be made by dissolving the GSK1940029 in the solvent of choice, which should be purged with an inert gas. GSK1940029 is soluble in DMSO.

Description

GSK1940029 is a stearyl-CoA desaturase (SCD) inhibitor.¹ It increases palmitate-induced phosphorylation of p70 ribosomal S6 kinase 1 (p70S6K1) and lactate dehydrogenase (LDH) release from AML12 hepatocytes when used at a concentration of 20 μM.² GSK1940029 (10 μM) also increases palmitate-induced increases in extracellular and intracellular triglyceride levels in AML12 hepatocytes, an effect that can be reversed by the mammalian target of rapamycin (mTOR) inhibitor torin 1 (Item No. 10997).

References

1. Bouillot, A.M.J., Laroze, A., and Trottet, L. 1,2,3-triazole derivatives for use as stearyl-coa desaturase inhibitors. *GlaxoSmithKline LLC WO 2009/060053* (2009).
2. Wang, J., Chen, Y., Song, Q., *et al.* mTORC1-IRE1α pathway activation contributes to palmitate-elicited triglyceride secretion and cell death in hepatocytes. *Exp. Biol. Med. (Maywood)* **245(14)**, 1268-1279 (2020).

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, 06/19/2023

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