

PRODUCT INFORMATION

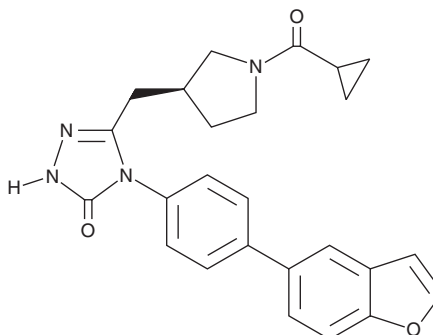


GSK2194069

Item No. 20022

CAS Registry No.: 1332331-08-4
Formal Name: 4-[4-(5-benzofuranyl)phenyl]-5-[[[(3S)-1-(cyclopropylcarbonyl)-3-pyrrolidinyl]methyl]-2,4-dihydro-3H-1,2,4-triazol-3-one

MF: C₂₅H₂₄N₄O₃
FW: 428.5
Purity: ≥98%
UV/Vis.: λ_{max}: 239 nm
Supplied as: A crystalline 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

GSK2194069 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK2194069 in the solvent of choice, which should be purged with an inert gas. GSK2194069 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GSK2194069 in these solvents is approximately 1.25, 20, and 30 mg/ml, respectively.

Description

GSK2194069 is an inhibitor of fatty acid synthase (FASN) with an IC₅₀ value of 7.7 nM in an assay detecting released CoA.¹ It is a competitive inhibitor that is selective for the KR domain over the KS domain of FASN, with IC₅₀ values of 29 and >10,000 nM, respectively, *in vitro*. GSK2194069 decreases incorporation of ¹³C-labeled acetate into cellular lipids in a dose-dependent manner in KATO-III, MKN-45, and SNU-1 gastric, and A549 non-small cell lung cancer cell lines. It also reduces cell growth of A549 cells (EC₅₀ = 15 nM) and inhibits lipid synthesis without decreasing FASN protein expression.

Reference

1. Hardwicke, M.A., Rendina, A.R., Williams, S.P., *et al.* A human fatty acid synthase inhibitor binds β-ketoacyl reductase in the keto-substrate site. *Nat. Chem. Biol.* **10**(9), 774-779 (2014).

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, 11/04/2022

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