

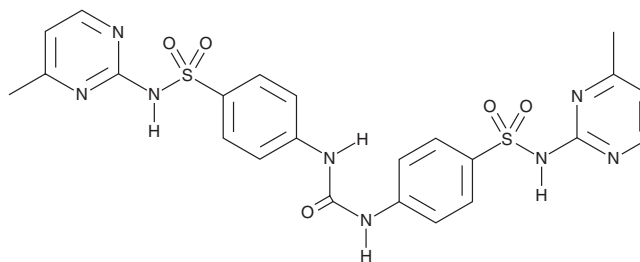
PRODUCT INFORMATION



GSK837149A

Item No. 19098

CAS Registry No.: 13616-29-0
Formal Name: 4,4'-(carbonyldiimino)bis[N-(4-methyl-2-pyrimidinyl)-benzenesulfonamide]
MF: C₂₃H₂₂N₈O₅S₂
FW: 554.6
Purity: ≥95%
UV/Vis.: λ_{max}: 288 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

GSK837149A is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK837149A in the solvent of choice. GSK837149A is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of GSK837149A in these solvents is approximately 5 mg/ml.

GSK837149A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK837149A should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GSK837149A has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A number of tumor cells and cell lines have been observed to have highly upregulated expression and activity of fatty acid synthase (FASN). GSK837149A is a selective, reversible inhibitor of FASN (IC₅₀ = 15.8 nM).¹ It acts by inhibiting the β-ketoacyl reductase activity of FASN.

Reference

1. Vázquez, M.J., Leavens, W., Liu, R., *et al.* Discovery of GSK837149A, an inhibitor of human fatty acid synthase targeting the β-ketoacyl reductase reaction. *FEBS J.* **275(7)**, 1556-1567 (2008).

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, 12/18/2023

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