

PRODUCT INFORMATION



GSK3326595

Item No. 27305

CAS Registry No.: 1616392-22-3

Formal Name: 6-[(1-acetyl-4-piperidinyl)amino]-
N-[(2S)-3-(3,4-dihydro-2(1H)-
isoquinolinyl)-2-hydroxypropyl]-4-
pyrimidinecarboxamide

MF: C₂₄H₃₂N₆O₃

FW: 452.6

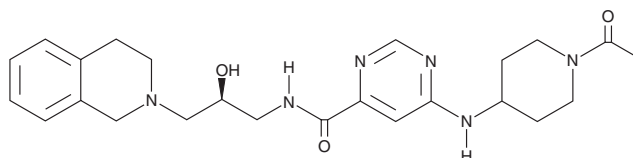
Purity: ≥98%

UV/Vis.: λ_{max}: 249 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

GSK3326595 is supplied as a solid. A stock solution may be made by dissolving the GSK3326595 in the solvent of choice, which should be purged with an inert gas. GSK3326595 is soluble in DMSO.

Description

GSK3326595 is an inhibitor of protein arginine methyltransferase 5 (PRMT5).¹ It inhibits methylation of peptides derived from the PRMT5 substrates histone H4 and H2A, SmD3, FUBP1, and HNRNPH1 by the PRMT5/MEP50 complex with IC₅₀ values ranging from 5.9 to 19.7 nM and is more than 4,000-fold selective for PRMT5/MEP50 over a panel of 20 methyltransferases. GSK3326595 (200 nM) induces alternative splicing of MDM4 in Z-138, MCF-7, JM1, and DOHH-2 cells and increases p53 and p21 protein levels in Z-138 cells in a concentration-dependent manner. GSK3326595 inhibits growth and symmetric dimethyl arginine (SDMA) methylation in several lymphoma and breast cancer cell lines. It also reduces tumor growth in a Z-138 mouse xenograft model when administered at doses of 25, 50, and 100 mg/kg twice per day.

Reference

1. Gerhart, S.V., Kellner, W.A., Thompson, C., *et al.* Activation of the p53-MDM4 regulatory axis defines the anti-tumour response to PRMT5 inhibition through its role in regulating cellular splicing. *Sci. Rep.* **8**(1), 9711 (2018).

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, 12/20/2022

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