

PRODUCT INFORMATION

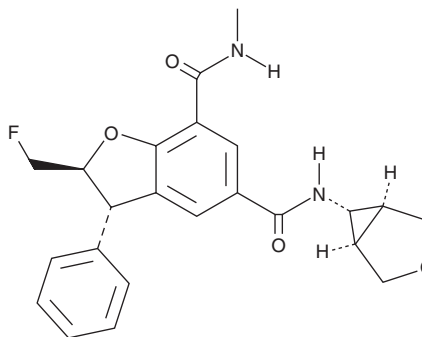


GSK973

Item No. 42057

CAS Registry No.: 2138473-38-6
Formal Name: (2S,3S)-2-(fluoromethyl)-2,3-dihydro-N⁷-methyl-N⁵-(1 α ,5 β ,6 β)-3-oxabicyclo[3.1.0]hex-6-yl-3-phenyl-5,7-benzofurandicarboxamide

MF: C₂₃H₂₃FN₂O₄
FW: 410.4
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

GSK973 is supplied as a solid. A stock solution may be made by dissolving the GSK973 in the solvent of choice, which should be purged with an inert gas. GSK973 is sparingly soluble (1-10 mg/ml) in methanol and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

GSK620 is a bromodomain and extra terminal domain family of proteins (BET) bromodomain 2 (BD2) inhibitor (IC₅₀s = 31.62, 15.85, 15.85, and 39.81 nM for BRD2, BRD3, BRD4, and BRDT_{BD2}, respectively).¹ It is selective for BD2 in bromodomain-containing protein 2 (BRD2), BRD3, BRD4, and BRDT over BD1 in the same BRDs (IC₅₀s = >10,000 nM for all).

Reference

1. Preston, A., Atkinson, S.J., Bamborough, P., *et al.* GSK973 Is an inhibitor of the second bromodomains (BD2s) of the bromodomain and extra-terminal (BET) family. *ACS Med. Chem. Lett.* **11**(8), 1581-1587 (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, 11/11/2024

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