

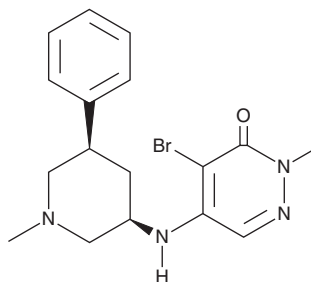
PRODUCT INFORMATION



GSK4027

Item No. 23421

CAS Registry No.: 2079896-25-4
Formal Name: 4-bromo-2-methyl-5-[[[(3R,5R)-1-methyl-5-phenyl-3-piperidyl]amino]-3(2H)-pyridazinone
MF: C₁₇H₂₁BrN₄O
FW: 377.3
Purity: ≥95%
UV/Vis.: λ_{max}: 228, 292 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

GSK4027 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK4027 in the solvent of choice, which should be purged with an inert gas. GSK4027 is soluble in DMSO.

Description

GSK4027 is a p300/CBP-associated factor (PCAF)/general control nonderepressible 5 (GCN5) bromodomain inhibitor ($K_i = 1.4$ nM; $IC_{50} = 60$ nM in a chromatin engagement assay).¹ It is greater than 18,000-fold selective for PCAF and GCN5 over BET family bromodomains and greater than 70-fold selective over other bromodomain families in a BROMOscan assay.

Reference

1. Humphreys, P.G., Bamborough, P., Chung, C.W., *et al.* Discovery of a potent, cell penetrant, and selective p300/CBP-associated factor (PCAF)/general control nonderepressible 5 (GCN5) bromodomain chemical probe. *J. Med. Chem.* **60**(2), 695-709 (2017).

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, 11/16/2022

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