

PRODUCT INFORMATION



BRD0639

Item No. 35625

CAS Registry No.: 2760881-74-9

Formal Name: (αS)-4-chloro-α-methyl-N-[4-methyl-3-[[[2-(2-pyridinyl)ethyl]amino]sulfonyl]phenyl]-6-oxo-1(6H)-pyridazineacetamide

MF: C₂₁H₂₂ClN₅O₄S

FW: 475.9

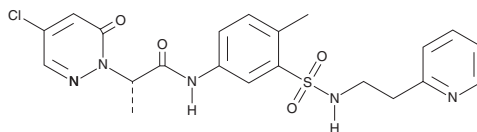
Purity: ≥98%

UV/Vis.: λ_{max}: 211, 252 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

BRD0639 is supplied as a solid. A stock solution may be made by dissolving the BRD0639 in the solvent of choice, which should be purged with an inert gas. BRD0639 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BRD0639 in these solvents is approximately 5 mg/ml.

Description

BRD0639 is an inhibitor of the protein-peptide interaction between protein arginine methyltransferase 5 (PRMT5) and PRMT5 binding motif (PBM; IC₅₀ = 13.8 μM), a conserved peptide motif contained in PRMT5 substrate adaptor proteins (SAPs) that mediates SAP binding to PRMT5.¹ It induces the formation of PRMT5 adducts in cells (EC₅₀ = 3 μM) and disrupts complex formation between PRMT5 and the SAP RIO kinase 1 (RIOK1) in permeabilized cells.

Reference

1. McKinney, D.C., McMillan, B.J., Ranaghan, M.J., *et al.* Discovery of a first-in-class inhibitor of the PRMT5–substrate adaptor interaction. *J. Med. Chem.* **64**(15), 11148-11168 (2022).

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, 03/29/2024

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