

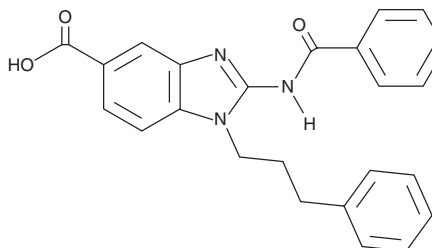
PRODUCT INFORMATION



BRD9539

Item No. 11788

CAS Registry No.: 1374601-41-8
Formal Name: 2-(benzoylamino)-1-(3-phenylpropyl)-1H-benzimidazole-5-carboxylic acid
MF: C₂₄H₂₁N₃O₃
FW: 399.4
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 230, 270, 329 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

BRD9539 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD9539 in the solvent of choice. BRD9539 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of BRD9539 in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

BRD9539 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD9539 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BRD9539 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD9539 is an inhibitor of euchromatin histone methyltransferase 2 (EHMT2), also known as G9a, with an IC₅₀ value of 6.3 μM.¹ It inhibits polycomb repressive complex 2 (PRC2) to a similar extent with 54 and 43% activity remaining for G9a and PRC2, respectively, when used at a concentration of 10 μM. It is selective for G9a and PRC2 over SU39H1 and NDMT1 up to a concentration of 40 μM. It is more potent than BRD4770 (Item No. 11787) in enzyme assays but has no activity in cell-based assays when used at concentrations of 5 and 10 μM.

Reference

1. Yuan, Y., Wang, Q., Paulk, J., *et al.* A small-molecule probe of the histone methyltransferase G9a induces cellular senescence in pancreatic adenocarcinoma. *ACS Chem. Biol.* **7**(7), 1152-1157 (2012).

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, 12/05/2022

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