

PRODUCT INFORMATION

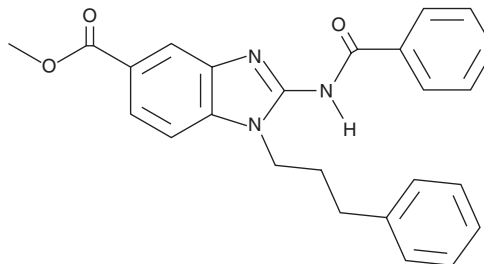


BRD4770

Item No. 11787

CAS Registry No.: 1374601-40-7
Formal Name: 2-(benzoylamino)-1-(3-phenylpropyl)-1H-benzimidazole-5-carboxylic acid, methyl ester

Synonym: HMTase Inhibitor VI
MF: C₂₅H₂₃N₃O₃
FW: 413.5
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 271, 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

BRD4770 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD4770 in the solvent of choice, which should be purged with an inert gas. BRD4770 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BRD4770 in these solvents is approximately 10 and 20 mg/ml, respectively.

BRD4770 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD4770 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD4770 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD4770 is an inhibitor of euchromatin histone methyltransferase 2 (EHMT2), also known as G9a, decreasing di- and trimethylation on lysine 9 of histone 3 (EC₅₀ = 5 μM).¹ Through its effects on EHMT2, BRD4770 induces senescence in the pancreatic cancer cell line PANC-1 without initiating apoptosis.¹ It also blocks both anchorage-dependent and -independent proliferation of PANC-1 cells.¹

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, 11/14/2022

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