

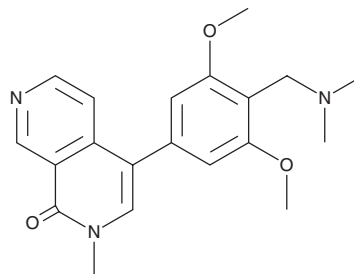
PRODUCT INFORMATION



BI-7273

Item No. 20311

CAS Registry No.: 1883429-21-7
Formal Name: 4-[4-[(dimethylamino)methyl]-3,5-dimethoxyphenyl]-2-methyl-2,7-naphthyridin-1(2H)-one
MF: C₂₀H₂₃N₃O₃
FW: 353.4
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 321 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

BI-7273 is supplied as a crystalline solid. A stock solution may be made by dissolving the BI-7273 in the solvent of choice. BI-7273 is soluble in organic solvents such as ethanol and DMSO. The solubility of BI-7273 in these solvents is approximately 0.2 and 0.1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of BI-7273 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of BI-7273 in PBS, pH 7.2, is approximately 0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BI-7273 is a potent BRD9 bromodomain inhibitor ($K_d = 15.4$ nM; $IC_{50} = 19$ nM) that less effectively inhibits the functionally and structurally related BRD7 bromodomain ($IC_{50} = 117$ nM).¹ It is without effect against the bromodomains of BRD2 and BRD4, as well as a panel of kinases. BI-7273 inhibits the growth of EOL-1 acute myeloid leukemia cells *in vitro* ($EC_{50} = 1.4$ μM).¹

Reference

1. Martin, L. J., Koegl, M., Bader, G., *et al.* Structure-based design of an *in vivo* active selective BRD9 inhibitor. *J. Med. Chem.* (2016).

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, 12/01/2022

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