

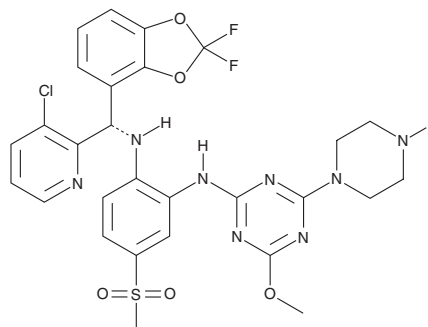
PRODUCT INFORMATION



DOT1L-IN-4

Item No. 36456

CAS Registry No.: 2565705-02-2
Formal Name: N¹-[(S)-(3-chloro-2-pyridinyl)(2,2-difluoro-1,3-benzodioxol-4-yl)methyl]-N²-[4-methoxy-6-(1-piperazinyl)-1,3,5-triazin-2-yl]-4-(methylsulfonyl)-1,2-benzenediamine
Synonym: DOT1-like Inhibitor 4
MF: C₂₈H₂₇ClF₂N₈O₅S
FW: 661.1
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 276 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

DOT1L-IN-4 is supplied as a solid. A stock solution may be made by dissolving the DOT1L-IN-4 in the solvent of choice, which should be purged with an inert gas. DOT1L-IN-4 is soluble in the organic solvent DMSO.

Description

DOT1L-IN-4 is an inhibitor of the non-SET domain-containing methyltransferase DOT1L (IC₅₀ = 0.11 nM).¹ It is selective for DOT1L over a panel of 11 additional histone methyltransferases at 5 μM.

Reference

1. Stauffer, F., Weiss, A., Scheufler, C., *et al.* New potent DOT1L inhibitors for *in vivo* evaluation in mouse. *ACS Med. Chem. Lett.* **10**(12), 1655-1660 (2019).

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, 12/19/2022

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