

PRODUCT INFORMATION

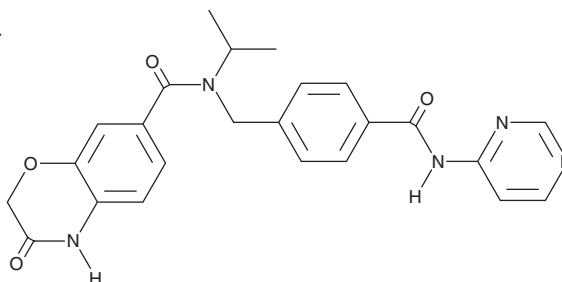


UNC7145

Item No. 40166

CAS Registry No.: 2561494-78-6
Formal Name: 3,4-dihydro-N-(1-methylethyl)-3-oxo-N-[[4-[[4-(pyrimidinylamino)carbonyl]phenyl)methyl]-2H-1,4-benzoxazine-7-carboxamide

MF: C₂₄H₂₃N₅O₄
FW: 445.5
Purity: ≥98%
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

UNC7145 is supplied as a solid. A stock solution may be made by dissolving the UNC7145 in the solvent of choice, which should be purged with an inert gas. UNC7145 is soluble in acetonitrile and DMSO.

Description

UNC7145 is an inactive derivative of the histone-lysine N-methyltransferase NSD2 inhibitor UNC6934 that has been used as a negative control for the histone methyltransferase activity of NSD2.¹

Reference

1. Dilworth, D., Hanley, R.P., Ferreira de Freitas, R., et al. A chemical probe targeting the PWWP domain alters NSD2 nucleolar localization. *Nat. Chem. Biol.* **18(1)**, 56-63 (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, 12/14/2023

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