

PRODUCT INFORMATION

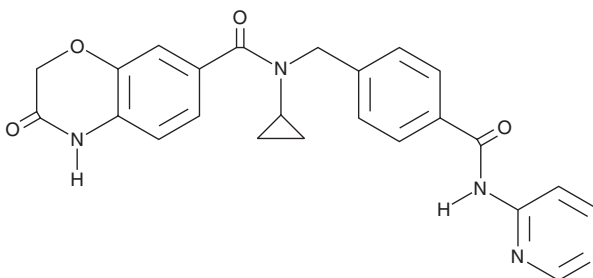


UNC6934

Item No. 28814

CAS Registry No.: 2561494-77-5
Formal Name: N-cyclopropyl-3,4-dihydro-3-oxo-N-[[4-[[4-(pyrimidinylamino)carbonyl]phenyl)methyl]-2H-1,4-benzoxazine-7-carboxamide

MF: C₂₄H₂₁N₅O₄
FW: 443.5
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

UNC6934 is supplied as a solid. A stock solution may be made by dissolving the UNC6934 in the solvent of choice, which should be purged with an inert gas. UNC6934 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

UNC6934 is an inhibitor of the protein-protein interaction between histone-lysine N-methyltransferase NSD2 and dimethylated lysine 36 histone 3 (H3K36me2) nucleosomes (IC₅₀ = 104 nM).¹ It binds to NSD2 (K_d = 92 nM). UNC6934 also inhibits the serotonin transporter (SERT) and μ-opioid receptor (K_{i,s} = 1.4 and 8 μM, respectively). It induces NSD2 nucleolus accumulation in U2OS cells when used at a concentration of 5 μM. See the Structural Genomics Consortium (SGC) website for more information.

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, 08/05/2024

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