

PRODUCT INFORMATION



NV03

Item No. 45383

CAS Registry No.: 2448341-58-8
Formal Name: N-[3-(diethylamino)propyl]-5,9-dimethyl-8-oxo-thieno[2',3':4,5]pyrrolo[1,2-d][1,2,4]triazine-7(8H)-acetamide

MF: C₁₉H₂₇N₅O₂S

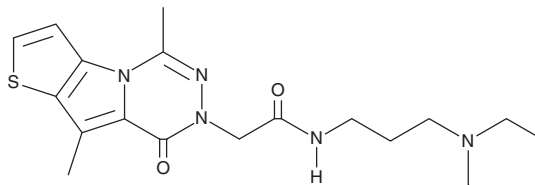
FW: 389.5

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

NV03 is supplied as a solid. A stock solution may be made by dissolving the NV03 in the solvent of choice, which should be purged with an inert gas. NV03 is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

Description

NV03 is an inhibitor of the protein-protein interaction between E3 ubiquitin-protein ligase UHRF1 (UHRF1) and histone H3 trimethylated lysine 9 (H3K9Me3).¹ It binds the tandem Tudor domain (TTD) of UHRF1 ($K_d = 2.4 \mu\text{M}$) and is selective for the UHRF1-TTD over a panel of 36 demethylases at 10 μM .

Reference

1. Senisterran, G., Zhu, H.Y., Luo, X., *et al.* Discovery of small-molecule antagonists of the H3K9me3 binding to UHRF1 tandem tudor domain. *SLAS Discov.* **23(9)**, 930-940 (2018).

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, 06/10/2026

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