

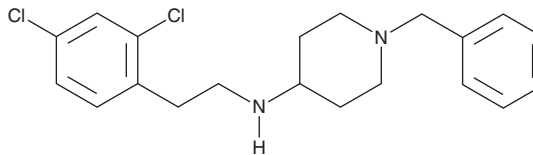
PRODUCT INFORMATION



NEDD8 Inhibitor M22

Item No. 37756

CAS Registry No.: 864420-54-2
Formal Name: N-[2-(2,4-dichlorophenyl)ethyl]-1-(phenylmethyl)-4-piperidinamine
Synonyms: NAE-IN-M22, NEDD8-activating Enzyme Inhibitor M22
MF: C₂₀H₂₄Cl₂N₂
FW: 363.3
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

NEDD8 inhibitor M22 is supplied as a solid. A stock solution may be made by dissolving the NEDD8 inhibitor M22 in the solvent of choice, which should be purged with an inert gas. NEDD8 inhibitor M22 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of NEDD8 inhibitor M22 in these solvents is approximately 5 and 10 mg/ml, respectively. NEDD8 inhibitor M22 is slightly soluble in DMSO.

Description

NEDD8 inhibitor M22 is an inhibitor of NEDD8-activating enzyme (NAE).¹ It inhibits NEDDylation of ubiquitin-conjugating enzyme 12 (UBC12) in a cell-free assay using recombinant human NAE in a concentration-dependent manner. NEDD8 inhibitor M22 selectively inhibits NEDDylation over SUMOylation and ubiquitination in A549 non-small cell lung cancer (NSCLC) cells and inhibits the growth of A549 cells (GI₅₀ = 5.5 μM). It inhibits tumor growth in an AGS gastric adenocarcinoma mouse xenograft model when administered at a dose of 60 mg/kg.

Reference

1. Lu, P., Liu, X., Yuan, X., *et al.* Discovery of a novel NEDD8 activating enzyme inhibitor with piperidin-4-amine scaffold by structure-based virtual screening. *ACS Chem. Biol.* **11**(7), 1901-1907 (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, 04/25/2023

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