

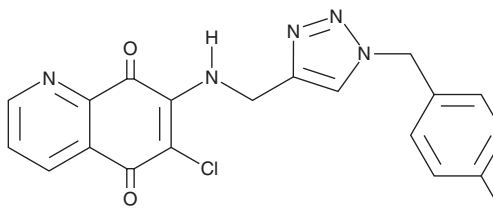
PRODUCT INFORMATION



M2N12

Item No. 37391

CAS Registry No.: 2376577-06-7
Formal Name: 6-chloro-7-[[[1-[(4-methylphenyl)methyl]-1H-1,2,3-triazol-4-yl]methyl]amino]-5,8-quinolinedione
MF: C₂₀H₁₆ClN₅O₂
FW: 393.8
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

M2N12 is supplied as a solid. A stock solution may be made by dissolving the M2N12 in the solvent of choice, which should be purged with an inert gas. M2N12 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of M2N12 in these solvents is approximately 10 mg/ml.

Description

M2N12 is a clickable inhibitor of cell division cycle 25 (Cdc25; IC₅₀s = 0.53, 1.39, and 0.09 μM for Cdc25A, Cdc25B, and Cdc25C, respectively).¹ It is composed of a naphthoquinone derivative functionalized with a clickable cyclic azide group. M2N12 inhibits the proliferation of A549 lung cancer and MDA-MB-231 and MCF-7 breast cancer cells (IC₅₀s = 3.92, 4.63, and 4.71 μM, respectively).

Reference

1. Jing, L., Wu, G., Hao, X., *et al.* Identification of highly potent and selective Cdc25 protein phosphatases inhibitors from miniaturization click-chemistry-based combinatorial libraries. *Eur. J. Med. Chem.* **183**, 111696 (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, 11/02/2023

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