

PRODUCT INFORMATION

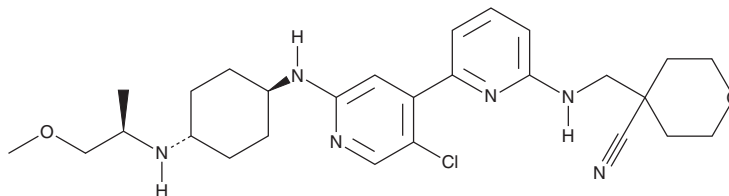


NVP-2

Item No. 34725

CAS Registry No.: 1263373-43-8
Formal Name: 4-[[[5'-chloro-2'-[[*trans*-4-[[[(1R)-2-methoxy-1-methylethyl]amino]cyclohexyl]amino][2,4'-bipyridin]-6-yl]amino]methyl]tetrahydro-2H-pyran-4-carbonitrile

MF: C₂₇H₃₇ClN₆O₂
FW: 513.1
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 255, 324 nm
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

NVP-2 is supplied as a solid. A stock solution may be made by dissolving the NVP-2 in the solvent of choice, which should be purged with an inert gas. NVP-2 is soluble in ethanol.

Description

NVP-2 is an inhibitor of the Cdk9/cyclin T1 complex (IC₅₀ = <0.514 nM).¹ It is selective for the Cdk9/cyclin T1 complex over nine additional Cdk complexes (IC₅₀s = 0.584->10 μM), as well as over a panel of 468 kinases at 1 μM. NVP-2 (250 nM) inhibits the proliferation of, and induces apoptosis in, MOLT-4 human acute lymphoblastic leukemia cells.

Reference

1. Olson, C.M., Jiang, B., Erb, M.A., *et al.* Pharmacological perturbation of CDK9 using selective CDK9 inhibition or degradation. *Nat. Chem. Biol.* **14**(2), 163-170 (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, 12/13/2022

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