

PRODUCT INFORMATION

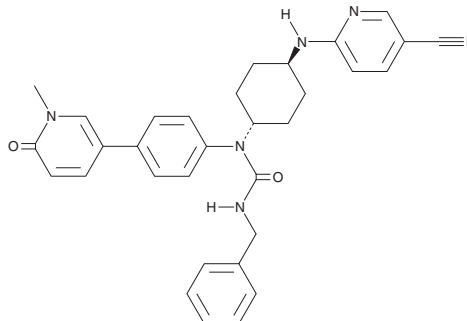


Cdk12-IN-2

Item No. 43703

CAS Registry No.: 2244987-03-7
Formal Name: N-[*trans*-4-[(5-cyano-2-pyridinyl)amino]cyclohexyl]-N-[4-(1,6-dihydro-1-methyl-6-oxo-3-pyridinyl)phenyl]-N'-(phenylmethyl)-urea
Synonyms: Cdk12 Inhibitor 2, Cyclin-dependent Kinase 12-IN-2, Cyclin-dependent Kinase 12 Inhibitor 2

MF: C₃₂H₃₂N₆O₂
FW: 532.6
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

Cdk12-IN-2 is supplied as a solid. A stock solution may be made by dissolving the Cdk12-IN-2 in the solvent of choice, which should be purged with an inert gas. Cdk12-IN-2 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Cdk12-IN-2 is an inhibitor of Cdk12 and Cdk13 (IC₅₀s = 0.052 and 0.01 μM, respectively).¹ It is selective for Cdk12 over Cdk2, -7, -8, and -9 (IC₅₀s = >100, >10, >10, and 16 μM, respectively). Cdk12-IN-2 inhibits RNA polymerase II phosphorylation in SK-BR-3 cells (IC₅₀ = 0.195 μM).

References

1. Ito, M., Tanaka, T., Toita, A., *et al.* Discovery of 3-benzyl-1-(*trans*-4-((5-cyanopyridin-2-yl)amino)cyclohexyl)-1-arylurea derivatives as novel and selective cyclin-dependent kinase 12 (CDK12) inhibitors. *J. Med. Chem.* **61**(17), 7710-7728 (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, 07/31/2025

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