

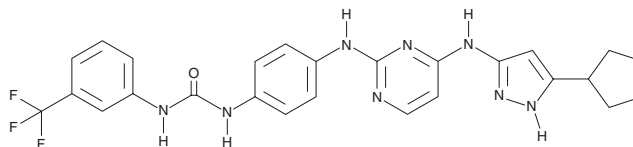
PRODUCT INFORMATION



CD532

Item No. 17054

CAS Registry No.: 1639009-81-6
Formal Name: N-[4-[[4-[(5-cyclopentyl-1H-pyrazol-3-yl)amino]-2-pyrimidinyl]amino]phenyl]-N'-[3-(trifluoromethyl)phenyl]-urea
MF: C₂₆H₂₅F₃N₈O
FW: 522.5
Purity: ≥98%
UV/Vis.: λ_{max}: 274 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

CD532 is supplied as a solid. A stock solution may be made by dissolving the CD532 in the solvent of choice, which should be purged with an inert gas. CD532 is soluble in the organic solvent ethanol.

Description

CD532 is an inhibitor of Aurora A kinase activity (IC₅₀ = 48 nM) and the protein-protein interaction between N-Myc and Aurora A kinase.¹ It also inhibits several cyclin-dependent kinases (CDKs), FGFRs, MEKs, and PDGFRs, as well as FLT3, KIT, and RET at 10 μM.² CD532 induces degradation of N-Myc in SK-N-BE(2) neuroblastoma cells (EC₅₀ = 223 nM).¹ It prevents S-phase entry in SK-N-BE(2) cells when used at a concentration of 1 μM.¹ CD532 (25 mg/kg) reduces tumor growth and increases survival in a MYCN-amplified SMS-KCN neuroblastoma mouse xenograft model.

References

1. Gustafson, W.C., Meyerowitz, J.G., Nekritz, E.A., *et al.* Drugging MYCN through an allosteric transition in Aurora kinase A. *Cancer Cell* **26(3)**, 414-427 (2014).
2. Lee, J.K., Phillips, J.W., Smith, B.A., *et al.* N-Myc drives neuroendocrine prostate cancer initiated from human prostate epithelial cells. *Cancer Cell* **29(4)**, 536-547 (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, 11/30/2022

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