

PRODUCT INFORMATION



NMS-P715

Item No. 31613

CAS Registry No.: 1202055-32-0
Formal Name: N-(2,6-diethylphenyl)-4,5-dihydro-1-methyl-8-[[4-[[[(1-methyl-4-piperidiny)amino]carbonyl]-2-(trifluoromethoxy)phenyl]amino]-1H-pyrazolo[4,3-h]quinazoline-3-carboxamide

MF: C₃₅H₃₉F₃N₈O₃

FW: 676.7

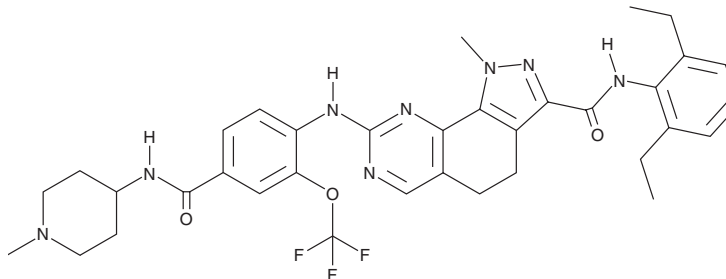
Purity: ≥95%

UV/Vis.: λ_{max}: 304 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

NMS-P715 is supplied as a solid. A stock solution may be made by dissolving the NMS-P715 in the solvent of choice, which should be purged with an inert gas. NMS-P715 is soluble in DMSO.

Description

NMS-P715 is an inhibitor of the spindle assembly checkpoint kinase Mps1/TTK (IC₅₀ = 0.182 μM).¹ It is selective for Mps1/TTK over a panel of 59 additional kinases (IC₅₀s = >5 μM for all). NMS-P715 (1 μM) accelerates mitosis, reducing the length of mitosis by approximately 3-fold, in U2OS cells. It induces aneuploidy and apoptosis in A2780 ovarian cancer cells when used at a concentration of 1 μM, as well as inhibits proliferation in a panel of 127 cancer cell lines, including colon, breast, renal, and melanoma cells, with IC₅₀ values ranging from 0.192 to 10 μM. NMS-P715 (90 mg/kg per day) reduces tumor growth in an A2780 mouse xenograft model.

Reference

1. Colombo, R., Caldarelli, M., Mennecozzi, M., *et al.* Targeting the mitotic checkpoint for cancer therapy with NMS-P715, an inhibitor of MPS1 kinase. *Cancer Res.* **70(24)**, 10255-10264 (2010).

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, 12/08/2022

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