

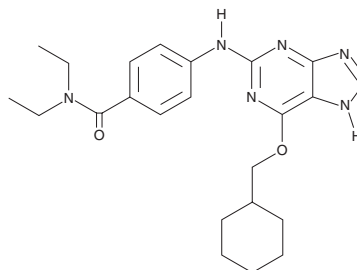
PRODUCT INFORMATION



NU 6140

Item No. 17271

CAS Registry No.: 444723-13-1
Formal Name: 4-[[6-(cyclohexylmethoxy)-9H-purin-2-yl]amino]-N,N-diethyl-benzamide
MF: C₂₃H₃₀N₆O₂
FW: 422.5
Purity: ≥98%
UV/Vis.: λ_{max}: 307 nm
Supplied as: A crystalline 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

NU 6140 is supplied as a crystalline solid. A stock solution may be made by dissolving the NU 6140 in the solvent of choice, which should be purged with an inert gas. NU 6140 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NU 6140 in these solvents is approximately 30 mg/ml.

NU 6140 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NU 6140 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. NU 6140 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NU 6140 is a cyclin-dependent kinase 2 (Cdk2) inhibitor (IC₅₀ = 0.41 μM) that demonstrates 10- to 36-fold selectivity against Cdk2-cyclin A compared to Cdk1-cyclin B, Cdk4-cyclin D, Cdk5-p25, or Cdk7-cyclin H.¹ NU 6140 has been found to induce cell-cycle arrest at the G₂-M phase and to potentiate the apoptotic effect of paclitaxel (Item No. 10461) in HeLa cells by inhibiting the anti-apoptotic protein, survivin.¹

Reference

1. Pennati, M., Campbell, A.J., Curto, M., *et al.* Potentiation of paclitaxel-induced apoptosis by the novel cyclin-dependent kinase inhibitor NU6140: A possible role for survivin down-regulation. *Mol. Cancer Ther.* 4(9), 1328-1337 (2005).

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, 11/14/2022

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