

PRODUCT INFORMATION



NVP-ADW742

Item No. 18390

CAS Registry No.: 475488-23-4
Formal Name: 5-[3-(phenylmethoxy)phenyl]-7-

[*trans*-3-(1-pyrrolidinylmethyl)
cyclobutyl]-7H-pyrrolo[2,3-d]
pyrimidin-4-amine

Synonyms: ADW742, GSK552602A

MF: C₂₈H₃₁N₅O

FW: 453.6

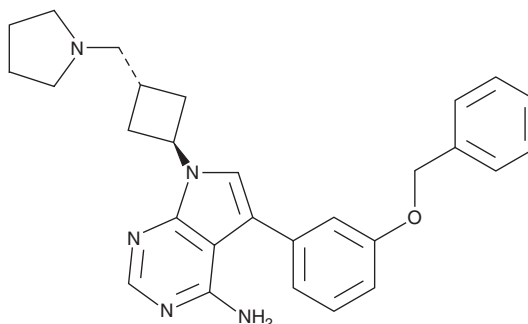
Purity: ≥98%

UV/Vis.: λ_{max}: 288 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

NVP-ADW742 is supplied as a crystalline solid. A stock solution may be made by dissolving the NVP-ADW742 in the solvent of choice, which should be purged with an inert gas. NVP-ADW742 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of NVP-ADW742 in these solvents is approximately 30 mg/ml.

NVP-ADW742 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NVP-ADW742 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. NVP-ADW742 has a solubility of approximately 0.09 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NVP-ADW742 is an insulin-like growth factor 1 receptor (IGF-1R) inhibitor with an IC₅₀ value of 0.17 μM.¹ It exhibits 6-fold greater selectivity for IGF-1R over the insulin receptor (InsR) kinase and minimal inhibitory activity against c-Kit, HER1, PDGFR, VEGFR2, or Bcr-Abl p210.¹ NVP-ADW742 has been shown to inhibit tumor proliferation both *in vitro* and in a mouse model of multiple myeloma.¹

Reference

1. Mitsiades, C. S., Mitsiades, N. S., McMullan, C. J., *et al.* Inhibition of the insulin-like growth factor receptor-1 tyrosine kinase activity as a therapeutic strategy for multiple myeloma, other hematologic malignancies, and solid tumors. *Cancer Cell* **5**(3), 221-230 (2004).

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, 12/02/2022

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