

PRODUCT INFORMATION



NVP-BHG712

Item No. 18489

CAS Registry No.: 940310-85-0
Formal Name: 4-methyl-3-[[1-methyl-6-(3-pyridinyl)-1H-pyrazolo[3,4-d]pyrimidin-4-yl]amino]-N-[3-(trifluoromethyl)phenyl]-benzamide

MF: C₂₆H₂₀F₃N₇O

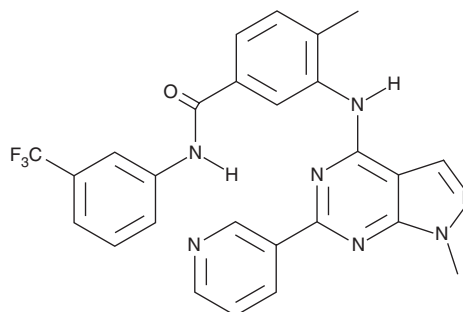
FW: 503.5

Purity: ≥98%

Stability: ≥2 years at -20°C

Supplied as: A crystalline solid

UV/Vis.: λ_{max}: 257 nm



Laboratory Procedures

For long term storage, we suggest that NVP-BHG712 be stored as supplied at -20°C. It should be stable for at least two years.

NVP-BHG712 is supplied as a crystalline solid. A stock solution may be made by dissolving the NVP-BHG712 in the solvent of choice. NVP-BHG712 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of NVP-BHG712 in ethanol is approximately 2 mg/ml and approximately 20 mg/ml in DMSO and DMF.

NVP-BHG712 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NVP-BHG712 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NVP-BHG712 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The EphB4 receptor tyrosine kinase (cloned from erythropoietin-producing hepatocellular carcinoma) and its ligand ephrinB2 play an important role in embryonic vessel development and vascular remodeling. Signaling from this pair is also involved in tumor angiogenesis. NVP-BHG712 is an orally bioavailable inhibitor of EphB4 kinase autophosphorylation (ED₅₀ = 25 nM).¹ It is selective for EphB4, demonstrating an ED₅₀ value of 4.2 μM at the related VEGFR2 receptor and ED₅₀ values >10 μM against a panel of 40 additional kinases.¹ At 3-30 mg/kg, NVP-BHG712 dose dependently inhibits VEGF-directed vessel formation in an *in vivo* mouse model of angiogenesis.¹

Reference

1. Martiny-Baron, G., Holzer, P., Billy, E., *et al.* The small molecule specific EphB4 kinase inhibitor NVP-BHG712 inhibits VEGF driven angiogenesis. *Angiogenesis* **13**, 259-267 (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, 10/30/2015

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