

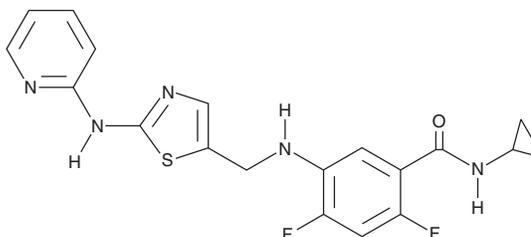
PRODUCT INFORMATION



BMS 605541

Item No. 43119

CAS Registry No.: 639858-32-5
Formal Name: N-cyclopropyl-2,4-difluoro-5-[[[2-(2-pyridinylamino)-5-thiazolyl]methyl]amino]-benzamide
MF: C₁₉H₁₇F₂N₅OS
FW: 401.4
Purity: ≥95%
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

BMS 605541 is supplied as a solid. A stock solution may be made by dissolving the BMS 605541 in the solvent of choice, which should be purged with an inert gas. BMS 605541 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

BMS 605541 is an inhibitor of VEGFR2 (IC₅₀ = 23 nM).¹ It is selective for VEGFR2 over 22 other kinases (IC₅₀s = 400->10,000 nM). BMS 605541 inhibits VEGF-induced proliferation in human umbilical vein endothelial cells (HUVECs; IC₅₀ = 25 nM). *In vivo*, BMS 605541 (90 or 180 mg/kg per day) decreases tumor volume in an L2987 lung adenocarcinoma mouse xenograft model.

Reference

1. Borzilleri, R.M., Bhide, R.S., Barrish, J.C., *et al.* Discovery and evaluation of N-cyclopropyl-2,4-difluoro-5-((2-(pyridin-2-ylamino)thiazol-5-ylmethyl)amino)benzamide (BMS-605541), a selective and orally efficacious inhibitor of vascular endothelial growth factor receptor-2. *J. Med. Chem.* **49(13)**, 3766-3769 (2006).

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, 06/25/2025

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