

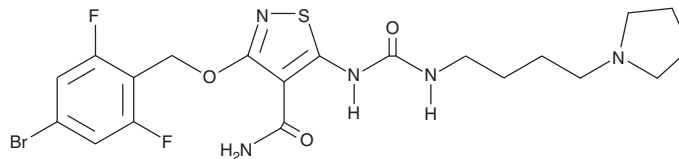
PRODUCT INFORMATION



CP 547,632

Item No. 21773

CAS Registry No.: 252003-65-9
Formal Name: 3-[[[4-(1-pyrrolidinyl)butyl]amino]carbonyl]-5-[[[4-(2,6-difluoro-3-bromophenyl)methoxy]amino]carbonyl]-4-isothiazolecarboxamide
MF: C₂₀H₂₄BrF₂N₅O₃S
FW: 532.4
Purity: ≥98%
UV/Vis.: λ_{max}: 258 nm
Supplied as: A solid
Storage: 4°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CP 547,632 is supplied as a solid. A stock solution may be made by dissolving the CP 547,632 in the solvent of choice. CP 547,632 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 94 mM.

Description

CP 547,632 is an orally bioavailable and potent inhibitor of the receptor tyrosine kinases VEGF receptor 2 (VEGFR2) and basic fibroblast growth factor (bFGF) with IC₅₀ values of 11 and 9 nM, respectively, in an enzyme assay.¹ It is selective for VEGFR2 and bFGF over EGFR, PDGF receptor β (PDGFRβ), and related tyrosine kinases. CP 547,632 inhibits VEGFR2 autophosphorylation induced by VEGF in porcine aortic endothelial cells transfected with VEGFR2 (IC₅₀ = 6 nM) and in a xenograft mouse model using NIH3T3/H-Ras cells (EC₅₀ = 590 nM). It decreases angiogenesis induced by VEGF or bFGF and suppresses tumor growth in athymic mice.

Reference

1. Beebe, J.S., Jani, J.P., Knauth, E., *et al.* Pharmacological characterization of CP-547,632, a novel vascular endothelial growth factor receptor-2 tyrosine kinase inhibitor for cancer therapy. *Cancer Res.* **63(21)**, 7301-7309 (2003).

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