

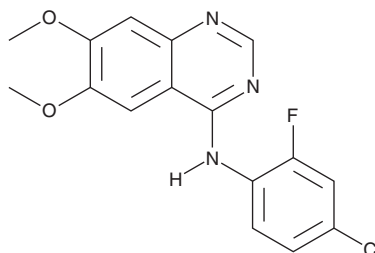
PRODUCT INFORMATION



ZM 306416

Item No. 16968

CAS Registry No.: 690206-97-4
Formal Name: N-(4-chloro-2-fluorophenyl)-6,7-dimethoxy-4-quinazolinamine
Synonym: CB-676475
MF: C₁₆H₁₃ClFN₃O₂
FW: 333.7
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 250, 331 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

ZM 306416 is supplied as a crystalline solid. A stock solution may be made by dissolving the ZM 306416 in the solvent of choice, which should be purged with an inert gas. ZM 306416 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ZM 306416 in ethanol is approximately 20 mg/ml and approximately 50 mg/ml in DMSO and DMF.

Description

ZM 306416 is an inhibitor of vascular endothelial growth factor (VEGF) receptor tyrosine kinase that antagonizes the activity of KDR and FLT VEGF receptors with IC₅₀ values of 100 nM and 2 μM, respectively.¹ It has also been reported to inhibit EGFR kinase with an IC₅₀ value of < 10 nM, reducing proliferation of EGFR-dependent cancer cells.²

References

1. Hennequin, L.F., Thomas, A.P., Johnstone, C., *et al.* Design and structure-activity relationship of a new class of potent VEGF receptor tyrosine kinase inhibitors. *J. Med. Chem.* **42(26)**, 5369-5389 (1999).
2. Antczak, C., Mahida, J.P., Bhinder, B., *et al.* A high-content biosensor-based screen identifies cell-permeable activators and inhibitors of EGFR function: Implications in drug discovery. *J. Biomol. Screen.* **17(7)**, 885-899 (2012).

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, 12/07/2022

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