

PRODUCT INFORMATION



FIIN-3

Item No. 20709

CAS Registry No.: 1637735-84-2
Formal Name: N-[4-[[[(2,6-dichloro-3,5-dimethoxyphenyl)amino]carbonyl][6-[[4-(4-methyl-1-piperazinyl)phenyl]amino]-4-pyrimidinyl]amino]methyl]phenyl]-2-propenamide

MF: C₃₄H₃₆Cl₂N₈O₄

FW: 691.6

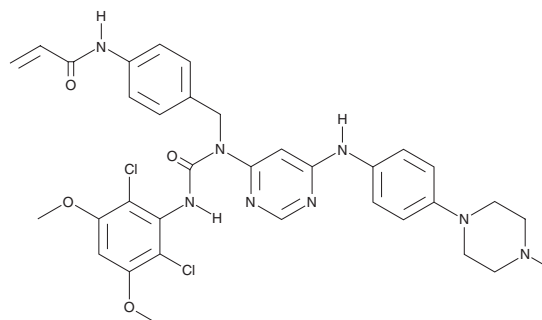
Purity: ≥98%

UV/Vis.: λ_{max}: 287 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

FIIN-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the FIIN-3 in the solvent of choice, which should be purged with an inert gas. FIIN-3 is sparingly soluble in methanol and acetonitrile.

Description

FIIN-3 is an inhibitor of FGF receptors (FGFRs; IC₅₀s = 13, 21, 31, and 35 nM for recombinant FGFR1-4, respectively).¹ It is selective for FGFRs over a panel of 456 kinases at a concentration of 1 μM, however, it does inhibit EGFR (IC₅₀ = 204 nM). FIIN-3 inhibits growth of Ba/F3 cells that are dependent on the kinase activity of wild-type FGFR1-4 as well as gatekeeper mutant FGFR2 and FGFR3 (EC₅₀s = <1-69 nM) and inhibits FGFR-dependent signaling in a concentration-dependent manner in FGFR2^{TEL/V564M}-dependent Ba/F3 cells. FIIN-3 also inhibits growth in a panel of cancer cell lines (EC₅₀s = 1.4-499 nM).

Reference

1. Tan, L., Wang, J., Tanizaki, J., *et al.* Development of covalent inhibitors that can overcome resistance to first-generation FGFR kinase inhibitors. *Proc. Natl. Acad. Sci. USA* **111**(45), E4869-E4877 (2014).

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, 11/10/2022

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