

PRODUCT INFORMATION

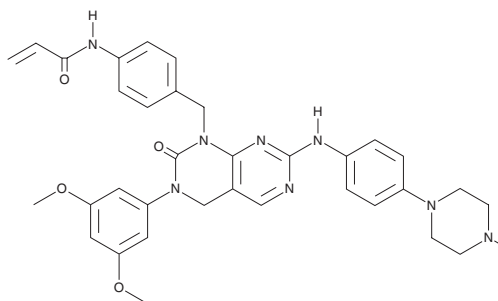


FIIN-2

Item No. 19837

CAS Registry No.: 1633044-56-0
Formal Name: N-[4-[[3-(3,5-dimethoxyphenyl)-3,4-dihydro-7-[[4-(4-methyl-1-piperazinyl)phenyl]amino]-2-oxopyrimido[4,5-d]pyrimidin-1(2H)-yl]methyl]phenyl]-2-propenamide

MF: C₃₅H₃₈N₈O₄
FW: 634.7
Purity: ≥98%
UV/Vis.: λ_{max}: 282 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-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the FIIN-2 in the solvent of choice. FIIN-2 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 10 mg/ml. FIIN-2 is also sparingly soluble in acetonitrile and methanol.

Description

FIIN-2 is an irreversible FGFR inhibitor with IC₅₀ values of 3.09, 4.3, 27, and 45.3 nM for FGFR1, FGFR2, FGFR3, and FGFR4, respectively.¹ It has been shown to inhibit cell proliferation of transformed Ba/F3 cell lines and demonstrates antiproliferative activity in cells dependent upon the gatekeeper mutants of FGFR1 or FGFR2.¹ By inhibiting FGFR in a zebrafish developmental model, FIIN-2 caused mild to severe alterations in tail morphogenesis similar to the phenotypes reported following genetic knockdown of FGFR.¹

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, 09/30/2022

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