

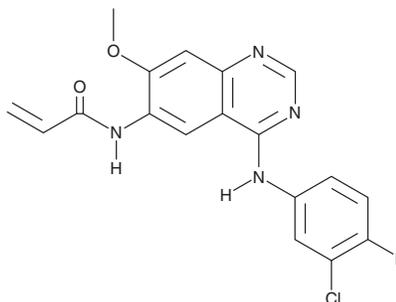
PRODUCT INFORMATION



PF-6274484

Item No. 22698

CAS Registry No.: 1035638-91-5
Formal Name: N-[4-[(3-chloro-4-fluorophenyl)amino]-7-methoxy-6-quinazoliny]-2-propenamide
MF: C₁₈H₁₄ClFN₄O₂
FW: 372.8
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 342 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

PF-6274484 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-6274484 in the solvent of choice, which should be purged with an inert gas. PF-6274484 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PF-6274484 in these solvents is approximately 25 and 33 mg/ml, respectively.

Description

PF-6274484 is an inhibitor of the EGF receptor (EGFR; IC₅₀s = 0.18 and 0.14 nM for wild-type EGFR and inhibitor-resistant EGFR^{L858R/T790M}, respectively).¹ It inhibits the autophosphorylation of wild-type EGFR in A549 cells and EGFR^{L858R/T790M} in H1975 cells (IC₅₀s = 5.8 and 6.6 nM, respectively).

Reference

1. Schwartz, P.A., Kuzmic, P., Solowiej, J., *et al.* Covalent EGFR inhibitor analysis reveals importance of reversible interactions to potency and mechanisms of drug resistance. *Proc. Nat. Acad. Sci. USA* **111**(1), 173-178 (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, 11/14/2022

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