

PRODUCT INFORMATION



PF-573228

Item No. 14924

CAS Registry No.: 869288-64-2
Formal Name: 3,4-dihydro-6-[[[4-[[[3-(methylsulfonyl)phenyl]methyl]amino]-5-(trifluoromethyl)-2-pyrimidinyl]amino]-2(1H)-quinolinone
Synonyms: FAK Inhibitor II, Focal Adhesion Kinase Inhibitor II

MF: C₂₂H₂₀F₃N₅O₃S

FW: 491.5

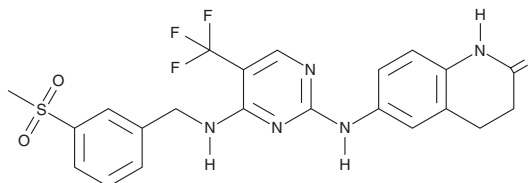
Purity: ≥95%

UV/Vis.: λ_{max}: 292 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-573228 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-573228 in the solvent of choice, which should be purged with an inert gas. PF-573228 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PF-573228 in these solvents is approximately 15 and 20 mg/ml, respectively.

Description

Focal adhesion kinases (FAKs) are non-receptor tyrosine kinases that play roles in regulating diverse processes, including cell adhesion, spreading, migration, proliferation, and apoptosis.¹ They are overexpressed in many types of cancer. PF-573228 inhibits FAK with IC₅₀ values of 4 and 30-100 nM for a purified recombinant catalytic fragment of FAK and in cultured cells, respectively.² It is 50-250-fold selective for FAK over other protein kinases.² PF-573228 can inhibit chemotactic and haptotactic migration of cells as well as prevent focal adhesion turnover.³

References

1. Parsons, J.T., Slack-Davis, J., Tilghman, R., *et al.* Focal adhesion kinase: Targeting adhesion signaling pathways for therapeutic intervention. *Clin. Cancer Res.* **14**(3), 627-632 (2008).
2. Medina, J.F., Haeggström, J., Kumlin, M., *et al.* Leukotriene A4: Metabolism in different rat tissues. *Biochim. Biophys. Acta* **961**, 203-212 (1988).
3. Slack-Davis, J.K., Martin, K.H., Tilghman, R.W., *et al.* Cellular characterization of a novel focal adhesion kinase inhibitor. *J. Biol. Chem.* **282**(20), 14845-14852 (2007).

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, 11/08/2022

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