

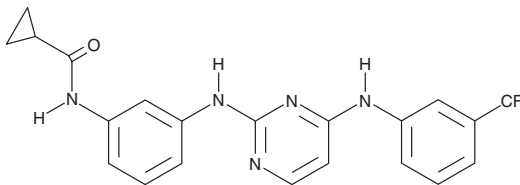
PRODUCT INFORMATION



Aurora Kinase Inhibitor III

Item No. 21851

CAS Registry No.: 879127-16-9
Formal Name: N-[3-[[4-[[3-(trifluoromethyl)phenyl]amino]-2-pyrimidinyl]amino]phenyl]-cyclopropanecarboxamide
MF: C₂₁H₁₈F₃N₅O
FW: 413.4
Purity: ≥98%
Supplied as: A 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

Aurora kinase inhibitor III is supplied as a solid. A stock solution may be made by dissolving the aurora kinase inhibitor III in the solvent of choice, which should be purged with an inert gas. Aurora kinase inhibitor III is soluble in organic solvents such as DMSO and ethanol. The solubility of aurora kinase inhibitor III in these solvents is approximately 20 and 50 mg/ml, respectively.

Description

Aurora kinase inhibitor III is a potent inhibitor of Aurora A kinase (IC₅₀ = 42 nM).¹ It is selective for Aurora A over BMX, BTK, IGF-1R, c-Src, TRKB, SYK, and EGFR (IC₅₀s = 386, 3,550, 591, 1,980, 2,510, 887, and >10,000 nM, respectively).

Reference

1. Zhang, Q., Liu, Y., Gao, F., *et al.* Discovery of EGFR selective 4,6-disubstituted pyrimidines from a combinatorial kinase-directed heterocycle library. *J. Am. Chem. Soc.* **128**(7), 2182-2183 (2006).

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, 04/05/2023

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