

PRODUCT INFORMATION

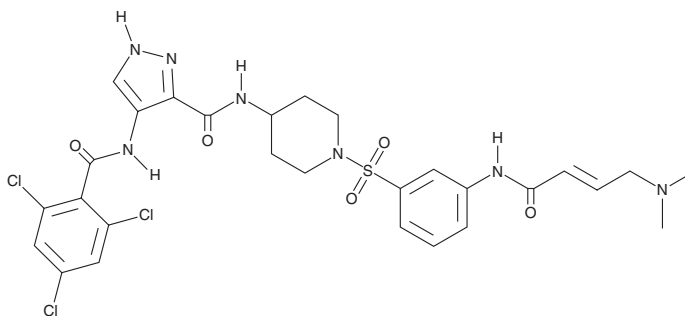


FMF-04-159-2

Item No. 39989

CAS Registry No.: 2364489-81-4
Formal Name: N-[1-[[[3-[[[(2E)-4-(dimethylamino)-1-oxo-2-buten-1-yl]amino]phenyl]sulfonyl]-4-piperidinyl]-4-[(2,4,6-trichlorobenzoyl)amino]-1H-pyrazole-3-carboxamide

MF: C₂₈H₃₀Cl₃N₇O₅S
FW: 683.0
Purity: ≥95%
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

FMF-04-159-2 is supplied as a solid. A stock solution may be made by dissolving the FMF-04-159-2 in the solvent of choice, which should be purged with an inert gas. FMF-04-159-2 is soluble in acetonitrile and DMSO.

Description

FMF-04-159-2 is an inhibitor of TAIRE family kinases.¹ It binds to Cdk14 and Cdk16 (IC₅₀s = 88 and 10.1 nM, respectively) and selectively inhibits Cdk14, Cdk16, Cdk17, and Cdk18 over a panel of 16 additional kinases at 1 μM but does not inhibit Cdk2 and Cdk10 by greater than 70% at 1 μM.

Reference

1. Ferguson, F.M., Doctor, Z.M., Ficarro, S.B., *et al.* Discovery of covalent CDK14 inhibitors with pan-TAIRE family specificity. *Cell. Chem. Biol.* **26**(6), 804-817 (2019).

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, 03/08/2024

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