

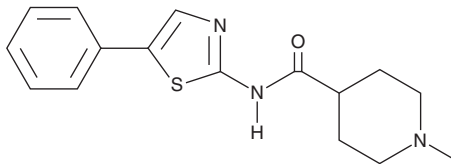
PRODUCT INFORMATION



DS12881479

Item No. 44121

CAS Registry No.: 2373065-59-7
Formal Name: 1-methyl-N-(5-phenyl-2-thiazolyl)-4-piperidinecarboxamide
MF: C₁₆H₁₉N₃OS
FW: 301.4
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

DS12881479 is supplied as a solid. A stock solution may be made by dissolving the DS12881479 in the solvent of choice, which should be purged with an inert gas. DS12881479 is soluble (≥10 mg/ml) in DMSO. DS12881479 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

DS12881479 is an inhibitor of MAP kinase-interacting serine/threonine protein kinase 1 (MNK1; IC₅₀s = 21 and 416 nM for the inactive and active forms, respectively).¹ It is selective for MNK1 over 46 other kinases but does inhibit FMS-related tyrosine kinase 3 (FLT3) and dual-specificity tyrosine phosphorylation-regulated kinase 1A (DYRK1A) at 5 μM.

Reference

1. Matsui, Y., Yasumatsu, I., Yoshida, K.I., *et al.* A novel inhibitor stabilizes the inactive conformation of MAPK-interacting kinase 1. *Acta Crystallogr. F Struct. Biol. Commun.* **74**(Pt. 3), 156-160 (2018).

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, 09/30/2025

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