

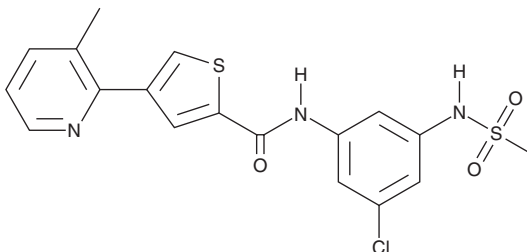
PRODUCT INFORMATION



ATX968

Item No. 44842

CAS Registry No.: 2973395-71-8
Formal Name: N-[3-chloro-5-[(methylsulfonyl)amino]phenyl]-4-(3-methyl-2-pyridinyl)-2-thiophenecarboxamide
Synonyms: DHX9-IN-2, DHX9 Inhibitor 2
MF: C₁₈H₁₆ClN₃O₃S₂
FW: 421.9
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

ATX968 is supplied as a solid. A stock solution may be made by dissolving the ATX968 in the solvent of choice, which should be purged with an inert gas. ATX968 is soluble (≥10 mg/ml) in DMSO.

Description

ATX968 is an inhibitor of DEAD-box helicase 9 (DDX9; IC₅₀ = 8 nM for helicase activity).¹ It is selective for DDX9 over DDX36, Brahma homolog (BRM), also known as SMARCA2, and Werner syndrome protein (WRN; IC₅₀s = >10, >10, and >100 μM, respectively), as well as a panel of 96 kinases at 10 μM but does not inhibit dual-specificity tyrosine phosphorylation-regulated kinase 1B (DYRK1B). ATX968 (1 μM) selectively induces unresolved nuclear R-loop accumulation in microsatellite instable-high with deficient mismatch repair (MSI-H/dMMR) LS411N colorectal carcinoma cells over microsatellite-stable with proficient mismatch repair (MSS/pMMR) NCI H747 colorectal carcinoma cells. It induces tumor regression in an LS411N mouse xenograft model when administered at a dose of 300 mg/kg.

Reference

1. Castro, J., Daniels, M.H., Brennan, D., *et al.* A potent, selective, small-molecule inhibitor of DHX9 abrogates proliferation of microsatellite instable cancers with deficient mismatch repair. *Cancer Res.* **85**(4), 758-776 (2025).

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, 01/22/2026

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