

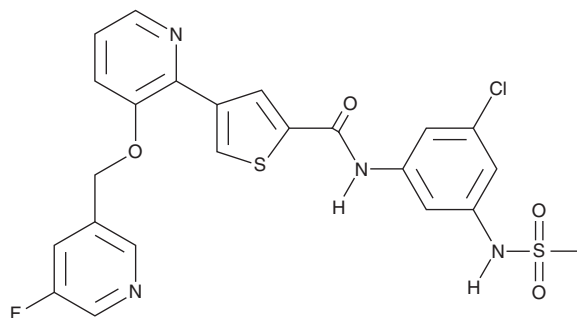
PRODUCT INFORMATION



DHX9-IN-6

Item No. 43559

CAS Registry No.: 2973401-56-6
Formal Name: N-[3-chloro-5-[(methylsulfonyl)amino]phenyl]-4-[3-[(5-fluoro-3-pyridinyl)methoxy]-2-pyridinyl]-2-thiophenecarboxamide
Synonyms: DEAD-box Helicase 9-IN-6, DEAD-box Helicase 9 Inhibitor 6, DHX9 Inhibitor 6
MF: C₂₃H₁₈ClFN₄O₄S₂
FW: 533.0
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

DHX9-IN-6 is supplied as a solid. A stock solution may be made by dissolving the DHX9-IN-6 in the solvent of choice, which should be purged with an inert gas. DHX9-IN-6 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DHX9-IN-6 is an inhibitor of DEAD-box helicase 9 (DDX9; IC₅₀ = 0.32 μM).¹ It reduces the proliferation of LS411N colorectal cancer cells (IC₅₀ = 0.0264 μM).

Reference

1. Daniels, M.H., Duncan, K.W., Sparling, B.A., *et al.* Inhibitors of RNA helicase DHX9 and uses thereof. WO2023/154519A1, Accent Therapeutics, Inc. (2023).

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, 08/07/2025

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