

PRODUCT INFORMATION

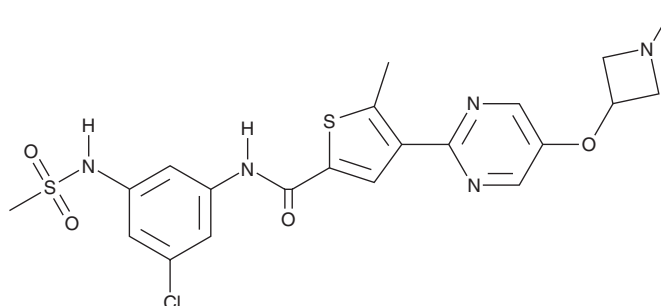


DHX9-IN-4

Item No. 45314

CAS Registry No.: 2973401-45-3
Formal Name: N-[3-chloro-5-[(methylsulfonyl)amino]phenyl]-5-methyl-4-[5-[(1-methyl-3-azetidinyloxy)-2-pyrimidinyl]-2-thiophenecarboxamide
Synonyms: DEAD-box Helicase 9-IN-4, DEAD-box Helicase 9 Inhibitor 4, DHX9 Inhibitor 4

MF: C₂₁H₂₂ClN₅O₄S₂
FW: 508.0
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-4 is supplied as a solid. A stock solution may be made by dissolving the DHX9-IN-4 in the solvent of choice, which should be purged with an inert gas. DHX9-IN-4 is slightly soluble (0.1-1 mg/ml) in acetonitrile and methanol.

Description

DHX9-IN-4 is an inhibitor of DEAD-box helicase 9 (DDX9; IC₅₀ = 3 nM).¹ It reduces the proliferation of LS411N cancer cells (IC₅₀ = 0.658 μM).

Reference

1. Daniels, M. H., Duncan, K. W., Sparling, B. A., *et al.* Inhibitors of RNA helicase DHX9 and uses thereof. (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, 03/17/2026

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