

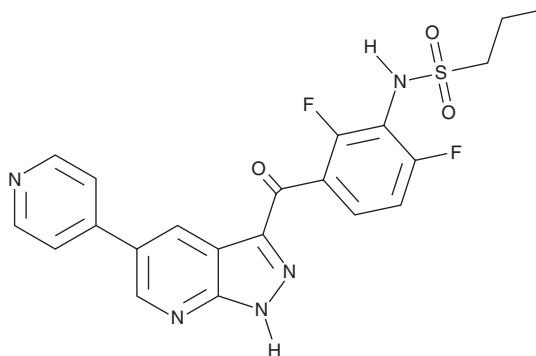
PRODUCT INFORMATION



HRX215

Item No. 41238

CAS Registry No.: 2369583-33-3
Formal Name: N-[2,6-difluoro-3-[[5-(4-pyridinyl)-1H-pyrazolo[3,4-b]pyridin-3-yl]carbonyl]phenyl]-1-propanesulfonamide
Synonyms: Darizmetinib, LN3348
MF: C₂₁H₁₇F₂N₅O₃S
FW: 457.5
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

HRX215 is supplied as a solid. A stock solution may be made by dissolving the HRX215 in the solvent of choice, which should be purged with an inert gas. HRX215 is soluble (≥10 mg/ml) in DMSO.

Description

HRX215 is an inhibitor of MAP kinase kinase 4 (MKK4; IC₅₀ = 0.02 μM).¹ It is selective for MKK4 over JNK1, B-RAF, and MKK7 (IC₅₀s = 7.07, 11.46, and 14.97 μM, respectively). HRX215 (10 mg/kg) decreases hepatocyte apoptosis in a mouse model of liver damage induced by carbon tetrachloride (CCl₄). It increases survival and hepatic size and regeneration, decreases intracranial pressure (ICP), and reduces blood ammonia, total bilirubin, and aspartate aminotransferase (AST) levels in a porcine model of partial hepatectomy-induced liver failure when administered at 5 mg/kg.

Reference

1. Zwirner, S., Rmilah, A.A.A., Klotz, S., et al. First-in-class MKK4 inhibitors enhance liver regeneration and prevent liver failure. *Cell* **187**(7), 1666-1684 (2024).

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, 04/23/2025

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