

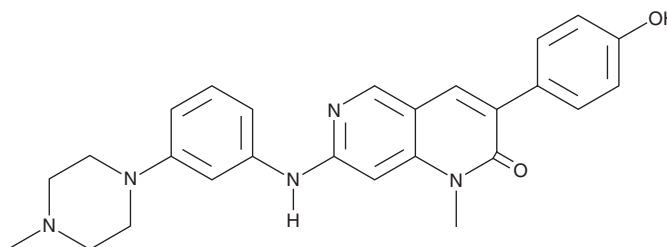
PRODUCT INFORMATION



UH15-38

Item No. 41056

CAS Registry No.: 2540881-21-6
Formal Name: 3-(3-hydroxyphenyl)-1-methyl-7-[[3-(4-methyl-1-piperazinyl)phenyl]amino]-1,6-naphthyridin-2(1H)-one
MF: $C_{26}H_{27}N_5O_2$
FW: 441.5
Purity: $\geq 95\%$
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

UH15-38 is supplied as a solid. A stock solution may be made by dissolving the UH15-38 in the solvent of choice, which should be purged with an inert gas. UH15-38 is soluble (≥ 10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Description

UH15-38 is an inhibitor of receptor-interacting serine/threonine kinase 3 (RIPK3; $IC_{50} = 0.02 \mu\text{M}$).¹ It is selective for RIPK3 over a panel of 50 receptors and enzymes at $1 \mu\text{M}$ and a panel of 77 kinases at $0.5 \mu\text{M}$ but does inhibit RIPK1 ($IC_{50} = 0.024 \mu\text{M}$) and 13 other kinases by greater than 90% at $0.5 \mu\text{M}$. UH15-38 inhibits LPS-induced necroptosis in RAW 264.7 macrophages and isolated mouse bone marrow-derived macrophages (BMDMs; IC_{50} s = 17 and 22 nM, respectively) and TNF- α -induced necroptosis in recombinant human RIPK3-expressing HeLa cells, Fas-associated death domain protein knockout (*FADD*^{-/-}) Jurkat cells, and U937 cells (IC_{50} s = 722, 205, and 329 nM, respectively). *In vivo*, UH15-38 (30 mg/kg) increases survival, decreases lung injury and immune cell bronchoalveolar fluid (BALF) infiltration, and inhibits decreases in arterial oxygen saturation in a mouse model of severe influenza A infection.

Reference

1. Gautam, A., Boyd, D.F., Nikhar, S., *et al.* Necroptosis blockade prevents lung injury in severe influenza. *Nature* **628**(8009), 835-843 (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, 03/19/2025

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