

PRODUCT INFORMATION



JH-X-119-01

Item No. 43041

CAS Registry No.: 2227368-54-7
Formal Name: N-[4-[[3-[(1-oxo-2-propen-1-yl)amino]benzoyl]amino]phenyl]-6-(1H-pyrazol-3-yl)-2-pyridinecarboxamide

MF: C₂₅H₂₀N₆O₃

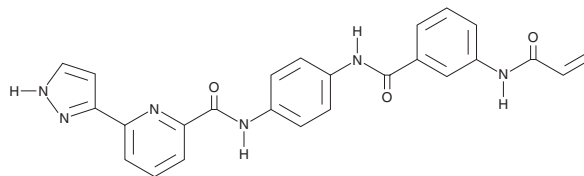
FW: 452.5

Purity: ≥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

JH-X-119-01 is supplied as a solid. A stock solution may be made by dissolving the JH-X-119-01 in the solvent of choice, which should be purged with an inert gas. JH-X-119-01 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

JH-X-119-01 is a covalent inhibitor of IL-1 receptor-associated kinase 1 (IRAK1; IC₅₀ = 9 nM).¹ It is selective for IRAK1 over IRAK4 (IC₅₀ = >10,000 nM). JH-X-119-01 is cytotoxic against a panel of mutant MyD88-expressing Waldenstrom's macroglobulinemia (WM), ABC subtype diffuse large B cell lymphoma (DLBCL), and lymphoma cells (EC₅₀s = 0.59-9.72 μM). It has a synergistic cytotoxic effect in MyD88-mutant-expressing BCWM.1 WM and TMD8 ABC DLBCL cells when used in combination with ibrutinib (Item No. 16274). JH-X-119-01 (5 and 10 mg/kg) also reduces lung injury and increases survival in a mouse model of LPS-induced sepsis.²

References

1. Hatcher, J.M., Yang, G., Wang, L., *et al.* Discovery of a selective, covalent IRAK1 inhibitor with antiproliferative activity in MYD88 mutated B-cell lymphoma. *ACS Med. Chem. Lett.* **11**(11), 2238-2243 (2020).
2. Pan, B., Gao, J., Chen, W., *et al.* Selective inhibition of interleukin-1 receptor-associated kinase 1 ameliorates lipopolysaccharide-induced sepsis in mice. *Int. Immunopharmacol.* **85**, 106597 (2020).

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, 03/21/2025

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