

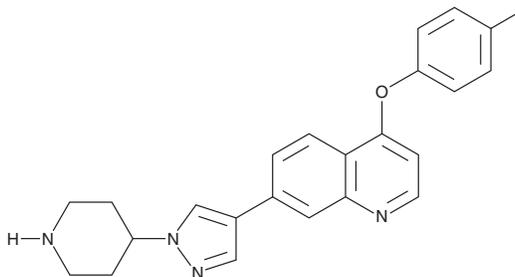
PRODUCT INFORMATION



HS-1371

Item No. 31033

CAS Registry No.: 2158197-70-5
Formal Name: 4-(4-methylphenoxy)-7-[1-(4-piperidinyl)-1H-pyrazol-4-yl]-quinoline
MF: C₂₄H₂₄N₄O
FW: 384.5
Purity: ≥98%
UV/Vis.: λ_{max}: 267 nm
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

HS-1371 is supplied as a solid. A stock solution may be made by dissolving the HS-1371 in the solvent of choice, which should be purged with an inert gas. HS-1371 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HS-1371 in DMF is approximately 30 mg/ml and approximately 5 mg/ml in ethanol and DMSO.

HS-1371 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HS-1371 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HS-1371 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HS-1371 is a receptor-interacting serine/threonine kinase 3 (RIPK3) inhibitor (IC₅₀ = 20.8 nM).¹ It inhibits basal RIPK3 autophosphorylation and TNF-α- or TRAIL-induced necroptosis in HT-29 cells when used at a concentration of 5 μM. It also inhibits necroptosis induced by a TNF-α Smac mimetic and the caspase inhibitor Z-VAD in HeLa cervical and NCI-H2009 lung cancer cells ectopically expressing RIPK3.

Reference

1. Park, H.-H., Park, S.-Y., Mah, S., *et al.* HS-1371, a novel kinase inhibitor of RIP3-mediated necroptosis. *Exp. Mol. Med.* **50(9)**, 1-15 (2018).

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, 12/14/2022

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