

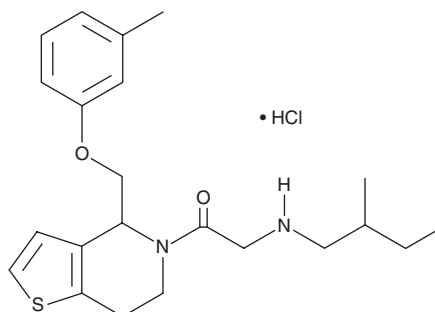
PRODUCT INFORMATION



RU-SKI 43 (hydrochloride)

Item No. 16972

CAS Registry No.: 1782573-67-4
Formal Name: 1-[6,7-dihydro-4-[(3-methylphenoxy)methyl]thieno[3,2-c]pyridin-5(4H)-yl]-2-[(2-methylbutyl)amino]-ethanone, monohydrochloride
MF: C₂₂H₃₀N₂O₂S • HCl
FW: 423.0
Purity: ≥98%
UV/Vis.: λ_{max}: 272 nm
Supplied as: A crystalline 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

RU-SKI 43 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the RU-SKI 43 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RU-SKI 43 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of RU-SKI 43 (hydrochloride) in these solvents is approximately 30 mg/ml.

RU-SKI 43 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RU-SKI 43 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. RU-SKI 43 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Hedgehog acyltransferase (Hhat) is an N-palmitoyltransferase that acylates Sonic hedgehog (Shh), which is critical for signaling through Shh. RU-SKI 43 is an inhibitor of Hhat (IC₅₀ = 0.85 μM).¹ It blocks palmitoylation of Shh without affecting palmitoylation of H-Ras or Fyn, myristoylation of c-Src, or acylation of Wnt3a.¹ RU-SKI 43 is cell-permeable and inhibits both autocrine and paracrine Shh-induced activation of Gli-mediated transcription.¹ In pancreatic cancer cells, RU-SKI 43 reduces both Gli1 activation and proliferation.²

References

1. Petrova, E., Rios-Esteves, J., Ourfelli, O., *et al.* Inhibitors of Hedgehog acyltransferase block Sonic Hedgehog signaling. *Nat. Chem. Biol.* **9(4)**, 247-249 (2013).
2. Petrova, E., Matevossian, A., and Resh, M.D. Hedgehog acyltransferase as a target in pancreatic ductal adenocarcinoma. *Oncogene* (2014).

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, 11/04/2022

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