

PRODUCT INFORMATION



RU-SKI 201 (hydrochloride)

Item No. 21101

CAS Registry No.: 2320262-09-5
Formal Name: 1-[6,7-dihydro-4-(6-methyl-2-pyridinyl)thieno[3,2-c]pyridin-5(4H)-yl]-2-[(2-methylbutyl)amino]-ethanone, dihydrochloride

MF: C₂₀H₂₇N₃OS • 2HCl

FW: 430.4

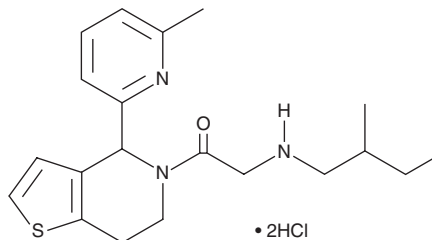
Purity: ≥95%

UV/Vis.: λ_{max}: 203, 266 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 201 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the RU-SKI 201 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RU-SKI 201 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of RU-SKI 201 (hydrochloride) in these solvents is approximately 30 mg/ml.

RU-SKI 201 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RU-SKI 201 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. RU-SKI 201 (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RU-SKI 201 is an inhibitor of hedgehog acyltransferase (Hhat; IC₅₀ = 0.2 μM for recombinant enzyme).¹ It inhibits palmitoylation of sonic hedgehog (Shh) in Shh-overexpressing HEK293 cells with a tagging IC₅₀ (TC₅₀) value of 0.87 μM. RU-SKI 201 inhibits hedgehog (Hh) signaling in H520, PANC-1, and MCF-7 cancer cells co-cultured with Shh-Light2 cells (IC₅₀s = 4.8, 7.8, and 8.5 μM, respectively).

Reference

1. Rodgers, U.R., Lanyon-Hogg, T., Masumoto, N., *et al.* Characterization of hedgehog acyltransferase inhibitors identifies a small molecule probe for hedgehog signaling by cancer cells. *ACS Chem. Biol.* **11**(12), 3256-3262 (2016).

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, 12/22/2022

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