

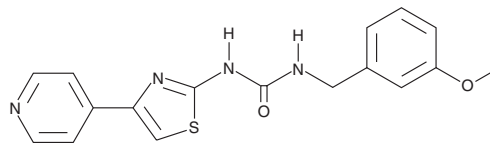
PRODUCT INFORMATION



RKI-1313

Item No. 22933

CAS Registry No.: 1342276-76-9
Formal Name: N-[(3-methoxyphenyl)methyl]-N'-[4-(4-pyridinyl)-2-thiazolyl]-urea
MF: C₁₇H₁₆N₄O₂S
FW: 340.4
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 265 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

RKI-1313 is supplied as a crystalline solid. A stock solution may be made by dissolving the RKI-1313 in the solvent of choice, which should be purged with an inert gas. RKI-1313 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of RKI-1313 in these solvents is approximately 30 mg/ml.

RKI-1313 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RKI-1313 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RKI-1313 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RKI-1313 is an inhibitor of Rho-associated kinase 1 (ROCK1) and 2 (IC₅₀s = 34 and 8 μM, respectively).¹ It has been used as a negative control for the ROCK inhibitor RKI-1447 (Item No. 16278).

Reference

1. Patel, R.A., Forinash, K.D., Pireddu, R., *et al.* RKI-1447 is a potent inhibitor of the Rho-associated ROCK kinases with anti-invasive and anti-tumor activities in breast cancer. *Cancer Res.* **72(19)**, 5025-5034 (2012).

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, 10/07/2022

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