

PRODUCT INFORMATION



8RK64

Item No. 41148

CAS Registry No.: 2705841-52-5

Formal Name: N-[5-(2-azidoacetyl)-4,5,6,7-tetrahydrothiazolo[5,4-c]pyridin-2-yl]-1-cyano-3S-pyrrolidinecarboxamide

MF: C₁₄H₁₆N₈O₂S

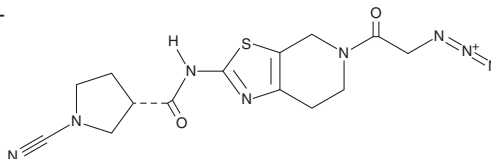
FW: 360.4

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

8RK64 is supplied as a solid. A stock solution may be made by dissolving the 8RK64 in the solvent of choice, which should be purged with an inert gas. 8RK64 is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly soluble (1-10 mg/ml) in DMSO.

Description

8RK64 is a covalent inhibitor of ubiquitin C-terminal hydrolase L1 (UCH-L1; IC₅₀ = 0.32 μM) and a click chemistry reagent.¹ It is selective for UCH-L1 over UCH-L3 and UCH-L5 (IC₅₀s = 216 and >1,000 μM, respectively). 8RK64 has been linked to fluorophores containing alkyne groups via copper(I)-catalyzed azide/alkyne cycloaddition (CuAAC) and used to visualize UCH-L1 activity in cells. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Kooij, R., Liu, S., Sapmaz, A., *et al.* Small-molecule activity-based probe for monitoring ubiquitin C-terminal hydrolase L1 (UCHL1) activity in live cells and zebrafish embryos. *J. Am. Chem. Soc.* **142**(39), 16825-16841 (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, 08/06/2024

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