

PRODUCT INFORMATION



KSQ-4279

Item No. 41862

CAS Registry No.: 2446480-97-1

Formal Name: 6-(4-cyclopropyl-6-methoxy-5-pyrimidinyl)-1-[[4-[1-(1-methylethyl)-4-(trifluoromethyl)-1H-imidazol-2-yl]phenyl]methyl]-1H-pyrazolo[3,4-d]pyrimidine

Synonyms: Ubiquitin-specific Protease 1 Inhibitor 1, USP1-IN-1, USP1 Inhibitor 1

MF: C₂₇H₂₅F₃N₈O

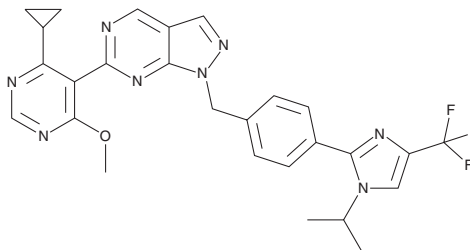
FW: 534.5

Purity: ≥98%

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

KSQ-4279 is supplied as a solid. A stock solution may be made by dissolving the KSQ-4279 in the solvent of choice, which should be purged with an inert gas. KSQ-4279 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

KSQ-4279 is an inhibitor of ubiquitin-specific protease 1 (USP1; IC₅₀ = 21.1 nM).¹ It reduces the viability of H1299 non-small cell lung cancer (NSCLC) cells when used at a concentration of 500 nM. KSQ-4279 (1 μM) induces cell cycle arrest at the S phase in MDA-MB-436 breast cancer cells.

Reference

1. Li, H., Liu, B.J., Xu, J., *et al.* Design, synthesis, and biological evaluation of pyrido[2,3-d]pyrimidin-7(8H)-one derivatives as potent USP1 inhibitors. *Eur. J. Med. Chem.* **275:116568**, (2024).

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, 09/26/2024

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