

PRODUCT INFORMATION



USP8-IN-3

Item No. 41600

CAS Registry No.: 2477651-10-6
Formal Name: N-(4-nitrophenyl)-N'-[1-[6-(trifluoromethyl)-2-pyridinyl]-4-piperidinyl]-thiourea

Synonyms: Ubiquitin-specific Protease 8-IN-3,
Ubiquitin-specific Protease 8 Inhibitor 3,
USP 8 Inhibitor 3

MF: C₁₈H₁₈F₃N₅O₂S

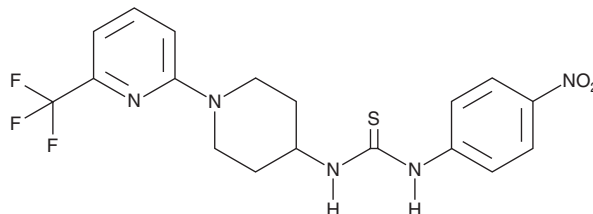
FW: 425.4

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

USP8-IN-3 is supplied as a solid. A stock solution may be made by dissolving the USP8-IN-3 in the solvent of choice, which should be purged with an inert gas. USP8-IN-3 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

USP8-IN-3 is an inhibitor of ubiquitin-specific protease 8 (USP8; IC₅₀ = 4 μM).¹ It inhibits the growth of H1975 non-small cell lung cancer (NSCLC) and GH3 rat pituitary cells (GI₅₀s = 6.01 and 37.03 μM).

Reference

1. Li, Z., Luo, C., Bian, J., et al. Preparing method and application of USP8 inhibitor. *China Pharmaceutical University, China*. CN111138358 (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, 09/20/2024

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