

PRODUCT INFORMATION



SJB2-043 analog (trifluoroacetate salt)

Item No. 38497

Formal Name: (R)-2-(2-chloro-6-fluoro-3-((1-methylpyrrolidin-3-yl)oxy)phenyl)naphtho[2,3-d]oxazole-4,9-dione, trifluoroacetate salt

Synonyms: OTUB1/Ubiquitin-specific Protease 8-IN-1, OTUB1/Ubiquitin-specific Protease 8 Inhibitor 1, OTUB1/USP8-IN-1, OTUB1/USP8 Inhibitor i

MF: C₂₂H₁₆ClFN₂O₄ • XCF₃COOH

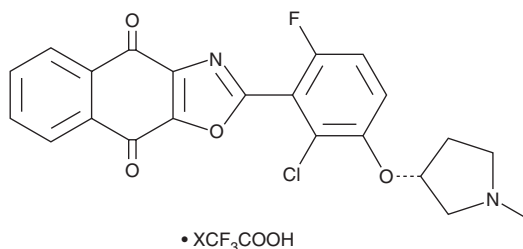
FW: 426.8

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

SJB2-043 analog (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SJB2-043 analog (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SJB2-043 analog (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

SJB2-043 analog is an inhibitor of OTUB1 and ubiquitin-specific protease 8 (USP8; IC₅₀s = 0.17 and 0.28 nM, respectively) and a derivative of the USP1/UAF1 complex and SARS-CoV-2 PL^{Pro} inhibitor SJB2-043 (Item No. 36605).¹ It is selective for OTUB1 and USP8 over USP1, USP2, USP7, USP36, and ataxin-3 (IC₅₀s = 1,370, 3,800, >10,000, 8,400, and >10,000 nM, respectively). SJB2-043 analog reduces viability in H1975, EBC-1, NCI H1703, H23, and A549 non-small cell lung cancer (NSCLC) cells (IC₅₀s = 118, 172, 145, 431, and 1,004 nM, respectively). *In vivo*, SJB2-043 analog (5 mg/kg twice daily) decreases tumor volume in an H1975 mouse xenograft model.

Reference

1. Tan, L., Shan, H., Han, C., *et al.* Discovery of potent OTUB1/USP8 dual inhibitors targeting proteostasis in non-small-cell lung cancer. *J. Med. Chem.* **65**(20), 13645-13659 (2022).

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, 08/30/2024

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