

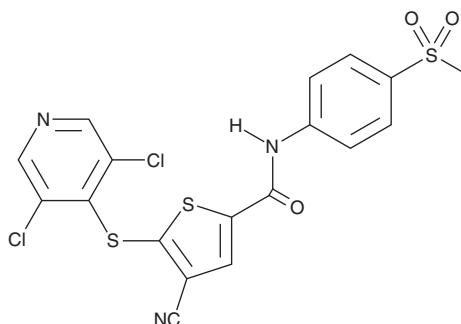
PRODUCT INFORMATION



USP7-USP47 Inhibitor

Item No. 25622

CAS Registry No.: 1247825-37-1
Formal Name: 4-cyano-5-[(3,5-dichloro-4-pyridinyl)thio]-N-[4-(methylsulfonyl)phenyl]-2-thiophenecarboxamide
MF: C₁₈H₁₁Cl₂N₃O₃S₃
FW: 484.4
Purity: ≥98%
UV/Vis.: λ_{max}: 237, 326 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

USP7-USP47 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the USP7-USP47 inhibitor in the solvent of choice. USP7-USP47 inhibitor is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 44 mg/ml.

Description

USP7-USP47 inhibitor is a dual inhibitor of ubiquitin-specific protease 7 (USP7) and USP47 (IC₅₀s = 0.42 and 1 μM, respectively).¹ It is selective for USP7 and USP47 over USP2, USP5, USP8, USP21, USP28, caspase-3, and cathepsin B (IC₅₀s = >31.6 μM). USP7-USP47 inhibitor inhibits the growth of HCT116 cells (EC₅₀ = 7.6 μM).

Reference

1. Weinstock, J., Wu, J., Cao, P., *et al.* Selective dual inhibitors of the cancer-related deubiquitylating proteases USP7 and USP47. *ACS Med. Chem. Lett.* **3**(10), 789-792 (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, 12/19/2022

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