

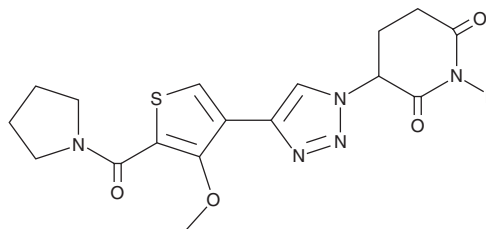
PRODUCT INFORMATION



TMX-4116

Item No. 43720

CAS Registry No.: 2766385-56-0
Formal Name: 3-[4-[4-methoxy-5-(1-pyrrolidinylcarbonyl)-3-thienyl]-1H-1,2,3-triazol-1-yl]-2,6-piperidinedione
MF: C₁₇H₁₉N₅O₄S
FW: 389.4
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

TMX-4116 is supplied as a solid. A stock solution may be made by dissolving the TMX-4116 in the solvent of choice, which should be purged with an inert gas. TMX-4116 is soluble in the organic solvent acetonitrile (≥10 mg/ml).

Description

TMX-4116 is a casein kinase 1α (CK1α) degrader.¹ It selectively induces E3 ubiquitin ligase-mediated degradation of CK1α over phosphodiesterase 6D (PDE6D), Ikaros family zinc finger protein 1 (IKZF1), and IKZF3 in MM.1S cells when used at concentrations ranging from 0.04 to 1 μM.

Reference

1. Teng, M., Lu, W., Donovan, K.A., *et al.* Development of PDE6D and CK1α degraders through chemical derivatization of FPFT-2216. *J. Med. Chem.* **65**(1), 747-756 (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, 07/28/2025

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