

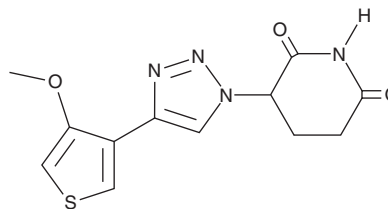
PRODUCT INFORMATION



FPFT-2216

Item No. 43551

CAS Registry No.: 2367619-87-0
Formal Name: 3-[4-(4-methoxy-3-thienyl)-1H-1,2,3-triazol-1-yl]-2,6-piperidinedione
MF: $C_{12}H_{12}N_4O_3S$
FW: 292.3
Purity: $\geq 98\%$
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

FPFT-2216 is supplied as a solid. A stock solution may be made by dissolving the FPFT-2216 in the solvent of choice, which should be purged with an inert gas. FPFT-2216 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

FPFT-2216 is a multi-protein degrader.¹ It degrades phosphodiesterase 6D (PDE6D), Ikaros family zinc finger protein 1 (IKZF1), IKZF3, and casein kinase 1 α (CK1 α) in MOLT-4 cells when used at a concentration of 1 μM .

Reference

1. Teng, M., Lu, W., Donovan, K.A., *et al.* Development of PDE6D and CK1 α 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, 08/04/2025

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