

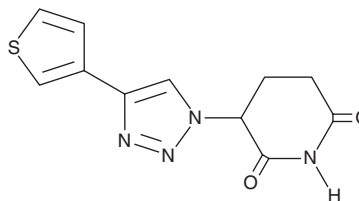
PRODUCT INFORMATION



TMX-4100

Item No. 44276

CAS Registry No.: 2367619-63-2
Formal Name: 3-[4-(3-thienyl)-1H-1,2,3-triazol-1-yl]-2,6-piperidinedione
MF: $C_{11}H_{10}N_4O_2S$
FW: 262.3
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

TMX-4100 is supplied as a solid. A stock solution may be made by dissolving the TMX-4100 in the solvent of choice, which should be purged with an inert gas. TMX-4100 is soluble (≥ 10 mg/ml) in DMSO.

Description

TMX-4100 is a molecular glue degrader of phosphodiesterase 6D (PDE6D).¹ It promotes the interaction between the E3 ubiquitin ligase cereblon and PDE6D, inducing the degradation of PDE6D in MOLT-4 T cell lymphoblasts, Jurkat leukemia cells, and MM.1S multiple myeloma cells with half-maximal degradation concentration (DC_{50}) values of <200 nM for all.

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, 10/07/2025

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