

PRODUCT INFORMATION



Dth

Item No. 38925

Formal Name: (Z)-4-((6-(4-(3,5-difluoro-4-hydroxybenzylidene)-2-methyl-5-oxo-4,5-dihydro-1H-imidazol-1-yl)hexyl)amino)-2-(2,6-dioxopiperidin-3-yl)isoindoline-1,3-dione

MF: $C_{30}H_{29}F_2N_5O_6$

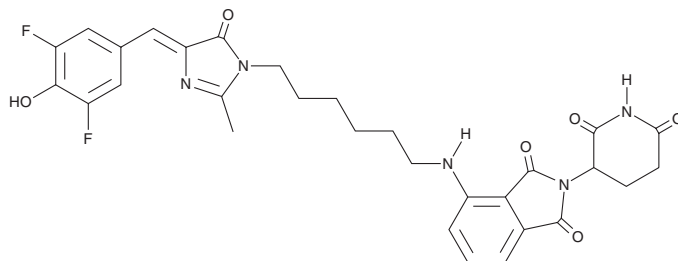
FW: 593.6

Purity: $\geq 95\%$

Supplied as: A 1 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Dth is a proteolysis-targeting chimera (PROTAC) composed of the fluorescent probe for RNA DFHBI (Item No. 34555) and the immunomodulatory compound thalidomide (Item No. 14610).¹ It binds to a genetically expressed RNA aptamer comprising Broccoli and a modifiable and selective protein targeting sequence and induces degradation of mCherry in HEK293T cells when used at a concentration of 20 μM . Dth (20 μM) induces degradation of the NF- κB subunit p50, as well as p50 and the E2F transcription factor subtype E2F1, in MCF-7 cells.

Reference

1. Xu, Y., Fu, D., Yuan, Y., *et al.* A heterobifunctional molecule recruits cereblon to an RNA scaffold and activates its PROTAC function. *Cell Rep. Phys. Sci.* **3**(10), 101064 (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, 01/26/2026

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