

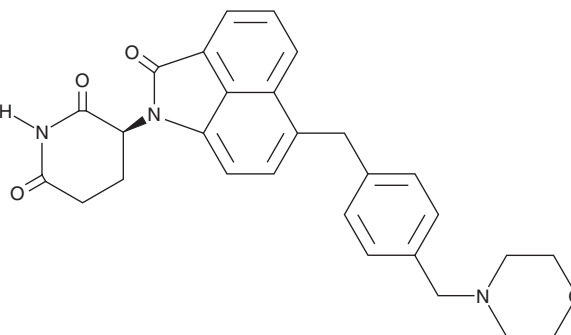
PRODUCT INFORMATION



CFT-7455

Item No. 42053

CAS Registry No.: 2504235-67-8
Formal Name: 3S-[6-[4-(4-morpholinylmethyl)phenyl]methyl]-2-oxobenz[cd]indol-1(2H)-yl]-2,6-piperidinedione
Synonym: Cemsidomide
MF: C₂₈H₂₇N₃O₄
FW: 469.5
Purity: ≥95%
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

CFT-7455 is supplied as a solid. A stock solution may be made by dissolving the CFT-7455 in the solvent of choice, which should be purged with an inert gas. CFT-7455 is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO.

Description

CFT-7455 is a cereblon ligand 1-containing degrader of Ikaros family zinc finger protein 1 (IKZF1) and IKZF3.¹ It reduces levels of IKZF1 and IKZF3 in NCI H929 B cell myeloma cells when used at a concentration of 0.1 nM. CFT-7455 decreases the growth of NCI H929 cells (GI₅₀ = 0.0743 nM). It reduces tumor volume in an RPMI-8226 plasmocytoma mouse xenograft model when administered at doses of 30 or 100 µg/kg per day.

Reference

1. Proia, D., Henderson, J.A., He, M., *et al.* Advantageous therapies for disorders mediated by karos or aiolos. (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, 12/16/2024

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