

PRODUCT INFORMATION



CPS2

Item No. 40457

CAS Registry No.: 2756741-90-7

Formal Name: 4-[[[5-amino-1-[(3-methyl-2-thienyl)carbonyl]-1H-1,2,4-triazol-3-yl]amino]-N-[2-[[1-[2-[2-[2-[[2-(2,6-dioxo-3-piperidiny)]-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]ethoxy]ethoxy]ethyl]-1H-1,2,3-triazol-4-yl]methoxy]ethyl]-benzenesulfonamide

MF: C₃₈H₄₂N₁₂O₁₀S₂

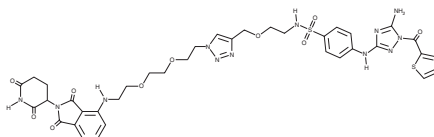
FW: 890.9

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

CPS2 is supplied as a solid. A stock solution may be made by dissolving the CPS2 in the solvent of choice, which should be purged with an inert gas. CPS2 is soluble in acetonitrile and methanol.

Description

CPS2 is a proteolysis-targeting chimera (PROTAC) composed of the Cdk2 binder J2, a derivative of the Cdk and Aurora kinase inhibitor JNJ-7706621 (Item No. 18494), connected to the cereblon inhibitor pomalidomide (Item No. 19877) by a linker region.¹ It selectively degrades Cdk2 (IC₅₀ = 24 nM) over seven additional CDKs at 500 nM but also degrades Aurora A kinase and Ikaros family zinc finger protein 1 (IKZF1) at 250 nM. CPS2 reduces Cdk2 levels and increases the percentage of cells expressing CD11b, a marker of differentiation, in primary acute myeloid lymphoma (AML) cells. It also decreases proliferation without reducing the viability of, as well as induces differentiation in, primary hematopoietic stem cells (HSCs).

Reference

1. Wang, L., Shao, X., Zhong, T., *et al.* Discovery of a first-in-class CDK2 selective degrader for AML differentiation therapy. *Nat. Chem. Biol.* **17**(5), 567-575 (2021).

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, 03/13/2024

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