

PRODUCT INFORMATION



BSJ-4-116

Item No. 39564

CAS Registry No.: 2519823-34-6

Formal Name: N-[7-[(3R)-3-[[5-chloro-4-[[2-[(1-methylethyl)sulfonyl]phenyl]amino]-2-pyrimidinyl]amino]-1-piperidinyl]heptyl]-2-[[2-(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]oxy]-acetamide

MF: C₄₀H₄₉ClN₈O₈S

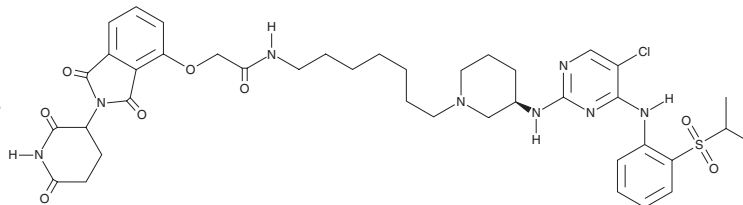
FW: 837.4

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

BSJ-4-116 is supplied as a solid. A stock solution may be made by dissolving the BSJ-4-116 in the solvent of choice, which should be purged with an inert gas. BSJ-4-116 is soluble in methanol.

Description

BSJ-4-116 is a proteolysis-targeting chimera (PROTAC) composed of a modified form of the multi-kinase degrader TL12-186 linked to the cereblon ligand and PROTAC synthetic intermediate E3 ligase ligand 2 (Item No. 36664).¹ It induces time- and concentration-dependent degradation of cyclin-dependent kinase 12 (Cdk12) in Jurkat T cells. BSJ-4-116 (50 nM) decreases expressions of various genes associated with DNA double-strand break repair in Jurkat cells. It inhibits the growth of MOLT-4 leukemia cells in a cereblon-dependent manner.

Reference

1. Jiang, B., Gao, Y., Che, J., *et al.* Discovery and resistance mechanism of a selective CDK12 degrader. *Nat. Chem. Biol.* **17**(6), 675-683 (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, 11/02/2023

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