

PRODUCT INFORMATION



BSJ-04-132

Item No. 41248

CAS Registry No.: 2349356-39-2
Formal Name: 7-cyclopentyl-2-[[[5-[4-[4-[[2-[(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]oxy]acetyl]amino]butyl]-1-piperazinyl]-2-pyridinyl]amino]-N,N-dimethyl-7H-pyrrolo[2,3-d]pyrimidine-6-carboxamide

MF: C₄₂H₄₉N₁₁O₇

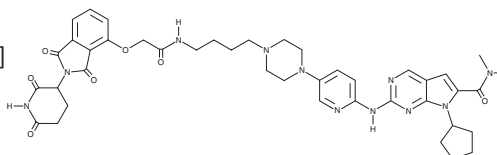
FW: 819.9

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

BSJ-04-132 is supplied as a solid. A stock solution may be made by dissolving the BSJ-04-132 in the solvent of choice, which should be purged with an inert gas. BSJ-04-132 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

BSJ-04-132 is a proteolysis-targeting chimera (PROTAC) containing the cyclin-dependent kinase 4 (Cdk4) and Cdk6 inhibitor ribociclib (LEE011; Item Nos. 17666 | 36414) conjugated to (±)-thalidomide (Item No. 14610).¹ It selectively induces the degradation of Cdk4 over Cdk6, as well as Ikaros family zinc finger protein 1 (IKZF1) and IKZF3, which are targets of other imide-based degraders, in a cereblon-dependent manner in Jurkat cells when used at a concentration of 1 μM.

Reference

1. Jiang, B., Wang, E.S., Donovan, K.A., *et al.* Development of dual and selective degraders of cyclin-dependent kinases 4 and 6. *Angew. Chem. Int. Ed. Engl.* **58(19)**, 6321-6326 (2019).

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, 11/15/2024

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