

PRODUCT INFORMATION



dBRD9

Item No. 45720

CAS Registry No.: 2170679-45-3

Formal Name: 2-[[[4-(1,2-dihydro-2-methyl-1-oxo-2,7-naphthyridin-4-yl)-2,6-dimethoxyphenyl]methyl]methylamino]-N-[2-[2-[2-[(2,6-dioxo-3-piperidinyl)-2,3-dihydro-1,3-dioxo-1H-isoindol-4-yl]amino]ethoxy]ethoxy]ethyl]-acetamide

MF: $C_{40}H_{45}N_7O_{10}$

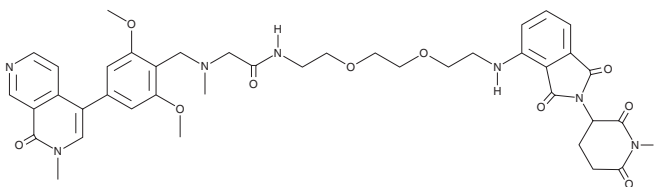
FW: 783.8

Purity: $\geq 98\%$

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

dBRD9 is supplied as a solid. A stock solution may be made by dissolving the dBRD9 in the solvent of choice, which should be purged with an inert gas. dBRD9 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

dBRD9 is a proteolysis-targeting chimera (PROTAC) composed of a bromodomain-containing protein 9 (BRD9) inhibitor conjugated to the cereblon (CRBN) inhibitor pomalidomide (Item No. 19877) via an alkyl ether linker.¹ It selectively degrades BRD9 over BRD4 and BRD7 at 100 nM and reduces the proliferation of MOLM-13 leukemia cells in a concentration-dependent manner.

Reference

1. Remillard, D., Buckley, D.L., Paulk, J., *et al.* Degradation of the BAF complex factor BRD9 by heterobifunctional ligands. *Angew. Chem. Int. Ed. Engl.* **56**(21), 5738-5743 (2017).

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, 06/01/2026

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