

# PRODUCT INFORMATION

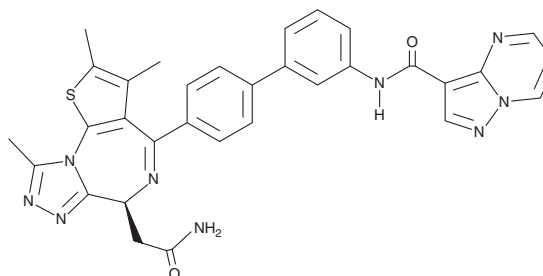


## PLX3618

Item No. 45519

**CAS Registry No.:** 2851986-76-8  
**Formal Name:** (6S)-2,3,9-trimethyl-4-[3'-[(pyrazolo[1,5-a]pyrimidin-3-ylcarbonyl)amino][1,1'-biphenyl]-4-yl]-6H-thieno[3,2-f][1,2,4]triazolo[4,3-a][1,4]diazepine-6-acetamide

**MF:** C<sub>32</sub>H<sub>27</sub>N<sub>9</sub>O<sub>2</sub>S  
**FW:** 601.7  
**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

PLX3618 is supplied as a solid. A stock solution may be made by dissolving the PLX3618 in the solvent of choice, which should be purged with an inert gas. PLX3618 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

PLX3618 is a monovalent direct degrader of bromodomain-containing protein 4 (BRD4).<sup>1</sup> It binds to both bromodomain 1 (BD1) and BD2 of BRD4 and mediates its degradation *via* recruitment of the E3 ubiquitin ligase DDB1- and CUL4-associated factor 11 (DCAF11) with a half-maximal degradation concentration (DC<sub>50</sub>) value of 12.2 nM. PLX3618 inhibits the proliferation of a variety of hematopoietic cancer cells and induces apoptosis in MV4-11 acute myeloid leukemia (AML) cells. It induces tumor regression and intratumoral BRD4 degradation in an MV4-11 mouse xenograft model when administered at a dose of 5 mg/kg.

### Reference

1. Parker, G.S., Toth, J.I., Fish, S., *et al.* Discovery of monovalent direct degraders of BRD4 that act via the recruitment of DCAF11. *Mol. Cancer Ther.* **23(10)**, 1446-1458 (2024).

#### 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, 04/21/2026

#### 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