

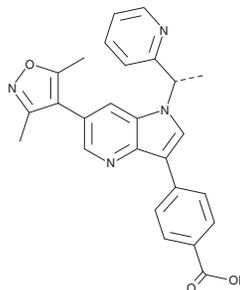
# PRODUCT INFORMATION



## PLX51107

Item No. 30742

**CAS Registry No.:** 1627929-55-8  
**Formal Name:** 4-[6-(3,5-dimethyl-4-isoxazolyl)-1-[(1S)-1-(2-pyridinyl)ethyl]-1H-pyrrolo[3,2-b]pyridin-3-yl]-benzoic acid  
**MF:** C<sub>26</sub>H<sub>22</sub>N<sub>4</sub>O<sub>3</sub>  
**FW:** 438.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 227, 251, 305 nm  
**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

PLX51107 is supplied as a solid. A stock solution may be made by dissolving the PLX51107 in the solvent of choice, which should be purged with an inert gas. PLX51107 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

### Description

PLX51107 is a bromodomain and extra terminal domain (BET) family protein inhibitor.<sup>1</sup> It binds to bromodomain 1 (BD1) in bromodomain-containing protein 2 (BRD2), BRD3, BRD4, and BRDT (K<sub>d</sub>s = 1.6, 2.1, 1.7, and 5 nM, respectively), as well as BD2 (K<sub>d</sub>s = 5.9, 6.2, 6.1, and 120 nM, respectively). It also binds to the bromodomains of CBP and p300 (K<sub>d</sub>s = ~100 nM for both). PLX51107 (0.1-10 μM) inhibits CpG-induced proliferation of primary chronic lymphocytic leukemia (CLL) cells. *In vivo*, PLX51107 (0.5-10 mg/kg) inhibits Ba/F3 cell-induced splenomegaly in mice. It also reduces tumor volume and increases survival in YUMM3.3 and D4M3.A mouse syngeneic B-RAF<sup>V600E</sup> mutant melanoma models.<sup>2</sup>

### References

- Ozer, H.G., El-Gamal, D., Powell, B., *et al.* BRD4 profiling identifies critical chronic lymphocytic leukemia oncogenic circuits and reveals sensitivity to PLX51107, a novel structurally distinct BET Inhibitor. *Cancer Discov.* **8**(4), 458-477 (2018).
- Erkes, D.A., Field, C.O., Capparelli, C., *et al.* The next-generation BET inhibitor, PLX51107, delays melanoma growth in a CD8-mediated manner. *Pigm. Cell Melanoma R.* **32**(5), 687-696 (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, 12/12/2022

#### 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