

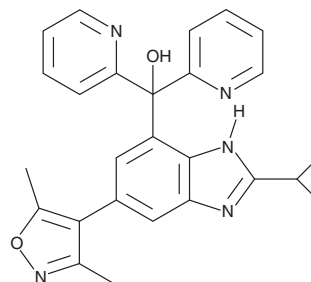
# PRODUCT INFORMATION



## Alobresib

Item No. 30726

**CAS Registry No.:** 1637771-14-2  
**Formal Name:** 2-cyclopropyl-5-(3,5-dimethyl-4-isoxazolyl)- $\alpha,\alpha$ -di-2-pyridinyl-1H-benzimidazole-7-methanol  
**Synonym:** GS-5829  
**MF:** C<sub>26</sub>H<sub>23</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 437.5  
**Purity:**  $\geq 95\%$   
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Alobresib is supplied as a solid. A stock solution may be made by dissolving the alobresib in the solvent of choice, which should be purged with an inert gas. Alobresib is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

Alobresib is an inhibitor of bromodomain and extra-terminal domain (BET) proteins.<sup>1,2</sup> It inhibits proliferation of MEC-1 chronic lymphocytic leukemia cells (CLL; IC<sub>50</sub> = 46.4 nM).<sup>1</sup> Alobresib (400 nM) decreases the levels of Blk and Myc and increases the levels of I $\kappa$ B $\alpha$  in patient-derived CLL cells co-cultured with nurse-like cells (NLC). It inhibits proliferation of USPC-ARK-1 and USPC-ARK-2 uterine serous carcinoma (USC) cells with IC<sub>50</sub> values of 31 and 27 nM, respectively.<sup>2</sup> Alobresib (10 and 20 mg/kg, p.o.) reduces tumor growth in a USPC-ARK-2 USC mouse xenograft model.

### References

1. Kim, E., Hacken, E.T., Sivina, M., *et al.* The BET inhibitor GS-5829 targets chronic lymphocytic leukemia cells and their supportive microenvironment. *Leukemia* **34(6)**, 1588-1598 (2020).
2. Bonazzoli, E., Predolini, F., Cocco, E., *et al.* Inhibition of BET bromodomain proteins with GS-5829 and GS-626510 in uterine serous carcinoma, a biologically aggressive variant of endometrial cancer. *Clin. Cancer Res.* **24(19)**, 4845-4853 (2018).

#### 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, 04/17/2024

#### 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