

# PRODUCT INFORMATION



## Valemetostat

Item No. 31674

**CAS Registry No.:** 1809336-39-7  
**Formal Name:** (2R)-7-chloro-N-[(1,2-dihydro-4,6-dimethyl-2-oxo-3-pyridinyl)methyl]-2-[trans-4-(dimethylamino)cyclohexyl]-2,4-dimethyl-1,3-benzodioxole-5-carboxamide

**Synonyms:** DS-3201b, (R)-OR-S2

**MF:** C<sub>26</sub>H<sub>34</sub>ClN<sub>3</sub>O<sub>4</sub>

**FW:** 488.0

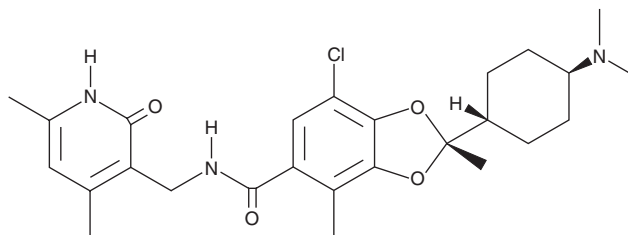
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 216 nm

**Supplied as:** A crystalline 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

Valemetostat is supplied as a crystalline solid. A stock solution may be made by dissolving the valemetostat in the solvent of choice, which should be purged with an inert gas. Valemetostat is soluble in DMSO.

### Description

Valemetostat is a dual inhibitor of enhancer of zeste homolog 1 (EZH1) and EZH2.<sup>1</sup> It inhibits polycomb repressive complex 2 (PRC2) containing EZH1 or EZH2 as the catalytic subunit in cell-free assays (IC<sub>50</sub>s = 8.4 and 2.5 nM, respectively). Valemetostat inhibits trimethylation of histone H3 lysine 27 in HCT116 cells (IC<sub>50</sub> = 0.44 nM). It inhibits the growth of Karpas-422 diffuse large B cell lymphoma (DLBCL) cells (GI<sub>50</sub> = 4.8 nM).

### Reference

1. Honma, D., Kanno, O., Watanabe, J., *et al.* Novel orally bioavailable EZH1/2 dual inhibitors with greater antitumor efficacy than an EZH2 selective inhibitor. *Cancer Sci.* **108**(10), 2069-2078 (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, 10/06/2022

#### 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