

# Tazemetostat

## EZH2 inhibitor

Tazemetostat is a potent and selective SAM-competitive inhibitor of the lysine methyltransferase EZH2. Tazemetostat displayed strong antiproliferative effects against SMARCB1-deleted malignant rhabdoid tumor (MRT) cell lines *in vitro*. This antitumor activity was also observed in SMARTCB1 mutant mouse xenografts. It displayed potent antitumor activity in various cancer models including non-Hodgkins lymphoma, pediatric glioma, small-cell carcinoma of the ovary, and synovial sarcomas. Tazemetostat has also been shown to control inflammatory genes by modulating IRF1, IRF8, and STAT1 levels suggesting therapeutic potential for the treatment of neuroinflammatory diseases associated with microglial activation.

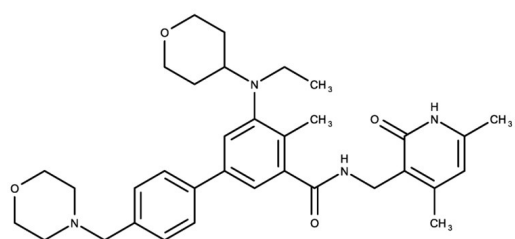
## Ordering Information

[Order Online »](#)

ENZ-CHM495-0025	25mg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-((4,6-Dimethyl-2-oxo-1,2-dihydropyridin-3-yl)methyl)-5-(ethyl(tetrahydro-2H-pyran-4-yl)amino)-4-methyl-4'-(morpholinomethyl)-[1,1'-biphenyl]-3-carboxamide

**Appearance** Off-white solid.

**CAS** 1403254-99-8

**Couple Target** Histone methyltransferase

**Couple Type** Inhibitor

**Formula**  $C_{34}H_{44}N_4O_4$

**Identity** Determined by NMR.

**MW** 572.75

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (>25 mg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)