

PRODUCT INFORMATION



Methylstat (hydrate)

Item No. 11091

Formal Name: 4-[hydroxy[4-[[[4-[[[(1-naphthalenylamino) carbonyl]oxy]methyl]phenyl]methyl]amino]butyl] amino]-4-oxo-2E-butenoic acid, methyl ester, hydrate

MF: C₂₈H₃₁N₃O₆ • XH₂O

FW: 505.6

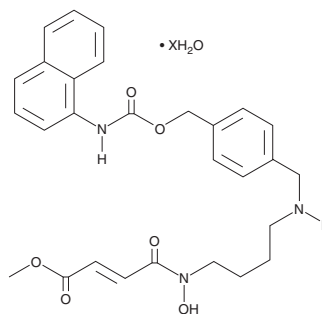
Purity: ≥98%

UV/Vis.: λ_{max}: 222 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

Methylstat (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the methylstat (hydrate) in the solvent of choice, which should be purged with an inert gas. Methylstat (hydrate) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of methylstat (hydrate) in these solvents is approximately 25 mg/ml.

Methylstat (hydrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methylstat (hydrate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Methylstat (hydrate) has a solubility of approximately 0.02 mg/ml in a 1:40 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Methylstat is a methyl ester prodrug of a Jumonji C domain-containing histone demethylase (JMJD) inhibitor that has favorable cell permeability.¹ The free acid of methylstat inhibits JMJD2A, JMJD2C, JMJD2E, PHF8, and JMJD3 with IC₅₀ values of approximately 4.3, 3.4, 5.9, 10, and 43 μM, respectively, in an *in vitro* assay. Methylstat inhibits growth of the JMJD2C-sensitive esophageal cancer cell line KYSE150 with a GI₅₀ of 5.1 μM, whereas the free acid of methylstat did not inhibit cell growth up to 100 μM. Methylstat induces histone hypermethylation at multiple sites in a concentration-dependent manner (EC₅₀ for H3K4me3 and H3K9me3 = 10.3 and 8.6 μM in KYSE150 cells and 6.7 and 6.3 μM in MCF-7 cells, respectively).

Reference

1. Luo, X., Liu, Y., Kubicek, S., *et al.* A selective inhibitor and probe of the cellular functions of Jumonji C domain-containing histone demethylases. *J. Am. Chem. Soc.* **133**(24), 9451-9456 (2011).

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, 12/16/2022

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