

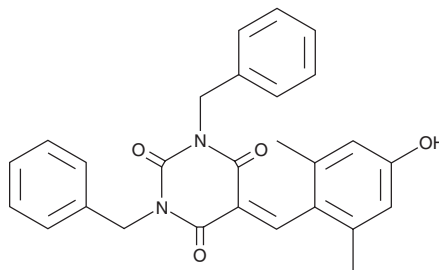
PRODUCT INFORMATION



EML 425

Item No. 30685

CAS Registry No.: 1675821-32-5
Formal Name: 5-[(4-hydroxy-2,6-dimethylphenyl)methylene]-1,3-bis(phenylmethyl)-2,4,6(1H,3H,5H)-pyrimidinetrione
MF: C₂₇H₂₄N₂O₄
FW: 440.5
Purity: ≥98%
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

EML 425 is supplied as a solid. A stock solution may be made by dissolving the EML 425 in the solvent of choice, which should be purged with an inert gas. EML 425 is soluble in the organic solvent DMSO at a concentration of approximately 40 mg/ml.

Description

EML 425 is a non-competitive inhibitor of the lysine acetyltransferase type 3 (KAT3) histone acetyltransferase p300 and CREB-binding protein (CBP; IC₅₀s = 2.9 and 1.1 μM, respectively).¹ It prevents glucose-induced cataract formation in isolated rat lens when used at a concentration of 200 μM.²

References

1. Milite, C., Feoli, A., Sasaki, K., *et al.* A novel cell-permeable, selective, and noncompetitive inhibitor of KAT3 histone acetyltransferases from a combined molecular pruning/classical isosterism approach. *J. Med. Chem.* **58**(6), 2779-2798 (2015).
2. Kanada, F., Takamura, Y., Miyake, S., *et al.* Histone acetyltransferase and Polo-like kinase 3 inhibitors prevent rat galactose-induced cataract. *Sci. Rep.* **9**(1), 20085 (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, 11/11/2022

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