

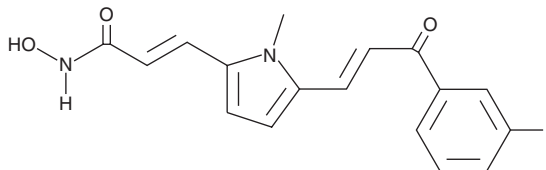
PRODUCT INFORMATION



MC 1568

Item No. 16265

CAS Registry No.: 852475-26-4
Formal Name: (2E)-3-[5-[(1E)-3-(3-fluorophenyl)-3-oxo-1-propen-1-yl]-1-methyl-1H-pyrrol-2-yl]-N-hydroxy-2-propenamamide
MF: C₁₇H₁₅FN₂O₃
FW: 314.3
Purity: ≥95%
UV/Vis.: λ_{max}: 435 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

MC 1568 is supplied as a crystalline solid. A stock solution may be made by dissolving the MC 1568 in the solvent of choice. MC 1568 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of MC 1568 in these solvents is approximately 1 and 0.5 mg/ml, respectively.

MC 1568 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MC 1568 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MC 1568 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

While class I HDACs are localized predominantly within the nucleus, class II HDACs shuttle into and out of the nucleus in response to intracellular signaling.¹ Class IIa HDACs, which includes HDAC4, 5, 7, and 9, commonly act as corepressors and play diverse roles in cell biology.² MC 1568 is a selective inhibitor of class IIa HDACs, with greater than 170-fold selectivity over class I HDACs, including HDAC1.³⁻⁵ It has been used in cells (1-10 μM) and in mice to elucidate the roles of class IIa HDACs in cell proliferation and differentiation, apoptosis, myogenesis, adipogenesis, and fibrosis.⁵⁻¹⁰

References

1. de Ruijter, A.J., van Gennip, A.H., Caron, H.N., et al. *Biochem. J.* **370**(Pt 3), 737-749 (2003).
2. Yang, X.-J. and Grégoire, S. *Mol. Cell. Biol.* **25**(8), 2873-2884 (2005).
3. Mai, A., Massa, S., Pezzi, R., et al. *J. Med. Chem.* **48**(9), 3344-3353 (2005).
4. Mai, A., Jelacic, K., Rotili, D., et al. *Mol. Pharmacol.* **72**(5), 1111-1123 (2007).
5. Nebbioso, A., Manzo, F., Miceli, M., et al. *EMBO Rep.* **10**(7), 776-782 (2009).
6. Duong, V., Bret, C., Altucci, L., et al. *Mol. Cancer Res.* **6**(12), 1908-1919 (2008).
7. Nebbioso, A., Dell'Aversana, C., Bugge, A., et al. *J. Mol. Endocrinol.* **45**(4), 219-228 (2010).
8. Wang, G., He, J., Zhao, J., et al. *PLoS One* **7**(12), 1-10 (2012).
9. Mannaerts, I., Eysackers, N., Onyema, O.O., et al. *PLoS One* **8**(1), 1-9 (2013).
10. Spallotta, F., Tardivo, S., Nanni, S., et al. *J. Biol. Chem.* **288**(32), 22915-22929 (2013).

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, 11/16/2022

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