

MS-275 (Entinostat)

Epigenetic Enzyme Inhibitor

Catalog # H83-904E

Lot # P3320-24

CAS # 209783-80-2

Product Description

Molecular Formula: C₂₁H₂₀N₄O₃

Appearance: Pale yellow powder

Melting Point: 159-160°C

Molecular Weight: 376.41

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (25 mg/ml, warm)

Alias

Pyridin-3-ylmethyl 4-((2-aminophenyl)carbamoyl)benzylcarbamate; Entinostat

Specific Activity

HDAC Inhibitor

Storage and Stability

Store desiccated as supplied at -20°C for up to 1 year.

Store solutions at -20°C for up to 1 week.

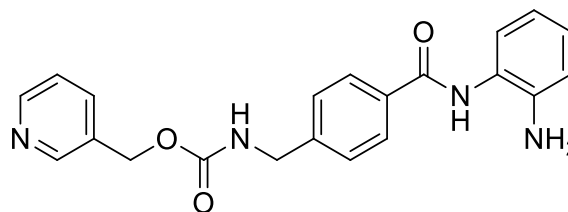
Scientific Background

A potent and selective non-hydroxamate HDAC inhibitor. Considered to be a class I selective inhibitor but displays some selectivity amongst class I members: HDAC1 (IC₅₀ = 0.3 μM); HDAC3 (IC₅₀ = 8 μM) with no inhibition of HDAC8 (at >100 μM)¹. Induces hyperacetylation of nuclear histones in tumor cells resulting in inhibition of proliferation of a variety of human tumor cell lines². Protects p53-deficient mice against ischemic injury³. Suppresses the p38 MAP kinase pathway in rheumatoid arthritic synovial fibroblasts⁴. Active in vivo.

References

1. E Hu et al. J. Pharmacol. Exp. Ther. 2003 307:720
2. A Saito et al. Proc. Natl. Acad. Sci. USA 1999 96:4592
3. SP Murphy et al. J. Neurochem. 2013 Oct 21 epub
4. QY Choo et al. Molecules 2013 18:14085

Molecular Structure



MS-275 (Entinostat)

Epigenetic Enzyme Inhibitor

Catalog #	H83-904E
CAS #	209783-80-2
Lot #	P3320-24
Purity	>98% (TLC)
Format	Pale yellow powder
Solubilization	May be dissolved in DMSO (25 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 1 week.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.