

Apicidin

Catalog Number P011-1MG

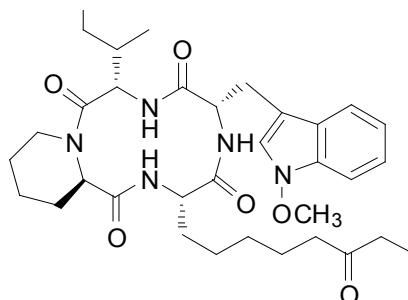
Catalog Number P011-5MG

FEATURES

- Potent HDAC inhibitor
- Blocks cell cycle progression at G1
- Induces activation of HIV-1 promoter

INTRODUCTION

Potent inhibitor of histone deacetylase (HDAC) activity and proliferation. Induces cell cycle arrest at the G1 phase. At 100nM it induces a long lasting hyperacetylation of histone H4 while that induced by trichostatin is transient. Stimulates apoptosis. Apoptosis is induced via induction of Fas/Fas ligand. Displays potent antiangiogenic effects and dramatically decreases HIF-1 α protein levels and transcriptional activity in human and mouse tumor cell lines. Apicidin is a fungal toxin that has broad spectrum activity against Apicomplexan parasites



FORM: White solid

MOLECULAR WEIGHT: 623.78

STORAGE: -20°C

FORMULA: C₃₄H₄₉N₅O₆

CAS NUMBER: 183506-66-3

OTHER NAMES: Cyclo[(2S)-2-amino-8-oxodecanoyl-1-methoxy-L-tryptophyl-L-isoleucyl-(2R)-2-piperidinexcarbonyl]

USES: Soluble in ethanol and DMSO to 10 mg/mL

REFERENCES:

Cao, H., and Stamatoyannopoulos, G. Histone deacetylase inhibitor FK228 is a potent inducer of human fetal hemoglobin. Am. J. Hematol. 81, 981-983 (2006).

Mazitschek, R., Patel, V., Wirth, DF., et al. Development of a fluorescence polarization based assay for histone deacetylase ligand discovery. Bioorg. Med. Chem. Lett. 18, 2809-2812 (2008).

Abbas, HK., Gronwald, JW., Plaisance, KL., et al. Histone deacetylase activity and phytotoxic effects following exposure of duckweed (*Lemna pausicostata* L.) to apicidin and HC-toxin. Phytopathology 91:12, 1141-1148 (2001).