
APHA Compound 8

(Catalog No. M45003)

Alternative Names

3-(1-Methyl-4-phenylacetyl-1H-2-pyrrolyl)-N-hydroxy-2-propenamide

Description

APHA Compound 8 is a potent novel hydroxamate HDAC inhibitor, is in the same structural class as SAHA, an investigational drug standard currently in clinical trial. The IC₅₀ for mouse HDAC1 is 0.5 μM. In a separate study, significant inhibition was found in the proliferation of MEL cells at 24 μM and differentiation, assayed as the accumulation of hemoglobin, at 5 μM.

CAS Number

676599-90-9

Chemical Formula

C₁₆H₁₆N₂O₃

Molecular Weight

284.31

Size

1 mg

Purity

>98%

Solubility

Soluble in DMSO: >10 mg/mL

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
APHA Compound 8	1 mg	M45003 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.