

## Vorinostat (SAHA)

Epigenetic Enzyme Inhibitor

### Catalog # H83-904M

Lot # P3320-43

CAS # 149647-78-9

### Product Description

Molecular Formula: C<sub>14</sub>H<sub>20</sub>N<sub>2</sub>O<sub>3</sub>

Appearance: White powder

Molecular Weight: 264.3

Purity: >98%; NMR (Conforms)

Solubilization: May be dissolved in DMSO (50 mg/ml); or Ethanol (2 mg/ml)

### Alias

N-Hydroxy-N'-phenyloctanediamide

### Specific Activity

HDAC Inhibitor

### Storage and Stability

Store desiccated as supplied at ambient temperature for up to 5 years. Store solutions at -20°C for up to 1 month.

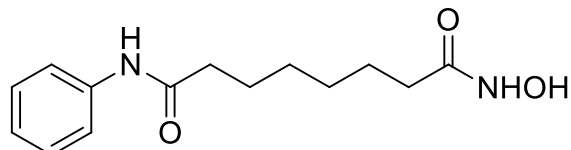
### Scientific Background

Potent and selective histone deacetylase inhibitor<sup>1</sup>.  
Induces apoptosis in a variety of tumor cell lines<sup>2</sup>.  
Synergizes with kinase inhibitors to kill CNS tumor cells<sup>3</sup>.  
Prevents diabetes in a mouse model by immunoregulation and β-cell protection<sup>4</sup>. Cell permeable.

### References

1. JA Vrana et al. Oncogene 1999 18:7016
2. LM Butler et al. Proc. Natl. Acad. Sci. USA 2002 99:11700
3. Y Tang et al. Cancer Biol. Ther. 2012 13:567
4. DP Christensen et al. Proc. Natl. Acad. Sci. USA 2014 111:1055

### Molecular Structure



## Vorinostat (SAHA)

Epigenetic Enzyme Inhibitor r

|                    |  |
|--------------------|--|
| Catalog #          | H83-904M   |
| CAS #              | 149647-78-9  |
| Lot #              | P3320-43   |
| Purity             | >98%   |
| Format             | White powder   |
| Solubilization     | May be dissolved in DMSO (50 mg/ml); or Ethanol (2 mg/ml)  |
| Storage & Shipping | Store desiccated as supplied at ambient temperature for up to 5 years. Store solutions at -20°C for up to 1 month. |

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

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