

## Apicidin

**Catalog No:** 14040, 14041

**Format:** 5 mg, 1 mg

**Chemical Properties:**

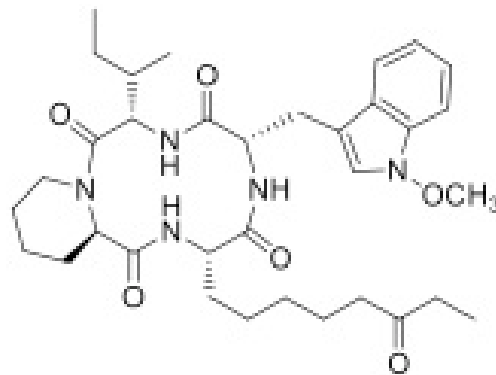
MW = 623.8

C<sub>34</sub>H<sub>49</sub>N<sub>5</sub>O<sub>6</sub>

CAS 183506-66-3

Physical Properties: White solid

**Names:** Apicidin, cyclo[(2S)-2-amino-8-oxodecanoyl-1-methoxy-L-tryptophyl-L-isoleucyl-(2R)-2-piperidinexcarbonyl]



**Pharmacology:** Cyclopeptide inhibitor of histone deacetylases (HDACs) (ref 1). Inhibits HeLa cell proliferation (IC<sub>50</sub> = 50-100 nM) and induces reversible transcriptional activation of p21 (WAF1) (ref 2). Decreases HIF1α protein levels and transcriptional activity in human and mouse tumor cell lines (ref 3). Cell permeable.

**Solubilization:** May be dissolved in DMSO (10 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

≥98% (HPLC); NMR (Conforms)

**References:**

1. SJ Darkin-Rattray *et al. Proc. Natl. Acad. Sci. USA* 1996, 93:13143
2. JW Han *et al. Cancer Res.* 2000, 60:6068
3. SH Kim *et al. Oncol. Rep.* 2007, 17:647

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.