

HPA (Hexyl-4-pentynoic acid)

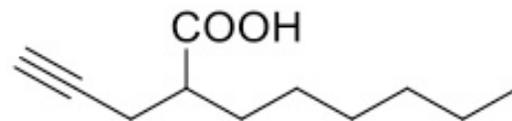
Catalog No: 14034, 14035**Format:** 50 mg, 10 mg**Chemical Properties:**

MW = 182.3

 $C_{11}H_{18}O_2$

CAS 96017-59-3

Physical Properties: Colorless oil

Names: HPA (Hexyl-4-pentynoic acid), 2-(2-propynyl)octanoic acid (racemic)

Pharmacology: Valproic acid analog. Histone deacetylase (HDAC) inhibitor ($IC_{50} = 13 \text{ M}$)(ref 1). More potent and robust than valproic acid at inducing histone hyperacetylation (600% at 50 M), HSP70 induction and protection against glutamate excitotoxicity in cultured neurons (ref 2). Cell permeable.

Solubilization: May be dissolved in DMSO (25 mg/ml); Ethanol (25 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>97% (TLC); NMR (Conforms)

References:

1. D Eikel *et al. Chem. Res. Toxicol.* 2006, 19:272
2. Y Leng *et al. Neurosci. Lett.* 2010, 476:127

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.