

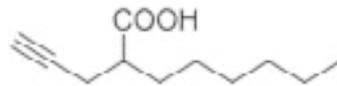
# Hexyl-4-pentynoic Acid (HPA)

Catalog Number P008-10MG

Catalog Number P008-50MG

## FEATURES

- Induces histone hyperacetylation
- 30 times lower IC50 than VPA
- Protects against glutamate-induced excitotoxicity



ARBOR  
ASSAYS

## INTRODUCTION

HPA causes histone hyperacetylation and protect against glutamate-induced excitotoxicity in cultured neurons. HPA is one of the most effective valproic acid derivatives in terms of teratogenic potential in ice, hyperacetylation of core histone 4 in treated F9 Cells, and one of the lowest concentrations of half-maximum effect in HDAC enzyme inhibition assays. Inhibits HDAC activity with an IC50 of 13  $\mu$ M (VPA= 398  $\mu$ M).

<b>FORM:</b>	Colorless Oil
<b>MOLECULAR WEIGHT:</b>	182.26
<b>STORAGE:</b>	Room temperature
<b>FORMULA:</b>	$C_{11}H_{18}O_2$
<b>CAS NUMBER:</b>	96017-59-3
<b>OTHER NAMES:</b>	$\pm$ 2-Hexyl-4-pentynoic acid, 2-(2-propynyl)octanoic acid (racemic), HPA
<b>USES:</b>	Soluble to 25 mg/mL in DMSO and Ethanol.

## REFERENCES:

Eikel D, Lampen A, Nau H. Teratogenic effects mediated by inhibition of histone deacetylases: evidence from quantitative structure activity relationships of 20 valproic acid derivatives. *Chem.Res. Toxicol.* 19:272–278 (2006).

Leng Y, Marinova Z, Reis-Fernandes MA, Nau H, Chuang DM. Potent neuroprotective effects of novel structural derivatives of valproic acid: potential roles of HDAC inhibition and HSP70 induction. *Neurosci Lett.* 476:3, 127-32. (2010).