Valproic Acid, sodium salt

Catalog No: 14021 Format: 5 g

Chemical Properties:

$$\begin{split} &\mathsf{MW} = 166.2\\ &\mathsf{C}_8\mathsf{H}_{15}\mathsf{O}_2\mathsf{Na}\\ &\mathsf{CAS}\ 1069\text{-}66\text{-}5\\ &\mathsf{Physical Properties:}\ &\mathsf{White powder}\\ &\mathsf{Names:}\ &\mathsf{Valproic \ acid, \ sodium \ salt, \ Sodium \ valproate, \ 2\text{-}Propylpentanoic \ acid}\\ &\mathsf{Na} \end{split}$$

Pharmacology: Histone deacetylase inhibitor ($IC_{50} = 400 \ \mu$ M). Demonstrates neuroprotective, anti-cancer and anti-inflammatory activity. Inhibits A β production and neuritic plaque formation in Alzheimer's mouse models. Clinically useful anti-convulsant.

Solubilization: May be dissolved in water (50 mg/ml)

Fluorescent Properties: N/A

Quality Control: >98% (TLC); NMR (Conforms)

References:

- 1. CJ Phiel et al. J. Biol. Chem. 2001 276:36734
- 2. HJ Kim et al. J. Pharmacol. Exp. Ther. 2007, 321:892
- 3. H Qing et al. J. Exp. Med. 2008, 205:2781

Storage and Guarantee: Store desiccated as supplied at 4°C 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.



