## Sodium phenylbutyrate



**Catalog No: 14032, 14033** Format: 5 g, 1 g

## **Chemical Properties:**

$$\begin{split} \text{MW} &= 186.2\\ \text{C}_{10}\text{H}_{11}\text{O}_2\text{ Na}\\ \text{CAS 1716-12-7}\\ \text{Physical Properties: White powder}\\ \textbf{Names: Phenylbutyrate Na, 4-Phenylbutyric acid sodium salt, 4-PB, PBNa} \end{split}$$

**Pharmacology:** Histone deacetylase (HDAC) inhibitor. Inhibits proliferation, migration and invasion of a number of cancer cell lines. Induces differentiation, and apoptosis. Active *in vivo*.

Solubilization: May be dissolved in DMSO (4 mg/ml); Water (25 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC; 5% CH<sub>3</sub>OH in CH<sub>2</sub>Cl<sub>2</sub>, R<sub>f</sub> = 0.36); NMR (Conforms)

References:

1. HH Engelhard et al. J. Neuro-Oncol. 1998, 37:97

2. IB Appelskog et al. Int. J. Oncol. 2004, 24:1419

**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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

