

Phenylbutyrate Na

Catalog No: 14033

Format: 1 g

Chemical Properties:

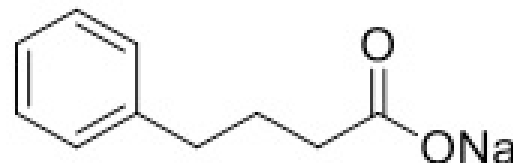
MW = 186.2

C₁₀H₁₁O₂ Na

CAS 1716-12-7

Physical Properties: White powder

Names: Phenylbutyrate Na, 4-Phenylbutyric acid sodium salt, 4-PB, PBNa



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% CH₃OH in CH₂Cl₂, R_f = 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 guaranteed for 6 months from date of arrival.