

Phenylbutyrate NA

Epigenetic Enzyme Inhibitor

Catalog # H83-904H

Lot # P3320-28
CAS # 1716-12-7

Product Description

Molecular Formula: C₁₀H₁₁NaO₂
Appearance: white powder
Molecular Weight: 186.2
Purity: >98% (TLC; 5% CH₃OH in CH₂Cl₂, R_f=0.36); NMR (Conforms)
Solubilization: May be dissolved in DMSO (4 mg/ml); or water (25 mg/ml)

Alias

4-Phenylbutyric acid, sodium salt

Specific Activity

HDAC Inhibitor

Storage and Stability

Store as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 months.

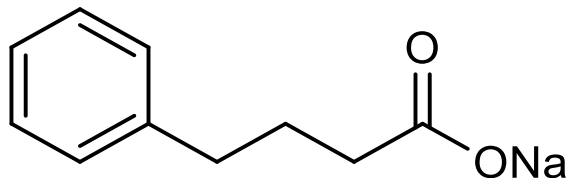
Scientific Background

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

References

1. HH Engelhard et al. J. Neuro-Oncol. 1998 37:97
2. IB Appelskog et al. Int. J. Oncol. 2004 24:1419

Molecular Structure



Phenylbutyrate NA

Epigenetic Enzyme Inhibitor

Catalog #	H83-904H
CAS #	1716-12-7
Lot #	P3320-28
Purity	white powder
Format	white powder
Solubilization	May be dissolved in DMSO (4 mg/ml); or water (25 mg/ml)
Storage & Shipping	Store as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.