

# Phenylbutyrate sodium

## HDAC inhibitor

Weak histone deacetylase (HDAC) inhibitor. Induces differentiation, growth arrest and apoptosis in a number of cell lines. Protects against cerebral ischemic injury and displays neuroprotective effect in a mouse model of Huntington's disease. Anti-neoplastic agent and transcriptional regulator. Also acts as an inducer of tumor cytostasis and differentiation. Shows chemical chaperone activity.

**Citations:** 11

[View Online »](#)

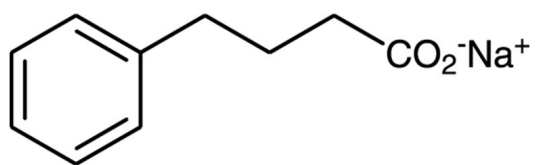
## Ordering Information

[Order Online »](#)

BML-EI320-0001	1g
----------------	----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Hygroscopic.
Long Term Storage	Ambient
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Sodium 4-phenylbutyrate, 4-Phenylbutyric acid sodium salt, SBP, 4PBA, PBNa
Appearance	White to yellowish solid.
CAS	1716-12-7
Couple Target	HDAC
Couple Type	Inhibitor
Formula	$C_{10}H_{11}O_2 \cdot Na$
MW	186.2
Purity	≥98% (UHPLC)
Solubility	Soluble in DMSO or water (25mg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)