

# Phenylbutyrate Sodium

HDAC inhibitor.

An aromatic fatty acid that induces cytostasis, differentiation and apoptosis in primary myeloid leukemic cells [1], human prostate cancer cell lines [2] and human colon carcinoma cells [3]. It also has chemopreventive effects against chemically induced colon carcinogenesis in experimental animals [4].

## Ordering Information

[Order Online »](#)

LKT-P2815-G005	5g
LKT-P2815-G025	25g

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      Ambient

Shipping      Ambient Temperature

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      Sodium 4-phenylbutyrate, 4-Phenylbutyric acid . Na, 4PBA

Appearance      White crystalline powder.

CAS      1716-12-7

Formula       $C_{10}H_{11}O_2 \cdot Na$

MW      186.18

Purity       $\geq 98\%$

Solubility      Soluble in water, ethanol, DMSO.



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)