

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-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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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