

# Sodium butyrate

**HDAC inhibitor. Apoptosis inducer.**

Short-chain fatty acid shown to be an inhibitor of histone deacetylases (HDACs). Induces apoptosis. Restores contextual memory in transgenic mouse model of Alzheimer's. Induces pancreatic progenitor formation. Directs differentiation of mouse embryonic stem cells to hepatocytes with Cytokine Activin A. Sodium butyrate interferes with signal transduction processes, as well as, inhibits c-myc splicing.

**Citations:** 7

[View Online »](#)

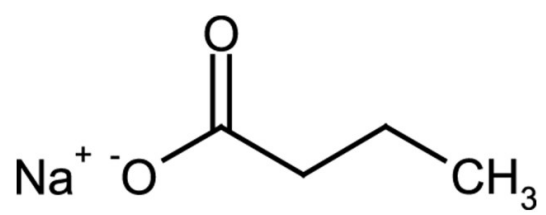
## Ordering Information

[Order Online »](#)

|                  |    |
|------------------|----|
| ALX-270-301-G001 | 1g |
|------------------|----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                  |
|-------------------|------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Hygroscopic.                                                     |
| Long Term Storage | -20°C                                                            |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|               |                              |
|---------------|------------------------------|
| Appearance    | White to off-white solid.    |
| CAS           | 156-54-7                     |
| Couple Target | HDAC                         |
| Couple Type   | Inhibitor                    |
| Formula       | $C_4H_7O_2Na$                |
| MW            | 110.1                        |
| Purity        | ≥95% (HPLC)                  |
| RTECS         | ET6400000                    |
| Solubility    | Soluble in water (100mg/ml). |

Last modified: May 29, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)