
Sodium Butyrate

(Catalog No. M41001)

Alternative Names

Butyric acid, Sodium salt, NaB

Description

A short chain fatty acid shown to be an inhibitor of histone deacetylase (HDAC). Induce apoptosis in colon carcinoma cell lines in a p53-independent pathway. Inhibits c-myc splicing and interferes with signal transduction processes, including the release of Ca²⁺ from intracellular stores. Downregulates the activity of serine/threonine protease kinase casein kinase II in HT29 cells.

CAS Number

156-54-7

Chemical Formula

C₄H₇O₂.Na

Molecular Weight

110.1

Size

1 g

Purity

≥99% by HPLC

Solubility

H₂O (100 mM)

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Sodium Butyrate	1 g	M41001 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.