
Valproic Acid

(Catalog No. M41002)

Alternative Names

Valproic acid sodium, Sodium Valproate, VPA

Description

A cell-permeable histone deacetylase 1 (HDAC1) inhibitor (IC₅₀ = 400 µM) that exhibits anticancer, anti-inflammatory and neuroprotective properties. Recently it has been reported that Valproic acid enables reprogramming of primary human fibroblasts with only two factors, Oct4 and Sox2, without the need for the oncogenes c-Myc or Klf4.

CAS Number

1069-66-5

Chemical Formula

C₈H₁₅O₂.Na

Molecular Weight

166.19

Size

500 mg

Purity

≥98% by titration

Solubility

H₂O (100 mM)

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Valproic Acid	500 mg	M41002-2

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