

# Sodium Valproate

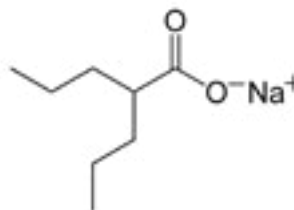
Catalog Number P001-5GM



ARBOR  
ASSAYS

## FEATURES

- Histone deacetylase inhibitor
- Anti-cancer, -inflammatory, neuroprotective effects
- Decreases A $\beta$  production



## INTRODUCTION

An anticonvulsant used in the treatment of epilepsy, anorexia nervosa, panic attack, anxiety disorder, post traumatic stress disorder, migraine and bipolar disorder. Valproate is also a histone deacetylase inhibitor (IC<sub>50</sub> = 400  $\mu$ M) that exhibits anticancer, anti-inflammatory and neuroprotective effects. Displays anticonvulsive activity via an increase in GABA levels and decreases A $\beta$  production in animal models of Alzheimer's disease. Also attenuates NMDA-mediated excitation, blocks voltage-gated Na<sup>+</sup> channels and modulates firing of neurons. Enables induction of pluripotent stem cells from somatic cells by Oct4 and Sox2.

Valproic acid also inhibits glycogen synthase kinase 3 (GSK3) and depletes cellular inositol-1,4,5-trisphosphate (1,4,5-IP<sub>3</sub>). Valproic acid shows promise in combination therapy for cancer and in treating Alzheimer's disease. Valproic acid (1 mM) also has pronounced effects on stem cell differentiation and self-renewal. Inhibits Class I HDACs with an IC<sub>50</sub> value of ~2 mM.

<b>FORM:</b>	White Powder
<b>MOLECULAR WEIGHT:</b>	166.2
<b>STORAGE:</b>	4°C, desiccated
<b>FORMULA:</b>	C <sub>8</sub> H <sub>15</sub> O <sub>2</sub> Na
<b>CAS NUMBER:</b>	1069-66-5
<b>OTHER NAMES:</b>	2-propyl-pentanoic acid, monosodium salt, 2-propylvaleric acid, monosodium salt, sodium 2-propylpentanoate
<b>USES:</b>	Soluble to 100 mM in water and 5 mg/ml in DMSO and DMF

## REFERENCES:

- Jung, G., Yoon, J., Moon, B., et al. Valproic acid induces differentiation and inhibition of proliferation in neural progenitor cells via the beta-catenin-Ras-ERK-p21Cip/WAF1 pathway. *BMC Cell Biol* 9(66) (2008).
- Bug, G., Gül, H., Schwarz, K., et al. Valproic acid stimulates proliferation and self-renewal of hematopoietic stem cells. *Cancer Res* 65(7) 2537-2541 (2005).
- Göttlicher, M., Minucci, S., Zhu, P., et al. Valproic acid defines a novel class of HDAC inhibitors inducing differentiation of transformed cells. *EMBO J* 20(24) 6969-6978 (2001).

**FOR RESEARCH USE ONLY**

info@ArborAssays.com • www.ArborAssays.com

10Aug12