

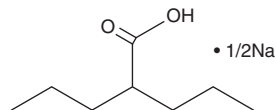
# PRODUCT INFORMATION



## Divalproex (sodium salt)

Item No. 22960

**CAS Registry No.:** 76584-70-8  
**Formal Name:** 2-propyl-pentanoic acid, sodium salt (2:1)  
**MF:**  $C_8H_{16}O_2 \cdot 1/2Na$   
**FW:** 155.7  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Divalproex (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the divalproex (sodium salt) in the solvent of choice. Divalproex (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of divalproex (sodium salt) in ethanol is approximately 30 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of divalproex (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of divalproex (sodium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Divalproex is a prodrug form of valproic acid (Item Nos. 35739 | 13033) that contains valproic acid and valproate sodium.<sup>1</sup> Formulations containing divalproex exhibit delayed gastrointestinal absorption and are converted to valproic acid in the intestine, which reduces gastric irritation and nervous system side effects. Formulations containing divalproex are widely used for the treatment of seizures and bipolar disorder.

### Reference

1. McGraw, D. Therapeutic drug monitoring with valproate—why product selection is an important factor. *Mental Health Clinician* 4(1), 31-34 (2014).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/13/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM