

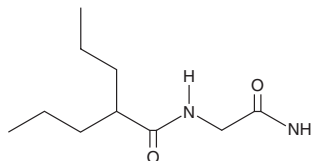
# PRODUCT INFORMATION



## Valrocemide

Item No. 30909

CAS Registry No.: 92262-58-3  
Formal Name: N-(2-amino-2-oxoethyl)-2-propyl-pentanamide  
Synonyms: N-Valproyl glycineamide, TV1901, VGD  
MF:  $C_{10}H_{20}N_2O_2$   
FW: 200.3  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Valrocemide is supplied as a solid. A stock solution may be made by dissolving the valrocemide in the solvent of choice, which should be purged with an inert gas. Valrocemide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of valrocemide in these solvents is approximately 25 mg/ml in ethanol and 30 mg/ml in DMSO and DMF.

Valrocemide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, valrocemide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Valrocemide has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Valrocemide is an anticonvulsant and a derivative of valproic acid (Item No. 13033).<sup>1</sup> *In vivo*, valrocemide protects against seizures induced by maximal electroshock (MES), pentylenetetrazole (PTZ; Item No. 18682), picrotoxin (Item No. 20771), or bicuculline (Item No. 11727) in mice ( $ED_{50}$ s = 151, 132, 275, and 248 mg/kg, respectively). It prevents sound-induced seizures in Frings audiogenic-sensitive mice ( $ED_{50}$  = 52 mg/kg) and MES-induced seizures in rats ( $ED_{50}$  = 73 mg/kg). Valrocemide also prevents seizures in corneally kindled rats ( $ED_{50}$  = 161 mg/kg).

### Reference

1. Isoherranen, N., Woodhead, J.H., White, H.S., *et al.* Anticonvulsant profile of valrocemide (TV1901): A new antiepileptic drug. *Epilepsia* 42(7), 831-836 (2001).

#### 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, 12/09/2022

#### 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