

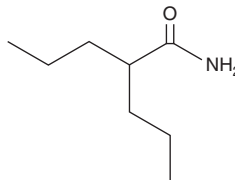
PRODUCT INFORMATION



Valpromide

Item No. 46269

CAS Registry No.: 2430-27-5
Formal Name: 2-propyl-pentanamide
Synonym: VPD
MF: $C_8H_{17}NO$
FW: 143.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Valpromide is supplied as a solid. A stock solution may be made by dissolving the valpromide in the solvent of choice, which should be purged with an inert gas. Valpromide is soluble (≥ 10 mg/ml) in ethanol and DMSO.

Description

Valpromide is a prodrug form of valproic acid (Item Nos. 35739 | 13033).^{1,2} It is metabolized to valproic acid via hydrolysis in the liver. Formulations containing valpromide have been used in the treatment of epilepsy and bipolar disorder.

References

1. Bialer, M. and Rubinstein, A. Pharmacokinetics of valpromide in dogs after various modes of administration. *Biopharm. Drug Dispos.* **5(2)**, 177-183 (1984).
2. Billig, H., Ziv, E., Bar-On, H., et al. The disposition of valpromide in rats and the isolated perfused rat liver. *Drug Metab. Dispos.* **18(2)**, 238-244 (1990).

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, 08/28/2026

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