

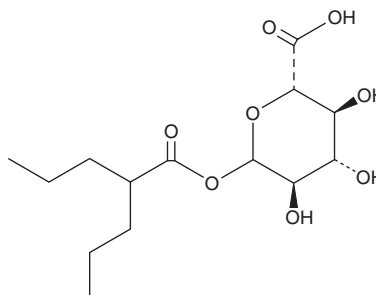
PRODUCT INFORMATION



Valproic Acid Acyl-D-Glucuronide

Item No. 22459

Formal Name: (2S)-3S,4S,5R-trihydroxy-6-((2-propylpentanoyl)oxy)tetrahydro-2H-pyran-2-carboxylic acid
Synonym: Valproic Acid Gluconide
MF: C₁₄H₂₄O₈
FW: 320.3
Purity: ≥95%
UV/Vis.: λ_{max}: 216 nm
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Valproic acid acyl-D-glucuronide is supplied as a lyophilized powder. A stock solution may be made by dissolving the valproic acid acyl-D-glucuronide in the solvent of choice, which should be purged with an inert gas. Valproic acid acyl-D-glucuronide is soluble in organic solvents such as methanol and DMSO.

Description

Valproic acid acyl-D-glucuronide is the major urinary metabolite of valproic acid (Item Nos. 35739 | 13033).¹

Reference

1. Dickinson, R. G., Hooper, W.D., and Eadie, M. pH-dependent rearrangement of the biosynthetic ester glucuronide of valproic acid to beta-glucuronidase-resistant forms. *Drug Metab. Dispos.* **12(2)**, 247-252 (1984).

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, 03/03/2023

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