

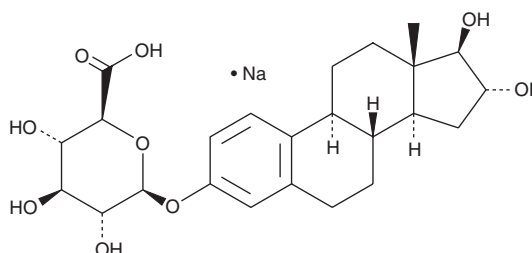
PRODUCT INFORMATION



Estriol 3-β-D-Glucuronide (sodium salt)

Item No. 14827

CAS Registry No.: 15087-06-6
Formal Name: 16α,17β-dihydroxyestra-1,3,5(10)-trien-3-yl, β-D-glucopyranosiduronic acid, monosodium salt
Synonym: Estriol sodium 3-glucuronate
MF: C₂₄H₃₂O₉ • Na
FW: 487.5
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Estriol 3-β-D-glucuronide (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the estriol 3-β-D-glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. Estriol 3-β-D-glucuronide (sodium salt) is slightly soluble in DMSO and methanol.

Description

Estriol 3-β-D-glucuronide is a metabolite of estriol (Item No. 10006484).¹ It is formed from estriol by the UDP-glucuronosyltransferase (UGT) isoform UGT1A10. Estriol 3-β-D-glucuronide binds to basolateral and canalicular liver plasma membranes with K_d values of 85 and 164 μM, respectively.² It competitively inhibits the hydrolysis of 4-methylumbelliferyl-β-D-glucuronide (4Mu-GlcU; Item No. 17203) and is a substrate for hydrolysis by Klotho-human IgG1 Fc protein (KLFc).³

References

1. Sneitz, N., Vanhermo, M., Mosorin, J., *et al.* Regiospecificity and stereospecificity of human UDP-glucuronosyltransferases in the glucuronidation of estriol, 16-epiestriol, 17-epiestriol, and 13-epiestradiol. *Drug Metab. Dispos.* **41**(3), 582-591 (2013).
2. Changchit, A., Durham, S.K., and Vore, M. Characterization of [³H]estradiol-17β-(β-D-glucuronide) binding sites in basolateral and canalicular liver plasma membranes. *Biochem. Pharmacol.* **40**(6), 1219-1225 (1990).
3. Tohyama, O., Imura, A., Freund, J.N., *et al.* Klotho is a novel β-glucuronidase capable of hydrolyzing steroid β-glucuronides. *J. Biol. Chem.* **279**(11), 9777-9784 (2004).

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, 12/05/2022

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