

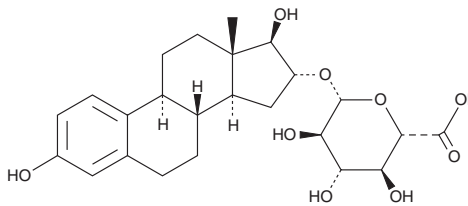
PRODUCT INFORMATION



Estriol 16 α -(β -D-Glucuronide)

Item No. 39140

CAS Registry No.: 1852-50-2
Formal Name: 3,17 β -dihydroxyestra-1,3,5(10)-trien-16 α -yl, β -D-glucoopyranosiduronic acid
Synonyms: Estriol 16-Glucuronidate, NSC 93823
MF: C₂₄H₃₂O₉
FW: 464.5
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Estriol 16 α -(β -D-glucuronide) is supplied as a solid. A stock solution may be made by dissolving the estriol 16 α -(β -D-glucuronide) in the solvent of choice, which should be purged with an inert gas. Estriol 16 α -(β -D-glucuronide) is soluble in ethanol:water (1:1) at a concentration of approximately 10 mg/ml.

Description

Estriol 16 α -(β -D-glucuronide) is a metabolite of the endogenous pregnancy estrogen estriol (Item No. 10006484).^{1,2} Plasma levels of estriol 16 α -(β -D-glucuronide) have been used as biomarkers of placental estriol production during pregnancy.²

References

- DiPietro, D.L. Assay of estriol-16 α -(β -D-glucuronide) in pregnancy urine with a specific antiserum. *Am. J. Obstet. Gynecol.* **125(6)**, 841-845 (1976).
- Ilgisonis, E.V., Shalina, R., Kasum-Zade, N., et al. Metabolomic markers for predicting preeclampsia in the first trimester of pregnancy: A retrospective study. *Molecules* **27(8)**, 2475 (2022).

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, 08/17/2023

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