

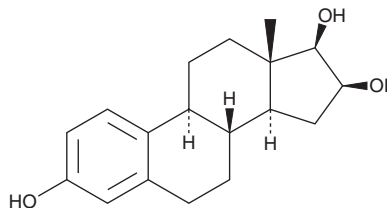
PRODUCT INFORMATION



16-Epiestriol

Item No. 33455

CAS Registry No.: 547-81-9
Formal Name: (16 β ,17 β)-estra-1,3,5(10)-triene-3,16,17-triol
Synonyms: 16-*epi* Estriol, 16-EpiE₃,
16 β -hydroxy-17 α -Estradiol, NSC 26646
MF: C₁₈H₂₄O₃
FW: 288.4
Purity: \geq 90%
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

16-Epiestriol is supplied as a solid. A stock solution may be made by dissolving the 16-epiestriol in the solvent of choice, which should be purged with an inert gas. 16-Epiestriol is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml.

Description

16-Epiestriol is a metabolite of the endogenous estrogen estrone (Item No. 10006485).¹ It is formed from estrone via a 16 β -hydroxy estrone intermediate by reduction of the C-17 ketone. 16-Epiestriol (200 μ g/ml) inhibits the growth of carbapenem-resistant *A. baumannii*.² It inhibits carrageenan-induced paw edema in rats when administered at a dose of 20 mg/kg.³ Unlike hydrocortisone, 16-epiestriol (240 μ g/animal) does not increase plasma or liver glucose levels in adrenalectomized rats.

References

1. Brinton, L.A., Trabert, B., Anderson, G.L., *et al.* Serum estrogens and estrogen metabolites and endometrial cancer risk among postmenopausal women. *Cancer Epidemiol. Biomarkers Prev.* **25(7)**, 1081-1089 (2016).
2. Skariyachan, S., Muddebihalkar, A.G., Badrinath, V., *et al.* Natural epiestriol-16 act as potential lead molecule against prospective molecular targets of multidrug resistant *Acinetobacter baumannii*-Insight from *in silico* modelling and *in vitro* investigations. *Infect. Genet. Evol.* **82**, 104314 (2020).
3. Latman, N.S., Kishore, V., and Bruot, B.C. 16-Epiestriol: An anti-inflammatory steroid without glycogenic activity. *J. Pharm. Sci.* **83(6)**, 874-877 (1994).

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, 02/06/2024

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