

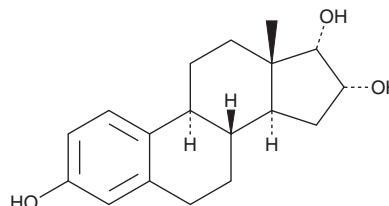
# PRODUCT INFORMATION



## 17-Epiestriol

Item No. 33601

**CAS Registry No.:** 1228-72-4  
**Formal Name:** estra-1,3,5(10)-triene-3,16 $\alpha$ ,17 $\alpha$ -triol  
**Synonyms:** 16 $\alpha$ -hydroxy-17 $\alpha$ -Estradiol,  
16 $\alpha$ ,17 $\alpha$ -Estriol, 17 $\alpha$ -Estriol,  
17-*epi*-Estriol, NSC 84051  
**MF:** C<sub>18</sub>H<sub>24</sub>O<sub>3</sub>  
**FW:** 288.4  
**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

17-Epiestriol is supplied as a solid. A stock solution may be made by dissolving the 17-epiestriol in the solvent of choice, which should be purged with an inert gas. 17-Epiestriol is slightly soluble in methanol.

### Description

17-Epiestriol is a metabolite of estrone (Item No. 10006485).<sup>1</sup> It is formed from estrone via a 16 $\alpha$ -hydroxy estrone (Item No. 15208) intermediate by reduction of the C-17 ketone. 17-Epiestriol binds to estrogen receptor  $\alpha$  (ER $\alpha$ ) and ER $\beta$  with relative binding affinities of 29 and 80 compared with 17 $\beta$ -estradiol (Item No. 10006315).

### 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. Kuiper, G.G.J.M., Lemmen, J.G., Carlsson, B., *et al.* Interaction of estrogenic chemicals and phytoestrogens with estrogen receptor  $\beta$ . *Endocrinology* **139(10)**, 4252-4263 (1998).

**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, 12/08/2022

### 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