

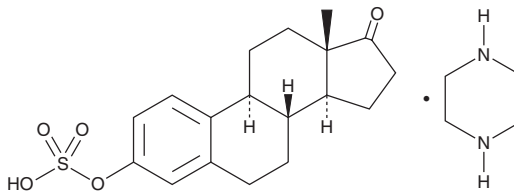
# PRODUCT INFORMATION



## Estropiate

Item No. 23806

**CAS Registry No.:** 7280-37-7  
**Formal Name:** 3-(sulfooxy)-estra-1,3,5(10)-trien-17-one, compd. with piperazine  
**Synonym:** Piperazine Estrone Sulfate  
**MF:**  $C_{18}H_{22}O_5S \cdot C_4H_{10}N_2$   
**FW:** 436.6  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Estropiate is supplied as a solid. A stock solution may be made by dissolving the estropiate in the solvent of choice. Estropiate is slightly soluble in the organic solvents DMSO and methanol, which should be purged with an inert gas.

### Description

Estropiate is a piperazine salt form of estrone 3-sulfate (Item No. 36989), which is an endogenous steroid that is converted into estrone (Item No. 10006485) *in vivo*.<sup>1</sup> Estropiate is also an inhibitor of organic anion transport polypeptide 1B1 (OATP1B1) that is selective for OATP1B1 over OATP1B3 ( $IC_{50}$ s = 0.06 and 19.3  $\mu M$ , respectively).<sup>2</sup> Formulations containing estropiate have been used in the treatment of menopausal symptoms and for the prevention of osteoporosis in high risk populations.

### References

1. Banerjee, N., Fonge, H., Mikahil, A., *et al.* Estrone-3-sulphate, a potential novel ligand for targeting breast cancers. *PLoS One* **8(5)**, e64069 (2013).
2. Gui, C., Obaidat, A., Chaguturu, R., *et al.* Development of a cell-based high-throughput assay to screen for inhibitors of organic anion transporting polypeptides 1B1 and 1B3. *Curr. Chem. Genomics* **4**, 1-8 (2010).

#### 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, 12/13/2022

#### 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