

# PRODUCT INFORMATION



## Estrone-3-Glucuronide (sodium salt hydrate)

Item No. 34417

**Formal Name:** 17-oxoestra-1,3,5(10)-trien-3-yl,  
β-D-glucopyranosiduronic acid,  
monosodium salt, hydrate

**Synonyms:** E<sub>1</sub>G, Estrone Glucuronide,  
Estrone β-D-Glucuronide

**MF:** C<sub>24</sub>H<sub>29</sub>O<sub>8</sub> • Na [XH<sub>2</sub>O]

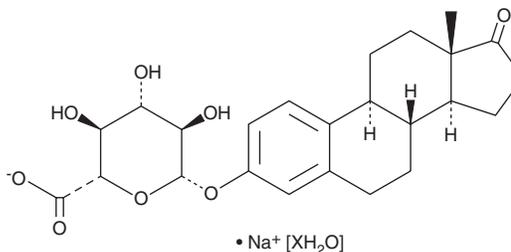
**FW:** 468.5

**Purity:** ≥98%

**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

Estrone-3-Glucuronide (sodium salt hydrate) is supplied as a solid. A stock solution may be made by dissolving the estrone-3-Glucuronide (sodium salt hydrate) in the solvent of choice, which should be purged with an inert gas. Estrone-3-Glucuronide (sodium salt hydrate) is soluble in ethanol and DMSO.

### Description

Estrone-3-Glucuronide is a metabolite of the primary premenopausal estrogen 17β-estradiol (Item Nos. ISO60155 | 10006315).<sup>1,2</sup> Urinary levels of estrone-3-Glucuronide peak approximately 36 hours prior to ovulation and have been used as a marker of fertility in women.

### References

1. Blackwell, L.F., Brown, J.B., and Cooke, D. Definition of the potentially fertile period from urinary steroid excretion rates. Part II. A threshold value for pregnanediol glucuronide as a marker for the end of the potentially fertile period in the human menstrual cycle. *Steroids* **63**(1), 5-13 (1998).
2. Adlercreutz, H., Brown, J., Collins, W., *et al.* The measurement of urinary steroid glucuronides as indices of the fertile period in women. *J. Steroid Biochem.* **17**(6), 695-702 (1982).

#### 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