

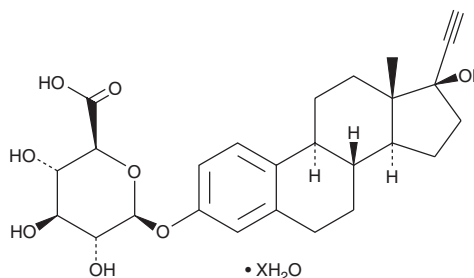
PRODUCT INFORMATION



Ethinyl Estradiol 3-β-D-Glucuronide (hydrate)

Item No. 22170

Formal Name:	(17α)-17-hydroxy-19-norpregna-1,3,5(10)-trien-20-yn-3-yl β-D-glucopyranosiduronic acid, hydrate
Synonym:	17α-Ethinylestradiol 3-Glucuronide
MF:	C ₂₆ H ₃₂ O ₈ • XH ₂ O
FW:	472.5
Purity:	≥96%
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

Ethinyl estradiol 3-β-D-glucuronide (hydrate) is supplied as a solid. A stock solution may be made by dissolving the ethinyl estradiol 3-β-D-glucuronide (hydrate) in the solvent of choice, which should be purged with an inert gas. Ethinyl estradiol 3-β-D-glucuronide (hydrate) is slightly soluble in organic solvents such as methanol and DMSO.

Description

Ethinyl estradiol 3-β-D-glucuronide is a metabolite of ethinyl estradiol (Item No. 10006486) that can be detected in urine.^{1,2} It is used as a marker of estrogen contamination in fresh and saltwater.²

References

1. Hadd, H. E., Slikker, W., Jr., and Helton, E.D. The synthesis and characterization of the glucopyranosiduronic acids of 17α-ethinylestradiol-17β. *J. Steroid Biochem.* **13(9)**, 1107-1114 (1980).
2. Griffith, D. R., Soule, M.C.K., Eglinton, T.I. *et al.* Steroidal estrogen sources in a sewage-impacted coastal ocean. *Environ. Sci.: Processes Impacts* **18(8)**, 931-1106 (2016).

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, 12/16/2022

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