

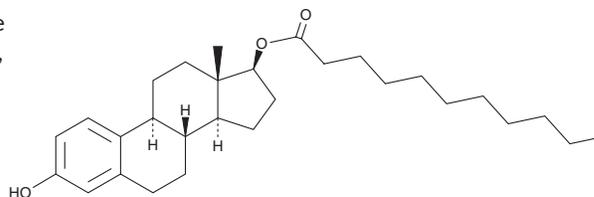
PRODUCT INFORMATION



Estradiol Undecylate

Item No. 9004927

CAS Registry No.: 3571-53-7
Formal Name: estra-1,3,5(10)-triene-3,17-diol, 17 β -undecanoate
Synonyms: β -Estradiol Undecylate, 17 β -Estradiol Undecylate, 17 β -Oestradiol Undecylate, E2 Undecylate
MF: C₂₉H₄₄O₃
FW: 440.7
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

Estradiol undecylate is supplied as a solid. A stock solution may be made by dissolving the estradiol undecylate in the solvent of choice, which should be purged with an inert gas. Estradiol undecylate is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Estradiol undecylate is a synthetic estrogen.^{1,2} It induces vaginal keratinization in ovariectomized rats when administered subcutaneously at a dose of 100 μ g/animal.¹ Estradiol undecylate (40 μ g/animal) induces vaginal estrus for 80 days in ovariectomized rats.²

References

1. Van Wauwe, J., Van Nyen, G., Coene, M.C., *et al.* Liarozole, an inhibitor of retinoic acid metabolism, exerts retinoid-mimetic effects *in vivo*. *J. Pharmacol. Exp. Ther.* **261**(2), 773-779 (1992).
2. Kuhl, H. and Taubert, H.D. A new class of long-acting hormonal steroid preparation: Synthesis of oligomeric estradiol derivatives. *Steroids* **22**(1), 73-87 (1973).

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, 12/04/2025

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