

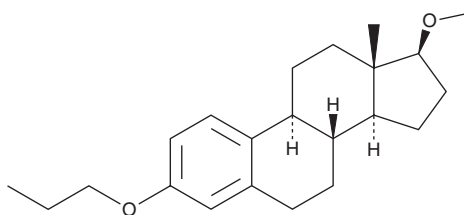
PRODUCT INFORMATION



Promestriene

Item No. 33426

CAS Registry No.: 39219-28-8
Formal Name: 17 β -methoxy-3-propoxy-estra-1,3,5(10)-triene
MF: C₂₂H₃₂O₂
FW: 328.5
Purity: \geq 95%
Supplied as: A crystalline 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

Promestriene is supplied as a crystalline solid. A stock solution may be made by dissolving the promestriene in the solvent of choice, which should be purged with an inert gas. Promestriene is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of promestriene in these solvents is approximately 5 and 15 mg/ml, respectively.

Description

Promestriene is a synthetic estrogen and derivative of estradiol (Item No. 10006315).¹ Topical application of a gel containing promestriene (1%) reduces vaginal atrophy and decreases in endometrial and myometrial wall thickness in ovariectomized rats.² Formulations containing promestriene have been used in the treatment of vaginal atrophy.

References

1. Del Pup, L., Di Francia, R., Cavaliere, C., *et al.* Promestriene, a specific topic estrogen. Review of 40 years of vaginal atrophy treatment: Is it safe even in cancer patients? *Anticancer Drugs* **24(10)**, 989-998 (2013).
2. López-Belmonte, J., Nieto, C., Estevez, J., *et al.* Comparative uterine effects on ovariectomized rats after repeated treatment with different vaginal estrogen formulations. *Maturitas* **72(4)**, 353-358 (2012).

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, 11/14/2022

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