

PRODUCT INFORMATION



Quinestrol

Item No. 10006320

CAS Registry No.: 152-43-2
Formal Name: 3-(cyclopentyloxy)-19-norpregna-1,3,5(10)-trien-20-yn-17 α -ol
Synonyms: Ethynyl Estradiol-3-cyclopentyl ether, W 3566
MF: C₂₅H₃₂O₂
FW: 3645
Purity: \geq 98%
UV/Vis.: λ_{max} : 202, 281 nm
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

Quinestrol is supplied as a crystalline solid. A stock solution may be made by dissolving the quinestrol in the solvent of choice, which should be purged with an inert gas. Quinestrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of quinestrol in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Quinestrol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Quinestrol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Quinestrol is a synthetic estrogen that is effective in hormone replacement therapy.^{1,2} It is a 3-cyclopentyl ether of ethynyl estradiol. After gastrointestinal absorption, it is stored in adipose tissue, where it is slowly released and metabolized in the liver to its active form, ethynyl estradiol. Quinestrol has found limited use in suppressing lactation in postpartum women and, in combination with synthetic progestogens, as contraceptive therapy, although additional studies are needed for both applications.

References

1. Baumgardner, S.B., Condrea, H., Daane, T.A., *et al.* Replacement estrogen therapy for menopausal vasomotor flushes. Comparison of Quinestrol and conjugated estrogens. *Obstet. Gynecol.* **51(4)**, 445-452 (1978).
2. Skouby, S.O. Criteria for the selection of an optimal estrogen replacement. *Gynecol. Endocrinol.* **15(1)**, 60-67 (2001).

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, 10/28/2022

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