

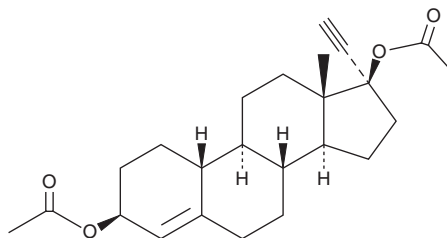
PRODUCT INFORMATION



Ethinodiol Diacetate

Item No. 29854

CAS Registry No.: 297-76-7
Formal Name: (3 β ,17 α)-19-norpregn-4-en-20-yne-3,17-diol, 3,17-diacetate
Synonym: SC-11800
MF: C₂₄H₃₂O₄
FW: 384.5
Purity: \geq 98%
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

Ethinodiol diacetate is supplied as a solid. A stock solution may be made by dissolving the ethynodiol diacetate in the solvent of choice, which should be purged with an inert gas. Ethynodiol diacetate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ethynodiol diacetate in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

Ethinodiol diacetate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ethynodiol diacetate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ethynodiol diacetate has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ethinodiol diacetate is a synthetic progestogen.¹ It prevents the development of osteoporosis in ovariectomized rats when administered at a dose of 9 μ g/animal per day. Ethynodiol diacetate inhibits copulatory ovulation in rabbits.² Formulations containing ethynodiol diacetate have been used as oral contraceptives.

References

1. Aitken, J.M., Armstrong, E., and Anderson, J.B. Osteoporosis after oophorectomy in the mature female rat and the effect of oestrogen and-or progestogen replacement therapy in its prevention. *J. Endocrinol.* **55(1)**, 79-87 (1972).
2. Suzukki, M., Hiroi, M., and Takahashi, K. Effects of ethynodiol diacetate (SC 11800) on the pregnancy and mating behavior in female rats. *Tohoku J. Exp. Med.* **107(4)**, 317-322 (1972).

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, 10/13/2022

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