

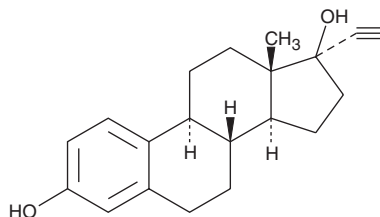
Product Information



Ethynyl Estradiol

Catalog No. 10006486

CAS Registry No.: 57-63-6
Formal Name: 19-norpregna-1,3,5(10)-trien-20-yne-3,17 α -diol
MF: C₂₀H₂₄O₂
FW: 296.4
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that ethynyl estradiol be stored as supplied at -20°C. It will be stable for at least two years.

Ethynyl estradiol is supplied as a crystalline solid. A stock solution may be made by dissolving the ethynyl estradiol in an organic solvent purged with an inert gas. Ethynyl estradiol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ethynyl estradiol in these solvents is at least 20 mg/ml.

Ethynyl estradiol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ethynyl estradiol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Ethynyl estradiol has a solubility of 0.125 mg/ml in a 1:7 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Estrogens direct the development of the female genotype in embryogenesis and at puberty. Estradiol is the major estrogen secreted by the premenopausal ovary.¹ Ethynyl estradiol is a synthetic analog of estradiol used commonly as an oral contraceptive, often in combination with a progestin such as norgestrel/levonorgestrel or desogestrel. Efficacy of oral administration of ethynyl estradiol is facilitated by the ethynyl substitution at the C-17 position, which inhibits first pass hepatic metabolism. Ethynyl estradiol is also rapidly and almost completely absorbed from the gastrointestinal tract.²

References

1. Gruber, C.J., Tschugguel, W., Schneeberger, C., *et al.* Production and actions of estrogens. *N. Engl. J. Med.* **346**(5), 340-352 (2002).
2. Coelingh Bennink, H.J.T. Are all estrogens the same? *Maturitas* **47**, 269-275 (2003).

Related Products

Estradiol EIA Kit - Cat. No. 582251 • Estradiol - Cat. No. 10006315 • Mifepristone - Cat. No. 10006317 • Levonorgestrel - Cat. No. 10006318 • Norgestrel - Cat. No. 10006319 • Tibolone - Cat. No. 10006321 • Estriol - Cat. No. 10006484 • Estrone - Cat. No. 10006485 • Estradiol Benzoate - Cat. No. 10006487 • Gestrinone - Cat. No. 10006488

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/17/2008

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com