

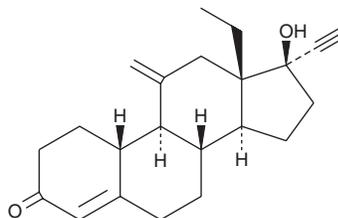
PRODUCT INFORMATION



Etonogestrel

Item No. 21062

CAS Registry No.: 54048-10-1
Formal Name: (17 α)-13-ethyl-17-hydroxy-11-methylene-18,19-dinorpregn-4-en-20-yn-3-one
Synonyms: ORG 3236, 3-keto Desogestrel, 3-oxo Desogestrel
MF: C₂₂H₂₈O₂
FW: 324.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 239 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

Etonogestrel is supplied as a crystalline solid. A stock solution may be made by dissolving the etonogestrel in the solvent of choice, which should be purged with an inert gas. Etonogestrel is soluble in organic solvents such as acetonitrile, methanol, and ethanol. The solubility of etonogestrel in these solvents is approximately 1 mg/ml.

Description

Etonogestrel is a progesterone and androgen receptor antagonist (IC₅₀s = 0.51 and 0.12 nM for isolated rabbit uterus and rat prostate cells, respectively) and an active metabolite of the synthetic progestogen desogestrel (Item No. 23651).^{1,2} It increases endometrial and myometrial vascularization, uterine weight, and uterine artery length, as well as decreases the number of hormone-secreting corpus luteal glands in non-ovariectomized guinea pigs when administered subcutaneously as a 1 mm sustained-release pellet.³ Formulations containing etonogestrel have been used in contraceptives to prevent pregnancy.

References

1. Stanczyk, F.Z. All progestins are not created equal. *Steroids* **68(10-13)**, 879-890 (2003).
2. Phillips, A., Demarest, K., Hahn, D.W., *et al.* Progestational and androgenic receptor binding affinities and in vivo activities of norgestimate and other progestins. *Contraception* **41(4)**, 399-410 (1990).
3. Krikun, G., Booth, C.J., Buchwalder, L., *et al.* Effects of etonogestrel treatment in the reproductive organs and uterine arteries of nonophorectomized guinea pigs. *Reprod. Sci.* **19(4)**, 400-406 (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, 03/21/2023

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