

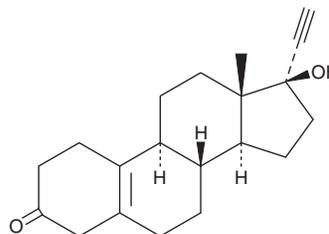
PRODUCT INFORMATION



Norethynodrel

Item No. 33230

CAS Registry No.: 68-23-5
Formal Name: 17 α -hydroxy-19-norpregn-5(10)-en-20-yn-3-one
Synonyms: Norethinodrel, NSC 15432, SC-4642
MF: C₂₀H₂₆O₂
FW: 298.4
Purity: \geq 90%
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

Norethynodrel is supplied as a solid. A stock solution may be made by dissolving the norethynodrel in the solvent of choice, which should be purged with an inert gas. Norethynodrel is slightly soluble in chloroform and methanol.

Description

Norethynodrel is a synthetic progestin.^{1,2} *In vivo*, norethynodrel (10 mg/animal per day) reduces fetal implantation and induces fetal resorption in mated female rats.² Formulations containing norethynodrel have been used as oral contraceptives.

References

1. Kubli-Garfias, C., González-Mariscal, G., Vásquez-Ramírez, R., *et al.* Bimodal binding and free energy of the progesterone receptor in the induction of female sexual receptivity by progesterone and synthetic progestins. *J. Steroid Biochem. Mol. Biol.* **133**, 43-50 (2013).
2. Davis, B.K. Termination of pregnancy in the rat with norethynodrel. *Nature* **197**, 308-309 (1963).

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, 03/13/2026

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