

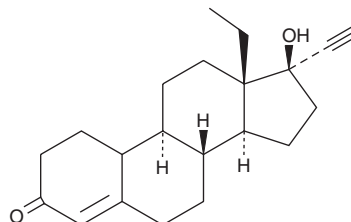
PRODUCT INFORMATION



Norgestrel

Item No. 10006319

CAS Registry No.: 6533-00-2
Formal Name: (±)-13-ethyl-17 α -hydroxy-18,19-dinorpregn-4-en-20-yn-3-one
Synonyms: (±)-Norgestrel, DL-Norgestrel, α -Norgestrel, SH 70850
MF: C₂₁H₂₈O₂
FW: 312.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 240 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Norgestrel is supplied as a crystalline solid. A stock solution may be made by dissolving the norgestrel in the solvent of choice. Norgestrel is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of norgestrel in these solvents is approximately 0.25, 5, and 10 mg/ml, respectively.

Norgestrel is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, norgestrel should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Norgestrel has a solubility of approximately 0.25 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Norgestrel is a synthetic progestin and a racemic mixture of dextronorgestrel and levonorgestrel (Item No. 10006318), of which levonorgestrel is the biologically active component.¹ *In vivo*, norgestrel administered *via* intrauterine device (IUD) prevents pregnancy in rats.² Formulations containing norgestrel have been used as contraceptives.

References

1. Jones, R.C., Singer, A.C., and Edgren, R.A. The biological activities of norgestrel and its enantiomers. *Int. J. Fertil.* **24(1)**, 39-43 (1979).
2. Jones, R.C. The effects of an intrauterine device containing norgestrel in the rat. *Fertil. Steril.* **25(9)**, 767-771 (1974).

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, 08/13/2019

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