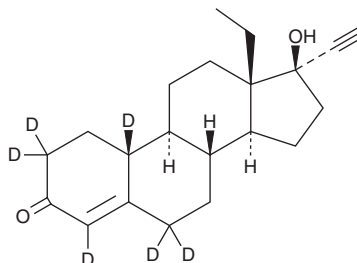


PRODUCT INFORMATION



Norgestrel-d₆ Item No. 28477

Formal Name:	(±)-13-ethyl-17 α -hydroxy-18,19-Dinorpregn-4-en-20-yn-3-one-2,2,4,6,6,10-d ₆
Synonyms:	(±)-Norgestrel-d ₆ , DL-Norgestrel-d ₆ , α -Norgestrel-d ₆
MF:	C ₂₁ H ₂₂ D ₆ O ₂
FW:	318.5
Chemical Purity:	≥98% (Norgestrel)
Deuterium Incorporation:	≥99% deuterated forms (d ₁ -d ₆); ≤1% d ₀
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years



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

Laboratory Procedures

Norgestrel-d₆ is intended for use as an internal standard for the quantification of norgestrel (Item No. 10006319) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Norgestrel-d₆ is supplied as a solid. A stock solution may be made by dissolving the norgestrel-d₆ in the solvent of choice, which should be purged with an inert gas. Norgestrel-d₆ is slightly soluble in chloroform (warmed) and methanol (sonicated).

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, 01/03/2023

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