

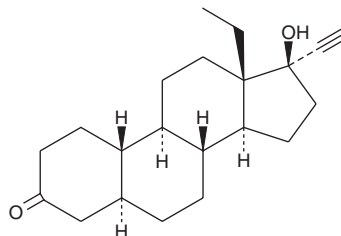
PRODUCT INFORMATION



5 α -dihydro Levonorgestrel

Item No. 27854

CAS Registry No.: 78088-19-4
Formal Name: (5 α ,17 α)-13-ethyl-17-hydroxy-18,19-dinorpregn-20-yn-3-one
Synonym: 5 α -DHLNG
MF: C₂₁H₃₀O₂
FW: 314.5
Purity: \geq 95%
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

5 α -dihydro Levonorgestrel is supplied as a solid. A stock solution may be made by dissolving the 5 α -dihydro levonorgestrel in the solvent of choice, which should be purged with an inert gas. 5 α -dihydro Levonorgestrel is slightly soluble in chloroform and methanol.

Description

5 α -dihydro Levonorgestrel is a metabolite of the synthetic progestin levonorgestrel (Item No. 10006318).^{1,2} It lacks *in vitro* and *in vivo* estrogenic activity, but induces copulatory behavior in male rats when administered at a dose of 1 mg per animal per day.

References

1. Santillán, R., Pérez-Palacios, G., Reyes, M., *et al.* Assessment of the oestrogenic activity of the contraceptive progestin levonorgestrel and its non-phenolic metabolites. *Eur. J. Pharmacol.* **427(2)**, 167-174 (2001).
2. Morali, G., Lemus, A.E., Munguía, R., *et al.* Hormone-like behavioral effects of levonorgestrel and its metabolites in the male rat. *Pharmacol. Biochem. Behav.* **73(4)**, 951-961 (2002).

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, 12/02/2022

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