

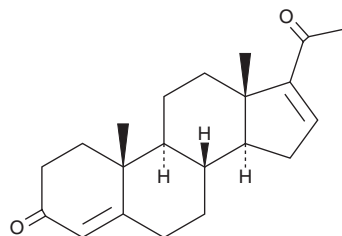
PRODUCT INFORMATION



16-Dehydroprogesterone

Item No. 43057

CAS Registry No.: 1096-38-4
Formal Name: pregna-4,16-diene-3,20-dione
Synonyms: NSC 9785, NSC 11037, Δ^{16} -Progesterone
MF: $C_{21}H_{28}O_2$
FW: 312.5
Purity: $\geq 98\%$
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

16-Dehydroprogesterone is supplied as a solid. A stock solution may be made by dissolving the 16-dehydroprogesterone in the solvent of choice, which should be purged with an inert gas. 16-Dehydroprogesterone is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

16-Dehydroprogesterone is a progestin and an active metabolite of pregnenolone (Item No. 19864).¹⁻³ It inhibits the proliferation of HL-60 leukemia and U937 lymphoma cells (IC_{50} s = 10.9 and 3.6 μM , respectively).² 16-Dehydroprogesterone (2 mg/kg), in combination with gonadotropin, increases ovary and uterine weights in prepuberal rhesus monkeys.³

References

1. Kwan, T.K., Taylor, and N.F., Gower, D.B. The use of steroid profiling in the resolution of pregnenolone metabolites from porcine testicular preparations. *J. Chromatogr.* **301**(1), 189-197 (1984).
2. Samudio, I., Konopleva, M., Safe, S., *et al.* Guggulsterones induce apoptosis and differentiation in acute myeloid leukemia: Identification of isomer-specific antileukemic activities of the pregnadienedione structure. *Mol. Cancer Ther.* **4**(12), 1982-1992 (2005).
3. Kar, A.B. and Chandra, H. Effect of some progestational steroids on the response of the ovary of prepuberal rhesus monkeys to exogenous gonadotrophin. *Steroids.* **6**(4), 463-472 (1965).

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, 04/08/2025

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