

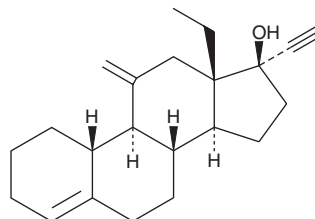
PRODUCT INFORMATION



Desogestrel

Item No. 23651

CAS Registry No.: 54024-22-5
Formal Name: (17 α)-13-ethyl-11-methylene-18,19-dinorpregn-4-en-20-yn-17-ol
Synonym: ORG 2969
MF: C₂₂H₃₀O
FW: 310.5
Purity: $\geq 95\%$
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

Desogestrel is supplied as a solid. A stock solution may be made by dissolving the desogestrel in the solvent of choice, which should be purged with an inert gas. Desogestrel is soluble in the organic solvent DMSO at a concentration of approximately 13.7 mg/ml. Desogestrel is soluble in methanol.

Description

Desogestrel is a synthetic progestogen.¹ It inhibits ovulation and prevents fertilization in rabbits and mice. Desogestrel inhibits rhodamine 123 (Item No. 16672) efflux, a measure of P-glycoprotein activity, *ex vivo* in human lymphocytes and CD8⁺ T cells in a dose-dependent manner.² Formulations containing desogestrel have been used as oral contraceptives and for the treatment of polycystic ovary syndrome.

References

1. van der Vies, J., and de Visser, J. Endocrinological studies with desogestrel. *Arzneimittelforschung* **33**(2), 231-236 (1983).
2. Fröhlich, M., Albermann, N., Sauer, A., *et al.* *In vitro* and *ex vivo* evidence for modulation of P-glycoprotein activity by progestins. *Biochem. Pharmacol.* **68**(12), 2409-2416 (2004).

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, 12/12/2022

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