

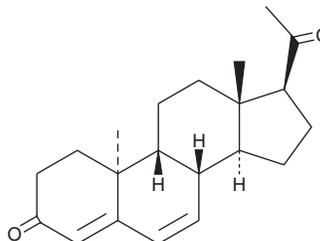
PRODUCT INFORMATION



Dydrogesterone

Item No. 20514

CAS Registry No.: 152-62-5
Formal Name: (9 β ,10 α)-pregna-4,6-diene-3,20-dione
Synonym: NSC 92336
MF: C₂₁H₂₈O₂
FW: 312.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 287 nm
Supplied as: A crystalline 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

Dydrogesterone is supplied as a crystalline solid. A stock solution may be made by dissolving the dydrogesterone in the solvent of choice. Dydrogesterone is soluble in organic solvents such as ethanol, acetonitrile, and methanol, which should be purged with an inert gas. The solubility of dydrogesterone in these solvents is approximately 1 mg/ml.

Description

Dydrogesterone is a synthetic progestogen that has no estrogenic, androgenic, glucocorticoid, or anabolic effects, although it is an agonist of progesterone receptors ($EC_{50} = 12.3$ nM) and an antagonist of androgen, mineralocorticoid, and glucocorticoid receptors (IC_{50} s = 28.6, 82.7, and 363 nM, respectively).^{1,2} Dydrogesterone protects against sound stress-induced abortion in mice when administered at a dose of 1.25 mg per animal prior to the stressor on day five of pregnancy.³ Formulations containing dydrogesterone have been used in hormone replacement therapy, in the treatment of menstrual disorders, and to prevent miscarriage.

References

1. Coelingh Bennink, H.J. and Boerrigter, P.J. Use of dydrogesterone as a progestogen for oral contraception. *Steroids* **68(10-13)**, 927-929 (2003).
2. Rižner, T.L., Brožič, P., Doucette, C., *et al.* Selectivity and potency of the retroprogesterone dydrogesterone *in vitro*. *Steroids* **76(6)**, 607-615 (2011).
3. Joachim, R., Zenclussen, A.C., Polgar, B., *et al.* The progesterone derivative dydrogesterone abrogates murine stress-triggered abortion by inducing a Th2 biased local immune response. *Steroids* **68(10-13)**, 931-940 (2003).

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, 12/19/2022

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