

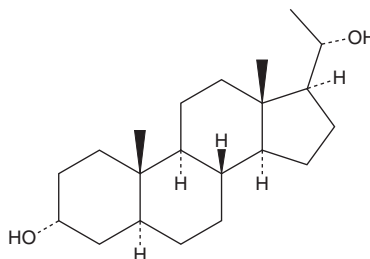
PRODUCT INFORMATION



5 α -Pregnane-3 α ,20 α -diol

Item No. 38890

CAS Registry No.: 566-58-5
Formal Name: (5 α)-pregnane-3 α ,20S-diol
Synonym: Allopregnanediol
MF: C₂₁H₃₆O₂
FW: 320.5
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

5 α -Pregnane-3 α ,20 α -diol is supplied as a solid. A stock solution may be made by dissolving the 5 α -pregnane-3 α ,20 α -diol in the solvent of choice, which should be purged with an inert gas. 5 α -Pregnane-3 α ,20 α -diol is slightly soluble in chloroform and methanol.

Description

5 α -Pregnane-3 α ,20 α -diol is an active metabolite of progesterone (Item No. 15876).¹ It is formed from progesterone via 5 α -pregnane-20 α -ol-3-one and 5 α -pregnane-3 α -ol-20-one intermediates. 5 α -Pregnane-3 α ,20 α -diol (10 μ M) increases intracellular calcium levels in washed isolated human platelets.²

References

1. Zhang, Y., Nadeau, M., Faucher, F., et al. Progesterone metabolism in adipose cells. *Mol. Cell. Endocrinol.* **298(1-2)**, 76-83 (2009).
2. Blackmore, P.F. Progesterone metabolites rapidly stimulate calcium influx in human platelets by a src-dependent pathway. *Steroids* **73(7)**, 738-750 (2008).

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, 06/13/2023

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