

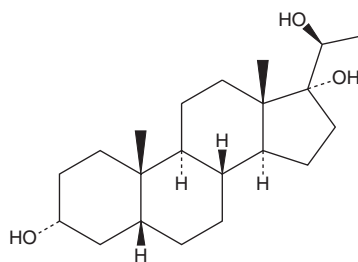
PRODUCT INFORMATION



Pregnanetriol

Item No. 33312

CAS Registry No.: 1098-45-9
Formal Name: (5 β)-pregnane-3 α ,17,20S-triol
Synonyms: NSC 52760, 5 β -Pregnane-3 α ,17 α ,20 α -triol
MF: C₂₁H₃₆O₃
FW: 336.5
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

Pregnanetriol is supplied as a solid. A stock solution may be made by dissolving the pregnanetriol in the solvent of choice, which should be purged with an inert gas. Pregnanetriol is soluble in methanol.

Description

Pregnanetriol is a metabolite of 17 α -hydroxyprogesterone (Item No. 33154).^{1,2} It is formed from 17 α -hydroxyprogesterone by reduction of the C-20 ketone.² Urinary levels of pregnanetriol are elevated in patients with 21-hydroxylase deficiency and congenital adrenal hyperplasia.^{3,4}

References

1. Kamrath, C., Hartmann, M.F., Boettcher, C., *et al.* Diagnosis of 21-hydroxylase deficiency by urinary metabolite ratios using gas chromatography-mass spectrometry analysis: Reference values for neonates and infants. *J. Steroid Biochem. Mol. Biol.* **156**, 10-16 (2016).
2. Schiffer, L., Barnard, L., Baranowski, E.S., *et al.* Human steroid biosynthesis, metabolism and excretion are differentially reflected by serum and urine steroid metabolomes: A comprehensive review. *J. Steroid Biochem. Mol. Biol.* **194**, 105439 (2019).
3. *Disorders of steroidogenesis guide to steroid profiling and biochemical diagnosis*. 1st ed., Springer International Publishing AG, Cham, Switzerland (2019).
4. Shackleton, C.H.L. Role of a disordered steroid metabolome in the elucidation of sterol and steroid biosynthesis. *Lipids* **47(1)**, 1-12 (2012).

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, 12/07/2022

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