

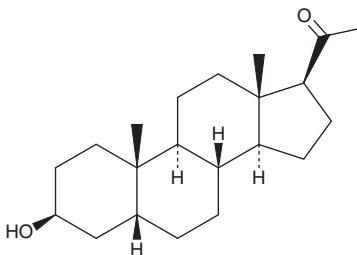
PRODUCT INFORMATION



Epipregnanolone

Item No. 34295

CAS Registry No.: 128-21-2
Formal Name: (5 β)-3 β -hydroxy-pregnan-20-one
Synonyms: NSC 21450, 5 β -Pregnan-3 β -ol-20-one
MF: C₂₁H₃₄O₂
FW: 318.5
Purity: \geq 98%
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

Epipregnanolone is supplied as a solid. A stock solution may be made by dissolving the epipregnanolone in the solvent of choice, which should be purged with an inert gas. Epipregnanolone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of epipregnanolone in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Epipregnanolone is a neurosteroid and an active metabolite of the steroid hormone pregnenolone (Item No. 19864).¹ It is enzymatically formed from pregnenolone via the intermediates progesterone (Item No. 15876) and 5 β -dihydroprogesterone in the placenta.² Epipregnanolone inhibits spontaneous contractions in myometrial strips isolated from at-term pregnant women (IC₅₀ = 156 μ M).³ Epipregnanolone (10 and 20 mg/kg) decreases operant alcohol self-administration in rats.⁴ Maternal plasma levels of epipregnanolone increase over the duration of pregnancy.²

References

1. Prince, R.J. and Simmonds, M.A. 5 β -pregnan-3 β -ol-20-one, a specific antagonist at the neurosteroid site of the GABA_A receptor-complex. *Neurosci. Lett.* **135**(2), 273-275 (1992).
2. Hill, M., Cibula, D., Havlíková, H., *et al.* Circulating levels of pregnanolone isomers during the third trimester of human pregnancy. *J. Steroid Biochem. Mol. Biol.* **105**(1-5), 166-175 (2007).
3. Perusquía, M. and Jasso-Kamel, J. Influence of 5 α - and 5 β -reduced progestins on the contractility of isolated human myometrium at term. *Life Sci.* **68**(26), 2933-2944 (2001).
4. O'Dell, L.E., Purdy, R.H., Covey, D.F., *et al.* Epipregnanolone and a novel synthetic neuroactive steroid reduce alcohol self-administration in rats. *Pharmacol. Biochem. Behav.* **81**(3), 543-550 (2005).

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, 07/07/2025

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