

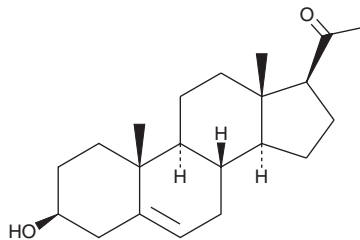
PRODUCT INFORMATION



Pregnenolone

Item No. 19864

CAS Registry No.: 145-13-1
Formal Name: 3 β -hydroxy-pregn-5-en-20-one
Synonyms: NSC 1616, NSC 18158,
3 β -hydroxy-5-Pregnen-20-one,
 Δ^5 -Pregnenolone
MF: C₂₁H₃₂O₂
FW: 316.5
Purity: $\geq 98\%$
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Pregnenolone is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Pregnenolone is a natural steroid hormone that serves as a precursor for a wide range of steroids, including mineralocorticoids, glucocorticoids, androgens, and estrogens.¹ Pregnenolone sulfate modulates NMDA receptor responses to exogenously applied glutamate and stimulates transient receptor potential melastatin 3 (TRPM3).²⁻⁴

References

1. Akhtar, M. K., Kelly, S. L. and Kaderbhai, M. A. Cytochrome b5 modulation of 17 α hydroxylase and 17-20 lyase (CYP17) activities in steroidogenesis. *J. Endocrinol.* **187**(2), 267-274 (2005).
2. Sedláček, M., Kořínek, M., Petrovič, M., et al. Neurosteroid modulation of ionotropic glutamate receptors and excitatory synaptic transmission. *Physiol. Res.* **57**(Suppl 3), S49-S57 (2008).
3. Smith, C. C., Gibbs, T. T., and Farb, D. H. Pregnenolone sulfate as a modulator of synaptic plasticity. *Psychopharmacology (Berl)* **231**(17), (2014).
4. Thiel, G., Müller, I., and Rössler, O. G. Signal transduction via TRPM3 channels in pancreatic β -cells. *J. Mol. Endocrinol.* **50**(3), R75-R83 (2013).

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, 12/15/2022

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