

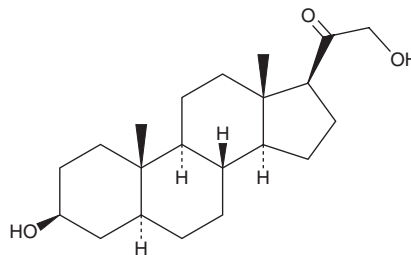
PRODUCT INFORMATION



3 β ,5 α -THDOC

Item No. 18431

CAS Registry No.: 567-01-1
Formal Name: (3 β ,5 α)-3,21-dihydroxy-pregnan-20-one
Synonyms: 3 β ,21-dihydroxy-5 α -Pregnan-20-one,
3 β ,5 α -Tetrahydrodeoxy Corticosterone
MF: C₂₁H₃₄O₃
FW: 334.5
Purity: \geq 95%
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

3 β ,5 α -THDOC is supplied as a crystalline solid. A stock solution may be made by dissolving the 3 β ,5 α -THDOC in the solvent of choice. 3 β ,5 α -THDOC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 3 β ,5 α -THDOC in these solvents is approximately 25 mg/ml.

3 β ,5 α -THDOC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3 β ,5 α -THDOC should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 3 β ,5 α -THDOC has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3 β ,5 α -THDOC is a neurosteroid.¹ It inhibits seizures induced by pilocarpine, but not pentylenetetrazol (PTZ; Item No. 18682), in mice (ED₅₀s = 45.9 and >100 mg/kg, respectively). 3 β ,5 α -THDOC has been found in the urine of women during the third trimester of pregnancy.²

References

1. Kokate, T.G., Cohen, A.L., Karp, E., *et al.* Neuroactive steroids protect against pilocarpine- and kainic acid-induced limbic seizures and status epilepticus in mice. *Neuropharmacology* **35(8)**, 1049-1056 (1996).
2. Pasqualini, J.R. and Bedrak, E. Identification of a new steroid: 3 β ,21-dihydroxy-5 α -pregnan-20-one in the urine of human pregnancy. *J. Clin. Endocrinol. Metab.* **38(2)**, 298-304 (1974).

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, 10/06/2022

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