

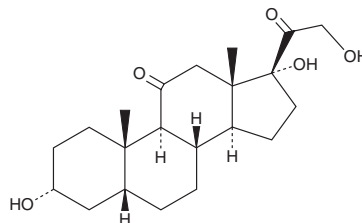
PRODUCT INFORMATION



Tetrahydrocortisone

Item No. 34344

CAS Registry No.: 53-05-4
Formal Name: (3 α ,5 β)-3,17,21-trihydroxy-pregnane-11,20-dione
Synonyms: NSC 76984, 5 β -Tetrahydrocortisone, THE, Urocortisone
MF: C₂₁H₃₂O₅
FW: 364.5
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

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

Description

Tetrahydrocortisone is a metabolite of cortisol (hydrocortisone; Item No. 20739).^{1,2} It is formed from cortisol via cortisone (Item No. 30763) as an intermediate.

References

1. Abel, S.M., Maggs, J.L., Back, D.J., *et al.* Cortisol metabolism by human liver *in vitro*—I. Metabolite identification and inter-individual variability. *J. Steroid. Biochem. Mol. Biol.* **43**(7), 713-719 (1992).
2. White, P.C., Mune, T., and Agarwal, A.K. 11 β -hydroxysteroid dehydrogenase and the syndrome of apparent mineralocorticoid excess. *Endocr. Rev.* **18**(1), 135-156 (1997).

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, 10/24/2022

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