

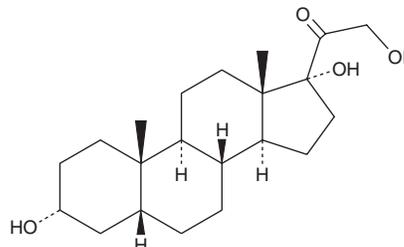
PRODUCT INFORMATION



Tetrahydro-11-deoxy Cortisol

Item No. 26501

CAS Registry No.: 68-60-0
Formal Name: 5 β -3 α ,17,21-trihydroxy-pregnan-20-one
Synonyms: NSC 53901, 5 β -Pregnane-3 α ,17 α ,21-triol-20-one, Tetrahydrodeoxycortisol, THS
MF: C₂₁H₃₄O₄
FW: 350.5
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

Tetrahydro-11-deoxy cortisol (THS) is supplied as a crystalline solid. A stock solution may be made by dissolving the THS in the solvent of choice. THS is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of THS in these solvents is approximately 30 mg/ml.

THS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, THS should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. THS has a solubility of approximately 0.2 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

THS is the primary urinary metabolite of 11-deoxycortisol.^{1,2} Urinary excretion of THS is elevated in patients with 11 β -hydroxylase deficiency, a condition resulting from mutations in the cytochrome P450 (CYP) isoform CYP11B1. Urinary levels of THS are also elevated in patients with adrenocortical carcinoma (ACC) and adrenocortical adenoma (ACA) but are higher in patients with ACC compared to ACA.^{2,3}

References

1. Keavney, B., Mayosi, B., Gaukrodger, N., *et al.* Genetic variation at the locus encompassing 11- β hydroxylase and aldosterone synthase accounts for heritability in cortisol precursor (11-deoxycortisol) urinary metabolite excretion. *J. Clin. Endocrinol. Metab.* **90**(2), 1072-1077 (2005).
2. Nguyen, H.-H., Eiden-Plach, A., Hannemann, F., *et al.* Phenotypic, metabolic, and molecular genetic characterization of six patients with congenital adrenal hyperplasia caused by novel mutations in the CYP11B1 gene. *J. Steroid Biochem. Mol. Biol.* **155**(Pt A), 126-134 (2016).
3. Arlt, W., Biehl, M., Taylor, A.E., *et al.* Urine steroid metabolomics as a biomarker tool for detecting malignancy in adrenal tumors. *J. Clin. Endocrinol. Metab.* **96**(12), 3775-3784 (2011).
4. Velikanova, L.I., Shafigullina, Z.R., Lisitsin, A.A., *et al.* Different types of urinary steroid profiling obtained by high-performance liquid chromatography and gas chromatography-mass spectrometry in patients with adrenocortical carcinoma. *Horm. Cancer* **7**(5-6), 327-335 (2016).

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, 12/19/2022

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