

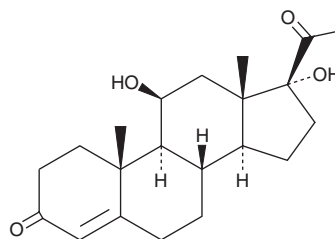
PRODUCT INFORMATION



21-Deoxycortisol

Item No. 30873

CAS Registry No.: 641-77-0
Formal Name: 11 β ,17-dihydroxy-pregn-4-ene-3,20-dione
Synonym: 21-Dehydrohydrocortisone
MF: C₂₁H₃₀O₄
FW: 346.5
Purity: \geq 95%
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

21-Deoxycortisol is supplied as a solid. A stock solution may be made by dissolving the 21-deoxycortisol in the solvent of choice, which should be purged with an inert gas. 21-Deoxycortisol is slightly soluble in chloroform and methanol.

Description

21-Deoxycortisol is a corticosteroid metabolite of 17-hydroxyprogesterone produced in the adrenal gland via 11-hydroxylation by 11 β -hydroxylase.^{1,2} Serum levels of 21-deoxycortisol are elevated in patients with congenital adrenal hyperplasia that are heterozygous for mutations in CYP2A21, the gene encoding steroid 21-hydroxylase, and have been used as a biomarker for the detection of 21-hydroxylase deficiencies.

References

1. Fiet, J., Villette, J.-M., Galons, H., *et al.* The application of a new highly-sensitive radioimmunoassay for plasma 21-deoxycortisol to the detection of steroid-21-hydroxylase deficiency. *Ann. Clin. Biochem.* **31**(Pt. 1), 56-64 (1994).
2. Cristoni, S., Cuccato, D., Sciannamblo, M., *et al.* Analysis of 21-deoxycortisol, a marker of congenital adrenal hyperplasia, in blood by atmospheric pressure chemical ionization and electrospray ionization using multiple reaction monitoring. *Rapid Commun. Mass Spectrom.* **18**(1), 77-82 (2004).

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, 11/04/2022

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