

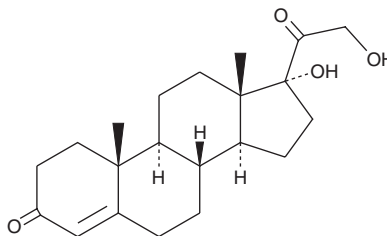
PRODUCT INFORMATION



11-deoxy Cortisol

Item No. 20775

CAS Registry No.: 152-58-9
Formal Name: 17,21-dihydroxy-pregn-4-ene-3,20-dione
Synonyms: Compound S, Cortexolone, Cortodoxone, NSC 18317, Reichstein's Substance S
MF: $C_{21}H_{30}O_4$
FW: 346.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

11-deoxy Cortisol is supplied as a solid. A stock solution may be made by dissolving the 11-deoxy cortisol in the solvent of choice, which should be purged with an inert gas. 11-deoxy Cortisol is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

11-deoxy Cortisol is a precursor in the biosynthesis of cortisol (Item No. 20739).¹ It is converted into cortisol by the cytochrome P450 (CYP) isoform CYP11B1, also known as 11 β -hydroxylase. Urinary levels of 11-deoxy cortisol and the ratio of 11-deoxy cortisol to cortisol are elevated in patients with 11 β -hydroxylase deficiency, a type of congenital adrenal hyperplasia (CAH) characterized by cortisol deficiency and hypertension.

Reference

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)

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, 06/15/2026

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