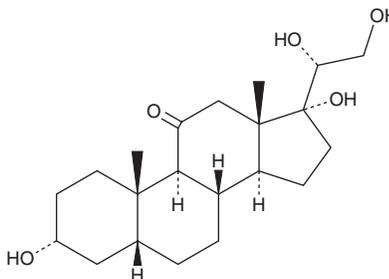


PRODUCT INFORMATION



β -Cortolone Item No. 36224

CAS Registry No.: 667-66-3
Formal Name: (5 β)-3 α ,17,20R,21-tetrahydroxy-pregnan-11-one
Synonyms: 20 β -Cortolone, NSC 56997
MF: C₂₁H₃₄O₅
FW: 366.5
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

β -Cortolone is supplied as a solid. A stock solution may be made by dissolving the β -cortolone in the solvent of choice, which should be purged with an inert gas. β -Cortolone is slightly soluble in chloroform (warmed) and methanol (warmed).

Description

β -Cortolone is a metabolite of the glucocorticoid cortisol (hydrocortisone; Item No. 20739).¹ It is formed from cortisol by 20 β -hydroxysteroid dehydrogenase (20 β -HSD) via a 5 β -tetrahydrocortisol (Item No. 34195) intermediate. Urinary levels of β -cortolone are decreased in postmenopausal women with endometrial adenocarcinoma compared to postmenopausal women without endometrial adenocarcinoma.²

References

1. Espiard, S., McQueen, J., Sherlock, M., *et al.* Improved urinary cortisol metabolome in Addison disease: A prospective trial of dual-release hydrocortisone. *J. Clin. Endocrinol. Metab.* **106(3)**, 814-825 (2021).
2. Bufa, A., Bíró, I., Poór, V., *et al.* Altered urinary profiles of endogenous steroids in postmenopausal women with adenocarcinoma endometrii. *Gynecol. Endocrinol.* **26(1)**, 10-15 (2010).

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, 02/23/2023

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