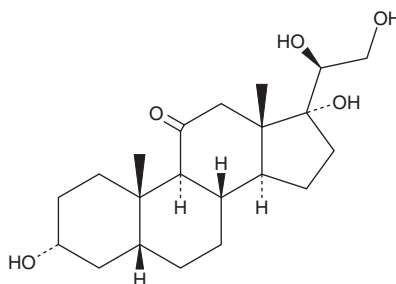


# PRODUCT INFORMATION



## $\alpha$ -Cortolone Item No. 35609

**CAS Registry No.:** 516-42-7  
**Formal Name:** (3 $\alpha$ ,5 $\beta$ ,20S)-3,17,20,21-tetrahydroxy-pregnan-11-one  
**Synonyms:** 20 $\alpha$ -Cortolone, NSC 59872  
**MF:** C<sub>21</sub>H<sub>34</sub>O<sub>5</sub>  
**FW:** 366.5  
**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

$\alpha$ -Cortolone is supplied as a solid. A stock solution may be made by dissolving the  $\alpha$ -cortolone in the solvent of choice, which should be purged with an inert gas.  $\alpha$ -Cortolone is soluble in ethanol and DMSO.

### Description

$\alpha$ -Cortolone is a metabolite of the glucocorticoid cortisol (hydrocortisone; Item No. 20739).<sup>1</sup> It is formed from cortisol by 20 $\alpha$ -hydroxysteroid dehydrogenase (20 $\alpha$ -HSD) via a tetrahydrocortisone (THE; Item No. 34344) intermediate. Urinary levels of  $\alpha$ -cortolone are decreased in patients with severe anorexia nervosa and in postmenopausal women with endometrial adenocarcinoma compared to postmenopausal women without endometrial adenocarcinoma.<sup>2,3</sup>

### 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. Wassif, W.S., McLoughlin, D.M., Vincent, R.P., *et al.* Steroid metabolism and excretion in severe anorexia nervosa: Effects of refeeding. *Am. J. Clin. Nutr.* **93(5)**, 911-917 (2011).
3. 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, 12/12/2022

#### 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