

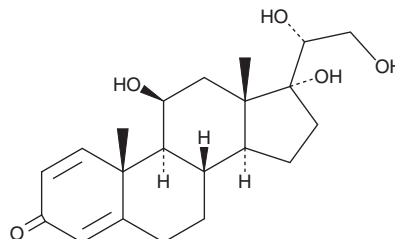
# PRODUCT INFORMATION



## 20 $\alpha$ -dihydro Prednisolone

Item No. 26502

**CAS Registry No.:** 15847-24-2  
**Formal Name:** 11 $\beta$ ,17,20R,21-tetrahydroxy-pregna-1,4-dien-3-one  
**Synonyms:** 20 $\alpha$ -DHP, 20 $\alpha$ -dihydro-PRED, 20(R)-hydroxy Prednisolone, 20 $\alpha$ -hydroxy Prednisolone  
**MF:** C<sub>21</sub>H<sub>30</sub>O<sub>5</sub>  
**FW:** 362.5  
**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

20 $\alpha$ -dihydro Prednisolone is supplied as a solid. A stock solution may be made by dissolving the 20 $\alpha$ -dihydro prednisolone in the solvent of choice, which should be purged with an inert gas. 20 $\alpha$ -dihydro Prednisolone is slightly soluble in DMSO and methanol.

### Description

20 $\alpha$ -dihydro Prednisolone is a metabolite of prednisolone (Item No. 20866).<sup>1,2</sup> It has been found in the urine of cattle following adrenocorticotrophic hormone (ACTH) administration but not in naturally stressed or non-stressed cattle.<sup>3</sup>

### References

1. Rodchenkov, G.M., Vedenin, A.N., Uralets, V.P., *et al.* Characterization of prednisone, prednisolone and their metabolites by gas chromatography-mass spectrometry. *J. Chromatogr.* **565(1-2)**, 45-51 (1991).
2. Matabosch, X., Pozo, O.J., Pérez-Mañá, C., *et al.* Detection and characterization of prednisolone metabolites in human urine by LC-MS/MS. *J. Mass. Spectrom.* **50(3)**, 633-642 (2015).
3. De Clercq, N., Van Meulebroek, L., Vanden Bussche, J., *et al.* The impact of stress on the prevalence of prednisolone in bovine urine: A metabolic fingerprinting approach. *J. Steroid. Biochem. Mol. Biol.* **154**, 206-216 (2015).

#### 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, 11/14/2022

#### 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