

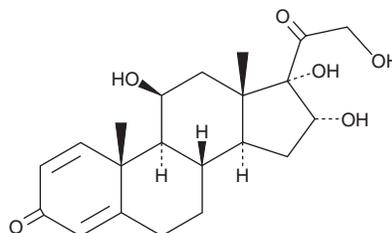
# PRODUCT INFORMATION



## 16 $\alpha$ -hydroxy Prednisolone

Item No. 27169

**CAS Registry No.:** 13951-70-7  
**Formal Name:** (11 $\beta$ ,16 $\alpha$ )-11,16,17,21-tetrahydroxy-pregna-1,4-diene-3,20-dione  
**MF:** C<sub>21</sub>H<sub>28</sub>O<sub>6</sub>  
**FW:** 376.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 243 nm  
**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

16 $\alpha$ -hydroxy Prednisolone is supplied as a solid. A stock solution may be made by dissolving the 16 $\alpha$ -hydroxy prednisolone in the solvent of choice. 16 $\alpha$ -hydroxy Prednisolone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of 16 $\alpha$ -hydroxy prednisolone in these solvents is approximately 30 mg/ml.

16 $\alpha$ -hydroxy Prednisolone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 16 $\alpha$ -hydroxy prednisolone should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 16 $\alpha$ -hydroxy Prednisolone has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

16 $\alpha$ -hydroxy Prednisolone is a metabolite of 22(R)-budesonide.<sup>1</sup> It is formed from 22(R)-budesonide by the cytochrome P450 (CYP) isoform CYP3A in human liver microsomes

### Reference

- Jönsson, G., Aström, A., and Andersson, P. Budesonide is metabolized by cytochrome P450 3A (CYP3A) enzymes in human liver. *Drug Metab. Dispos.* **23(1)**, 137-142 (1995).

#### 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, 03/29/2023

#### 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