

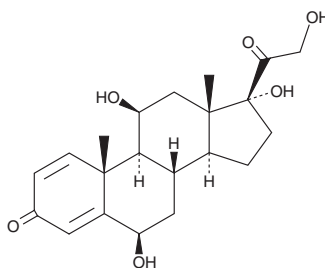
PRODUCT INFORMATION



6 β -hydroxy Prednisolone

Item No. 35720

CAS Registry No.: 16355-29-6
Formal Name: 6 β ,11 β ,17,21-tetrahydroxy-
pregna-1,4-diene-3,20-dione
MF: C₂₁H₂₈O₆
FW: 376.4
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

6 β -hydroxy Prednisolone is supplied as a solid. A stock solution may be made by dissolving the 6 β -hydroxy prednisolone in the solvent of choice, which should be purged with an inert gas. 6 β -hydroxy Prednisolone is soluble in methanol and DMSO.

Description

6 β -hydroxy Prednisolone is a metabolite of prednisolone (Item No. 20866), the active metabolite of prednisone (Item No. 20677).¹ It is formed from prednisolone primarily by the cytochrome P450 (CYP) isoform CYP3A4 and to a lesser extent by CYP3A5

Reference

1. Skauby, R.H., Bergan, S., Andersen, A.M., *et al.* In vitro assessments predict that CYP3A4 contributes to a greater extent than CYP3A5 to prednisolone clearance. *Basic Clin. Pharmacol. Toxicol.* **129(6)**, 427-436 (2021).

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, 12/13/2022

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