

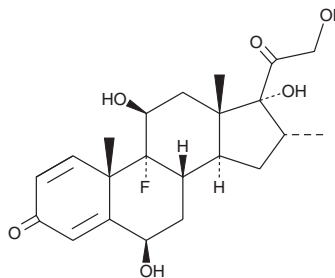
PRODUCT INFORMATION



6 β -hydroxy Dexamethasone

Item No. 22099

CAS Registry No.: 55879-47-5
Formal Name: 9-fluoro-6 β ,11 β ,17,21-tetrahydroxy-16 α -methyl-pregna-1,4-diene-3,20-dione
MF: C₂₂H₂₉FO₆
FW: 408.5
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

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

Description

6 β -hydroxy Dexamethasone is an active metabolite of the synthetic glucocorticoid dexamethasone (Item No. 11015).¹ It is formed from dexamethasone by the cytochrome P450 (CYP) isoform CYP3A4. 6 β -hydroxy Dexamethasone (1 mg/kg) reduces cortisol (hydrocortisone; Item No. 20739) production in rats.²

References

1. Gentile, D.M., Tomlinson, E.S., Maggs, J.L., *et al.* Dexamethasone metabolism by human liver in vitro. Metabolite identification and inhibition of 6-hydroxylation. *J. Pharmacol. Exp. Ther.* **277**(1), 105-112 (1996).
2. Ritchie, J.C., Owens, M.J., Mayer, H., *et al.* Preliminary studies of 6 β -hydroxydexamethasone and its importance in the DST. *Biol. Psychiatry* **32**(9), 825-833 (1992).

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, 03/28/2023

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