

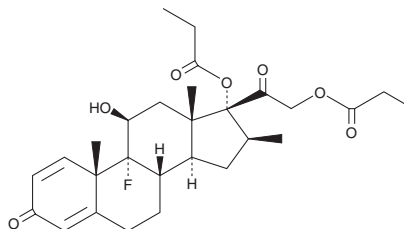
PRODUCT INFORMATION



Betamethasone Dipropionate

Item No. 39499

CAS Registry No.: 5593-20-4
Formal Name: 9-fluoro-11 β -hydroxy-16 β -methyl-17,21-bis(1-oxopropoxy)-pregna-1,4-diene-3,20-dione
Synonyms: Betamethasone 17,21-Dipropionate, β -Methasone Dipropionate, SCH 11460
MF: C₂₈H₃₇FO₇
FW: 504.6
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

Betamethasone dipropionate is supplied as a solid. A stock solution may be made by dissolving the betamethasone dipropionate in the solvent of choice, which should be purged with an inert gas. Betamethasone dipropionate is soluble in acetonitrile and DMSO.

Description

Betamethasone dipropionate is a synthetic corticosteroid and prodrug form of betamethasone (Item No. 20363).^{1,2} It binds to the glucocorticoid receptor (IC₅₀ = 5.25 nM).¹ Betamethasone dipropionate decreases the number of circulating eosinophils in dogs and reduces croton oil-induced exudate formation in a granuloma pouch assay in rats.³ Formulations containing betamethasone dipropionate have been used in the treatment of dermatoses.

References

1. Ueno, H., Maruyama, A., Miyake, M., *et al.* Synthesis and evaluation of antiinflammatory activities of a series of corticosteroid 17 α -esters containing a functional group. *J. Med. Chem.* **34(8)**, 2468-2473 (1991).
2. Li, L., Zhang, M., P., M., *et al.* Development of a simple method for simultaneous determination of tazarotene and betamethasone dipropionate and their metabolites using LC-MS method and its application to dermatopharmacokinetic study. *Biomed. Chromatogr.* **33(8)**, e4557 (2019).
3. Collins, E.J., Aschenbrenner, J., and Nakahama, M. Biologic action of betamethasone 17,21-dipropionate in rat and dog. *Steroids* **20(5)**, 543-554 (1972).

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, 03/05/2024

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