

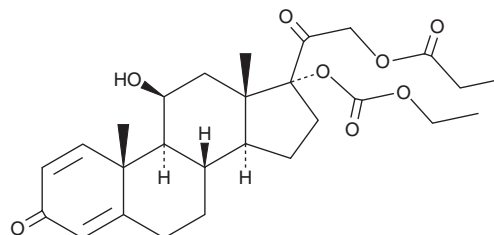
PRODUCT INFORMATION



Prednicarbate

Item No. 23668

CAS Registry No.: 73771-04-7
Formal Name: (11 β)-17-[(ethoxycarbonyl)oxy]-11-hydroxy-21-(1-oxopropoxy)-pregna-1,4-diene-3,20-dione
Synonyms: Prednisolone 17-(ethyl carbonate) 21-propionate, S 77-0777
MF: C₂₇H₃₆O₈
FW: 488.6
Purity: \geq 98%
Supplied as: A solid
Storage: 4°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Prednicarbate is supplied as a solid. A stock solution may be made by dissolving the prednicarbate in the solvent of choice, which should be purged with an inert gas. Prednicarbate is slightly soluble in organic solvents such as chloroform, ethanol, and methanol.

Description

Prednicarbate is a synthetic corticosteroid and a derivative of prednisolone (Item No. 20866).¹ It has anti-inflammatory activity comparable to that of desoximetasone in various murine models when administered topically but lacks activity when administered subcutaneously. Topical administration of prednicarbate (2 drops of a 0.1% solution per day) induces wound healing in hairless mice.² Formulations containing prednicarbate have been used to treat chronic hand eczema.³

References

1. Alpermann, H.G., Sandow, J., and Vogel, H.G. Animal experiment studies on the topical and systemic effectiveness of prednisolone-17-ethylcarbonate-21-propionate. *Z. Hautkr.* **61(1)**, 7-17 (1986).
2. Bodor, N.S., Kiss-Buris, S.T., and Buris, L. Novel soft steroids: Effects on cell growth *in vitro* and on wound healing in the mouse. *Steroids* **56(8)**, 434-439 (1991).
3. Soost, S., Abdollahnia, M., Kostev, K., *et al.* Topical therapy of hand eczema - analysis of the prescription profile from dermatologists in private practice. *J. Dtsch. Dermatol. Ges.* **10(3)**, 180-184 (2012).

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, 10/12/2022

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