

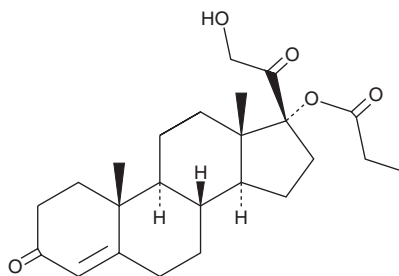
PRODUCT INFORMATION



Cortisolone 17 α -propionate

Item No. 28392

CAS Registry No.: 19608-29-8
Formal Name: 21-hydroxy-17-(1-oxopropoxy)-pregn-4-ene-3,20-dione
Synonym: CB-03-01
MF: C₂₄H₃₄O₅
FW: 402.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 240, 295 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

Cortisolone 17 α -propionate is supplied as a solid. A stock solution may be made by dissolving the cortisolone 17 α -propionate in the solvent of choice, which should be purged with an inert gas. Cortisolone 17 α -propionate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cortisolone 17 α -propionate in DMSO is approximately 10 mg/ml and approximately 20 mg/ml in ethanol and DMF.

Cortisolone 17 α -propionate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cortisolone 17 α -propionate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Cortisolone 17 α -propionate has a solubility of approximately 0.1 mg/ml in a 1:9 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cortisolone 17 α -propionate is a peripherally selective androgen receptor antagonist.¹ Cortisolone 17 α -propionate binds to the androgen receptor (AR) and inhibits AR-regulated transcription in a reporter assay.² It also inhibits androgen-regulated lipid synthesis in primary human sebocytes. Cortisolone 17 α -propionate inhibits inflammatory cytokine production from primary human sebocytes and scalp dermal papilla cells.^{2,3} It inhibits testosterone propionate-induced flank organ enlargement in Syrian golden hamsters by 84% when administered topically at a dose of 400 μ g/animal.

References

1. Celasco, G., Moro, L., Bozzella, R., *et al.* Biological profile of cortisolone 17 α -propionate (CB-03-01), a new topical and peripherally selective androgen antagonist. *Arzneimittelforschung* **54(12)**, 881-886 (2004).
2. Rosette, C., Agan, F.J., Mazzetti, A., *et al.* Cortisolone 17 α -propionate (clascoterone) is a novel androgen receptor antagonist that inhibits production of lipids and inflammatory cytokines from sebocytes in vitro. *J. Drugs Dermatol.* **18(5)**, 412-418 (2019).
3. Rosette, C., Rosette, N., Mazezzetti, A., *et al.* Cortisolone 17 α -propionate (clascoterone) is an androgen receptor antagonist in dermal papilla cells in vitro. *J. Drugs Dermatol.* **18(2)**, 197-201 (2019).

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, 11/21/2022

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