

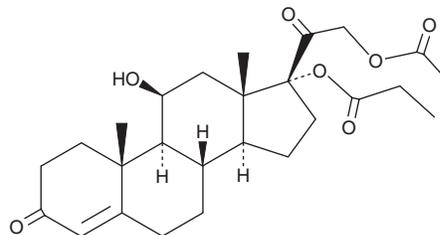
PRODUCT INFORMATION



Hydrocortisone Aceponate

Item No. 35359

CAS Registry No.: 74050-20-7
Formal Name: (11 β)-21-(acetyloxy)-11-hydroxy-17-(1-oxopropoxy)-pregn-4-ene-3,20-dione
Synonym: Hydrocortisone 17-Propionate 21-Acetate
MF: C₂₆H₃₆O₇
FW: 460.6
Purity: \geq 98%
UV/Vis.: λ_{max} : 242 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

Hydrocortisone aceponate is supplied as a solid. A stock solution may be made by dissolving the hydrocortisone aceponate in the solvent of choice, which should be purged with an inert gas. Hydrocortisone aceponate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of hydrocortisone aceponate in DMF is approximately 3 mg/ml and approximately 10 mg/ml in ethanol and DMSO.

Description

Hydrocortisone aceponate is a corticosteroid.^{1,2} Topical application of hydrocortisone aceponate (0.0584%) reduces lesion size and pruritis in cats with allergic dermatitis.¹ It also induces remission and increases time to relapse in dogs with spontaneous atopic dermatitis.² Formulations containing hydrocortisone aceponate have been used in the treatment of veterinary dermatoses.

References

- Schmidt, V., Buckley, L.M., McEwan, N.A., *et al.* Efficacy of a 0.0584% hydrocortisone aceponate spray in presumed feline allergic dermatitis: An open label pilot study. *Vet. Dermatol.* **23(1)**, 11-16 (2012).
- Lourenço, A.M., Schmidt, V., Braz, B.S., *et al.* Efficacy of proactive long-term maintenance therapy of canine atopic dermatitis with 0.0584% hydrocortisone aceponate spray: A double-blind placebo controlled pilot study. *Vet. Dermatol.* **27(2)**, 88e25-92e25 (2016).

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, 12/12/2022

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