

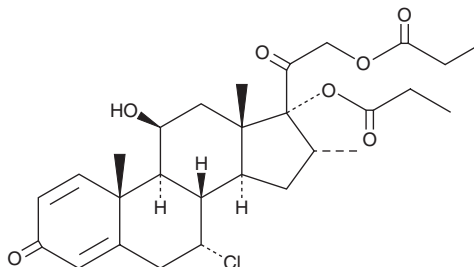
PRODUCT INFORMATION



Alclometasone Dipropionate

Item No. 23901

CAS Registry No.: 66734-13-2
Formal Name: 7 α -chloro-11 β -hydroxy-16 α -methyl-17,21-bis(1-oxopropoxy)-pregna-1,4-diene-3,20-dione
Synonyms: ACM, Alclometasone
17,21-Dipropionate, SCH 22219
MF: C₂₈H₃₇ClO₇
FW: 521.0
Purity: \geq 98%
UV/Vis.: λ_{max} : 244 nm
Supplied as: A crystalline 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

Alclometasone dipropionate (ACM) is supplied as a crystalline solid. A stock solution may be made by dissolving the ACM in the solvent of choice. ACM is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ACM in these solvents is approximately 30 mg/ml.

ACM is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ACM should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. ACM has a solubility of approximately 0.2 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ACM is a corticosteroid.¹ It reduces cutaneous anaphylaxis reactions induced by tuberculin or albumin in mice when administered topically (20 μ l of a 0.1% solution). ACM also inhibits androgen-dependent cytochrome P450 activity and the O-depropylation activity of 7-alkoxy-coumarin O-dealkylase in a dose-dependent manner in male rats but has no effect on hepatic drug metabolism in female rats or mice of both sexes.² Formulations containing ACM have been used to treat radiation and allergic contact dermatitis.

References

1. Sengoku, T., Morita, K., Sakuma, S., *et al.* Possible inhibitory mechanism of FK506 (tacrolimus hydrate) ointment for atopic dermatitis based on animal models. *Eur. J. Pharmacol.* **379(2-3)**, 183-189 (1999).
2. Nishibe, Y., Hamataka, M., Hasegawa, T., *et al.* Species and sex differences in the inhibitory action of the corticosteroid alclometasone dipropionate on the hepatic drug-metabolizing system. *Japan J. Pharmacol.* **50(4)**, 435-443 (1989).

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, 12/22/2022

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