

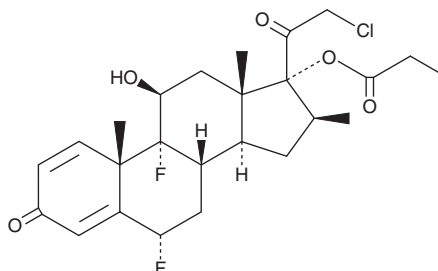
PRODUCT INFORMATION



Halobetasol Propionate

Item No. 23781

CAS Registry No.: 66852-54-8
Formal Name: 21-chloro-6 α ,9-difluoro-11 β -hydroxy-16 β -methyl-17-(1-oxopropoxy)-pregna-1,4-diene-3,20-dione
Synonyms: BMY 30056, CGP 14458
MF: C₂₅H₃₁ClF₂O₅
FW: 485.0
Purity: \geq 95%
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

Halobetasol propionate is supplied as a solid. A stock solution may be made by dissolving the halobetasol propionate in the solvent of choice, which should be purged with an inert gas. Halobetasol propionate is slightly soluble in methanol and DMSO.

Description

Halobetasol propionate is a synthetic glucocorticoid.¹ It inhibits dermatitis induced by ultraviolet light in guinea pigs (ED₅₀ = 1.45 μ g/animal) and by oxazolone in mice (ED₅₀ = 0.8 μ g/animal).² Halobetasol propionate also has anti-inflammatory effects on croton oil-induced ear edema in rats (ED₅₀s = 10 μ g/ml and 0.5 μ g/ear) and mice (ED₅₀s = 1.2 and <0.32 μ g/ear after 6 and 24 hours of treatment, respectively). Topical formulations containing halobetasol propionate have been used for inflammatory and pruritic dermatitis.

References

1. Avery, A.M. and Chittiboyina, A.G. Anti-inflammatory glucocorticoids. In *Burger's medicinal chemistry, drug discovery and development*. 35-152, 7th edition, Wiley (2010).
2. Yawalkar, S., Wiesenberg-Boettcher, I., Gibson, J.R., et al. Dermatopharmacologic investigations of halobetasol propionate in comparison with clobetasol 17-propionate. *J. Am. Acad. Dermatol.* **25**(6), 1137-1144 (1991).

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, 12/19/2022

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