

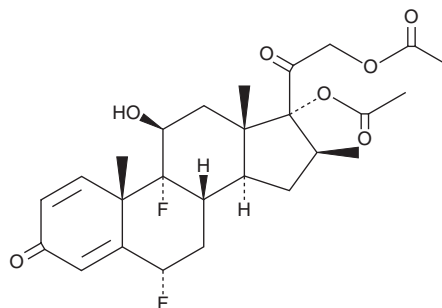
PRODUCT INFORMATION



Diflorasone Diacetate

Item No. 23808

CAS Registry No.: 33564-31-7
Formal Name: 17,21-bis(acetyloxy)-6 α ,9-difluoro-11 β -hydroxy-16 β -methyl-pregna-1,4-diene-3,20-dione
Synonym: U-34865
MF: C₂₆H₃₂F₂O₇
FW: 494.5
Purity: \geq 98%
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

Diflorasone diacetate is supplied as a solid. A stock solution may be made by dissolving the diflorasone diacetate in the solvent of choice, which should be purged with an inert gas. Diflorasone diacetate is slightly soluble in methanol and chloroform.

Description

Diflorasone is a corticosteroid with anti-inflammatory activity.¹ It reduces production of the inflammatory cytokines Il-6, Il-4, and macrophage inflammatory protein-2 (Mip-2) and completely inhibits swelling in the ears of toluene-2,4-diisocyanate-sensitized mice following topical administration of 20 μ L of a 0.05% solution. Diflorasone inhibits RANTES and thymus and activation regulated chemokine (TARC) in a dinitrochlorobenzene-induced mouse model of allergic contact dermatitis.² It also inhibits proliferation of human keratinocytes *in vitro* and reduces epidermal thickening in a mouse model of psoriasis induced by imiquimod (Item No. 14956).³

References

1. Bäumer, W., Gorr, G., Hoppmann, J., *et al.* AWD 12-281, a highly selective phosphodiesterase 4 inhibitor, is effective in the prevention and treatment of inflammatory reactions in a model of allergic dermatitis. *J. Pharm. Pharmacol.* **55(8)**, 1107-1114 (2003).
2. Bäumer, W., Seegers, U., Braun, M., *et al.* TARC and RANTES, but not CTACK, are induced in two models of allergic contact dermatitis. Effects of cilmilast and diflorasone diacetate on T-cell-attracting chemokines. *Br. J. Dermatol.* **151(4)**, 823-830 (2004).
3. Schaper, K., Dickhaut, J., Japtok, L., *et al.* Sphingosine-1-phosphate exhibits anti-proliferative and anti-inflammatory effects in mouse models of psoriasis. *J. Dermatol. Sci.* **71(1)**, 29-36 (2013).

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