

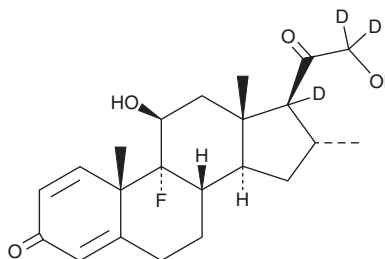
PRODUCT INFORMATION



Desoximetasone-d₃

Item No. 28159

Formal Name: 9-fluoro-11β,21-dihydroxy-16α-methylpregna-1,4-diene-3,20-dione-d₃
Synonym: 17-deoxy Dexamethasone-d₃
MF: C₂₂H₂₆D₃FO₄
FW: 379.5
Chemical Purity: ≥98% (Desoximetasone)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Desoximetasone-d₃ is intended for use as an internal standard for the quantification of desoximetasone (Item No. 23945) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Desoximetasone-d₃ is supplied as a solid. A stock solution may be made by dissolving the desoximetasone-d₃ in the solvent of choice, which should be purged with an inert gas. Desoximetasone-d₃ is soluble in organic solvents such as methanol and DMSO.

Description

Desoximetasone is a glucocorticoid.¹ *In vivo*, desoximetasone (0.001-0.1 mg) reverses increases in epidermal cell proliferation induced by phorbol 12-myristate 13-acetate (TPA; Item No. 10008014) in SENCAR mice when administered topically. It also reduces TPA-induced expression of IL-6, COX-2, iNOS, and c-Jun mRNA levels in the skin of SENCAR mice, indicating anti-inflammatory action. Formulations containing desoximetasone have been used in the treatment of psoriasis and atopic dermatitis.

Reference

1. Kowalczyk, P., Junco, J.J., Kowalczyk, M.C., *et al.* The effects of dissociated glucocorticoids RU24858 and RU24782 on TPA-induced skin tumor promotion biomarkers in SENCAR mice. *Mol. Carcinog.* **53(6)**, 488-497 (2014).

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, 12/05/2022

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