

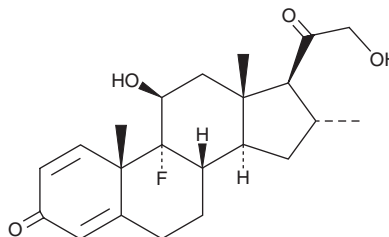
# PRODUCT INFORMATION



## Desoximetasone

Item No. 23945

**CAS Registry No.:** 382-67-2  
**Formal Name:** 9-fluoro-11 $\beta$ ,21-dihydroxy-16 $\alpha$ -methyl-pregna-1,4-diene-3,20-dione  
**Synonyms:** A 41-304,  
17-deoxy Dexamethasone,  
HOE 304  
**MF:** C<sub>22</sub>H<sub>29</sub>FO<sub>4</sub>  
**FW:** 376.5  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 237 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

Desoximetasone is supplied as a crystalline solid. A stock solution may be made by dissolving the desoximetasone in the solvent of choice. Desoximetasone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of desoximetasone in these solvents is approximately 3, 30, and 25 mg/ml, respectively.

Desoximetasone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, desoximetasone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Desoximetasone has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Desoximetasone is a glucocorticoid.<sup>1</sup> *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, 12/22/2022

#### 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