

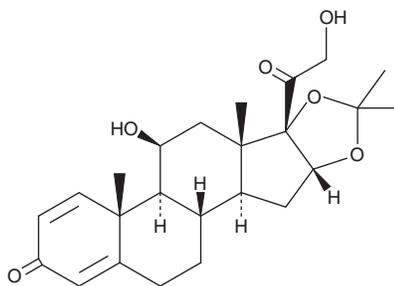
PRODUCT INFORMATION



Desonide

Item No. 21368

CAS Registry No.: 638-94-8
Formal Name: (11 β ,16 α)-11,21-dihydroxy-16,17-
[[1-methylethylidene]bis(oxy)]-
pregna-1,4-diene-3,20-dione
Synonym: Budesonide acetonide
MF: C₂₄H₃₂O₆
FW: 416.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 243 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

Desonide is supplied as a crystalline solid. A stock solution may be made by dissolving the desonide in the solvent of choice, which should be purged with an inert gas. Desonide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of desonide in these solvents is approximately 10, 25, and 20 mg/ml, respectively.

Desonide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, desonide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Desonide has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Desonide is a synthetic corticosteroid.¹ It is a nonfluorinated analog of triamcinolone acetonide (Item No. 18026). Formulations containing desonide are used topically to reduce inflammation in conditions that include allergic reactions and psoriasis.^{1,2}

References

1. Stewart, W.D., Runikis, J.O., Verma, S.C., *et al.* Problems in selection of topical anti-inflammatory corticosteroids. *Can. Med. Assoc. J.* **108(1)**, 33-38 (1973).
2. Iqbal, J., Husain, A., and Gupta, A. Photochemistry of desonide, a non-fluorinated steroidal anti-inflammatory drug. *Chem. Pharm. Bull. (Tokyo)* **54(6)**, 836-838 (2006).

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, 11/30/2022

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