

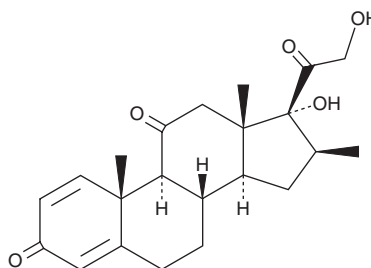
# PRODUCT INFORMATION



## Meprednisone

Item No. 30289

**CAS Registry No.:** 1247-42-3  
**Formal Name:** (16 $\beta$ )-17,21-dihydroxy-16-methyl-pregna-1,4-diene-3,11,20-trione  
**Synonyms:** Methylprednisone, NSC 527579, SCH 4358  
**MF:** C<sub>22</sub>H<sub>28</sub>O<sub>5</sub>  
**FW:** 372.5  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 238 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

Meprednisone is supplied as a crystalline solid. A stock solution may be made by dissolving the meprednisone in the solvent of choice, which should be purged with an inert gas. Meprednisone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of meprednisone in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Meprednisone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, meprednisone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Meprednisone has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Meprednisone is a corticosteroid.<sup>1</sup> It is a methylated form of prednisone (Item No. 20677) and a metabolite of methylprednisolone (Item No. 15013).<sup>1,2</sup> Formulations containing meprednisone have been used in the treatment of drug-induced or seasonal allergies.

### References

1. Jacob, S.E. and Steele, T. Corticosteroid classes: A quick reference guide including patch test substances and cross-reactivity. *J. Am. Acad. Dermatol.* **54**(4), 723-727 (2006).
2. Haughey, D.B. and Jusko, W.J. Effect of ketoconazole on methylprednisolone pharmacokinetics and receptor/gene-mediated pharmacodynamics. *J. Pharmacol. Exp. Ther.* **258**(2), 826-832 (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, 10/10/2022

#### 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