

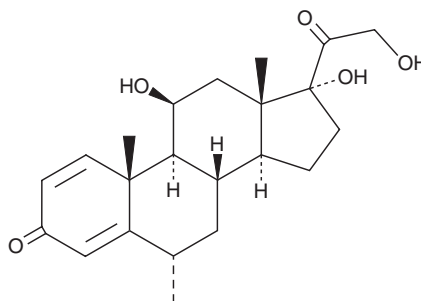
PRODUCT INFORMATION



Methylprednisolone

Item No. 15013

CAS Registry No.: 83-43-2
Formal Name: 11 β ,17,21-trihydroxy-6 α -methyl-pregna-1,4-diene-3,20-dione
Synonyms: 6 α -Methylprednisolone, NSC 19987, U-7532
MF: C₂₂H₃₀O₅
FW: 374.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 243 nm
Supplied as: A crystalline 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

Methylprednisolone is supplied as a crystalline solid. A stock solution may be made by dissolving the methylprednisolone in the solvent of choice. Methylprednisolone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of methylprednisolone in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Methylprednisolone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methylprednisolone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Methylprednisolone 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

Methylprednisolone is a synthetic glucocorticoid, an agonist of glucocorticoid and mineralocorticoid receptors, and an active metabolite of the prodrug 6 α -methylprednisolone 21-hemisuccinate (MPS; Item Nos. 11800 | 35742).¹ It transactivates glucocorticoid and mineralocorticoid receptors in a reporter assay using CV-1 cells (EC₅₀s = 2.92 and 2.31 nM for the human receptors, respectively). Methylprednisolone (40 mg/kg) reduces endotoxin-induced complement activation in rabbits.² It induces T cell apoptosis and reduces the number of T cell spinal cord infiltrates in a rat model of experimental autoimmune encephalomyelitis (EAE) when administered at a dose of 50 mg/kg.³ It decreases the number of reactive spinal astrocytes, improves proprioceptive limb placing and coordination, and reduces spinal edema and the number of spinal necrotic cells in a rat model of spinal crush injury.⁴ Formulations containing methylprednisolone have been used in the treatment of immune and autoimmune disorders, such as rheumatoid arthritis, systemic lupus erythematosus, psoriasis, corneal ulcers, ulcerative colitis, and allergies, among others.

References

1. Grossmann, C., Scholz, T., Rochel, M., et al. *Eur. J. Endocrinol.* **151**(3), 397-406 (2004).
2. Bult, H., Herman, A.G., and Rampart, M. *Br. J. Pharmacol.* **84**(2), 317-327 (1985).
3. Schmidt, J., Gold, R., Schönrock, L., et al. *Brain* **123**(Pt 7), 1431-1441 (2000).
4. Jiang, S., Khan, M.I., Middlemiss, P.J., et al. *Int. J. Immunopathol. Pharmacol.* **17**(3), 353-366 (2004).

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, 09/20/2023

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