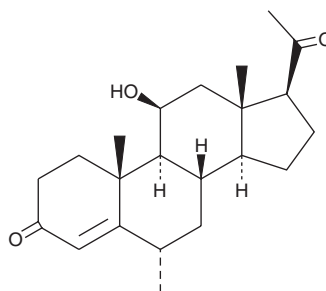


PRODUCT INFORMATION

6 α -methyl-11 β -hydroxy Progesterone

Item No. 19533

CAS Registry No.: 2668-66-8
Formal Name: 11 β -hydroxy-6 α -methyl-pregn-4-ene-3,20-dione
Synonyms: Medrysone, NSC 63278, Visudrisone
MF: C₂₂H₃₂O₃
FW: 344.5
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 242 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

6 α -methyl-11 β -hydroxy Progesterone is supplied as a crystalline solid. A stock solution may be made by dissolving the 6 α -methyl-11 β -hydroxy progesterone in the solvent of choice, which should be purged with an inert gas. 6 α -methyl-11 β -hydroxy Progesterone is soluble in organic solvents such as ethanol, methanol, and acetonitrile. The solubility of 6 α -methyl-11 β -hydroxy progesterone in these solvents is approximately 1 mg/ml.

Description

6 α -methyl-11 β -hydroxy Progesterone is an anti-inflammatory steroid.¹ It is a hydroxylated progesterone that lacks progestational and androgenic activity. It inhibits phytohemagglutinin-activated T cell proliferation *in vitro* with an IC₅₀ value of 2.9 $\mu\text{g/ml}$.² 6 α -methyl-11 β -hydroxy Progesterone also inhibits TNF- α production in LPS-stimulated THP-1 cells (IC₅₀ = 30.54 $\mu\text{g/ml}$). Topical formulations containing 6 α -methyl-11 β -hydroxyprogesterone have been used to treat inflammatory ocular diseases.¹

References

- Schwartz, A., and Leopold, I.H. Topical anti-inflammatory steroids and intraocular pressure: The place of medrysone. *Drugs* **2**(1), 1-4 (1971).
- Bano, S., Wahab, A.-t., Yousuf, S., *et al.* New anti-inflammatory metabolites by microbial transformation of medrysone. *PLoS One* **11**(4), e0153951 (2016).

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, 11/29/2022

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