

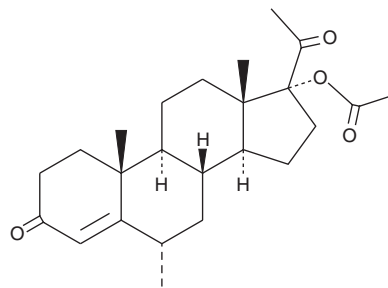
PRODUCT INFORMATION



Medroxyprogesterone 17-acetate

Item No. 23664

CAS Registry No.: 71-58-9
Formal Name: 17-(acetyloxy)-6 α -methyl-pregn-4-ene-3,20-dione
Synonyms: NSC 21171, NSC 26386, U-8839
MF: C₂₄H₃₄O₄
FW: 386.5
Purity: \geq 98%
Supplied as: A 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

Medroxyprogesterone 17-acetate is supplied as a solid. A stock solution may be made by dissolving the medroxyprogesterone 17-acetate in the solvent of choice. Medroxyprogesterone 17-acetate is slightly soluble in methanol and chloroform.

Description

Medroxyprogesterone 17-acetate is a synthetic progestogen.¹⁻³ It prevents fertilization and increases the rate of transport of eggs from the fallopian tubes to the uterus in female ferrets when administered prior to ovulation.¹ Medroxyprogesterone 17-acetate reversibly blocks ovulation in rats when injected on the last day of diestrus.² It also has anti-androgenic activity in rats, decreasing plasma testosterone (Item Nos. 15645 | ISO60154) levels via induction of hepatic testosterone reductase activity.³ Medroxyprogesterone 17-acetate exhibits immunosuppressive effects *in vitro* and *in vivo*, inhibiting the production of IFN- γ by CD2/CD3/CD28-stimulated peripheral blood mononuclear cells (PBMCs) at concentrations \geq 10 nM and extending the survival of rabbit skin allografts.^{4,5} Injectable formulations containing medroxyprogesterone 17-acetate have been used as contraceptives.⁶

References

1. Chang, M.C. Effects of medroxyprogesterone acetate and of ethinyl oestradiol on the fertilization and transportation of ferret eggs. *J. Reprod. Fertil.* **13(1)**, 173-174 (1967).
2. Dickmann, Z. Short- and long-term effects of a single injection of depo-medroxyprogesterone acetate (provera) on the vaginal smear, ovulation and mating in the rat. *J. Reprod. Fertil.* **32(3)**, 447-451 (1973).
3. Albin, J., Vittek, J., Gordon, G.G., *et al.* On the mechanism of the anti-androgenic effect of medroxyprogesterone acetate. *Endocrinology* **93(2)** (1973).
4. Huijbregts, R.P., Michel, K.G., and Hel, Z. Effect of progestins on immunity: Medroxyprogesterone but not norethisterone or levonorgestrel suppresses the function of T cells and pDCs. *Contraception* **90(2)**, 123-129 (2014).
5. Turcotte, J.G., Haines, R.F., Brody, G.L., *et al.* Immunosuppression with medroxyprogesterone acetate. *Transplantation* **6(2)**, 248-260 (1968).
6. Hofmeyr, G.J., Singata-Madliki, M., Lawrie, T.A., *et al.* Effects of the copper intrauterine device versus injectable progestin contraception on pregnancy rates and method discontinuation among women attending termination of pregnancy services in South Africa: A pragmatic randomized controlled trial. *Reprod. Health* **13**, 42 (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, 10/25/2022

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