

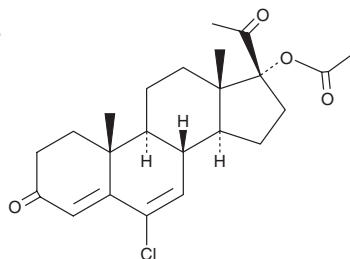
PRODUCT INFORMATION



Chlormadinone Acetate

Item No. 29392

CAS Registry No.: 302-22-7
Formal Name: 17-(acetyloxy)-6-chloro-pregna-4,6-diene-3,20-dione
Synonym: NSC 92338
MF: C₂₃H₂₉ClO₄
FW: 404.9
Purity: ≥98%
UV/Vis.: λ_{max}: 283 nm
Supplied as: A 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

Chlormadinone acetate is supplied as a solid. A stock solution may be made by dissolving the chlormadinone acetate in the solvent of choice, which should be purged with an inert gas. Chlormadinone acetate is soluble in the organic solvent chloroform at a concentration of approximately 50 mg/ml.

Description

Chlormadinone acetate is a synthetic progestin.¹ It binds to progesterone, androgen, and glucocorticoid receptors *in vitro* (K_s = 2.5, 3.8, and 16 nM, respectively, for the human receptors). Chlormadinone acetate increases the number of endometrial glands and uterine weight in β-estradiol-primed rabbits when administered at a dose of 45 μg/kg per day for five days. Chlormadinone acetate reduces testosterone-induced increases in the seminal vesicle weight of castrated male rats when administered at doses of 4.6 and 21.5 mg/kg per day for eight days.

Reference

1. Schneider, J., Kneip, C., and Jahnel, U. Comparative effects of chlormadinone acetate and its 3α- and 3β-hydroxy metabolites on progesterone, androgen and glucocorticoid receptors. *Pharmacology* **84**(2), 74-81 (2009).

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, 10/05/2022

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