

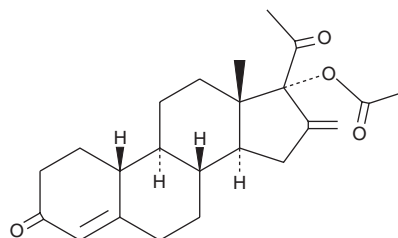
PRODUCT INFORMATION



Segesterone Acetate

Item No. 26441

CAS Registry No.: 7759-35-5
Formal Name: 17-(acetyloxy)-16-methylene-19-norpregn-4-ene-3,20-dione
Synonym: ST-1435
MF: $C_{23}H_{30}O_4$
FW: 370.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Segesterone acetate is supplied as a solid. A stock solution may be made by dissolving the segesterone acetate in the solvent of choice, which should be purged with an inert gas. Segesterone acetate is soluble in DMSO.

Description

Segesterone acetate is a synthetic progestin that binds to the progesterone receptor ($EC_{50} = 50.3$ nM).¹ It increases uterine weight in rabbits ($ED_{50} = \sim 1$ μ g), maintains pregnancy in 40-50% of rats when administered at doses ranging from 0.03 to 0.1 mg per day, and completely inhibits spontaneous ovulation in rats at a dose of 10 μ g per day. It also binds to the glucocorticoid receptor ($EC_{50} = 56$ nM) but does not have *in vivo* glucocorticoid activity. Segesterone acetate (8 μ g, i.c.v., *via* osmotic minipump over four weeks) increases cell proliferation in the adult female mouse brain, increasing the production of new neurons and oligodendrocytes.² Sustained-release formulations containing segesterone have been used as contraceptives.

References

1. Kumar, N., Koide, S.S., Tsong, Y.-Y., *et al.* Nestorone®: A progestin with a unique pharmacological profile. *Steroids* **65(10-11)**, 629-636 (2000).
2. Chen, S., Kumar, N., Mao, Z., *et al.* Therapeutic progestin segesterone acetate promotes neurogenesis: Implications for sustaining regeneration in female brain. *Menopause* **25(10)**, 1138-1151 (2018).

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, 10/11/2022

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