

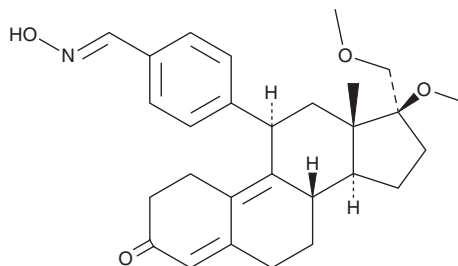
PRODUCT INFORMATION



Asoprisnil

Item No. 35170

CAS Registry No.: 199396-76-4
Formal Name: [C(E)]-4-[(11 β ,17 β)-17-methoxy-17-(methoxymethyl)-3-oxoestra-4,9-dien-11-yl]-benzaldehyde, 1-oxime
Synonym: J867
MF: C₂₈H₃₅NO₄
FW: 449.6
Purity: \geq 95%
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

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

Description

Asoprisnil is a selective progesterone receptor modulator (SPRM).¹ It selectively binds to the progesterone receptor over the glucocorticoid receptor. Asoprisnil inhibits the proliferation of vaginal epithelium and induces mucification of the vaginal epithelium in guinea pigs. It also inhibits ovulation in guinea pigs when administered at doses greater than 1 mg/kg.²

References

- Schubert, G., Elger, W., Kaufmann, G., *et al.* Discovery, chemistry, and reproductive pharmacology of asoprisnil and related 11 β -benzaloxime substituted selective progesterone receptor modulators (SPRMs). *Semin. Reprod. Med.* **23(1)**, 58-73 (2005).
- DeManno, D., Elger, W., Garg, R., *et al.* Asoprisnil (J867): A selective progesterone receptor modulator for gynecological therapy. *Steroids* **68(10-13)**, 1019-1032 (2003).

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, 12/12/2022

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