

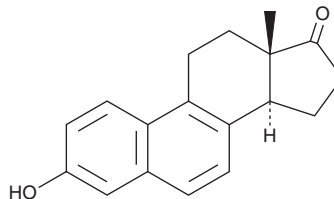
PRODUCT INFORMATION



Equilenin

Item No. 23322

CAS Registry No.: 517-09-9
Formal Name: 3-hydroxy-estra-1,3,5,7,9-pentaen-17-one
Synonyms: (+)-Equilenin, d-Equilenin, NSC 9901
MF: C₁₈H₁₈O₂
FW: 266.3
Purity: ≥98%
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

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

Description

Equilenin is an equine estrogen found in the urine of pregnant mares.¹ Unlike naturally occurring human estrogens, equilenin has no effect on nitric oxide synthase (eNOS) activity or expression and does not increase production of nitric oxide (NO). It is metabolized to 4-hydroxyequilenin, a carcinogenic compound that induces transformation of C3H 10T½ murine fibroblasts.² Formulations containing equilenin conjugates have been widely used as hormone replacement therapies for post-menopausal women.

References

1. Novensa, L., Selent, J., Pastor, M., *et al.* Equine estrogens impair nitric oxide production and endothelial nitric oxide synthase transcription in human endothelial cells compared with the natural 17β-estradiol. *Hypertension* **56**(3), 405-411 (2010).
2. Spink, D.C., Zhang, F., Hussain, M.M., *et al.* Metabolism of equilenin in MCF-7 and MDA-MB-231 human breast cancer cells. *Chem. Res. Toxicol.* **14**(5), 572-581 (2001).

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, 10/04/2022

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