

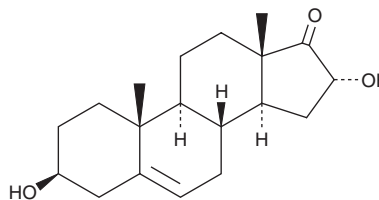
PRODUCT INFORMATION



16 α -hydroxy Dehydroepiandrosterone

Item No. 34116

CAS Registry No.: 1232-73-1
Formal Name: 3,16-dihydroxy-(3 β ,16 α)-androst-5-en-17-one
Synonym: 16 α -hydroxy DHEA
MF: C₁₉H₂₈O₃
FW: 304.4
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

16 α -hydroxy Dehydroepiandrosterone is supplied as a solid. A stock solution may be made by dissolving the 16 α -hydroxy dehydroepiandrosterone in the solvent of choice, which should be purged with an inert gas. 16 α -hydroxy Dehydroepiandrosterone is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

16 α -hydroxy Dehydroepiandrosterone is a metabolite of the endogenous steroid hormone dehydroepiandrosterone (Item No. 15728).¹ 16 α -hydroxy Dehydroepiandrosterone is formed from dehydroepiandrosterone via 16-hydroxylation by the cytochrome P450 (CYP) isoforms CYP3A4 and CYP3A5 in adult human liver microsomes, as well as by fetal recombinant CYP3A7. It is a precursor to fetal estrogens, including estriol (Item No. 10006484).²

References

1. Miller, K.K.M., Cai, J., Ripp, S.L., *et al.* Stereo- and regioselectivity account for the diversity of dehydroepiandrosterone (DHEA) metabolites produced by liver microsomal cytochromes P450. *Drug Metab. Dispos.* **32(3)**, 305-313 (2004).
2. Hampl, R. and Starka, L. Minireview: 16 α -Hydroxylated metabolites of dehydroepiandrosterone and their biological significance. *Endocr. Regul.* **34(3)**, 161-163 (2000).

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, 02/06/2024

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