

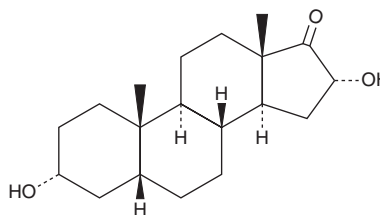
PRODUCT INFORMATION



16 α -hydroxy Etiocholanolone

Item No. 27683

CAS Registry No.: 14167-50-1
Formal Name: (3 α ,5 β ,16 α)-3,16-dihydroxy-androstan-17-one
MF: C₁₉H₃₀O₃
FW: 306.4
Purity: \geq 95%
Supplied as: A crystalline 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 Etiocholanolone is supplied as a crystalline solid. A stock solution may be made by dissolving the 16 α -hydroxy etiocholanolone in the solvent of choice, which should be purged with an inert gas. 16 α -hydroxy Etiocholanolone is soluble in organic solvents such as acetonitrile, ethanol, and methanol. The solubility of 16 α -hydroxy etiocholanolone in these solvents is approximately 1 mg/ml.

Description

16 α -hydroxy Etiocholanolone is a metabolite of 16 α -hydroxydehydroisoandrosterone (16 α -DHEA) and androstenedione.^{1,2}

References

1. Lai, E.Y. and Solomon, S. The *in vivo* metabolism of 16 α -hydroxydehydroisoandrosterone in man. *Biochemistry* **6**(7), 2040-2052 (1967).
2. Christakoudi, S., Cowan, D.A., Christakudis, G., et al. 21-hydroxylase deficiency in the neonate - trends in steroid anabolism and catabolism during the first weeks of life. *J. Steroid Biochem. Mol. Biol.* **138**, 334-347 (2013).

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, 12/22/2022

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