

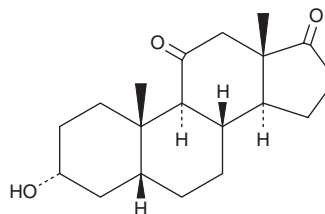
PRODUCT INFORMATION



11-oxo Etiocholanolone

Item No. 35009

CAS Registry No.: 739-27-5
Formal Name: (5 β)-3 α -hydroxy-androstane-11,17-dione
Synonyms: 5 β -Androstane-3 α -ol-11,17-dione,
11-keto Etiocholanolone, NSC 53896
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

11-oxo Etiocholanolone is supplied as a solid. A stock solution may be made by dissolving the 11-oxo etiocholanolone in the solvent of choice, which should be purged with an inert gas. 11-oxo Etiocholanolone is soluble in ethanol.

Description

11-oxo Etiocholanolone is a metabolite of cortisol (hydrocortisone; Item No. 20739) and 11-keto testosterone.^{1,2} It is formed from cortisol *via* cortisone (Item No. 30763) and tetrahydrocortistone (Item No. 34344) intermediates, and it is formed from 11-keto testosterone *via* 5 β -11-keto dihydrotestosterone and 3 α -11-keto etiocholanediol intermediates. Urinary levels of 11-oxo etiocholanolone increase during pregnancy and are elevated in patients with uterine leiomyomas.^{3,4} Fecal levels of 11-oxo etiocholanolone have been used as a marker of stress in domestic livestock and free-range ruminants.⁵

References

- Schiffer, L., Barnard, L., Baranowski, E.S., *et al.* Human steroid biosynthesis, metabolism and excretion are differentially reflected by serum and urine steroid metabolomes: A comprehensive review. *J. Steroid Biochem. Mol. Biol.* **194**, 105439 (2019).
- Barnard, L., Nikolaou, N., Louw, C., *et al.* The A-ring reduction of 11-ketotestosterone is efficiently catalysed by AKR1D1 and SRD5A2 but not SRD5A1. *J. Steroid Biochem. Mol. Biol.* **202**, 105724 (2020).
- Birke, G., Gemzell, C.A., Plantin, L.O., *et al.* Plasma levels of 17-hydroxycorticosteroids and urinary excretion pattern of keto-steroids in normal pregnancy. *Acta Endocrinol. (Copenh.)* **27(4)**, 389-402 (1958).
- Jung, B.H., Bai, S.W., and Chung, B.C. Endogenous urinary steroids in premenopausal women with uterine leiomyomas. *Int. J. Gynaecol. Obstet.* **84(1)**, 55-60 (2004).
- Molina-García, L., Pérez, J.M., Sarasa, M., *et al.* HPLC-QTOF method for quantifying 11-ketoetiocholanolone, a cortisol metabolite, in ruminants' feces: Optimization and validation. *Ecol. Evol.* **8(18)**, 9218-9228 (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, 11/03/2022

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