

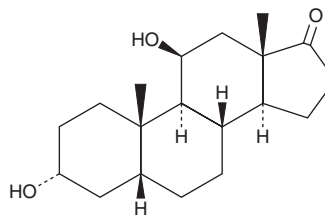
# PRODUCT INFORMATION



## 11 $\beta$ -hydroxy Etiocholanolone

Item No. 33190

CAS Registry No.: 739-26-4  
Formal Name: (5 $\beta$ )-3 $\alpha$ ,11 $\beta$ -dihydroxy-androstan-17-one  
Synonym: NSC 53897  
MF: C<sub>19</sub>H<sub>30</sub>O<sub>3</sub>  
FW: 306.4  
Purity:  $\geq 98\%$   
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

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

### Description

11 $\beta$ -hydroxy Etiocholanolone is a metabolite of hydrocortisone (cortisol; Item No. 20739).<sup>1</sup> Fecal levels of 11 $\beta$ -hydroxy etiocholanolone are increased in response to environmental stress in giraffes (*G. camelopardalis*) and plains zebras (*E. quagga*).<sup>2,3</sup>

### References

1. Shackleton, C.H.L., Neres, M.S., Hughes, B.A., *et al.* 17-Hydroxylase/C17,20-lyase (CYP17) is not the enzyme responsible for side-chain cleavage of cortisol and its metabolites. *Steroids* **73**(6), 652-656 (2008).
2. Bashaw, M.J., Sicks, F., Palme, R., *et al.* Non-invasive assessment of adrenocortical activity as a measure of stress in giraffe (*Giraffa camelopardalis*). *BMC Vet. Res.* **12**(1), 235 (2016).
3. Seeber, P.A., Franz, M., Dehnhard, M., *et al.* Plains zebra (*Equus quagga*) adrenocortical activity increases during times of large aggregations in the Serengeti ecosystem. *Horm. Behav.* **102**, 1-9 (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, 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