

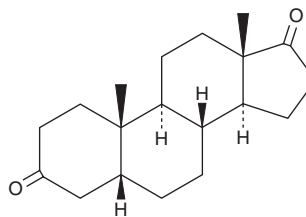
# PRODUCT INFORMATION



## Etiocholanedione

Item No. 34358

CAS Registry No.: 1229-12-5  
Formal Name: (5 $\beta$ )-androsterane-3,17-dione  
Synonyms: 5 $\beta$ -Androstanedione, Etiocholane-3,17-dione  
MF: C<sub>19</sub>H<sub>28</sub>O<sub>2</sub>  
FW: 288.4  
Purity:  $\geq$ 98%  
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

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

### Description

Etiocholanedione is a neurosteroid.<sup>1</sup> It inhibits glycine-induced chloride currents ( $I_{Gly}$ ) in isolated rat hippocampal pyramidal neurons when used at a concentration of 50  $\mu$ M. Etiocholanedione (0.1, 0.3, and 0.5 mM) also inhibits glucose exchange transport in isolated human erythrocytes.<sup>2</sup>

### References

1. Bukanova, J.V., Solntsera, I., and Kudova, E. Neurosteroids as selective inhibitors of glycine receptor activity: Structure-activity relationship study on endogenous androstanes and androstenes. *Front. Mol. Neurosci.* **13**, 44 (2020).
2. Lacko, L., Wittke, B., and Geck, P. Interaction of steroids with the transport system of glucose in human erythrocytes. *J. Cell. Physiol.* **86 Suppl 2(3 Pt 2)**, 673-680 (1975).

#### 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, 12/06/2022

#### 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