

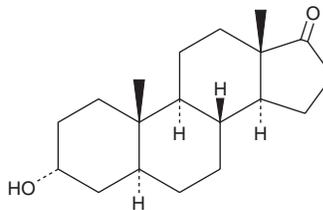
# PRODUCT INFORMATION



## Androsterone

Item No. 15872

**CAS Registry No.:** 53-41-8  
**Formal Name:** (5 $\alpha$ )-3 $\alpha$ -hydroxy-androstan-17-one  
**Synonyms:** Androkinin, NSC 9898  
**MF:** C<sub>19</sub>H<sub>30</sub>O<sub>2</sub>  
**FW:** 290.4  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 295 nm  
**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

Androsterone is supplied as a crystalline solid. A stock solution may be made by dissolving the androsterone in the solvent of choice, which should be purged with an inert gas. Androsterone is soluble in organic solvents such as ethanol, methanol, and acetonitrile. The solubility of androsterone in these solvents is approximately 1 mg/ml.

### Description

Androsterone is a steroid hormone produced in the body from 5 $\alpha$ -reduced metabolites of dehydroepiandrosterone and possesses 7-fold weaker androgen potency compared to that of testosterone (Item No. 15645).<sup>1</sup> It is the major androgen excreted in urine as a metabolite of testosterone and is used as the international reference standard for androgenic activity.<sup>1</sup>

### Reference

1. de Gruyter, W. Androsterone. *Concise encyclopedia biology*. Scott, T., editor, de Gruyter (1996).

**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, 11/16/2022

### 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