

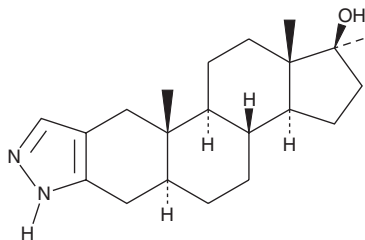
# PRODUCT INFORMATION



## Stanozolol

Item No. 21167

**CAS Registry No.:** 10418-03-8  
**Formal Name:** (5 $\alpha$ ,17 $\beta$ )-17-methyl-2'H-androst-2-eno[3,2-c]pyrazol-17-ol  
**Synonyms:** Androstanazol, NSC 43193, NSC 233046, WIN 14,833  
**MF:** C<sub>21</sub>H<sub>32</sub>N<sub>2</sub>O  
**FW:** 328.5  
**Purity:**  $\geq$ 98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 224 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

Stanozolol (Item No. 21167) is an analytical reference standard that is categorized as a synthetic anabolic androgenic steroid. Stanozolol has low urinary excretion; therefore, its metabolites are usually used for detection.<sup>1</sup> Formulations containing stanozolol are used as a treatment for hereditary angioedema.<sup>2</sup> It has also been abused in athletics to improve performance. Stanozolol is regulated as a Schedule III compound in the United States. This product is intended for research and forensic applications.

### References

1. Schanzer, W. and Donike, M. Metabolism of anabolic steroids in man: Synthesis and use of reference substances for identification of anabolic steroid metabolites. *Analytica Chim. Acta* **275**, 23-48 (1993).
2. Sloane, D.E., Lee, C.W., and Sheffer, A.L. Hereditary angioedema: Safety of long-term stanozolol therapy. *J. Allergy Clin. Immunol.* **120**, 654-658 (2007).

#### 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, 06/22/2023

#### 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