

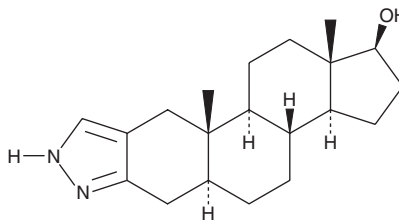
# PRODUCT INFORMATION



## [3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol

Item No. 17943

CAS Registry No.: 99996-65-3  
Formal Name: (5 $\alpha$ ,17 $\beta$ )-1'H-androstano[3,2-c]pyrazol-17-ol  
MF: C<sub>20</sub>H<sub>30</sub>N<sub>2</sub>O  
FW: 314.5  
Purity:  $\geq$ 98%  
Stability:  $\geq$ 2 years at -20°C  
Supplied as: A crystalline solid  
UV/Vis.:  $\lambda_{\text{max}}$ : 223 nm



### Laboratory Procedures

For long term storage, we suggest that [3,2-c]pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol be stored as supplied at -20°C. It should be stable for at least two years.

[3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol is supplied as a crystalline solid. A stock solution may be made by dissolving the [3,2-c]pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol in the solvent of choice. [3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of [3,2-c]pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

[3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, [3,2-c]pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. [3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol has a solubility of approximately 0.1 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

[3,2-c]Pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol is an anabolic androgenic steroid (AAS) that is closely related to stanozolol. Both steroids are regulated as Schedule III compounds by the US Drug Enforcement Agency. The physiological and toxicological properties of [3,2-c]pyrazole-5 $\alpha$ -androstan-17 $\beta$ -ol are not known, although the effects of AAS, including deleterious effects on the hypothalamic-pituitary-gonadal axis, are well known.<sup>1,2</sup> This product is intended for forensic and research applications.

### References

1. Cunningham, R.L., Lumia, A.R., and McGinnis, M.Y. Androgenic anabolic steroid exposure during adolescence: Ramifications for brain development and behavior. *Horm. Behav.* **64**(2), 350-356 (2013).
2. Nieschlag, E. and Vorona, E. Mechanisms in endocrinology: Medical consequences of doping with anabolic androgenic steroids: Effects on reproductive functions. *Eur. J. Endocrinol.* **173**, R47-R58 (2015).

#### 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, 10/28/2015

#### 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