

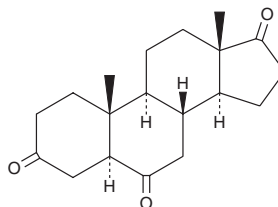
PRODUCT INFORMATION



5 α -Androstan-3,6,17-trione

Item No. 17945

CAS Registry No.: 2243-05-2
Formal Name: 5 α -androstan-3,6,17-trione
MF: C₁₉H₂₆O₃
FW: 302.4
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that 5 α -androstan-3,6,17-trione be stored as supplied at -20°C. It should be stable for at least two years.

5 α -Androstan-3,6,17-trione is supplied as a crystalline solid. A stock solution may be made by dissolving the 5 α -androstan-3,6,17-trione in the solvent of choice. 5 α -Androstan-3,6,17-trione is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 5 α -androstan-3,6,17-trione in these solvents is approximately 5, 30, and 50 mg/ml, respectively.

5 α -Androstan-3,6,17-trione is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 5 α -androstan-3,6,17-trione should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 5 α -Androstan-3,6,17-trione has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

5 α -Androstan-3,6,17-trione is an analog of 4-androstene-3,6,17-trione (4-AT, 6-OXO), an irreversible aromatase inhibitor that is used to increase serum androgen levels.^{1,2} 5 α -Androstan-3,6,17-trione is considered a designer drug and is regulated in the United States. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Numazawa, M., Tsuji, M., and Mutsumi, A. Studies on aromatase inhibition with 4-androstene-3,6,17-trione: Its 3 beta-reduction and time-dependent irreversible binding to aromatase with human placental microsomes. *J. Steroid Biochem.* **28(3)**, 337-344 (1987).
2. Rohle, D., Wilborn, C., Taylor, L., et al. Effects of eight weeks of an alleged aromatase inhibiting nutritional supplement 6-OXO (androst-4-ene-3,6,17-trione) on serum hormone profiles and clinical safety markers in resistance-trained, eugonadal males. *J. Int. Soc. Sports Nutr.* **4(13)**, (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, 10/29/2015

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