

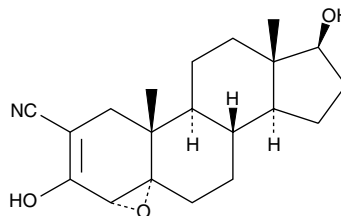
Product Information



Trilostane

Item No. 14164

CAS Registry No.: 13647-35-3
Formal Name: 4 α ,5 α -epoxy-3,17 β -dihydroxy-androst-2-ene-2-carbonitrile
Synonyms: Desopan, Modrefen, Vetoryl, WIN 24,540
MF: C₂₀H₂₇NO₃
FW: 329.4
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{\max} : 252 nm



Laboratory Procedures

For long term storage, we suggest that trilostane be stored as supplied at -20°C. It should be stable for at least two years.

Trilostane is supplied as a crystalline solid. A stock solution may be made by dissolving the trilostane in the solvent of choice. Trilostane is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of trilostane in these solvents is approximately 16 and 20 mg/ml, respectively.

Trilostane is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, trilostane should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Trilostane 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.

3 β -hydroxysteroid dehydrogenase (3 β -HSD) type 1 and type 2 isoforms are key enzymes for the biosynthesis of all active steroid hormones. 3 β -HSD1 (type I) is expressed in placenta and peripheral tissues including breast tumors, whereas 3 β -HSD2 (type 2) is expressed in the adrenal gland, ovary, and testis. Trilostane is an inhibitor of the 3 β -HSDs: 3 β -HSD1 and 3 β -HSD2 with K_i values of 0.10 and 1.60 μ M, respectively.¹ Trilostane has been approved for use in the treatment of Cushing's syndrome in dogs to reduce cortisol, aldosterone, and corticosterone levels.² Because human 3 β -HSD (type 1) is a critical enzyme in the conversion of DHEA to estradiol in breast tumors, trilostane is also of interest for the treatment of breast cancer in postmenopausal women.³

References

1. Thomas, J.L., Mack, V.L., Glow, J.A., *et al.* Structure/function of the inhibition of human 3 β -hydroxysteroid dehydrogenase type 1 and type 2 by trilostane. *J. Steroid Biochem. Mol. Biol.* **111**(1-2), 66-73 (2008).
2. Machida, T., Uchida, E., Matsuda, K., *et al.* Aldosterone-, corticosterone- and cortisol-secreting adrenocortical carcinoma in a dog: Case report. *J. Vet. Med. Sci.* **70**(3), 317-320 (2008).
3. Thomas, J.L., Bucholtz, K.M., and Kacsoh, B. Selective inhibition of human 3 β -hydroxysteroid dehydrogenase type 1 as a potential treatment for breast cancer. *J. Steroid Biochem. Mol. Biol.* **125**(1-2), 57-65 (2011).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/14164

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 05/21/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com