

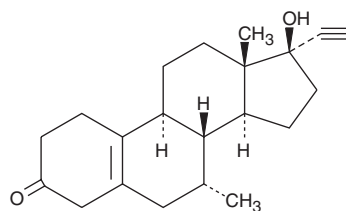
Product Information



Tibolone

Catalog No. 10006321

CAS Registry No.: 5630-53-5
Formal Name: 17 α -hydroxy-7 α -methyl-19-norpregn-5(10)-en-20-yn-3-one
Synonyms: Livial, Liviella
MF: C₂₁H₂₈O₂
FW: 312.5
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

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

Tibolone is supplied as a crystalline solid. A stock solution may be made by dissolving the tibolone in an organic solvent purged with an inert gas. Tibolone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of tibolone in these solvents is 2, 20, and 30 mg/ml, respectively.

Tibolone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tibolone should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Tibolone has a solubility of 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Tibolone is an estrogen-like compound used for the treatment of the symptoms associated with menopausal transition (i.e., climacteric symptoms) and also for the treatment of osteoporosis.¹ Three major metabolites of tibolone are responsible for its tissue selective mechanism of action. Conversion into 3 α - and 3 β -hydroxy-tibolone results in estrogenic effects in brain, vagina, and bone. The Δ^4 isomer has progestogenic and androgenic effects and does not cause estrogenic stimulation in the endometrium.¹ A two-year longitudinal study indicates that low doses (1.25-2.5 mg) of tibolone effectively relieve climacteric symptoms and prevent loss of bone mass in early postmenopausal women.²

References

1. Kloosterboer, H.J. Tissue-selectivity: The mechanism of action of tibolone. *Maturitas* **48**, S30-S40 (2004).
2. Gambacciani, M., Ciaponi, M., Cappagli, B., *et al.* A longitudinal evaluation of the effect of two doses of tibolone on bone density and metabolism in postmenopausal women. *Gynecol. Endocrinol.* **18**, 9-16 (2004).

Related Products

Estradiol EIA Kit - Cat. No. 582251 • Progesterone EIA Kit - Cat. No. 582601 • Estradiol - Cat. No. 10006315 • Levonorgestrel - Cat. No. 10006318 • Norgestrel - Cat. No. 10006319 • Tibolone - Cat. No. 10006321 • Estriol - Cat. No. 10006484 • Ethinyl Estradiol - Cat. No. 10006486 • Estradiol Benzoate - Cat. No. 10006487 • Gestrinone - Cat. No. 10006488

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

MATERIAL 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 Material 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, 11/17/2009

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