

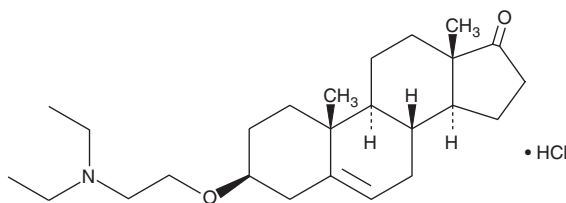
# Product Information



## U-18666A

Item No. 10009085

**Cas Registry No.:** 3039-71-2  
**Formal Name:** 3 $\beta$ -[2-(diethylamino)ethoxy]-androst-5-en-17-one, monohydrochloride  
**MF:** C<sub>25</sub>H<sub>41</sub>NO<sub>2</sub> • HCl  
**FW:** 424.1  
**Purity:**  $\geq$ 95%  
**Stability:**  $\geq$ 2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that U-18666A be stored as supplied at -20°C. It should be stable for at least two years.

U-18666A is supplied as a crystalline solid. A stock solution may be made by dissolving the U-18666A in an organic solvent purged with an inert gas. U-18666A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of U-18666A in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of U-18666A can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of U-18666A in PBS, pH 7.2, is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Impaired trafficking of cholesterol to or from the endoplasmic reticulum (ER) has been implicated in many disease states including Neimann-Pick disease type C (NPC), Alzheimers disease, and atherosclerosis. U-18666A is a cell permeable drug that inhibits cholesterol trafficking. It inhibits cholesterol transport from late endosomes/lysosomes to the ER, but not cholesterol transport to the plasma membrane as demonstrated in many cell types including macrophages, primary cortical neurons, and primary fibroblasts.<sup>1-3</sup> In macrophages, micromolar concentrations of U-18666A inhibit multiple pathways of cholesterol trafficking from late endosomes, whereas nanomolar concentrations impair cholesterol trafficking to the ER, a response similar to that found in NPC. U-18666A inhibits oxidosqualene cyclase at high (>0.5  $\mu$ M) concentrations and oral doses (10 mg/kg) induces cataracts in rats.<sup>1</sup>

### References

1. Koh, C.H.V. and Cheung, N.S. Cellular mechanism of U18666A-mediated apoptosis in cultured murine cortical neurons: Bridging Niemann-Pick disease type C and Alzheimer's disease. *Cell. Signal.* **18(11)**, 1844-1853 (2006).
2. Tabas, I. Apoptosis and plaque destabilization in atherosclerosis: The role of macrophage apoptosis induced by cholesterol. *Cell Death and Differentiation* **11**, S12-S16 (2004).
3. Feng, B., Yao, P.M., Li, Y., *et al.* The endoplasmic reticulum is the site of cholesterol-induced cytotoxicity in macrophages. *Nature Cell Biology* **5(9)**, 781-792 (2003).

### Related Products

NPC1L1 Polyclonal Antibody - Item No. 10005385 • Levonorgestrel - Item No. 10006318 • Norgestrel - Item No. 10006319 • PCSK9 (human) Polyclonal Antibody - Item No. 10007185 • SREBP-2 Polyclonal Antibody - Item No. 10007663 • LDL Receptor Polyclonal Antibody - Item No. 10007665

### 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

**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, 09/07/2011