

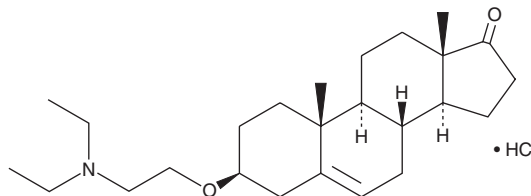
PRODUCT INFORMATION



U-18666A

Item No. 10009085

CAS Registry No.: 3039-71-2
Formal Name: 3 β -[2-(diethylamino)ethoxy]-
androst-5-en-17-one,
monohydrochloride
MF: C₂₅H₄₁NO₂ • 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.

Description

Impaired trafficking of cholesterol to or from the endoplasmic reticulum (ER) has been implicated in many disease states including Niemann-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.¹⁻³ 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 μ M) concentrations and oral doses (10 mg/kg) induces cataracts in rats.¹

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).

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, 09/09/2015

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