

# U-18666A

## Oxido-squalene cyclase inhibitor

A useful tool for inhibition of cholesterol transport from late endosomes/lysosomes to the endoplasmic reticulum. Inhibits cholesterol transport in a variety of cell types such as macrophages and sympathetic neurons. May be used to induce a Neimann-Pick disease type C (NPC) or NPC-like phenotype in a variety of cell types. Inhibits cholesterol biosynthesis at the level of oxido-squalene cyclase (OSC). Induction of cataracts is a common toxic feature of OSC inhibitors such as U18666A.

Citations: 6

[View Online »](#)

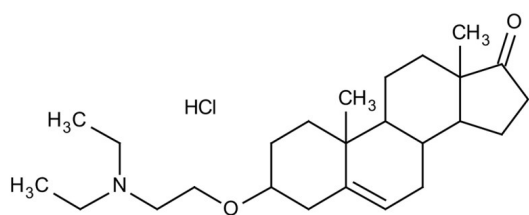
## Ordering Information

[Order Online »](#)

BML-S200-0010	10mg
BML-S200-0050	50mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** Ambient

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name**  $3\beta$ -(2-Diethylaminoethoxy)-Androstenone . HCl

**Appearance** White solid.

**CAS** 3039-71-2

**Couple Target** Oxido-squalene cyclase

**Couple Type** Inhibitor

**Formula**  $C_{25}H_{41}NO_2 \cdot HCl$

**MW** 424.1

**Purity**  $\geq 99\%$  (TLC)



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)