

# Pyripyropene A

ACAT2 inhibitor

Highly specific inhibitor of acyl-coenzyme A:cholesterol acetyltransferase 2 (ACAT2).

Citations: 15

[View Online »](#)

## Ordering Information

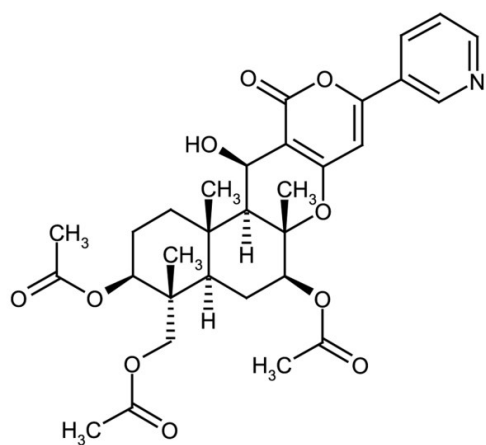
[Order Online »](#)

ALX-350-134-C500

500µg

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** (+)-Pyripyropene A, PPPA

**Appearance** White to brown powder.

**CAS** 147444-03-9

**Couple Target** ACAT

**Couple Type** Inhibitor

**Formula**  $C_{31}H_{37}NO_{10}$

**MW** 583.6

**Purity** ≥85% (HPLC)

**Solubility** Soluble in DMSO, chloroform, methanol or ethyl acetate; insoluble in water or hexane.

**Source** Isolated from *Aspergillus fumigatus* FO-1289.

**Last modified: May 29, 2024**



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-  
fr@enzolifesciences.com](mailto:info-fr@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)