

3-O-Acetyl-11-keto- β -boswellic acid

Inhibitor of lipoxygenase and topoisomerase

Shows antitumor and anti-inflammatory activities. Potent non-redox, noncompetitive 5-lipoxygenase and topoisomerase I and IIa inhibitor leading to apoptosis. 10-times more potent than 3-O-acetyl- β -boswellic acid (Prod. No. ALX-350-308). Exhibits *in vivo* efficacy in tumor growth and inhibition.

Citations: 7

[View Online »](#)

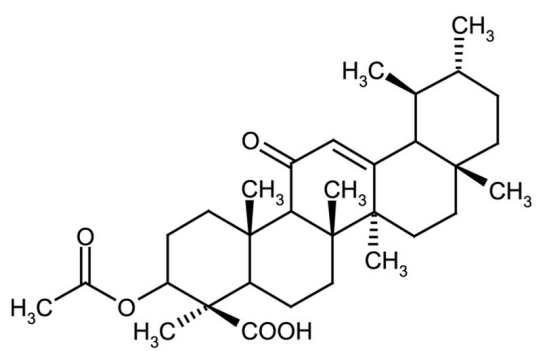
Ordering Information

[Order Online »](#)

ALX-350-310-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AKβBA
Appearance	White to off-white solid.
CAS	67416-61-9
Couple Target	Lipoxygenase, Topoisomerase
Couple Type	Inhibitor
Formula	$C_{32}H_{48}O_5$
MW	512.7
Purity	≥98% (HPLC)
Solubility	Soluble in acetone, dichloromethane, diethyl ether or DMSO.
Source	Isolated from <i>Boswellia serrata</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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