

# Bufotalin

Na<sup>+</sup>/K<sup>+</sup> ATPase inhibitor

Less potent Na<sup>+</sup>/K<sup>+</sup>-ATPase inhibitor than other bufadienolides. Potent immunosuppressor.

Citations: 2

[View Online »](#)

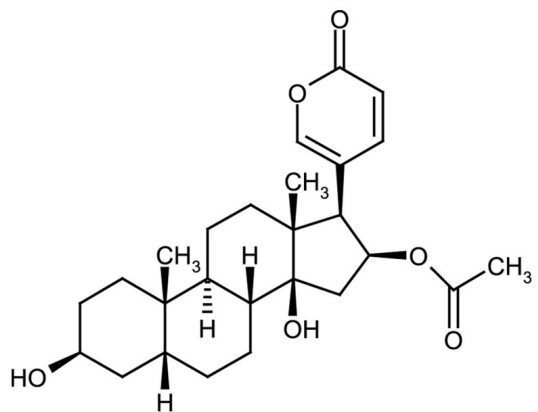
Ordering Information

[Order Online »](#)

ALX-350-282-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	(3β,5β,16β)-16-(Acetyloxy)-3-,14-dihydroxybufa-20,22-dienolide, 3β,14,16β-Trihydroxy-5β bufa-20,22-dienolide 16-acetate
Appearance	White to off-white solid.
CAS	471-95-4
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{26}H_{36}O_6$
MI	14: 1477
MW	444.6
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol or chloroform.
Source	Isolated from <i>Bufo</i> venom.

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-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)