

Aconitine

Na⁺ channel activator

A lipid-soluble neurotoxin that acts as a sodium channel modulator. Binds with high affinity to the open state of the voltage-sensitive sodium channels at site 2, thereby causing a persistent activation of the sodium channels, which become refractory to excitation.

Citations: 2

[View Online »](#)

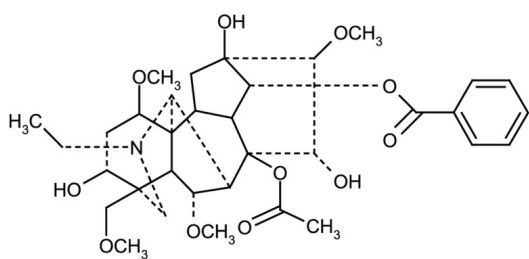
Ordering Information

[Order Online »](#)

ALX-550-232-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(1 α ,3 α ,6 α ,14 α ,15 α ,16 β)-20-Ethyl-1,6,16-trimethoxy-4-(methoxymethyl)aconitane-3,8,13,14,15-pentol 8-acetate 14-benzoate
Appearance	White to off-white powder.
CAS	302-27-2
Couple Target	Sodium
Couple Type	Activator
Formula	C ₃₄ H ₄₇ NO ₁₁
MI	14: 118
MW	645.7
Purity	≥98% (HPLC)
RTECS	AR5960000
Solubility	Soluble in 100% ethanol (35mg/ml); almost insoluble in water (0.3mg/ml).

Source



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-eu@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

Isolated from *Aconitum karacolum*.

Last modified: May 29, 2024