

Vindoline

Lacks physiological activity itself, but is contained as the pentacyclic moiety in the antineoplastic agents vinblastine and vincristine.

Citations: 4

[View Online »](#)

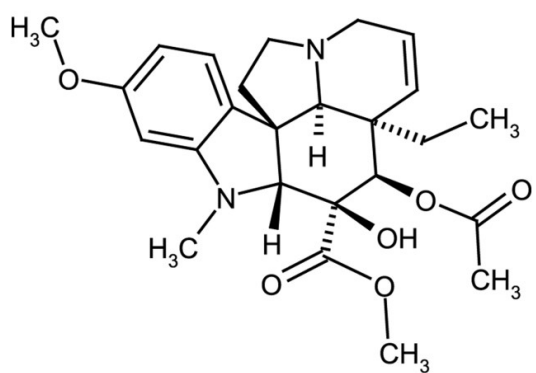
Ordering Information

[Order Online »](#)

ALX-350-102-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Very hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(2β,3β,4β,5α,12β,19α)-4-(acetyloxy)-6,7-didehydro-3-hydroxy-16-methoxy-1-methyl-Aspidospermidine-3-carboxylic acid methyl ester
Appearance	White to off-white crystalline powder.
CAS	2182-14-1
Formula	$C_{25}H_{32}N_2O_6$
MI	14: 9988
MW	456.5
Purity	≥90% (assay on dried basis)
RTECS	CJ0120000
Solubility	Soluble in 100% ethanol and water (30mg/ml).
Source	Isolated from <i>Catharanthus roseus</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-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