

# Catharanthine . tartrate

Starting material for the synthesis of the antitumor drugs vinblastine and vincristine. It is less active as an inhibitor of tubulin self-assembly into microtubules than the latter two compounds.

Citations: 2

[View Online »](#)

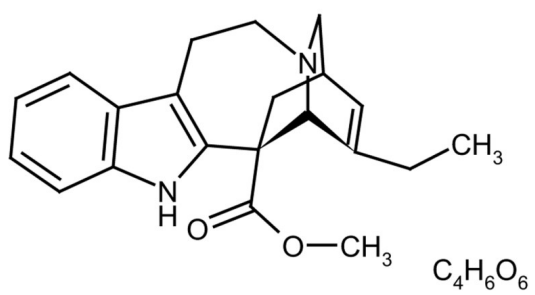
## Ordering Information

[Order Online »](#)

ALX-350-101-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. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	3,4-Didehydroibogamine-18-carboxylic acid methyl ester . tartrate, 7-Ethyl-9,10,12,13-tetrahydro-6,9-methano-5H-pyrido[1',2':1,2]azepino[4,5-b]indole-6(6aH)-carboxylic acid methyl ester . tartrate
Appearance	White to off-white crystalline powder.
CAS	4168-17-6
Formula	$C_{21}H_{24}N_2O_2 \cdot C_4H_6O_6$
MI	<b>14:</b> 1904
MW	336.4 . 150.1
Purity	≥90% (assay on dried basis)
Solubility	Soluble in 100% ethanol.
Source	Isolated from <i>Catharanthus roseus</i> .

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)