

Catharanthine . tartrate

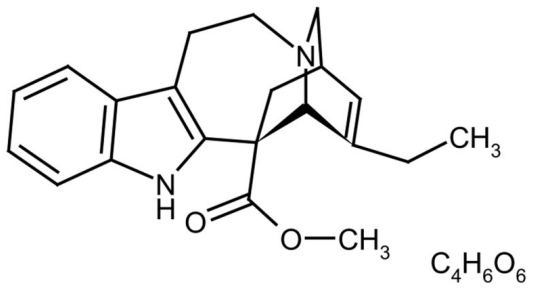
Starting material for the synthesis of the antitumor drugs vinblastine and vincristine. It is less active as an inhibitor of tubulin self-assembly into microtubules than the latter two compounds.

Citations: 2

[View Online »](#)

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	14: 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

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-be@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