

Vinblastine . sulfate

Cell cycle blocker

Plant alkaloid that arrests the cell cycle in G2/M phase by blocking mitotic spindle formation. Inhibits normal microtubule assembly and causes aberrant tubulin polymerization. Blocks binding of tubulin to microtubule proteins. Induces apoptosis. Triggers Raf-1 activation, phosphorylation of Bcl-2 family proteins and induction of p53 expression. Cancer chemotherapeutic agent.

Citations: 16

[View Online »](#)

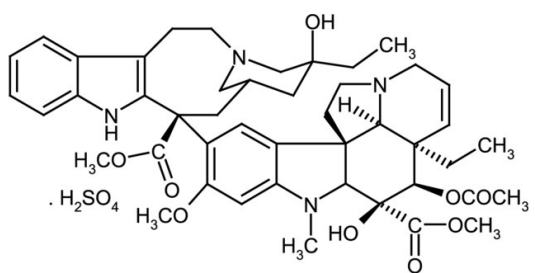
Ordering Information

[Order Online »](#)

ALX-350-257-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Vincaleukoblastine . sulfate

Appearance White to light yellow crystalline powder.

CAS 143-67-9

Formula $C_{46}H_{58}N_4O_9 \cdot H_2SO_4$

MI 14: 9982

MW 909.1

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, methanol or water (25mg/ml).

Source Semisynthetic from catharanthine or vindoline.

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)