

Vincristine . sulfate

Microtubule inhibitor

Plant alkaloid that arrests the cell cycle in G2/M phase by blocking mitotic spindle formation. Depolymerizes microtubules and blocks binding of tubulin to microtubule proteins. Causes aberrant tubulin polymerization. Induces apoptosis. Triggers Raf-1 activation, phosphorylation of Bcl-2 family proteins and induction of p53 expression. Cancer chemotherapeutic agent. Posttranscriptionally downregulates HIF-1 α expression.

Citations: 20

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-T117-0025	25mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	VCR . sulfate, Leukocristine . sulfate
Appearance	White to off-white powder.
CAS	2068-78-2
Couple Target	HIF, Microtubule
Couple Type	Inhibitor
Formula	$C_{46}H_{56}N_4O_{10} \cdot H_2SO_4$
MI	14: 9986
MW	825.0 . 98.1
Purity	≥98% TLC
RTECS	OH6340000
Solubility	Soluble in water (25mg/ml) or methanol.

Technical Info / Product Notes

Note: Product is not sterile.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com

Last modified: May 29, 2024