

Colchicine

Inhibitor of tubulin polymerization

Depolymerizes microtubules and inhibits tubulin polymerization. Induces apoptosis in a variety of cells.

Citations: 20

[View Online »](#)

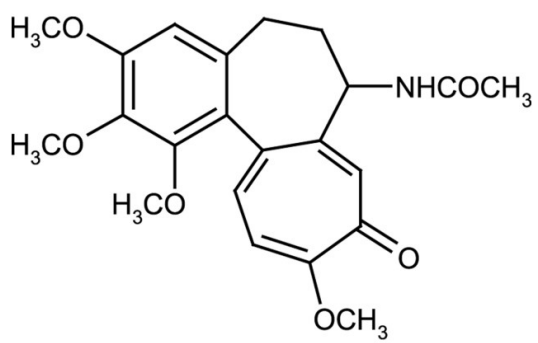
Ordering Information

[Order Online »](#)

ALX-380-033-G001	1g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	64-86-8
Formula	$C_{22}H_{25}NO_6$
MI	14: 2471
MW	399.4
Purity	≥94% (HPLC)
RTECS	GH0700000
Solubility	Soluble in DMSO (14mg/ml), 100% ethanol (7mg/ml), water (7mg/ml) or dimethylformamide (12mg/ml).
Source	Isolated from <i>Colchicum autumnale</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-
be@enzolifesciences.com](mailto:info-be@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)