

Colcemid

Depolymerizes microtubules and inhibits their formation. Induces apoptosis.

Citations: 15

[View Online »](#)

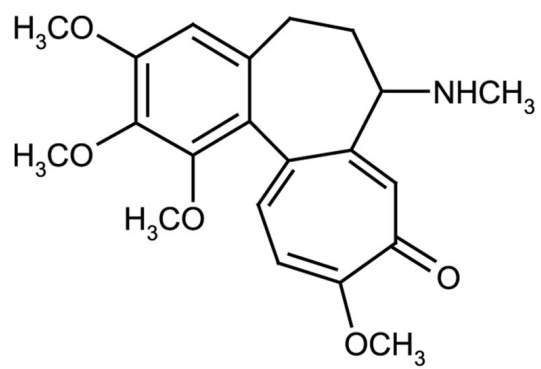
Ordering Information

[Order Online »](#)

ALX-430-033-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for at least 6 months when stored at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Demecolcine
Appearance	Light yellow powder.
CAS	477-30-5
Formula	$C_{21}H_{25}NO_5$
Identity	Determined by 1H-NMR and TLC.
MI	14: 2887
MW	371.4
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol, chloroform or DMSO.
Source	Semisynthetic from colchicine (Prod. No. ALX-380-033).

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)