

Malvidin chloride

Cell cycle inhibitor

Antioxidant flavonoid. Antitumor compound. Induces cell cycle arrest in the G2/M-phase.

Citations: 2

[View Online »](#)

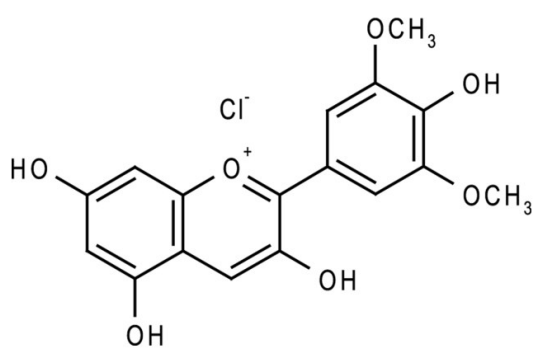
Ordering Information

[Order Online »](#)

ALX-385-013-M010	10mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3,4',5,7-Tetrahydroxy-3',5'-dimethoxyflavylium chloride
Appearance	Reddish brown to black powder.
CAS	643-84-5
Formula	$C_{17}H_{15}O_7Cl$
MI	14: 5715
MW	366.8
Purity	≥97% (HPLC)
RTECS	LK9900000
Solubility	Soluble in 100% ethanol; slightly soluble in water.

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)