

Delphinidin chloride (high purity)

Antioxidant

Anthocyanidin with antioxidant effect found in pigmented fruits and vegetables. Shown to inhibit angiogenesis and endothelial cell apoptosis by stimulating nitric oxide (NO) production. Inhibits solar radiation (UVB)-mediated oxidative stress, reducing DNA damage. Delphinidin induces necrosis in hepatocellular carcinoma cells with no caspase activation.

Citations: 8

[View Online »](#)

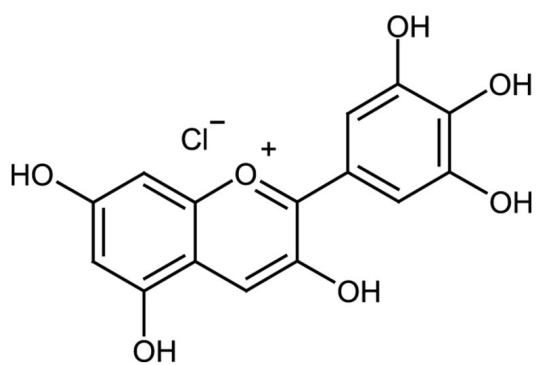
Ordering Information

[Order Online »](#)

ALX-385-028-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 and moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Red to dark brown powder.
CAS	528-53-0
Formula	$C_{15}H_{11}ClO_7$
MW	338.7
Purity	≥97% (HPLC)
RTECS	DK1310000
Solubility	Soluble in water or DMSO.

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)