

Cyanidin chloride

Nitric oxide (NO) scavenger

Antioxidant flavonoid. Nitric oxide (NO) scavenger.

Citations: 4

[View Online »](#)

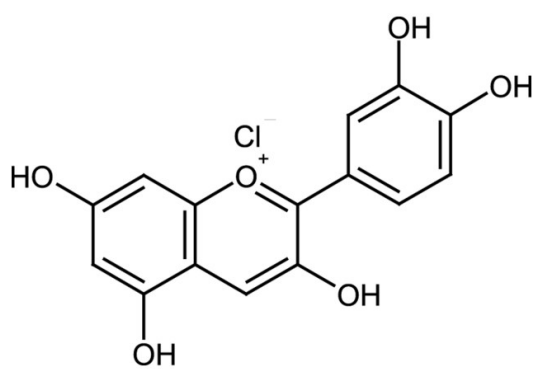
Ordering Information

[Order Online »](#)

ALX-385-003-M010	10mg
ALX-385-003-M050	50mg

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

Alternative Name 3,3',4',5,7-Pentahydroxy-flavylium chloride

Appearance Red to black solid.

CAS 528-58-5

Formula $C_{15}H_{11}ClO_6$

Identity Determined by 1H -NMR, ^{13}C -NMR and MS.

MI 14: 2687

MW 323.7

Purity $\geq 97\%$ (HPLC)

RTECS LK9824000

Solubility Soluble in methanol or 0.1% hydrochloric acid.

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)