

3,5-Di-O- caffeoylquinic acid

Antioxidant

Antioxidant. Shows antiproliferative activity.

Citations: 5

[View Online »](#)

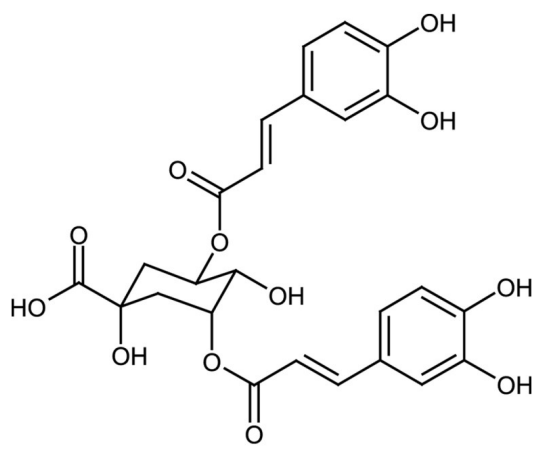
Ordering Information

[Order Online »](#)

ALX-350-320-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	3,5-CQA, Isochlorogenic acid
Appearance	Off-white to yellow/brown crystalline solid.
CAS	2450-53-5
Formula	$C_{25}H_{24}O_{12}$
Identity	Determined by 1H-NMR, 13C-NMR and MS.
MW	516.5
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.
Source	Isolated from <i>Cynara scolymus</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com