

(±)-Taxifolin

Antioxidant flavonoid

Antioxidant flavonoid.

Citations: 7

[View Online »](#)

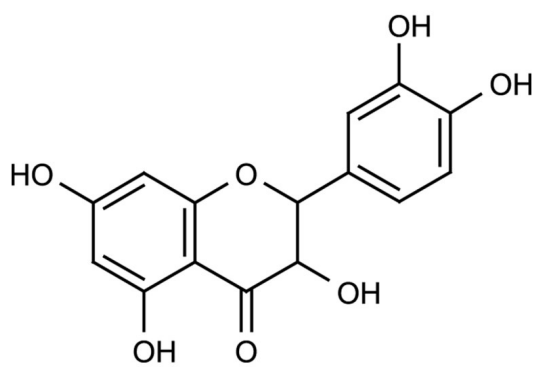
Ordering Information

[Order Online »](#)

ALX-385-006-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-Dihydroquercetin, (±)-3,3',4',5,7-Pentahydroxyflavanone, Taxifoliol, Distylin, Catechin hydrate
Appearance	White to slightly brown powder.
CAS	17654-26-1
Formula	$C_{15}H_{12}O_7$
Identity	Determined by NMR.
MW	304.3
Purity	≥95% (HPLC)
RTECS	LK6920000
Solubility	Soluble in 100% ethanol.

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