

Pseudohypericin

Inhibitor of PKC and dopamine β -hydroxylase

Inhibitor of protein kinase C (PKC) (IC_{50} =15 μ g/ml). Antiviral and antiretroviral activity. Potent inhibitor of dopamine- β -hydroxylase (IC_{50} =3 μ g/ml).

Citations: 8

[View Online »](#)

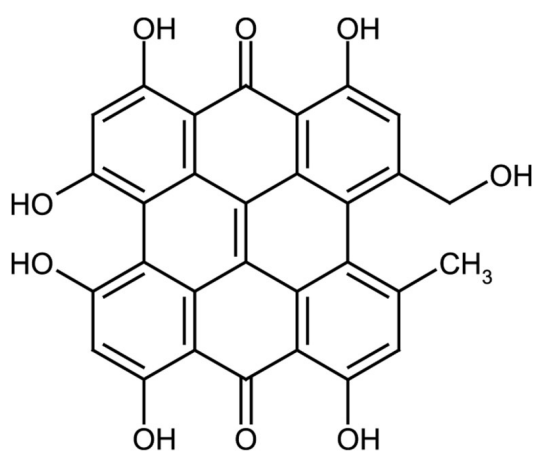
Ordering Information

[Order Online »](#)

ALX-350-276-M001	1mg
ALX-350-276-M005	5mg

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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Black to black-red powder.
CAS	55954-61-5
Couple Target	Dopamine beta-hydroxylase, PKC
Couple Type	Inhibitor
Formula	$C_{30}H_{16}O_9$
MW	520.5
Purity	≥98% (HPLC, TLC)
Solubility	Soluble in methanol, 100% ethanol, DMSO, acetone and other organic solvents.
Source	Isolated from <i>Hypericum perforatum</i> .

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)