

Hypericin (native)

PKC inhibitor

A potent and selective inhibitor of protein kinase C ($IC_{50}=3.4\mu M$). Anti-viral and anti-retroviral activity. Displays antineoplastic and antitumor activities. Most activities require light activation. Induces apoptosis in neuroblastoma cells at $1\mu M$ or above ($IC_{50}=0.1\mu M$ at 29 mW/cm^2 (tungsten light)). Inhibition of other kinases (e.g. EGFR-PTK, CKII and MAP kinase). Bright red fluorescence emissino and photostability. Antidepressant.

Citations: 17

[View Online »](#)

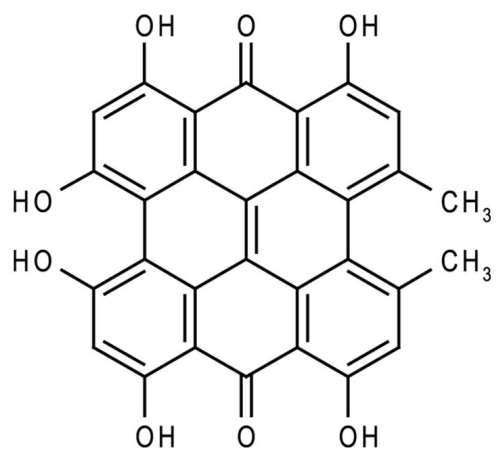
Ordering Information

[Order Online »](#)

ALX-350-030-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. High purity hypericin tautomerizes to 1,6-dioxo tautomer in the presence of water contained in solvents.
Handling	Protect from light and oxygen.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Dark brown to black-red powder.
CAS	548-04-9
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{30}H_{16}O_8$
MI	14: 4863
MW	504.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol, methanol, acetone, ethylmethylketone, pyridine or other organic solvents and bases.
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)