

# 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  $29\text{ mW/cm}^2$  (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 »](#)

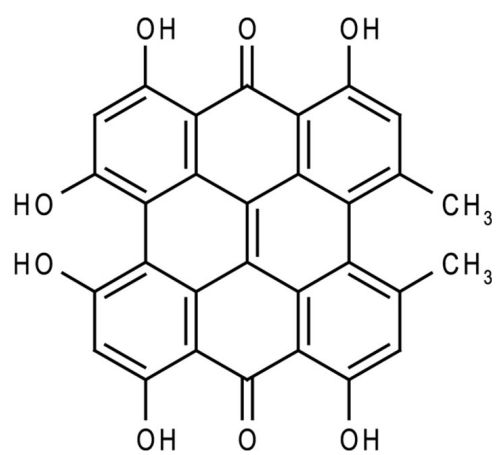
## Ordering Information

[Order Online »](#)

ALX-350-030-M001	1mg
ALX-350-030-M005	5mg
ALX-350-030-M010	10mg

## 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	<b>14:</b> 4863
MW	504.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol, methanol, acetone, ethylmethyleketone, pyridine or other organic solvents and bases.
Source	Isolated from <i>Hypericum perforatum</i> .



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)