Hypericin (native)

PKC inhibitor

A potent and selective inhibitor of protein kinase C (IC $_{50}$ =3.4µM). Anti-viral and anti-retroviral actitity. Displays antineoplastic and antitumor actitities. Most activities require light activation. Induces apoptosis in neuroblastoma cells at 1µM or above (IC50=0.1µM at 29 mW/cm² (tungsten light)). Inhibitition of other kinases (e.g. EGFR-PTK, CKII and MAP kinase). Bright red fluorescence emissino and photostability. Antidepressant.

Citations: 17

View Online »

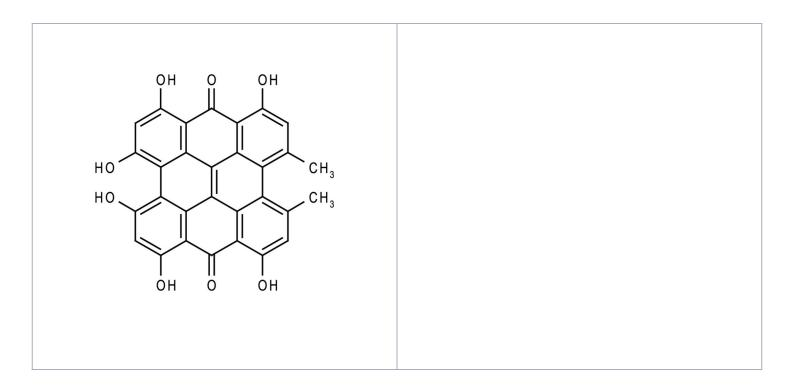
Ordering Information

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 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 *Hypericum perforatum*.

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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