

Ionomycin (free acid)

Calcium ionophore

Potent and highly selective calcium ionophore. Ca^{2+} /ionomycin complex forms between pH 7-9.5 which has intense UV absorption. Induces apoptotic neuronal degeneration in embryonic cortical neurons and cell cycle arrest at G_1 phase. Useful tool for increasing intracellular Ca^{2+} concentration. More effective than A23187. Non fluorescent. Has been used at 4nM and 2-4 μM (whole cells).

Citations: 32

[View Online »](#)

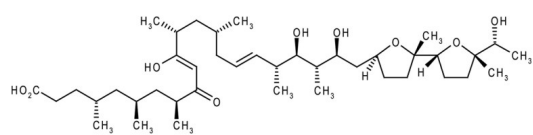
Ordering Information

[Order Online »](#)

ALX-450-006-M001	1mg
ALX-450-006-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for 1 year when stored at +4°C.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Waxy yellow to colorless material
CAS	56092-81-0
Formula	$C_{41}H_{72}O_9$
MW	709.0
Purity	≥90% (HPLC, TLC)
Solubility	Soluble in DMSO or methanol.
Source	Isolated from <i>Streptomyces conglobatus</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-
fr@enzolifesciences.com](mailto:info-fr@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)