

# Ionomycin . calcium salt

Calcium ionophore

For technical information and literature references see Ionomycin (free acid) (Prod. No. ALX-450-006).

Citations: 13

[View Online »](#)

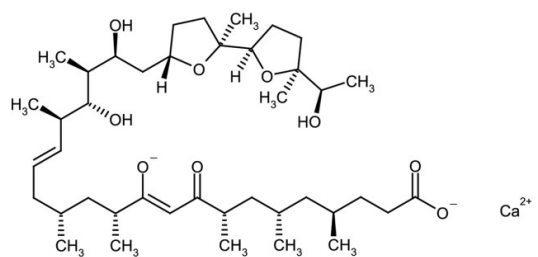
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-450-007-M001 | 1mg |
| ALX-450-007-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 | -20°C                                                            |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|            |                                                                      |
|------------|----------------------------------------------------------------------|
| Appearance | White to off-white solid.                                            |
| CAS        | 56092-82-1                                                           |
| Formula    | $C_{41}H_{70}O_9 \cdot Ca$                                           |
| MW         | 707.0 . 40.1                                                         |
| Purity     | ≥97% (HPLC, TLC)                                                     |
| Solubility | 10mg/ml soluble in DMSO or 100% ethanol, 5mg/ml soluble in methanol. |
| Source     | Isolated from <i>Streptomyces conglobatus</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)