

# Thapsigargin

## ATPase inhibitor

Thapsigargin is a potent inhibitor of sarco/endoplasmic reticulum  $\text{Ca}^{2+}$ -ATPases (SERCA), which are linked with endoplasmic reticulum stress, and induces the release of intracellular stored  $\text{Ca}^{2+}$  without hydrolysis of inositolphospholipid ( $\text{IC}_{50}=30\text{nM}$ ). Inhibition of SERCA by thapsigargin reveals a significant change in intracellular  $\text{Ca}^{2+}$  homeostasis and pH regulation in tumor cells, and may be used to distinguish between discrete intracellular  $\text{Ca}^{2+}$  pools. Thapsigargin increases  $\text{Ca}^{2+}$ -dependent  $\text{Na}^{+}$  influx in human platelets in a dose-dependent manner. Thapsigargin-induced tumor promotion and down-regulation of the EGF receptor is independent of PKC activation. Thapsigargin also stimulates nitric oxide (NO) production and as a result, may promote apoptosis.

Citations: 75

[View Online »](#)

## Ordering Information

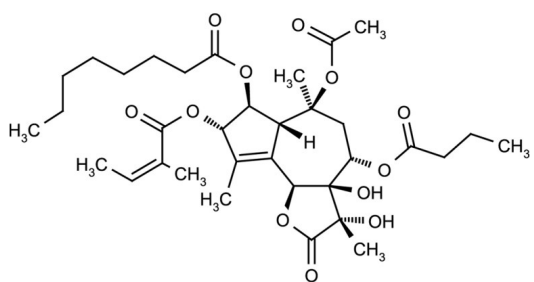
[Order Online »](#)

BML-PE180-0001	1mg
BML-PE180-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)

- Potent inhibitor of SERCA
- Cell-permeable tumor promoter



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C.
Handling	USE CAUTION – product is a tumor promoter. Solutions may lose activity after one week. For longer term storage, dissolve in acetone or ethanol, aliquot and evaporate solvent to dryness with a stream of nitrogen or argon. Store aliquots of neat compound at -20°C or colder. Protect from light and oxygen.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	Colorless wax or white solid.
CAS	67526-95-8
Couple Target	ATPase
Couple Type	Inhibitor
Formula	$C_{34}H_{50}O_{12}$
MI	14: 9272
MW	650.8
Purity	≥95% (HPLC, TLC)
Solubility	Soluble in acetone, DMSO (20mg/ml), or 100% ethanol (20mg/ml).
Source	Isolated from <i>Thapsia garganica</i> .

**Technical Info / Product Notes** Replacement for **ADI-908-298**.



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)