

Thapsigargin

ATPase inhibitor

Thapsigargin is a potent inhibitor of sarco/endoplasmic reticulum Ca^{2+} -ATPases (SERCA), which are linked with endoplasmic reticulum stress, and induces the release of intracellular stored Ca^{2+} without hydrolysis of inositolphospholipid ($\text{IC}_{50}=30\text{nM}$). Inhibition of SERCA by thapsigargin reveals a significant change in intracellular Ca^{2+} homeostasis and pH regulation in tumor cells, and may be used to distinguish between discrete intracellular Ca^{2+} pools. Thapsigargin increases Ca^{2+} -dependent 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

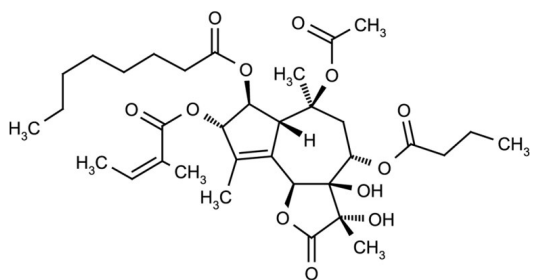
[Order Online »](#)

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024