

# BAPTA-AM

## PKC activator

BAPTA is a  $\text{Ca}^{2+}$  chelator with 105-fold greater affinity for  $\text{Ca}^{2+}$  than for  $\text{Mg}^{2+}$ . BAPTA-AM is a cell-permeable derivative which is widely used as an intracellular calcium sponge. Induces inactivation of protein kinase C. Also inhibits thapsigargin-induced apoptosis in rat thymocytes.

Citations: 23

[View Online »](#)

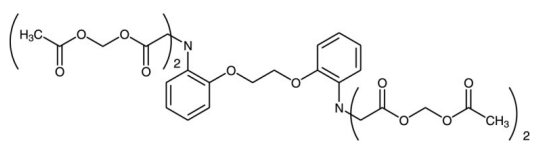
## Ordering Information

[Order Online »](#)

BML-CA411-0025	25mg
----------------	------

## 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

Alternative Name	1,2-bis-(o-Aminophenoxy)-ethane-N,N,N',N'-tetraacetic acid, tetraacetoxymethyl ester
Appearance	White to off-white powder.
CAS	126150-97-8
Couple Target	PKC
Couple Type	Activator
Formula	$C_{34}H_{40}N_2O_{18}$
Identity	Determined by NMR.
MW	764.7
Purity	≥90% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), dimethyl formamide (25mg/ml) or 100% ethanol (5mg/ml, warm)

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-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)