

# Xestospongin C

## IP3 receptor inhibitor

Xestonspongin C is a structurally novel marine alkaloid isolated from the Okinawan sponge *Xestospongia* sp. It is a potent, cell-permeable inhibitor of IP3 receptor-mediated  $\text{Ca}^{2+}$  release ( $\text{IC}_{50} = 358 \text{ nM}$ ). However, since xestospongin C also inhibits voltage-dependent  $\text{Ca}^{2+}$  and  $\text{K}^{+}$  currents at concentrations similar to those which inhibit the IP3 receptor, it can only be regarded as a selective blocker of the IP3 receptor in permeabilized cells and not in cells with intact plasma membranes. Has vasodilatory properties.

Citations: 20

[View Online »](#)

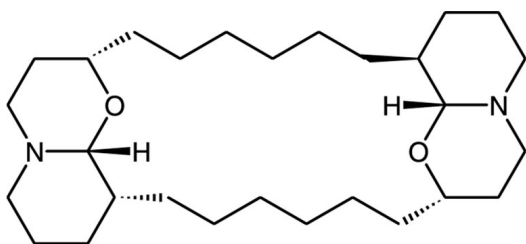
## Ordering Information

[Order Online »](#)

BML-CA409-0050	50µg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	White to off-white powder.
CAS	88903-69-9
Couple Target	IP3 receptor
Couple Type	Inhibitor
Formula	$C_{28}H_{50}N_2O_2$
MW	446.7
Purity	≥90% (TLC)
Solubility	Soluble in DMSO or 100% ethanol at 2mM.
Source	Synthetic.

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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)