

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 Ca^{2+} release ($\text{IC}_{50} = 358 \text{ nM}$). However, since xestospongin C also inhibits voltage-dependent Ca^{2+} and 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 »](#)

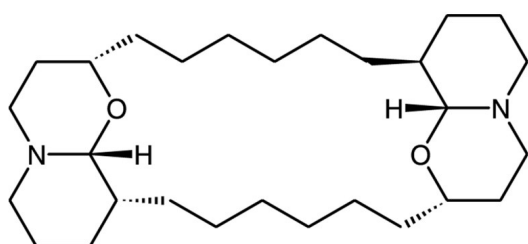
Ordering Information

[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.



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)