

Sphingosylphosphorylcholine

Induces release of Ca^{2+} from intracellular stores

Sphingosylphosphorylcholine induces release of Ca^{2+} from intracellular stores via an IP_3 independent pathway in permeabilized pancreatic acinar cells. It is an extremely potent mitogen in numerous cell types and induces neurite outgrowth in mouse neuroblastoma cells. It rapidly activates transcription factor AP-14 and transiently activates MAP kinase.

Citations: 6

[View Online »](#)

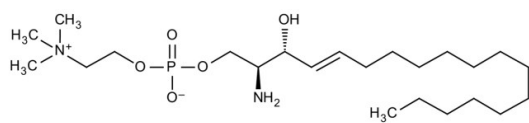
Ordering Information

[Order Online »](#)

BML-SL141-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Waxy solid.

CAS 82970-80-7

Couple Target MAPK

Couple Type Activator

Formula $C_{23}H_{49}N_2O_5P$

MW 464.6

Purity ≥98% (TLC)

Solubility Soluble in Chloroform: Methanol (2:1).

Last modified: August 19, 2025



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)