

Sphingomyelinase (Staphylococcus aureus)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE108-0010	10U
BML-SE108-0050	50U

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Sphingomyelin cholinephosphohydrolase
Application Notes	May be used exogenously to initiate the sphingomyelin signaling pathway. Has very little activity on phosphatidylcholine.
CAS	9031-54-3
Formulation	Liquid. Supplied in 50% glycerol/0.25 M phosphate buffer.
Source	Isolated from <i>Staphylococcus aureus</i> .
Specific Activity	Specific activity: 100-300 U/mg protein. One unit will hydrolyze 1 µmol of TNPAL-sphingomyelin per minute at pH 7.4 and 37°C
UniProt ID	P09978

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-
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)