

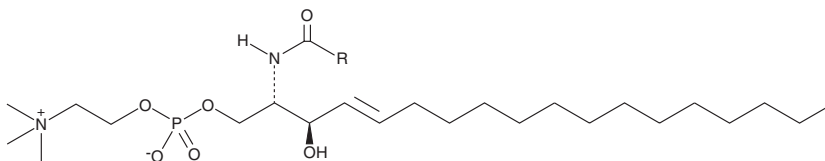
PRODUCT INFORMATION



Sphingomyelin (bovine buttermilk)

Item No. 31559

CAS Registry No.: 475662-40-9
Synonym: SMs (buttermilk)
MF: $C_{46}H_{93}N_2O_6P$ (for tricosanoyl)
FW: 801.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Sphingomyelin (bovine buttermilk) is supplied as a solid. A stock solution may be made by dissolving the sphingomyelin (bovine buttermilk) in the solvent of choice, which should be purged with an inert gas. Sphingomyelin (bovine buttermilk) is soluble in ethanol and chloroform.

Description

Sphingomyelins (SMs) are bioactive sphingolipids found in mammalian cell membranes.¹ SMs make up 2-15% of the total organ phospholipid population but are found at higher concentrations in the brain and myelin sheaths surrounding peripheral nerves. They interact with cholesterol to control its distribution within cellular membranes and maintain cholesterol homeostasis in cells. SMs undergo hydrolysis by sphingomyelinase to form ceramides, which are sphingolipid mediators of intracellular signaling.² This product contains SM molecular species with primarily C22:0, C23:0, and C24:0 fatty acyl chain lengths. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Slotte, J.P. and Ramstedt, B. The functional role of sphingomyelin in cell membranes. *Eur. J. Lipid Sci. Technol.* **109**(10), 977-981 (2007).
2. Shayman, J.A. Sphingolipids. *Kidney Int.* **58**(1), 11-26 (2000).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/20/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM