

PRODUCT INFORMATION



C20 Sphingomyelin (d18:1/20:0)

Item No. 24450

CAS Registry No.: 121999-68-6

Formal Name: 4-hydroxy-7S-[(1R,2E)-1-hydroxy-2-hexadecen-1-yl]-N,N,N-trimethyl-9-oxo-3,5-dioxo-8-aza-4-phosphaoctacosan-1-aminium, 4-oxide, inner salt

Synonyms: N-eicosanoyl-D-erythro-Sphingosylphosphorylcholine, SM(d18:1/20:0)

MF: $C_{43}H_{87}N_2O_6P$

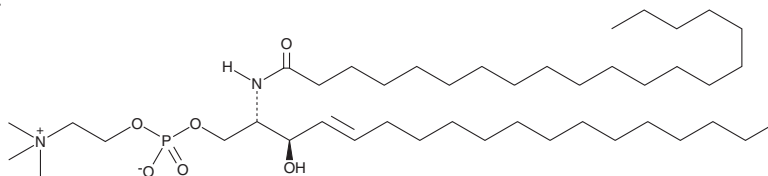
FW: 759.2

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

C20 Sphingomyelin (d18:1/20:0) is supplied as a solid. A stock solution may be made by dissolving the C20 sphingomyelin (d18:1/20:0) in the solvent of choice. C20 Sphingomyelin (d18:1/20:0) is soluble in ethanol, methanol, and a 14:1 solution of chloroform:methanol. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Description

C20 Sphingomyelin is a naturally occurring sphingolipid.¹⁻⁴ Levels of C20 sphingomyelin are upregulated in the hippocampus of rats with diabetes induced by streptozotocin (Item No. 13104) and in human plasma where it is positively correlated with insulin resistance in obese humans.² C20 sphingomyelin is also upregulated in the liver of a mouse model of Niemann-Pick type C1 disease, a neurodegenerative cholesterol-sphingolipid lysosomal storage disorder.¹ The plasma concentration of C20 sphingomyelin is decreased in men with prostate cancer.³

References

1. Praggastis, M., Tortelli, B., Zhang, J., *et al.* A murine Niemann-Pick C1 I1061T knock-in model recapitulates the pathological features of the most prevalent human disease allele. *J. Neurosci.* **35**(21), 8091-8106 (2015).
2. Hanamatsu, H., Ohnishi, S., Sakai, S., *et al.* Altered levels of serum sphingomyelin and ceramide containing distinct acyl chains in young obese adults. *Nutr. Diabetes* **4**(10), e141 (2014).
3. Awwad, H.M., Ohlmann, C.H., Stoeckle, M., *et al.* Choline-phospholipids inter-conversion is altered in elderly patients with prostate cancer. *Biochimie* **126**, 108-114 (2016).
4. Fiedorowicz, A., Prokopiuk, S., Zendzian-Piotrowska, M., *et al.* Sphingolipid profiles are altered in prefrontal cortex of rats under acute hyperglycemia. *Neuroscience* **256**, 282-291 (2014).

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, 12/01/2022

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