

PRODUCT INFORMATION



Globotetraosylceramide (porcine RBC)

Item No. 24881

CAS Registry No.: 11034-93-8

Formal Name: 1-O-[O-2-(acetylamino)-2-deoxy-β-D-galactopyranosyl-(1→3)-O-α-D-galactopyranosyl-(1→4)-O-β-D-galactopyranosyl-(1→4)-β-D-glucopyranosyl]-ceramide

Synonyms: Gb₄Cer (porcine RBC), Globoside (porcine RBC), Globotetrahexosylceramide (porcine RBC)

MF: C₆₈H₁₂₆N₂O₂₃ (for tetracosanoyl)

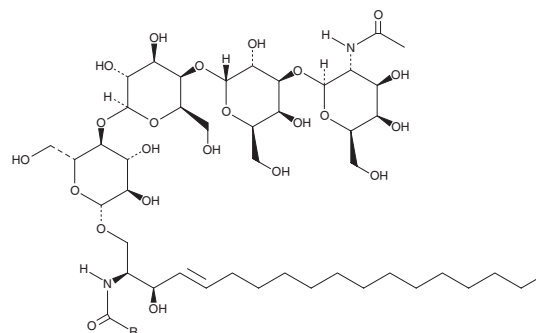
FW: 1,339.7

Purity: ≥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

Globotetraosylceramide (porcine RBC) is supplied as a solid. A stock solution may be made by dissolving the globotetraosylceramide (porcine RBC) in the solvent of choice. Globotetraosylceramide (porcine RBC) is soluble in DMSO, methanol (warmed), and a 2:1 solution of chloroform:methanol.

Description

Globotetraosylceramides are bioactive neutral glycosphingolipids. They are the major glycolipids in human erythrocytes.¹ They act as receptors for the Shiga toxins Stx1, Stx2, and Stx2e, the cytotoxic protein pierisin-1, and parvovirus B19.²⁻⁴ Globotetraosylceramides increase the expression of proteins responsible for enamel deposition, including ameloblastin, amelogenin, and enamelin, in dental epithelial cells and activate the ERK and p38 MAPK signaling pathways.⁵ Levels of globotetraosylceramides are elevated in fibroblasts from patients with salt and pepper syndrome, a neurocutaneous condition characterized by intellectual disability and hyper- and hypo-pigmented skin.⁶ This product contains globotetraosylceramides with primarily C24:0, C22:0, and 2-hydroxy C24:0 fatty acyl chains. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Yamakawa, T., Yokoyama, S., and Kiso, N. Structure of main globoside of human erythrocytes. *J. Biochem.* **52**(3), 228-229 (1962).
2. Müthing, J., Scheweppe, C.H., Karch, H., *et al.* Shiga toxins, glycosphingolipid diversity, and endothelial cell injury. *Thromb. Haemost.* **101**(2), 252-264 (2009).
3. Matsushima-Hibiya, Y., Watanabe, M., Hidari, J.I.-P.J., *et al.* Identification of glycosphingolipid receptors for pierisin-1, a guanine-specific ADP-ribosylating toxin from the cabbage butterfly. *J. Biol. Chem.* **278**(11), 9972-9978 (2003).
4. Nasir, W., Nilsson, J., Ologsson, S., *et al.* Parvovirus B19 VLP recognizes globoside in supported lipid bilayers. *Virology* **456-457**, 364-369 (2014).
5. Nakamura, T., Chiba, Y., Naruse, M., *et al.* Globoside accelerates the differentiation of dental epithelial cells into ameloblasts. *Int. J. Oral Sci.* **8**(4), 205-212 (2016).
6. Boccuto, L., Aoki, K., Flanagan-Steet, H., *et al.* A mutation in a ganglioside biosynthetic enzyme, ST3GAL5, results in salt & pepper syndrome, a neurocutaneous disorder with altered glycolipid and glycoprotein glycosylation. *Hum. Mol. Genet.* **23**(2), 418-433 (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, 11/04/2022

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