

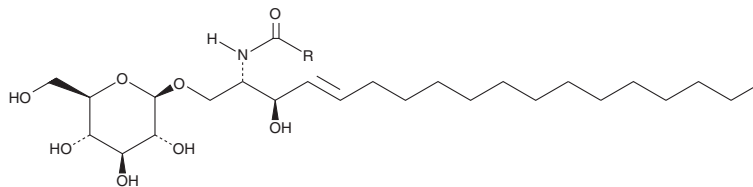
# PRODUCT INFORMATION



## Glucosylceramide (Gaucher's spleen)

Item No. 23207

**CAS Registry No.:** 85305-87-9  
**Formal Name:** 1-O-β-D-glucopyranosyl-ceramide  
**Synonyms:** Gaucher cerebroside, GL1a, GluCers (Gaucher's spleen), Glucocerebrosides, Glucosylceramides (gaucher's spleen)  
**MF:** C<sub>48</sub>H<sub>93</sub>NO<sub>8</sub> (for lignoceryl)  
**FW:** 812.3  
**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

Glucosylceramide (Gaucher's spleen) is supplied as a solid. A stock solution may be made by dissolving the glucosylceramide (Gaucher's spleen) in the solvent of choice, which should be purged with an inert gas. Glucosylceramide (Gaucher's spleen) is soluble in a 2:1 solution of chloroform:methanol.

### Description

Glucosylceramides are formed by the tethering of glucose to a ceramide by glucosylceramide synthase.<sup>1</sup> They are present in neuronal and non-neuronal mammalian tissues and are found at low quantities in a large number of plant species, where they comprise 5-30% of total lipids in the plant plasma membrane, and in fungi.<sup>1-3</sup> Glucosylceramide levels decrease during cold acclimatization in plants and glucosylceramides in fungi are involved in the regulation of virulence and act as elicitors in plants, stimulating plant defense mechanisms.<sup>3,4</sup> Glucosylceramides are precursors in the synthesis of lactosylceramides and gangliosides. Increased levels of glucosylceramides are associated with obesity-induced insulin resistance in mice and with neuronal deficits observed in neuronopathic Gaucher disease.<sup>1,5</sup> This product contains glucosylceramide molecular species with primarily C16:0, C22:0, 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

- Holland, W.L. and Summers, S.A. Sphingolipids, insulin resistance, and metabolic disease: New insights from *in vivo* manipulation of sphingolipid metabolism. *Endocr. Rev.* **29(4)**, 381-402 (2008).
- Cahoon, E.B. and Lynch, D.V. Analysis of Glucocerebrosides of Rye (*Secale cereale* L. cv Puma) Leaf and Plasma Membrane. *Plant Physiol.* **95(1)**, 56-68 (1991).
- Lynch, D.V. and Dunn, T.M. An introduction to plant sphingolipids and a review of recent advances in understanding their metabolism and function. *New Phytol.* **161(3)**, 667-702 (2004).
- Rollin-Pinheiro, R., Bernardino, M.C., and Barreto-Bergter, E. Sphingolipids: Functional and biological aspects in mammals, plants, and fungi. *Analysis of membrane lipids*. Prasad, R. and Singh, A., editors Springer (2020).
- Dai, M., Liou, B., Swope, B., *et al.* Progression of behavioral and CNS deficits in a viable murine model of chronic neuronopathic Gaucher disease. *PLoS One* **11(9)**, e0162367 (2016).

#### 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, 03/15/2024

#### 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