

# PRODUCT INFORMATION



## C18 Galactosylceramide-d<sub>35</sub> (d18:1/18:0-d<sub>35</sub>)

Item No. 24467

**CAS Registry No.:** 145176-92-7  
**Formal Name:** N-[(1S,2R,3E)-1-[(β-D-galactopyranosyloxy)methyl]-2-hydroxy-3-heptadecenyl]-octadecanamide-d<sub>35</sub>

**Synonyms:** Galactosylceramide (d18:1/18:0-d<sub>35</sub>), GalCer(d18:1/18:0-d<sub>35</sub>), N-octadecanoyl Psychosine-d<sub>35</sub>

**MF:** C<sub>42</sub>H<sub>46</sub>D<sub>35</sub>NO<sub>8</sub>

**FW:** 763.3

**Chemical Purity:** ≥98% (C18 Galactosylceramide)

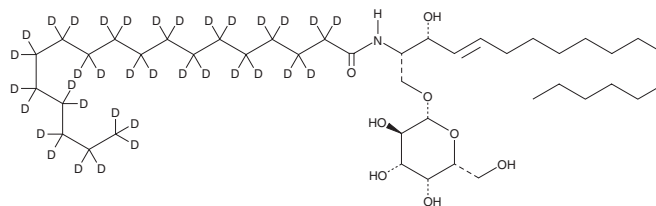
**Deuterium**

**Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>35</sub>); ≤1% d<sub>0</sub>

**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

C18 Galactosylceramide-d<sub>35</sub> (d18:1/18:0-d<sub>35</sub>) is intended for use as an internal standard for the quantification of C18 galactosylceramide by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

C18 Galactosylceramide-d<sub>35</sub> (d18:1/18:0-d<sub>35</sub>) is supplied as a solid. A stock solution may be made by dissolving the C18 galactosylceramide-d<sub>35</sub> (d18:1/18:0-d<sub>35</sub>) in the solvent of choice, which should be purged with an inert gas. C18 Galactosylceramide-d<sub>35</sub> (d18:1/18:0-d<sub>35</sub>) is soluble in ethanol (warmed), chloroform, and a 2:1 solution of chloroform:methanol.

### Description

C18 Galactosylceramide is a naturally occurring sphingolipid that has been found in the cerebellum of rats and humans.<sup>1,2</sup> It is selectively localized with cholesterol over potassium, sodium, or phosphocholine in rat cerebellum white matter.<sup>1</sup>

### References

1. Börner, K., Nygren, H., Hagenhoff, B., *et al.* Distribution of cholesterol and galactosylceramide in rat cerebellar white matter. *Biochem. Biophys. Acta* **1761**(3), 335-344 (2006).
2. Boutin, M., Sun, Y., Shacka, J.J., *et al.* Tandem mass spectrometry multiplex analysis of glucosylceramide and galactosylceramide isoforms in brain tissues at different stages of Parkinson Disease. *Anal. Chem.* **88**(3), 1856-1863 (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, 11/10/2022

#### 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