

PRODUCT INFORMATION



Neutral Glycosphingolipid Mixture

Item No. 29358

Supplied as: A solution in chloroform:methanol (2:1)

Storage: -20°C

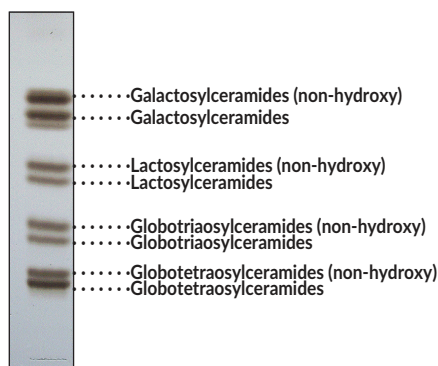
Stability: ≥2 years

Concentration: 1 mg/ml

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

Description and Data

Galactosylceramides, lactosylceramides, globotriaosylceramides, and globotetraosylceramides are neutral membrane glycosphingolipids that contain a saccharide moiety attached to a ceramide containing N-acyl hydroxy or non-hydroxy fatty acids. These glycosphingolipids have roles in synaptic signaling and immune activation and accumulate in a variety of lysosomal storage disorders, including Fabry and Krabbe diseases.¹⁻⁴ Neutral Glycosphingolipid Mixture contains hydroxy- and non-hydroxy galactosylceramides, lactosylceramides, globotriaosylceramides, and globotetraosylceramides. It is intended for use as a reference standard for the detection of these glycosphingolipids in various sample types by TLC or LC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation.



References

1. Boggs, J.M., Gao, W., Zhao, J., *et al.* Participation of galactosylceramide and sulfatide in glycosynapses between oligodendrocyte or myelin membranes. *FEBS Lett.* **584**(9), 1771-1778 (2010).
2. Birkholz, A.M., Howell, A.R., and Kronenberg, M. The α and Ω of galactosylceramides in T cell immune function. *J. Biol. Chem.* **290**(25), 15365-15370 (2015).
3. Wenger, D.A., Rafi, M.A., and Luzzi, P. Krabbe disease: One hundred years from the bedside to the bench to the bedside. *J. Neurosci. Res.* **94**(11), 982-989 (2016).
4. Feldt-Rasmussen, U., Rasmussen, A.K., Mersebach, H., *et al.* Fabry disease: A new challenge in endocrinology and metabolism? *Eur. J. Endocrinol.* **146**(6), 741-742 (2002).

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, 04/29/2029

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