PRODUCT INFORMATION



Monogalactosyldiacylglyceride (hydrogenated) (plant)

Item No. 17886

CAS Registry No.: Formal Name:	41670-62-6 2,3- <i>bis</i> [(1-oxooctadecyl)oxy]propyl-β-	
i offilar Name.	D-galactopyranoside	
Synonyms:	MGDG, Monogalactosyldiacylglycerol	
MF:	C ₄₅ H ₈₆ O ₁₀ (for stearoyl)	
FW:	787.2	
Purity:	≥95%	
UV/Vis.:	λ _{max} : 210 nm	но
Supplied as:	A solid	ОН
Storage:	-20°C	
Stability:	≥4 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

Monogalactosyldiacylglyceride (hydrogenated) (plant) is supplied as a solid. A stock solution may be made by dissolving the monogalactosyldiacylglyceride (hydrogenated) (plant) in the solvent of choice, which should be purged with an inert gas. The solubility of the monogalactosyldiacylglyceride (hydrogenated) (plant) in chloroform:methanol (1:1) is approximately 10 mg/ml.

Description

Monogalactosyldiacylglyceride is a minor galactolipid present in oligodendrocytes and myelin, as well as comprises a major fraction of the lipids in photosynthesizing plants and green microorganisms.^{1,2} It is used as a marker for myelination and has also been shown to stimulate PKC α activity in oligodendrocytes.¹ This product contains monogalactosyldiacylglyceride molecular species with C18:0 and C16:0 fatty acyl chain lengths.

References

- 1. Schmidt-Schultz, T. and Althaus, H.H. Monogalactosyl diglyceride, a marker for myelination, activates oligodendroglial protein kinase C. J. Neurochem. 62(4), 1578-1585 (1994).
- 2. Rosenberg, A., Gouaux, J., and Milch, P. Monogalactosyl and digalactosyl diglycerides from heterotrophic, hetero-autotrophic, and photobiotic Euglena gracilis. J. Lipid Res. 7(6), 733-738 (1996).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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.

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, 02/02/2023

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