PRODUCT INFORMATION



C22 Glucosylceramide-d₄ (d18:1/22:0-d₄)

Item No. 9003462

CAS Registry No.:	2738376-81-1	
Formal Name:	N-[(1S,2R,3E)-1-[(β-D-glucopyranosyloxy)methyl]-	
	2-hydroxy-3-heptadecen-1-yl]-docosanamide-	
	12,12,13,13-d ₄	
Synonyms:	N-Docosanoyl-glucopsychosine-d₄,	
	N-Docosanoyl- β -glucosylsphingosine-d ₄ ,	
	GlcCer(d18:1/22:0-d ₄), GluCer(18:1/22:0-d ₄),	
	N-C22:0-Glucocerebroside-d ₄ ,	
	C22 Glucosylceramide- d_4 ,	~
	Glucosylceramide-d ₄ (d18:1/22:0-d ₄)	но' 🔪
MF:	$C_{46}H_{85}D_4NO_8$	но.
FW:	788.2	
Chemical Purity:	≥98% (C22 Glucosylceramide)	
Deuterium		
Incorporation:	≥99% deuterated forms (d ₁ -d ₄); ≤1% d ₀	
Supplied as:	A crystalline 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

C22 Glucosylceramide- d_4 (d18:1/22:0- d_4) is intended for use as an internal standard for the quantification of C22 glucosylceramide (Item No. 23210) 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).

Description

C22 Glucosylceramide is an endogenous glucosylceramide.¹⁻³ Glucosylceramides are major constituents of skin lipid membranes where they play a role in maintaining the water permeability barrier. They are precursors in the synthesis of lactosylceramide (Item No. 16983), as well as oligoglycolipids and gangliosides. Phospholipase A_2 (PLA₂) type XIIA knockdown increases C22 glucosylceramide expression in rat brain.³ It is also increased in the brain, but not the liver or spinal cord, of mice fed a methionine-restricted diet.² In human athletes, plasma levels of C22 glucosylceramide increase during exercise and return to basal levels during recovery.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

- 1. Bergman, B.C., Brozinick, J.T., Strauss, A., et al. Serum sphingolipids: Relationships to insulin sensitivity and changes with exercise in humans. Am. J. Physiol. Endocrinol. Metab. 309(4), E398-E408 (2015).
- 2. Jové, M., Ayala, V., Ramírez-Núñez, O., et al. Specific lipidome signatures in central nervous system from methionine-restricted mice. J. Proteome Res. 12(6), 2679-2689 (2013).
- 3. Ee, S.-M., Lo, Y.-L., Shui, G., et al. Distribution of secretory phospholipase A₂ XIIA in the brain and its role in lipid metabolism and cognition. Mol. Neurobiol. 50(1), 60-75 (2014).

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.

WARRANTY AND LIMITATION OF REMEDY

uyer 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, 11/12/2024

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