

PRODUCT DATA SHEET

N-Hexanoyl-NBD-galactosylceramide

Catalog number: 1621

Common names: N-C6:0-NBD-*beta*-D-Galactosylsphingosine; N-C6:0-NBD-Cerebroside; N-C6:0-NBD-Galactosylceramide, fluorescent; N-(NBD-D-Aminocaproyl)-galactosylsphingosine

Source: semisynthetic, bovine

Solubility: chloroform/methanol, 5:1; methanol

CAS number: 170212-26-7

Molecular Formula: C₃₆H₅₉N₅O₁₁

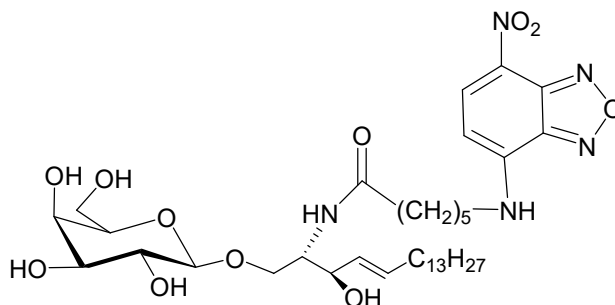
Molecular Weight: 738

Storage: -20°C

Purity: TLC >98%; identity confirmed by MS

TLC System: chloroform/methanol/DI water (65:25:3)

Appearance: solid



Application Notes:

This cerebroside product is a fluorescent labeled glycosphingolipid containing a galactose (galactocerebroside) attached to a ceramide acylated with a fluorescent NBD fatty acid. Galactocerebroside is found primarily in neuronal tissues and are the major glycosphingolipids in the central nervous system. They are the largest single component of the myelin sheath of nerves and seem to act, along with other molecules, to form part of the structural support of the myelin sheath.¹ Cerebroside is involved in a very wide range of biological activities such as cell agglutination, intracellular communication, cellular development, and antitumor/cytotoxic effects.² Galactocerebroside can be metabolized into sulfatide which is also abundant in the nervous system and myelin sheath. Due to the relatively high melting point of cerebroside (much greater than physiological body temperature) they have a para-crystalline structure. Krabbe's disease (globoid cell leukodystrophy) is characterized by a deficiency in the enzyme galactocerebrosidase, which is responsible for degrading galactocerebroside. This leads to an accumulation of cerebroside and psychosine (which is very cytotoxic and can result in demyelination of nerves and loss of axonal conductivity). This fluorescent standard from Matreya is excellent for use in the identification and isolation of cerebroside in the study of Krabbe's disease and other studies.³

Selected References:

1. M. Sheldon, D. Lyudmila, "Cycloserine-induced decrease of cerebroside in myelin" *Lipids*, Vol. 33:4 pp. 441-443, 1998
2. X. Zhou, L. Tang and Y. Liu "An Isomeric Mixture of Novel Cerebroside Isolated from *Impatiens pritzellii* Reduces Lipopolysaccharide-Induced Release of IL-18 from Human Peripheral Blood Mononuclear Cells" *Lipids*, Vol. 44:8 pp. 759-763, 2009
3. K. Zama et al. "Simultaneous quantification of glucosylceramide and galactosylceramide by normal-phase HPLC using *O*-phthalaldehyde derivatives prepared with sphingolipid ceramide *N*-deacylase" *Glycobiology*, vol. 19 pp. 767-775, 2009

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.