
Ceramide Monoclonal Antibody [MID15B4]

(Cat. No. A-0549)

Background

A ceramide is a waxy lipid molecule that is found in high concentration within the cell membranes of cells. They are a component of sphingomyelin, a major lipid in the lipid bilayer. Ceramides and other sphingolipids found in cell membranes can participate in a variety of cellular signaling, including proliferation and differentiation, as well as apoptosis or cell death and cell-cycle arrest.

Description

Mouse monoclonal antibody to Ceramide, Clone MID15B4.

Formulation

Liquid. In PBS, pH 7.2, containing 0.5M sodium chloride, 0.1% BSA and 0.09% sodium azide.

Immunogen

Ceramide (sphingosine-[trans-D-erythro-2-amino-4-octadecene-1,3-diol]) conjugated to BSA.

Specificity

Species Independent

Isotype

IgM

Purification

Serum

Storage

Store at 4°C. Avoid multiple freeze/thaw cycles.

Suggested amount:

20µl/test.

Application

ELISA (1:10), FC, ICC, IHC (1:10)

Ordering Information

Product	Size	Cat. No.
Ceramide Monoclonal Antibody [MID15B4]	1 ml	A-0549-001

This product is for research purposes only. Not intended for use in diagnostic procedures.