

Sphingomyelin Assay Kit ab133118

[2 References](#) [画像数 1](#)

製品の概要

製品名 Sphingomyelin Assay Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
plasma	16			= 3.9%
serum	16			= 2.9%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
plasma	16			= 2.5%
serum	16			= 2.4%

サンプルの種類 Serum, Plasma

アッセイタイプ Quantitative

検出範囲 5000 µg/dl - 50000 µg/dl

種交差性 **交差種:** Mammals, Other species

製品の概要 Sphingomyelin Assay Kit (ab133118) provides a specific, sensitive, and convenient method for quantifying sphingomyelin in plasma or serum. In this assay, sphingomyelinase is first used to hydrolyze sphingomyelin to phosphorylcholine and ceramide. Alkaline phosphatase then generates choline from the phosphorylcholine and the newly formed choline is used to generate hydrogen peroxide in a reaction catalyzed by choline oxidase. Finally, with peroxidase as a catalyst, hydrogen peroxide reacts with DAOS and 4-aminoantipyrine to generate a blue color with an optimal absorption at OD595 nm.

Under the standardized conditions of the assay described in this booklet, the dynamic range of the kit is 5-50 mg/dL sphingomyelin.

特記事項

Sphingomyelin (ceramide phosphorylcholine) is an important lipid component of cell membranes and lipoproteins. It consists of a ceramide moiety linked via a phosphodiester bond to phosphorylcholine. Sphingomyelinases are a family of enzymes that can hydrolyze sphingomyelin

into ceramide and phosphorylcholine. Ceramides have been implicated as key mediators in signaling pathways, with outcomes as diverse as cell proliferation, differentiation, growth arrest, and apoptosis. An inherited deficiency of acid sphingomyelinase activity results in the sphingomyelin storage disorder Niemann- Pick disease. This disease results in the accumulation of sphingomyelin in cells, tissues, and fluids. Since sphingomyelin has been implicated in the pathogenesis of several diseases, including atherosclerosis, sensitive and reliable techniques for its quantification are of considerable importance.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム

Microplate reader

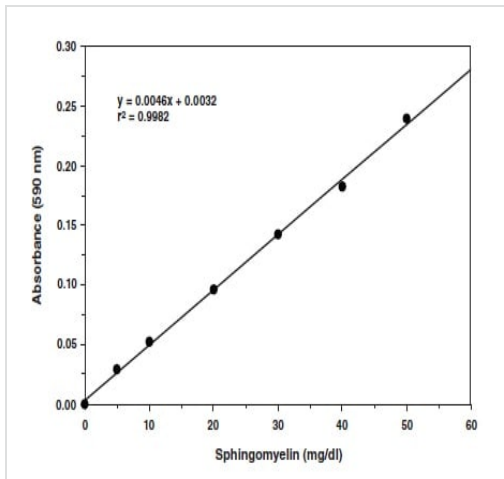
製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	96 tests
96-Well Plate (Colorimetric Assay)	1 unit
96-Well Plate Cover	1 unit
SM Alkaline Phosphatase	1 vial
SM Buffer (5X)	1 vial
SM Color Detector	2 vials
SM Enzyme Mixture	2 vials
Sphingomyelin Standard	1 vial
Sphingomyelinase	1 vial

画像



Typical standard curve obtained with sphingomyelin provided following the protocol.

Sphingomyelin standard curve.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors