

Sphingomyelin Assay Kit (Fluorometric) ab138877

6 References 1 Image

Overview

Product name	Sphingomyelin Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Cell culture extracts, Whole Blood
Assay type	Quantitative
Sensitivity	1 µM
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Abcam's Sphingomyelin Assay Kit (Fluorometric) (ab138877) provides the most sensitive method for detecting neutral SM activity or screening SM inhibitors. The kit uses AbRed Indicator as a fluorogenic probe to indirectly quantify the phosphocholine produced from the hydrolysis of sphingomyelin (SM) by sphingomyelinase (SMase). It can be used for measuring the amount of SM in blood, cell extracts or other solutions. The fluorescence intensity of AbRed Indicator is proportional to the formation of phosphocholine, therefore to the amount of SM. AbRed Indicator enables the assay readable by either a fluorescence reader or an absorbance reader.</p> <p>The kit is an optimized “mix and read” assay that can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step.</p>
Notes	<p>Sphingomyelin (SM) is largely found in the exoplasmic leaflet of the cell membrane, primarily in nervous tissue. It plays an important role in signal transduction. Sphingomyelin accumulates abnormally in Niemann-Pick disease and Abetalipoproteinemia.</p>
Platform	Microplate reader

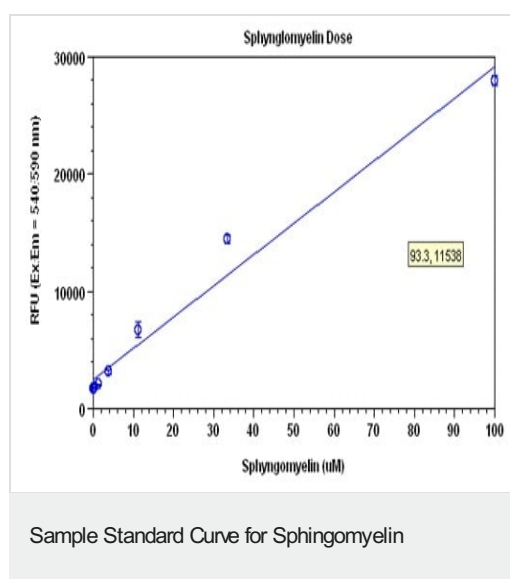
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
AbRed Indicator	1 vial
Assay Buffer	1 x 5ml
DMSO	1 x 200µl

Components	100 tests
Enzyme Mix	1 vial
SMase Reaction Buffer	1 x 20ml
50mM Sphingomyelin standard	1 x 20µl
Sphingomyelinase	1 vial

Images



Sphingomyelin dose response was measured on a solid black 96-well plate with ab138877 using a fluorescence microplate reader. As low as 1 µM sphingomyelin can be detected with 60 minutes incubation (n=3).

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.com/abpromise> or contact our technical team.

Terms and conditions

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