

Lipid Extraction Kit (Chloroform Free) ab211044

[13 References](#) [2 Images](#)

Overview

Product name	Lipid Extraction Kit (Chloroform Free)
Sample type	Serum, Plasma, Tissue, Adherent cells, Suspension cells
Assay time	0h 30m
Product overview	Abcam's Lipid Extraction Kit (Chloroform-Free) (ab211044) provides a chloroform-free method of lipid isolation. The recovered organic phase can be dried and resuspended for downstream lipid analysis. This protocol overcomes the difficulties of lower phase extraction and the dangers of using chloroform while resulting in highly efficient lipid extraction. Each kit provides sufficient reagents to isolate up to 50 lipid preparations based on a 25 µL sample size.

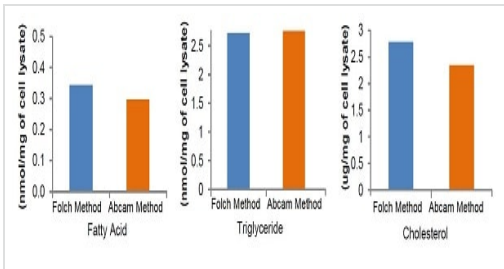
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K216 Lipid Extraction Kit (Chloroform Free). K216-50 is the same size as the 50 test size of ab211044.</p> <p>The term lipids describe a group of naturally occurring molecules that includes monoglycerides, diglycerides, triglycerides, phospholipids, fat-soluble vitamins, fats and sterols. Lipids define and preserve cellular membrane integrity and play an important role in cellular processes such as membrane trafficking, signal transduction and energy storage. Dysregulation of lipid metabolism is linked to many diseases like cancer, diabetes and heart disease. Traditional lipid extraction is performed using the Folch method, in which lipids are extracted into an organic phase. Isolating lipids from this organic phase is tricky and can result in contamination with material from other phases. These contaminants may have a negative impact on downstream lipid analysis. Furthermore, the Folch method uses chloroform, which is toxic, as the organic phase solvent.</p>
--------------	--

Properties

Storage instructions	Store at room temperature. Please refer to protocols.
-----------------------------	---

Components	50 tests
Lipid Extraction Buffer	1 x 25ml
Lysis Buffer II	1 x 5ml

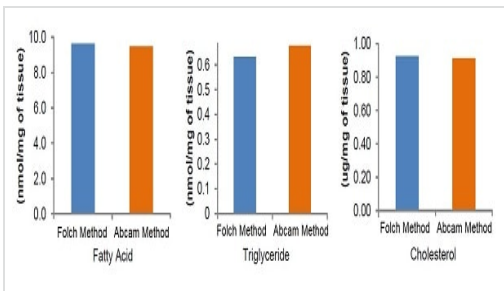
Images



Analysis of Total Free Fatty Acid, Triglyceride and Cholesterol on Extracted Lipids using Adipocytes Cell Lysate

Analysis of Total Free Fatty Acid, Triglyceride and Cholesterol on Extracted Lipids using Adipocytes Cell

Lysate. Comparison of Total Free Fatty Acid, Triglyceride and Cholesterol using Folch Method and Abcam's Lipid Extraction Method.



Analysis of Total Free Fatty Acid, Triglyceride and Cholesterol on Extracted Lipids using Rat Liver

Analysis of Total Free Fatty Acid, Triglyceride and Cholesterol on Extracted Lipids using Rat Liver Tissue.

Comparison of Total Free Fatty Acid, Triglyceride and Cholesterol using Folch Method and Abcam's Lipid Extraction Method.

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