

Adipogenesis Detection Assay Kit (Colorimetric/Fluorometric) ab102513

15 References **5 图像**

概述

产品名称	Adipogenesis Detection Assay试剂盒(Colorimetric/Fluorometric)
检测方法	Colorimetric/Fluorometric
样品类型	Tissue, Adherent cells, Suspension cells
检测类型	Quantitative
灵敏度	> 0.2 nM
范围	0.2 nM - 10 nM
检测时间	0h 40m
产品概述	Adipogenesis Detection Assay Kit (Colorimetric/Fluorometric) (ab102513) quantifies triglyceride accumulation in cells and tissues. In the assay, triglycerides are efficiently solubilized then hydrolyzed to glycerol which is subsequently oxidized to convert the probe to generate color (ODmax = 570 nm) and fluorescence (Ex/Em = 535/587 nm). The kit can detect 0.2 - 10 nmol of triglyceride in <1,000 differentiated 3T3-L1 cells. The high detection sensitivity and the convenient microplate assay format make the kit a convenient tool for studying the effect of adipogenesis inducers or to screen inhibitor compounds.

Visit our [FAQs page](#) for tips and troubleshooting.

说明 This product is manufactured by BioVision, an Abcam company and was previously called K610 Adipogenesis Colorimetric/Fluorometric Assay Kit. K610-100 is the same size as the 100 test size of ab102513.

Adipogenesis is the process of differentiation of different cell types into adipocytes, the primary fat storage cell type. The accumulation of adipocytes is the basis for obesity, a significant risk factor in many diseases, including diabetes, atherosclerosis, cancer and cardiovascular disease, etc. Adipocytes accumulate triglycerides, in the form of lipid droplets which can be measured.

平台 Microplate reader

性能

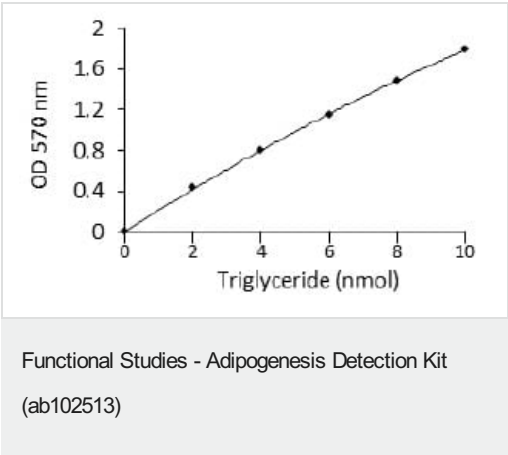
存放说明 Store at -20°C. Please refer to protocols.

组件	100 tests	2000 tests
Adipogenesis Lipid Extraction Solution	1 x 10ml	20 x 10ml
Assay Buffer V	1 x 25ml	20 x 25ml
Enzyme Mix VI	1 vial	20 vials
Lipase I	1 vial	20 vials
OxiRed Probe	1 x 200µl	20 x 200µl
Triglyceride Standard	1 x 0.3ml	20 x 0.3ml

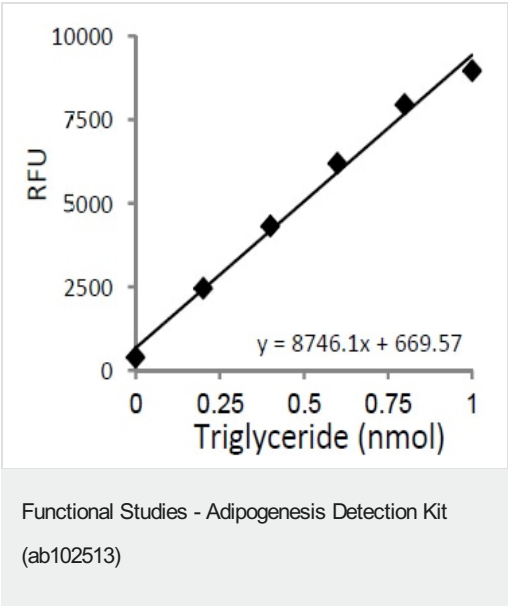
相关性

Adipogenesis is the process of differentiation of different cell types into adipocytes, the primary fat storage cell type. The accumulation of adipocytes is the basis for obesity, a significant risk factor in many diseases, including diabetes, atherosclerosis, cancer and cardiovascular disease, etc. Adipocytes accumulate triglycerides, in the form of lipid droplets which can be measured.

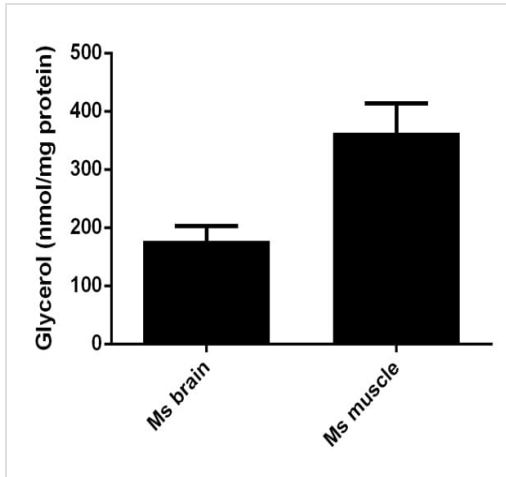
图片



Example of a standard curve obtained with ab102513.

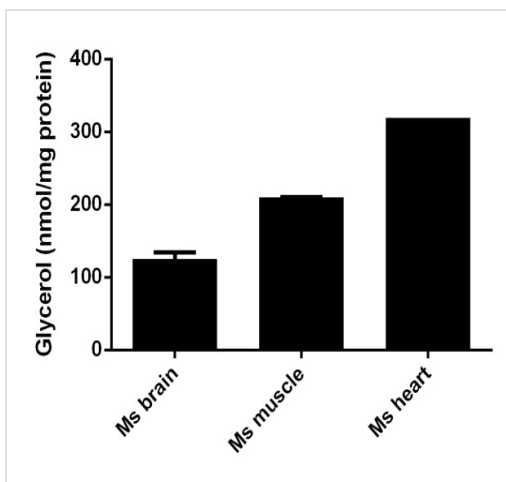


Example of a standard curve obtained with ab102513.



Glycerol measured fluorometrically in mouse tissue lysates showing quantity (nmol) per mg protein of tested sample

Functional Studies - Adipogenesis Detection Kit
(ab102513) - Fluorometric



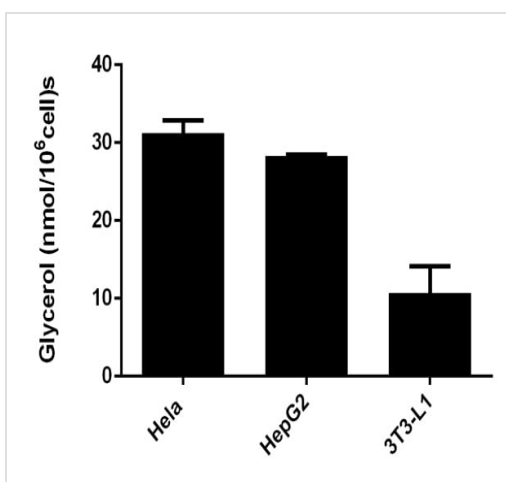
Glycerol measured colourimetrically in mouse tissue lysates showing quantity (nmol) per mg protein of tested sample.

Ms brain value retrieved from 1:9 dilution of original 12.4mg/mL sample.

Ms muscle value retrieved from 1:9 dilution of original 7.5mg/mL sample.

Ms heart value retrieved from 1:27 dilution of original 16.4mg/mL sample.

Functional Studies - Adipogenesis Detection Kit
(ab102513) - Colourimetric



Glycerol measured colourimetrically in cell lysates showing quantity (nmol) per 1 mln of tested cells

Functional Studies - Adipogenesis Detection Kit
(ab102513) - Colourimetric

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.cn/abpromise> or contact our technical team.

Terms and conditions

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