

Adipogenesis Detection Assay Kit (Colorimetric/Fluorometric) ab102513

15 References **画像数 5**

製品の概要

製品名	Adipogenesis Detection Assay Kit (Colorimetric/Fluorometric)
検出方法	Colorimetric/Fluorometric
サンプルの種類	Tissue, Adherent cells, Suspension cells
アッセイタイプ	Quantitative
検出感度	> 0.2 nM
検出範囲	0.2 nM - 10 nM
全工程の試験時間	0h 40m
製品の概要	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.

特記事項	<p>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.</p> <p>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.</p>
------	---

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

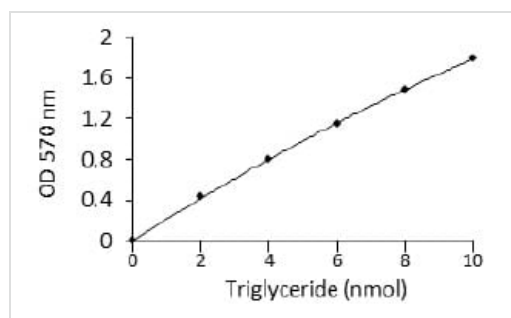
保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	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

関連性

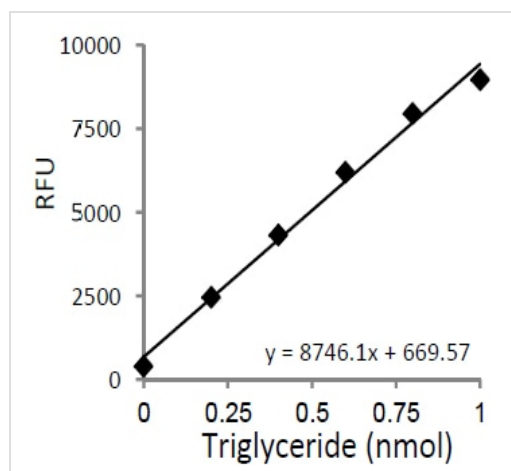
Adipogenesis is the process of differentiation of different cell types into adipocytes, the primary fat storage cell type. The accumulation of adipocytes is the basis for obesity, a significant risk factor in many diseases, including diabetes, atherosclerosis, cancer and cardiovascular disease, etc.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

画像



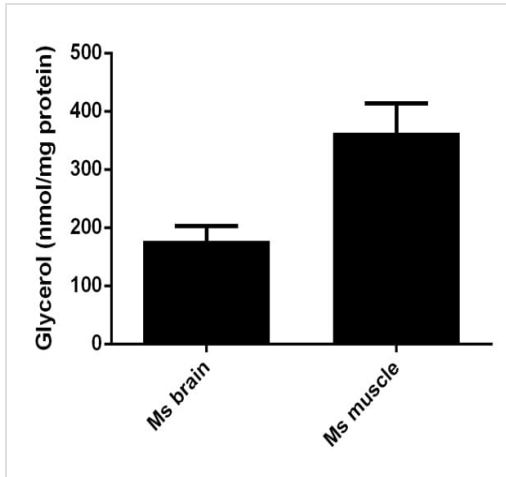
Example of a standard curve obtained with ab102513.

Functional Studies - Adipogenesis Detection Kit
(ab102513)



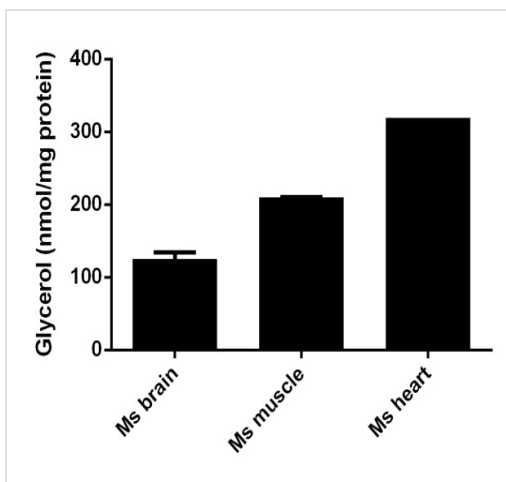
Example of a standard curve obtained with ab102513.

Functional Studies - Adipogenesis Detection Kit
(ab102513)



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

Functional Studies - Adipogenesis Detection Kit
(ab102513) - Fluorometric



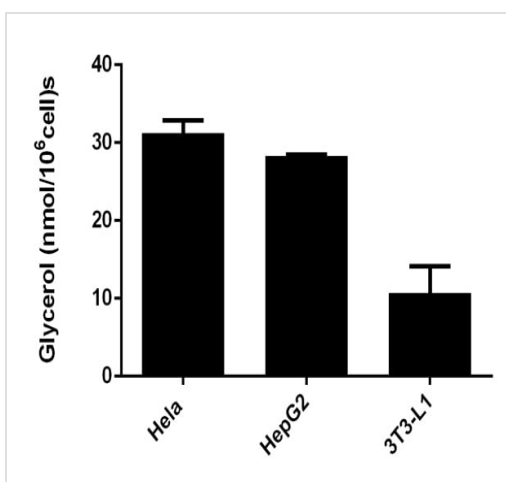
Glycerol measured colourimetrically in mouse tissue lysates showing quantity (nmol) per mg protein of tested sample.

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.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