

Adipolysis Assay Kit ab133115

[3 References](#) [3 Images](#)

Overview

Product name	Adipolysis Assay Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Adherent cells
Assay type	Quantitative
Species reactivity	Reacts with: Mammals, Other species
Product overview	Abcam's Adipolysis Assay Kit (ab133115) provides a convenient tool for studying the adipolysis of triglycerides in differentiated 3T3-L1 cells, using a common adipogenesis induction procedure. This kit will allow investigators to screen compounds involved in lipid storage and metabolism. Isoproterenol is included in the kit as a positive control for screening pharmaceuticals that regulate FFA release from adipocytes.

Triglycerides stored in these lipid droplets can be stimulated by known factors or compounds to release FFA and glycerol. The latter can be measured by incubation with glycerol kinase, glycerol phosphate oxidase, and horseradish peroxidase in the presence of a colorimetric substrate to generate a chromophore detectable at 540 nm. The amount of glycerol released into the medium is proportional to the level of triglyceride storage and/or degree of adipolysis.

Notes

Obesity is a risk factor in many of the major chronic diseases such as diabetes mellitus, and cancer. The ability to regulate the number of fat cells and/or the size of fat cells is a key in the development and physiology of obesity and also in the origin of chronic disease. Adipose tissue plays an important role in energy homeostasis. Mammals have both brown adipose tissue (BAT), which utilizes lipids to generate heat in a process known as thermogenesis, and white adipose tissue (WAT) which stores excess energy as triglycerides in lipid droplets. Triglycerides can be broken down into free fatty acids (FFA) and glycerol, through a process of adipolysis (also referred as lipolysis).

Adipolysis is a highly regulated process which allows appropriate delivery of FFA to meet energy needs. Appropriate FFA levels are important for optimal human health but excess FFAs are generally associated with development of insulin resistance, hypertriglyceridemia, reduced hepatic insulin clearance, and impaired β -cell insulin secretion. Adipolysis is under tight hormonal control.

3T3-L1 cells are a well-characterized model for studying lipid metabolism, including the uptake and release of fatty acids from fat cells. This model shares similar morphology, gene expression, and metabolism with adipocytes *in vivo*. During terminal differentiation, the fibroblast-like preadipocytes undergo a series of morphological and biochemical changes to eventually

accumulate lipid droplets.

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.

Platform

Microplate reader

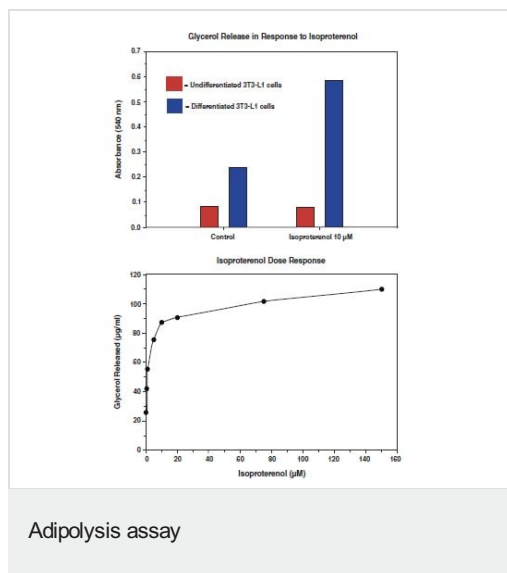
Properties

Storage instructions

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

Components	1 kit
Adipolysis Assay Dexamethasone Solution (1000X)	1 vial
Adipolysis Assay IBMX Solution (1000X)	1 vial
Adipolysis Assay Insulin Solution (1000X)	1 vial
Adipolysis Assay Isoproterenol Solution	1 vial
Free Glycerol Assay Reagent (10X)	2 vials
Glycerol Standard Solution	1 vial

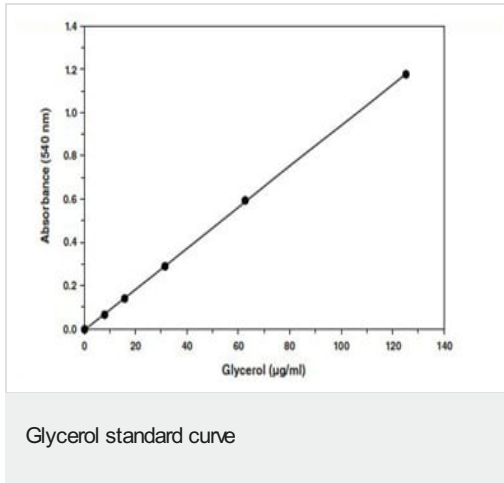
Images



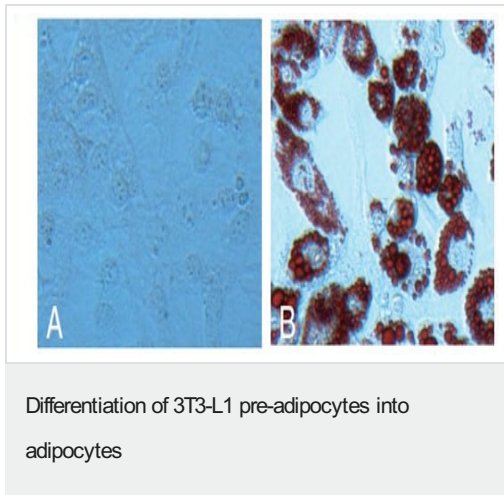
Differentiated 3T3-L1 cells were treated with increasing concentrations of isoproterenol for 48 hours and the culture supernatant was assayed for free glycerol.

Top panel: glycerol release from undifferentiated (*red*) and differentiated (*blue*) 3T3-L1 cells as determined using Free Glycerol Assay Reagent.

Bottom panel: isoproterenol dose response curve showing glycerol released after 24 hours *versus* isoproterenol concentration.



Example of standard curve of typical data produced with 3T3-L1 culture medium.



Upon induction, 3T3-L1 cells differentiate into adipocytes-like cells and they are ready to be used for adipolysis experiments.

Panel A: non-differentiated 3T3-L1 cells show no staining of Oil Red O Solution.

Panel B: increased lipid droplet accumulation indicates that more than 80% of preadipocytes were differentiated into adipocytes four days after weaning the cells from Induction Medium to Insulin Medium.

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