

Adipogenesis Assay Kit (Cell-Based) ab133102

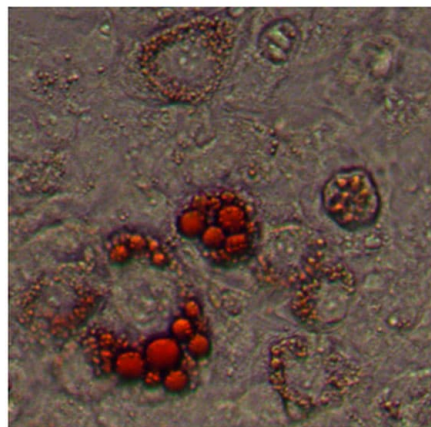
6 References 2 图像

概述	
产品名称	Adipogenesis Assay试剂盒(Cell-Based)
检测方法	Colorimetric
样品类型	Adherent cells
检测类型	Cell-based
种属反应性	与反应: Mammals, Other species
产品概述	Adipogenesis Assay Kit (ab133102) provides the reagents required for studying the induction and inhibition of adipogenesis in the established 3T3-L1 model using the adipogenesis induction procedure. This kit can also be used to screen drug candidates involved in adipogenesis. The classic Oil Red O staining for lipid droplets is used in this kit as an indicator of the degree of adipogenesis and can be quantified with a plate reader after the dye is conveniently extracted from the lipid droplet.
说明	<p>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. 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. Adipocytes are derived from multipotent mesenchymal precursor cells commit to preadipocytes and then either remain dormant or proceed to become differentiated adipocytes.</p> <p>3T3-L1 cells are a well-characterized cell line to study the differentiation of adipocytes, and other mechanisms such as insulin-induced glucose uptake and mechanisms of obesity development. This model system has greatly advanced the understanding of the molecular basis and signaling pathways of adipogenesis. During terminal differentiation, the fibroblast-like preadipocytes undergo a series of morphological and biochemical changes to eventually accumulate lipid droplets. This <i>in vitro</i> differentiated adipocytes share similar morphology with adipocytes <i>in vivo</i>.</p>
平台	Microplate reader
性能	
存放说明	Please refer to protocols.

组件	1 kit
Adipogenesis Assay Dexamethasone Solution	1 x 500μl
Adipogenesis Assay Insulin Solution (1,000X)	1 x 1.5ml
Cell Based Assay IBMX Solution (1,000X)	1 x 500μl
Fixative (10X)	1 x 10ml
Lipid Droplets Assay Dye Extraction Solution	1 x 30ml
Lipid Droplets Assay Oil Red O Solution	1 x 25ml
Lipid Droplets Assay Wash Solution	6 x 30ml

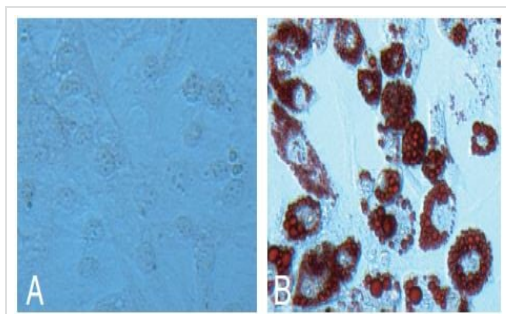
相关性 Adipogenesis is the process of differentiation of different cell types into adipocytes, the primary fat storage cell type. The accumulation of adipocytes is the basis for obesity, a significant risk factor in many diseases, including diabetes, atherosclerosis, cancer and cardiovascular disease, etc.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

图片



Lipid droplet accumulation in the differentiated 3T3-L1 cells visualized by Oil Red O Solution staining.

Adipogenesis Assay Kit (Cell-Based) (ab133102)



Functional Studies - Adipogenesis Assay Kit (Cell-Based) (ab133102)

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

Panel A: Non-differentiated 3T3-L1 cells were not stained by Oil Red O Solution.

Panel B: more than 80% of preadipocytes were differentiated four days after weaning the cells from induction medium to insulin medium. Lipid droplet accumulation in the differentiated cells can be visualized by Oil Red O Solution staining.

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