



## **Adipogenesis Assay Kit**

Catalog #: ENZ-KIT103-0005

5 x 24 Well or 225 cm<sup>2</sup> Tissue Culture Plate Surface Area

## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## TABLE OF CONTENTS

Description .....	1
Background .....	1
Safety Warning and Precautions .....	3
Materials Supplied .....	3
Storage .....	4
Other Materials Needed .....	4
Optional Material .....	4
Important Technical Notes .....	5
Workflow Diagram .....	6
Assay Procedure .....	6
Typical Results .....	9
References .....	10
Contact Information .....	13

## DESCRIPTION

The Adipogenesis Assay Kit is a complete kit for the qualitative determination of differentiation of 3T3L1 cells into adipocytes.

All of reagents needed to induce and detect differentiation are provided. This kit can be used to either test agonists that induce adipogenesis or antagonists that inhibit it from occurring. The protocol provided is for use with 24 well plates, but can be adapted for use with other plate sizes. The stain provided allows for visual confirmation and can also be easily extracted allowing measurement using a spectrophotometer or microplate reader. Please read the complete kit insert before performing this assay.

## BACKGROUND

Obesity is a serious health problem that leads to elevated triglyceride levels, which can lead to cardiovascular disease, diabetes, hypertension, and hypercholesterolemia<sup>5,13</sup>. It is characterized by the enlargement of adipose tissue to store excess triglyceride intake<sup>5</sup>. This can occur through two different mechanisms, known as hyperplasia (increase in the number of adipocytes) and hypertrophy (increase in the volume of adipocytes)<sup>3</sup>. Triglycerides accumulate in lipid droplets that are embedded in the cytoplasm of the adipocyte<sup>13</sup>. Lipid droplets are intracellular organelles that are involved in the regulation of the storage and hydrolysis of triglycerides and cholesterol esters, which are both neutral lipids<sup>13</sup>. Lipid droplets are found in two types of adipose tissue. White adipose tissue is made up of white adipocytes, which are composed of a single, large lipid droplet that takes up most of the cell volume and functions in neutral lipid storage<sup>9,10</sup>. Brown adipose tissue is made up of brown adipocytes, which are composed of several small lipid droplets with increased mitochondria in the cytoplasm<sup>9,10</sup>. This tissue functions in maintaining body temperature during long periods of cold exposure<sup>10</sup>.

Essential to understanding and treating disease, is the identification of drugs that modulate differentiation and adipogenesis. Agonists of PPAR $\gamma$  play an important role in decreasing insulin resistance<sup>6</sup>. Rosiglitazone, a thiazolidinedione that binds to PPAR $\gamma$  and sensitizes fat cells to insulin, is used clinically as an anti-diabetic drug in the treatment of type 2 diabetes<sup>6,12</sup>. It enhances adipogenesis and stimulates hyperplasia by speeding up differentiation and increasing the degree of differentiation<sup>7,8</sup>. Antagonists also play an important role in

treating obesity. Pro-inflammatory cytokines, such as TNF- $\alpha$  and IL-1 inhibit adipogenesis by decreasing the expression of PPAR $\gamma$  and CEBP $\alpha$ <sup>9,11</sup>. Inhibitors of PPAR $\gamma$  have been identified as a potential target of anti-obesity drugs<sup>2</sup>.

One of the most utilized models for the study of differentiation of fibroblast into adipocytes is the 3T3-L1 cell line<sup>1,5,9</sup>. Differentiation requires three components, which include insulin or insulin-like growth factor, dexamethasone (DXM), and 3-isobutyl-1-methylxanthine (IBMX)<sup>5,7,9</sup>. Insulin induces proliferation and differentiation of preadipocytes<sup>8,9</sup>. DXM is a glucocorticoid, which activates CEBP $\beta$  (CCAAT/enhancer binding protein- $\beta$ )<sup>5,8</sup>. IBMX is a competitive, nonselective phosphodiesterase inhibitor that increases intracellular cAMP levels and activates CEBP $\delta$ <sup>5,8</sup>. It also increases intracellular protein kinase A levels, which is important for the transcriptional activation of PPAR $\gamma$  (peroxisome proliferator-activated receptor- $\gamma$ )<sup>8</sup>. The combination of DXM and IBMX regulates PPAR $\gamma$ , which is a nuclear hormone receptor that is highly expressed in adipose tissue and promotes adipogenesis<sup>2,3,12</sup>. CEBP $\beta$  and CEBP $\delta$  induce transcription of CEBP $\alpha$  and PPAR, which in turn activates the expression of adipocyte-specific genes<sup>11</sup>. During the differentiation process, 3T3-L1 cells undergo a post-confluent mitosis, which occurs 24 hours after induction with insulin, DXM, and IBMX, followed by growth arrest<sup>5</sup>. After growth arrest occurs, the cells are committed to becoming adipocytes and express late markers of differentiation at day 3, after induction<sup>5</sup>. Growth arrest of the cells is a requirement for terminal adipocyte differentiation<sup>5,9</sup>. After 5-7 days post-induction, the cell morphology changes from the elongated, fibroblastic cells to rounded cells, and lipid droplets begin to accumulate<sup>4,5</sup>.

## SAFETY WARNINGS & PRECAUTIONS

*FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.*

- Refer to the products MSDS before use.
- Reagents should be treated as though they are potentially toxic. They should be handled with care and disposed of properly.
- Observe good laboratory practices. The use of protective equipment including, nitrile gloves, lab coat, and protective eyewear is recommended.
- Solutions in DMSO can be readily absorbed through the skin.
- Adipogenesis Dye will stain clothing and exposed skin.
- Isopropanol is flammable. Keep away from open flames.



Handle  
with care

## MATERIALS SUPPLIED

1. **Differentiation Solution, 65  $\mu$ L, Catalog No. 80-2634:**  
A 1000X solution of dexamethasone and 3-Isobutyl-1-methylxanthine (IBMX) in DMSO.
2. **Insulin Solution, 130  $\mu$ L, Catalog No. 80-2635:**  
A 1000X solution of insulin in 0.02N HCl.
3. **Enhancer Solution, 65  $\mu$ L, Catalog No. 80-2639:**  
A 1000X solution of rosiglitazone in DMSO.
4. **Cell Fixative, 28 mL, Catalog No. 80-2636:**  
A 1X solution of 3.7% formaldehyde in phosphate buffered saline.
5. **Adipogenesis Dye, 28 mL, Catalog No. 80-2637:**  
A 1X solution of Oil Red O lipid dye in 60% isopropanol.
6. **Extraction Solution, 28 mL, Catalog No. 80-2638:**  
A 1X solution containing 60% isopropanol.



Reagents require separate storage conditions.

## STORAGE

All components of this kit are shipped at  $-20^{\circ}\text{C}$ . Upon receipt the Differentiation Solution, Insulin Solution, and Enhancer Solution should be stored at or below  $-20^{\circ}\text{C}$ . The Cell Fixative, Adipogenesis Dye, and Extraction Solution can be stored at room temperature. If stored properly all reagents should be stable until the kits expiration date. Induction media is stable for 2 weeks at  $4^{\circ}\text{C}$ .

**Note:** The Cell Fixative and Adipogenesis Dye will form insoluble precipitates especially upon freezing. Removal of the precipitate from the Cell Fixative is recommended but not required for performance; however removal of the precipitate from the dye is required for assay performance. Filtering either reagent will have no effect on their performance in the assay.

## OTHER MATERIALS NEEDED

1. Preadipocyte cell line (Ex. 3T3-L1, ATCC # CL-173)
2. 24 well or other size tissue culture treated plate.
3. Tissue culture media for growing 3T3-L1 cells (DMEM +10% FBS)
4. Deionized or distilled water.
5. Precision pipets for volumes between  $5\ \mu\text{L}$  and  $1,000\ \mu\text{L}$ .
6. Spectrophotometer or microplate reader capable of reading 96 well plates at 490-520nm.
7.  $0.2\ \mu\text{M}$  PES syringe filter

## OPTIONAL MATERIAL

1. Collagen I, rat tail (thin layer coating), Enzo Catalog No. ALX-522-440-0050.

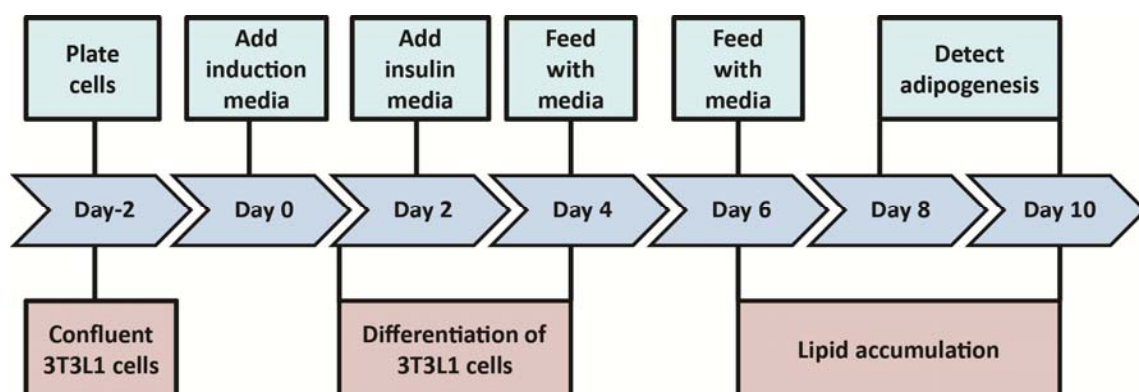
## IMPORTANT TECHNICAL NOTES

1. The 3T3-L1 cell line must be maintained as a subconfluent monolayer prior to use in the assay. However, actively growing cells will not undergo adipogenesis.
2. Low passage cells are highly recommended. Higher passage number cells ( $\geq 10$ ) have been shown to cause increased OD's in untreated cells and may dislodge from the plate when treated. Improperly maintained cells may adversely affect assay performance.
3. Other preadipocyte cell lines may be used but it is up to the user to validate the results.
4. This kit was validated using 24 well plates, however if the use of smaller plates is required the user must determine the appropriate culture conditions. Use tremendous care when changing the media. As stated above the cells have a tendency to pull away from the edges and form a clump of cells. This may lead to variability in the data. The use of collagen coated plates may help.
5. Fetal bovine serum lots may vary potentially causing increased background. The use of calf serum may help to eliminate background in the assay.
6. The components of the kit are not sterile. They should be sterile filtered through a 0.2  $\mu\text{M}$  PES filter before adding to cells to prevent contamination. The use of antibiotics in the media does not affect assay performance.
7. Absorbance values may vary due to plate reader differences and plate characteristics.
8. The Cell Fixative and Adipogenesis Dye will likely form insoluble precipitates which should be removed by centrifugation or filtering through a 0.2  $\mu\text{M}$  PES filter.
9. Adipogenesis Dye can be read between 490nm - 520nm. The Adipogenesis Dye will stick to the plastic so it is recommended that non-specific background be subtracted from a blank plate.



- It is recommended that only the amount of induction media or insulin media needed for each assay be made. However, if there is extra it may be stored at 4°C for up to two weeks without significant loss of activity. The stability of drugs that are being used as test conditions in the assay must be evaluated by the user.

## WORKFLOW DIAGRAM



Basic Workflow of Adipogenesis Assay. Procedure on top row, cell characterization on bottom row.

## ASSAY PROCEDURE

### Cell Culture (volumes are for 1 x 24 well plate):

**Note:** For different size plates adjust the volumes accordingly.

- Plate 1 mL of subconfluent 3T3L1 cells @  $0.25 - 0.5 \times 10^6$  / mL ( $\leq$  passage 10) onto the appropriate number of wells of a 24 well plate.
- Grow 2 - 4 days at 37°C, 5% CO<sub>2</sub> until they are 2 days past confluent (This is Day 0).
- On Day 0 prepare Induction Media by adding 12  $\mu$ L of differentiation solution, 12  $\mu$ L enhancer solution, and 12  $\mu$ L of insulin solution to 12 mL of DMEM + 10% FBS + 1X Penstrep.
- Aspirate media from wells.
- For induced wells: Add 0.5 mL Induction Media (from step 3) per well.

For uninduced wells: Add 0.5 mL DMEM + 10% FBS + 1X PenStrep per well.

6. Incubate for 2 days at 37°C, 5% CO<sub>2</sub>.
7. Prepare Insulin Media by adding 12 µL of insulin solution to 12 mL of DMEM + 10% FBS + 1X penstrep.
8. Carefully aspirate media from wells. **(Note: Differentiated cells have a tendency to detach from wells).**
9. For induced wells: Add 0.5 mL Insulin Media (from step 7) per well.  
  
For uninduced wells: Add 0.5 mL DMEM + 10% FBS + 1X PenStrep per well.
10. Incubate for 2 days at 37°C, 5% CO<sub>2</sub>.
11. Remove media from all wells and feed all cells with 1 mL of DMEM + 10% FBS + 1X penstrep.
12. Feed the cells every 2 days with 1mL of DMEM + 10% FBS + 1X penstrep for at least 4-6 more days. At 8-10 days post induction, full differentiation into adipocytes should be achieved. Fully differentiated cells have large oil droplets that are visible and often appear viscous when feeding. If cells do not appear to be fully differentiated, continue incubation and feed every 2 days with DMEM + 10 FBS + 1X penstrep.

**Staining:**

**The Adipogenesis Dye can stick to the plastic causing non-specific background to occur. To correct for this perform the staining procedure on several empty well from an empty 24 well plate. The average value can then be subtracted from both uninduced and induced samples to remove non-specific background staining.**

13. If needed centrifuge or filter the Adipogenesis Dye using a 0.2 µM filter if precipitate is observed. **Precipitated dye will lead to variability in the results.**
14. Aspirate media from cells.
15. Add 200 µL of Cell Fixative to all wells and allow the plate to sit at room temperature for 30 minutes. Remove cell fixative and add 200 µL of Adipogenesis Dye per well, and allow the plate to sit at room temperature for 30min.

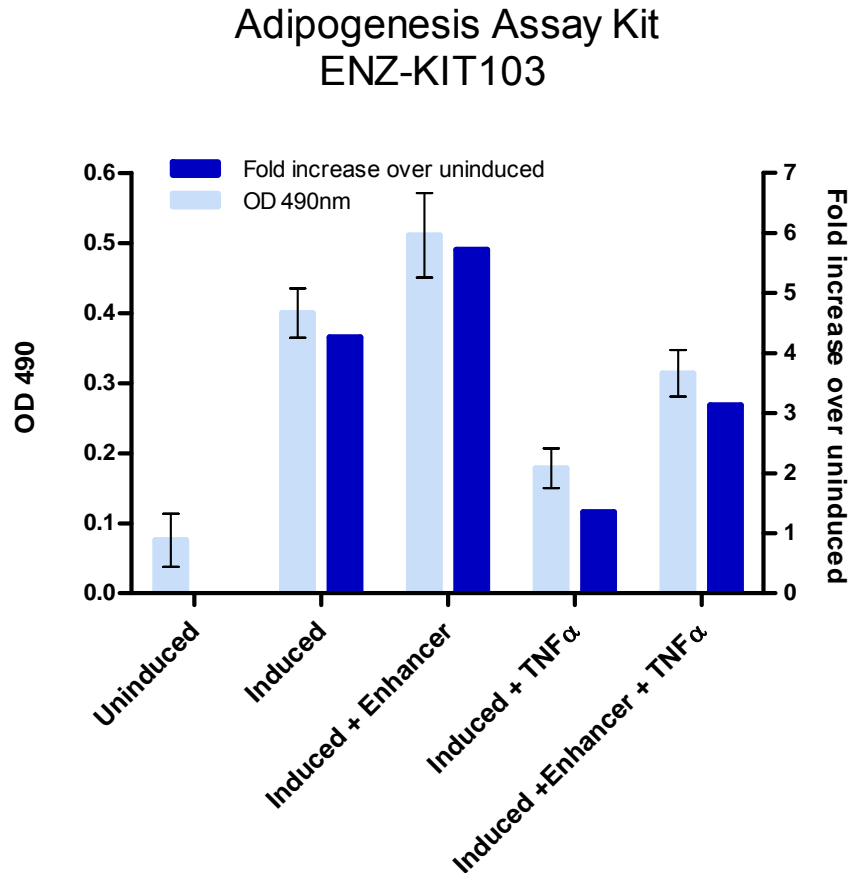


# Product Manual

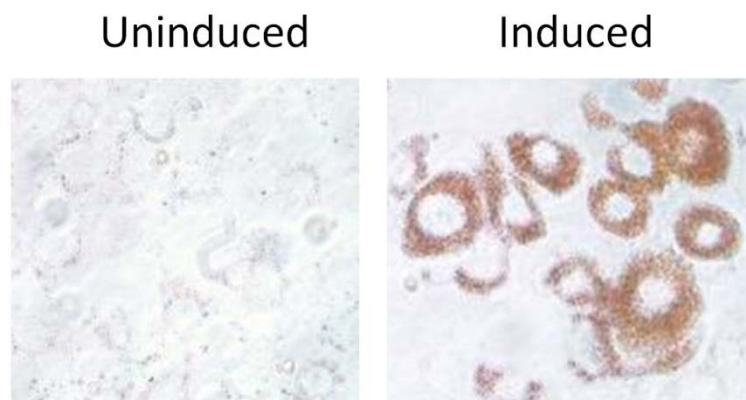
16. Remove the stain and rinse each well with 3X with 1 mL distilled water.
17. Allow the wells to dry completely.
18. If desired the stained cells can be visualized using microscopy.
19. Once the wells are dry add 200  $\mu$ L of Extraction Solution per well and incubate for 30min at room temperature with shaking (if possible).
20. Transfer contents of each well to a new 96 well plate and read at 490nm.
21. Remove non-specific background by subtracting the average OD 490nm readings obtained from empty wells from each sample well.

## TYPICAL RESULTS

The results shown below are for illustration only.



**Figure 1. Induction Levels:** OD 490nm values were obtained after extraction and the average background was subtracted from each well. The fold increase over the uninduced condition was determined using the following calculation:  $(\text{Induced OD 490nm} - \text{Uninduced OD 490nm}) / \text{Uninduced OD 490nm}$



**Figure 2. Microscopic Analysis of Induction:** 3T3-L1 cells were treated with and without induction media. After 10 days they were stained and imaged at 40X magnification.

**REFERENCES**

1. H. Green, et al., *Cell*, (1975) 5(1):19–27.
2. H.M. Wright, et al., *J. Biol. Chem.*, (2000) 275(3):1873-1877.
3. J.E.B. Reusch, et al., *Mol. Cell. Biol.*, (2000) 1008-1020.
4. J.L. Ramírez-Zacarías, et al., *Histochemistry.*, (1992) 97:493-497.
5. J.M. Ntambi, et al., *J.Nutr.*, (2000) 130: 3122S-3126S.
6. J.T. Crossno Jr., et al., *J. Clin. Invest.*, (2006) 116(12):3220-3228.
7. K. Zebisch, et al., *Anal. Biochem.*, (2012) 425:88-90.
8. M.A. Scott, et al., *Stem Cells Dev.*, (2011) 20(10):1793-1804.
9. M.E. Symonds, “*Adipose Tissue Biology*”, (2012) New York: Springer.
10. M. Rosenwald, et al., *Adipocyte.*, (2014) 3(1):4-9.
11. S. Kurebayashi, et al., *Endocr. J.*, (2001) 48(2):249-253.
12. T. Albrektsen, et al., *Diabetes.*, (2002) 51:1042-1051.
13. V. Rizzatti, et al., *Eur. J. Histochem.*, (2013) 57:e24.



# Product Manual

## NOTES

---



# Product Manual

## NOTES

---



# Product Manual

**[www.enzolifesciences.com](http://www.enzolifesciences.com)**

***Enabling Discovery in Life Science<sup>®</sup>***

**GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**EUROPE/ASIA**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
[info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

For local distributors and detailed product information visit us online:

**[www.enzolifesciences.com](http://www.enzolifesciences.com)**