

# Adipogenesis assay kit

**Convenient method for the enhanced induction and detection of adipogenesis in the classic 3T3L-1 model.**

The Adipogenesis Assay Kit provides all the reagents needed to induce and detect adipogenesis. Optimized Differentiation, Insulin, and Enhancer solutions ensure maximal adipocyte induction. The Adipogenesis Dye allows for visual confirmation of differentiation and can also be easily extracted allowing measurement using a spectrophotometer or microplate reader.

The ability to regulate the cell cycle and differentiation of adipocytes are key in the development and physiology of obesity. One of the most utilized models for the study of differentiation of fibroblast into adipocytes is the 3T3-L1 cell line. Differentiation of 3T3-L1 cells into adipocytes requires three primary components: insulin or insulin-like growth factor, dexamethasone (DXM), and 3-isobutyl-1-methylxanthine (IBMX). In addition, rosiglitazone, a thiazolidinedione that binds to PPAR $\gamma$  and sensitizes fat cells to insulin, enhances adipogenesis.

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. After growth arrest occurs, the cells are committed to becoming adipocytes and express late markers of differentiation at day 3, after induction. Growth arrest of the cells is a requirement for terminal adipocyte differentiation. After 5-7 days post-induction, the cell morphology changes from the elongated, fibroblastic cells to rounded cells, and lipid droplets begin to accumulate.

## Ordering Information

[Order Online »](#)

ENZ-KIT103-0005	5x24 wells
-----------------	------------

## Manuals, SDS & CofA

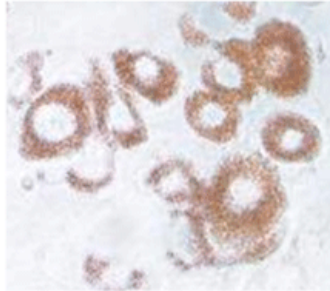
[View Online »](#)

- Provides optimized reagents for adipocyte differentiation and subsequent colorimetric lipid oil detection
- Contains enhancer solution for the optimal induction of adipogenesis
- For drug-screening and testing agonists/antagonists of adipogenesis

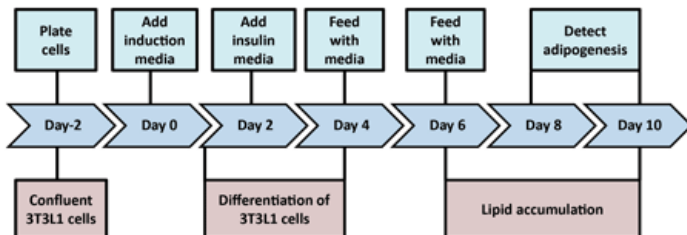
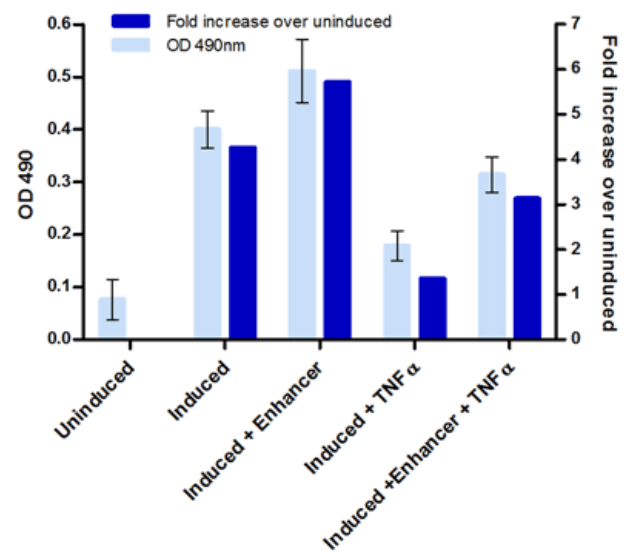
Uninduced



Induced



## Adipogenesis Assay Kit ENZ-KIT-103



## Handling & Storage

<b>Use/Stability</b>	Store Differentiation Solution, Insulin Solution, and Enhancer Solution at -20°C. All other reagents can be stored at room temp.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Application</b>	Colorimetric detection
<b>Application Notes</b>	For the induction and colorimetric detection of adipogenesis.
<b>Contents</b>	1000x Differentiation Solution (Dexamethasone and IBMX) 1000x Insulin Solution 1000x Enhancer Solution (Rosiglitazone) 1x Cell Fixative 1x Adipogenesis Dye (Oil Red O) 1x Extraction Solution
<b>Quantity</b>	Contains enough reagents to conduct experiments for: 5 x 24 well plates or 225 cm <sup>2</sup> tissue culture plate surface area



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)