



# SensoLyte<sup>®</sup> ONPG $\beta$ -Galactosidase Assay Kit

## *\*Colorimetric\**

Revision Number: 1.1	Last updated: October 2014
<b>Catalog #</b>	<b>AS-72134</b>
<b>Kit Size</b>	500 Assays (96-well)

- **Optimized Performance:** This kit is optimized to detect  $\beta$ -galactosidase activity.
- **Enhanced Value:** It provides ample reagents to perform 500 assays in a 96-well format.
- **High Speed:** The entire process can be completed within 30 minutes.
- **Assured Reliability:** Detailed protocol is provided.

---

## Kit Components, Storage and Handling

Component	Description	Quantity
Component A	Substrate solution	50 mL
Component B	$\beta$ -galactosidase enzyme	0.1 mg/mL, 20 $\mu$ L
Component C	Lysis buffer	100 mL
Component D	Stop solution	30 mL
Component E	DTT	1 M, 5 mL

### Other Materials Required (but not provided)

- 96-well microplate: Clear microplates provide better signal for absorbance reading.
- Microplate reader: Capable of detecting absorbance at 420 nm.

### Storage and Handling

- Store all components at -20°C.
- Keep component A away from light.
- Component D can be stored at room temperature for convenience.

---

## Introduction

Reporter enzymes are commonly used in cell biology to study transcriptional activity of genes.  $\beta$ -galactosidase, is one of the first and most popular reporter enzymes.  $\beta$ -galactosidase, encoded by the *lacZ* gene of *E. coli*, catalyzes the hydrolysis of  $\beta$ -galactosides into monosaccharides. It is widely used as a reporter enzyme to study gene expression, protein-protein interactions,<sup>1</sup> and normalization of transfection efficiency.<sup>2</sup>

The SensoLyte<sup>®</sup> ONPG  $\beta$ -Galactosidase Assay Kit utilizes the widely used chromogenic substrate, O-nitrophenyl  $\beta$ -D-galactopyranoside (ONPG) to determine the activity of  $\beta$ -galactosidase in transfected cells or tissues. The colorless compound is hydrolyzed into galactose and ortho-nitrophenol in the presence of  $\beta$ -galactosidase. Upon  $\beta$ -galactosidase protease cleavage, the o-nitrophenol produces a yellow color that can be detected at absorbance 420 nm. The intensity of the color produced is proportional to the concentration of  $\beta$ -galactosidase; thus can be used to measure  $\beta$ -galactosidase activity.

---

## Protocol

### 1. Prepare working solutions.

Note: Warm all kit components to room temperature before starting the experiment.

1.1 Prepare cell extracts: Aspirate culture medium from cells. Wash cells by using PBS or Hanks' Balanced Salts Solution (HBSS). Add 50  $\mu$ L of lysis buffer (Component C) to the microplate wells. Incubate at room temperature for 10 min to allow cell lysis.

1.2 Substrate solution: Dilute DTT (Component E) 1:10 in substrate solution (Component A). Use this DTT-containing substrate solution in all subsequent steps. Prepare fresh substrate solution for each experiment. Adjust volume as needed.

Table 1. ONPG substrate solution for one 96-well plate (100 assays)

Components	Volume
ONPG substrate solution (Component A)	45 mL
DTT (Component E)	5 mL
Total volume	50 mL

1.3 Prepare dilutions of  $\beta$ -galactosidase standard (optional): Dilute  $\beta$ -galactosidase (Component B) to 1000 ng/mL (1:100) in lysis buffer (Component C). Then make 2-fold serial dilutions to get concentration of 500, 250, 125, 62.5, 31.25, 15.63 ng/mL, include a blank control.

### 2. Set up the enzymatic reaction.

2.1 Add 10  $\mu$ L of cell extracts containing  $\beta$ -galactosidase to the wells.

Note: The amount of cell extract can be adjusted depending on the level of enzyme in the samples. Use lysis buffer to dilute test samples.

2.2 Set up  $\beta$ -galactosidase standard (optional): Add 10  $\mu\text{L}$  of serially diluted  $\beta$ -galactosidase reference solutions to the wells. The final amounts of  $\beta$ -galactosidase are 10, 5, 2.5, 1.25, 0.625, 0.3125, 0.156, and 0 nanogram/well.

2.3 Simultaneously establish the following control wells, as deemed necessary:

- Negative control contains 10  $\mu\text{L}$  of biological sample without  $\beta$ -galactosidase.

### 3. Detect $\beta$ -galactosidase activity.

3.1 Add 90  $\mu\text{L}$ /well of substrate solution into each well. Mix the reagents completely by shaking the plate gently for 30 sec.

3.2 Measure absorbance signal: Incubate reaction for 30 min at 37°C. Optional: Add 50  $\mu\text{L}$  of stop solution (Component D) to each well. Mix the reagents and measure absorbance at 420 nm.

3.3 Data Analysis:

- The absorbance reading from the blank control well is used as the background absorbance. Subtract the background reading from the readings of the other wells containing substrate.
- To evaluate the amount of  $\beta$ -galactosidase in the samples, use the enzyme standard curve as shown in Figure 1.

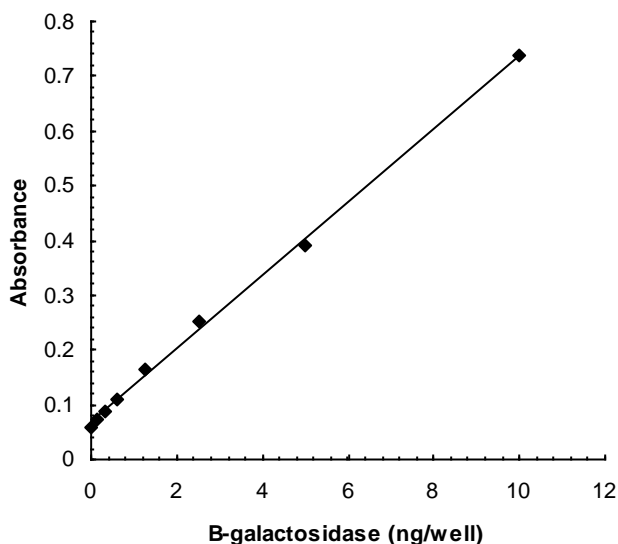


Figure 1. Detection of  $\beta$ -galactosidase with the SensoLyte<sup>®</sup> ONPG  $\beta$ -galactosidase Activity Assay Kit. The detection limit can reach as low as 0.1 ng.

---

## References

1. Rossi, F. et al. *PNAS* 94, 8405 (1997).
2. Thompson CD et al. *Biotechniques* 27, 824 (1999).