

# $\beta$ -Galactosidase Activity Assay Kit (Colorimetric)

Catalog #: MA-BGAL

ISO 13485:2016

## Introduction

$\beta$ -galactosidase is one of the first and most popular reporter enzymes which are commonly used in cell biology to study the transcriptional activity of genes.  $\beta$ -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, and normalization of transfection efficiency.

$\beta$ -Galactosidase Activity Assay Kit provides a simple, reproducible, and sensitive tool for measuring  $\beta$ -Galactosidase activity in cell lysates, plasma, serum, and other biological liquid samples. In this assay, O-nitrophenyl  $\beta$ -D-galactopyranoside (ONPG) 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 an absorbance of 420 nm. The optical density measured is directly proportional to the activity of  $\beta$ -Galactosidase present in the sample.

## Storage

The entire kit should be stored below -20 °C for up to 6 months from the date of shipment. For prepared reagent storage, see the table below.

Component	Size / Description	Storage After Preparation
Microplate	A 96-well (12 strips x 8 wells) plate	RT
Pre-Assay Buffer	15 ml	2–8 °C, 3 months
$\beta$ -Galactosidase Standard	1 vial (10 $\mu$ L)	-20 °C, 3 months
Substrate	3 vials	Use immediately
$\beta$ -Mercaptoethanol	1 vial (100 $\mu$ L)	2–8 °C, 3 months
Stop Solution	15 ml	2–8 °C, 3 months

RT = room temperature

## Additional Materials Required

1. Microplate reader capable of measuring absorbance at 420 nm
2. Incubator at 37 °C
3. Precision pipettes to deliver 2  $\mu$ l to 1 ml volumes
4. Tubes to prepare sample dilutions
5. 15 ml conical tubes

## Sample Tips and General Considerations

NOTE: Optimal methods of sample preparation will need to be determined by each researcher empirically based on researched literature and knowledge of the samples.

- If not using fresh samples, freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot samples prior to initial storage.
- It is strongly recommended to add a protease inhibitor cocktail to cell and tissue lysate samples.
- Avoid sonication of 1 ml or less as this can quickly heat and denature proteins.
- Most samples will not need to be concentrated. If concentration is required, a spin column concentrator with a chilled centrifuge is recommended.

### 1. Cell lysates can be prepared as follows:

For attached cells, remove supernatant from cell culture and wash cells twice with cold 1X PBS (for suspension cells, pellet the cells by spinning down the cells at 1,000 x g for 10 min), making sure to remove any remaining PBS before adding lysis buffer. Solubilize the cells at  $2 \times 10^7$  cells/ml in lysis buffer containing protease inhibitors. Pipette up and down to resuspend cells and rock the lysates gently at 2–8 °C for 30 minutes. Transfer extracts to microfuge tubes and centrifuge at 14,000 x g for 10 minutes.

It is recommended that sample protein concentrations should be determined using a total protein assay. Lysates should be used immediately or aliquot and stored at -70 °C. Thawed lysates should be kept on ice prior to use.

General tips for preparing lysate samples can be viewed on the online Resources page of the website:

<https://www.raybiotech.com/tips-on-sample-preparation/>

### 2. Plasma samples:

Collect blood with an anticoagulant such as citrate, EDTA, or oxalate and mix by inversion. Centrifuge the blood at 1000 x g at 4°C for 10 minutes. Collect plasma supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage. The sample may be assayed directly or diluted as necessary in Assay Buffer.

### 3. Serum samples:

Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage. The sample may be assayed directly or diluted as necessary in Assay Buffer.

***NOTE:*** Optimal experimental conditions for samples must be determined by the investigator. A set of serial dilutions is recommended for samples to achieve optimal assay results.

## Reagent Preparation

REAGENT	PREPARATION
Microplate	No preparation
Pre-Assay Buffer	
$\beta$ -Mercaptoethanol	
Stop Solution	
<b>Assay Buffer</b>	Immediately before using, add 17.5 $\mu$ L $\beta$ -Mercaptoethanol into 10 mL Pre-Assay Buffer to create Assay Buffer. Mix well. Prepare additional Assay Buffer as needed.
$\beta$ -Galactosidase Standard	See "Standard Preparation" section.
Substrate	Quickly spin down each vial <u>immediately before use</u> . Reconstitute with 1 ml Assay Buffer.
<b>Working Solution</b>	Add 1mL reconstituted Substrate solution into 1800 $\mu$ L Assay Buffer to make the <b>Working Solution</b> . This is enough to assay 38 wells. Prepare the additional vials as needed. <u>The Working Solution should be used immediately after preparation. Keep it in the dark, on ice.</u>

## Standard Preparation

To prepare a dilution series of  $\beta$ -Galactosidase Standard in the concentration range of 0 U/L – 160 U/L (see table below).

1. Label 8 microtubes #1 through 8 with the following concentrations: 160, 80, 40, 20, 10, 5, 2.5, 0 U/L.
2. Pipette 196  $\mu$ L Assay Buffer into labeled tube #1, and 100  $\mu$ L Assay Buffer into labeled tube #2 through #8.
3. Briefly spin down the  $\beta$ -Galactosidase standard tube provided, and pipette 2  $\mu$ L of the  $\beta$ -Galactosidase standard into 285.3  $\mu$ L Assay Buffer to make 8000 U/L stock. Mix well.
4. Pipette 4  $\mu$ L of 8000 U/L stock into tube#1, mix thoroughly, and quick spin to make a 160 U/L  $\beta$ -Galactosidase Standard.
5. To make the 80 U/L standard, pipette 100  $\mu$ L of tube #1 into the tube labeled #2. Mix thoroughly and quick spin.
6. Repeat this step with each successive concentration, preparing a dilution series as shown in the table below. Each time, use 100  $\mu$ L of the prior concentration until the 2.5 U/L is reached. Mix each tube thoroughly before the next transfer. Standards should be prepared fresh and used immediately.

Labeled Tubes	$\beta$ -Galactosidase Standard ( $\mu$ L)	Assay Buffer ( $\mu$ L)	Standard Activity (U/L)
1	4 $\mu$ L of 8000 U/L stock	196 $\mu$ L	160
2	100 $\mu$ L of Tube #1	100 $\mu$ L	80
3	100 $\mu$ L of Tube #2	100 $\mu$ L	40
4	100 $\mu$ L of Tube #3	100 $\mu$ L	20
5	100 $\mu$ L of Tube #4	100 $\mu$ L	10
6	100 $\mu$ L of Tube #5	100 $\mu$ L	5
7	100 $\mu$ L of Tube #6	100 $\mu$ L	2.5
8	0 $\mu$ L	100 $\mu$ L	0

## Assay Procedure

Each  $\beta$ -Galactosidase Standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

1. Add 20  $\mu$ L of the diluted  $\beta$ -Galactosidase Standards or samples to the 96-well microtiter plate.
2. Initiate the reaction by adding 70  $\mu$ L Working Solution (See "Reagent Preparation") to each well. Carefully shake the plate for a few seconds to mix.
3. Incubate the plate for 30 minutes at 37 °C in the dark with mixing.
4. Add 120  $\mu$ L of Stop Solution to each well.
5. Measure the absorbance at 420nm using a plate reader.

## Calculation of Results

### Subtract the blanks

Average the absorbance value of the blank wells (Standard 0 U/L) and subtract this from the absorbance values of all the other wells. These are the corrected absorbances.

### Plotting the standard curves

Make a plot of corrected absorbance at 420nm as a function of  $\beta$ -Galactosidase activity.

### Determination of sample $\beta$ -Galactosidase activity

$$\beta\text{-Galactosidase (U/L)} = \frac{OD_{\text{Sample}} - OD_{\text{Blank}}}{\text{slope}} \times DF$$

OD<sub>Sample</sub> = Optical density (OD) reading of the Sample

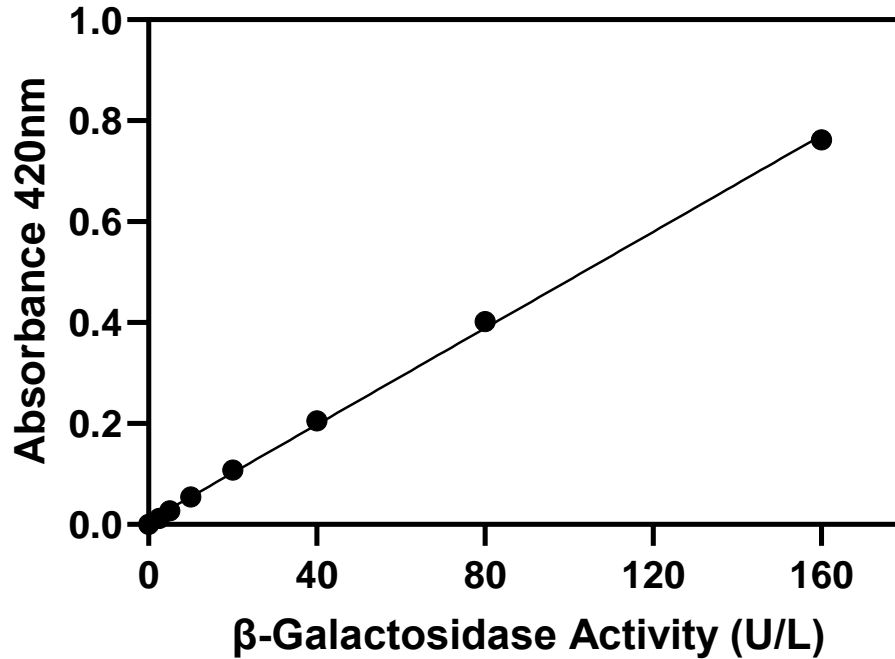
OD<sub>Blank</sub> = Optical density (OD) reading of the Blank (Standard 0 U/L)

Slope is from the plot of  $\beta$ -Galactosidase activity vs. Absorbance (as shown in Typical data below).

DF = Sample Dilution factor (DF = 1 for undiluted Samples)

*Note:* If the calculated  $\beta$ -Galactosidase activity of the sample is higher than 160 U/L, dilute the sample in assay buffer and repeat the assay.

## A. Typical Data



This standard curve is for demonstration only. A standard curve must be run with each assay.

## B. Reproducibility

Intra-assay Precision (Precision within an assay):

To assess intra-assay precision, 16 wells per sample (total of 4 samples) were tested on a single plate. The intra-assay coefficient of variation was found to be 2.3%.

Inter-assay Precision (Precision between assays):

To assess inter-assay precision, 4 samples were tested in separate assays (n=4). The inter-assay coefficient of variation was found to be 4.7%.

This product is for research use only.