

## Product Information

### **$\beta$ -GALACTOSIDASE (GAL) FLUORESCENT ACTIVITY DETECTION KIT**

Product No. **GAL-F**

Technical Bulletin MB-550  
November 1998

### TECHNICAL BULLETIN

#### **Product Description**

Reporter genes are "markers" widely used for analysis of mutationally altered genes as well as gene regulation and localization. The expressed reporter genes are detected by biochemical activity assay, immunological assay or by histochemical staining of tissue sections or cells<sup>1</sup>.

The *E. coli*  $\beta$ -Gal ( $\beta$ -galactosidase) gene is extensively used as a gene fusion marker for analysis of gene expression in eukaryotes. The *E. coli*  $\beta$ -galactosidase enzyme hydrolyzes D-galactose from various  $\beta$ -galactosides. The compound, 4-methylumbelliferyl-galactoside (4-MU-Gal) is a  $\beta$ -Gal substrate which upon hydrolysis produces the fluorescent 4-methylumbelliferone (4-MU). It is used in this fluorescent activity detection test designed for transfected cells or tissues expressing the *E. coli*  $\beta$ -Gal enzyme<sup>2</sup>.

The  $\beta$ -galactosidase hydrolyzes the 4-MU-Gal to galactose and the fluorescent compound 4-MU. The reaction is stopped with sodium carbonate solution; raising the pH also serves to enhance the fluorescence of the 4-MU. The 4-MU can be excited at 365 nm; its emission maximum is at 455 nm. *E. coli*  $\beta$ -galactosidase serves as a positive control enzyme.

This kit includes all buffers, substrate and reagents required for highly sensitive, easy to perform quantitative activity assays.

#### **Unit Definition**

One unit of  $\beta$ -galactosidase will release one pmole of 4-MU from 4-MU-Gal per minute at pH 7.0 and 37 °C. Note: This unit definition is different from Sigma's usual units for  $\beta$ -galactosidase.

#### **Reagents Provided**

This kit is sufficient for 200 tests.

- 5X Lysis Buffer, Product No. L5784 25 ml  
250 mM HEPES, pH 7.4, containing  
25 mM CHAPS
- 2X Assay Buffer, Product No. F4672 10 ml  
200 mM sodium phosphate buffer, pH 7.3,  
containing 2 mM MgCl<sub>2</sub> and 100 mM  
 $\beta$ -mercaptoethanol
- 4-Methylumbelliferyl  $\beta$ -D-Galacto-  
pyranoside, (4-MU-Gal substrate),  
Product No. M1633 10 mg
- 4-Methylumbelliferone Standard,  
(4-MU), Product No. M1508 25 mg
- $\beta$ -Galactosidase Positive Control 100  $\mu$ l  
From *E. coli*, Product No. G4538
- 5X Stop Solution, Product No. S5930 100 ml  
1 M sodium carbonate

#### Reagents and Equipment Required but Not Provided (Sigma product numbers are given where appropriate)

- Temperature controlled 37 °C water bath
- 10 mm x 75 mm glass test tubes
- Microcentrifuge tubes
- Fluorimeter and cuvetts
- Dimethyl sulfoxide (DMSO), Product No. D8418
- Reagents for determination of protein concentration

### Precautions and Disclaimer

Sigma's  $\beta$ -Galactosidase Fluorescent Activity Detection Kit is for laboratory use only. Not for drug, household or other uses.

### Storage

Store kit at  $-20\text{ }^{\circ}\text{C}$ .

### Preparation Instructions

Preparation of working solutions

1. Lysis buffer:  
Mix thawed 5X lysis buffer thoroughly before diluting 1:5 with deionized water.
2. Substrate solution:  
Dissolve 4-MU-Gal at a concentration of 11.4 mg/ml (34 mM) in DMSO. Store solution at  $-20\text{ }^{\circ}\text{C}$  up to one year.
3. 2X Assay buffer with substrate:  
Mix thawed 2X assay buffer thoroughly. To 0.5 ml 2X assay buffer add 25  $\mu\text{l}$  of substrate solution (34 mM 4-MU-Gal solution). Use freshly prepared solutions.
4. Calibration standard:  
Dissolve 2.0 mg 4-MU in 10 ml deionized water to make a 1 mM stock solution. Dilute 100  $\mu\text{l}$  of the stock solution to a volume of 10 ml with deionized water (1:100) to make a 10  $\mu\text{M}$  intermediate stock solution. Dilute 50  $\mu\text{l}$  of the intermediate stock solution to a volume of 5 ml with 1X stop solution to make the 100 nM working solution. Use freshly prepared solutions.
5. Positive control enzyme solution:  
Dilute 10  $\mu\text{l}$   $\beta$ -galactosidase positive control to a volume of 1 ml with 1X lysis buffer. Dilute a 10  $\mu\text{l}$  aliquot of the first dilution to a volume of 1 ml with 1X lysis buffer. Dilute a 10  $\mu\text{l}$  aliquot of the second dilution to a volume of 1 ml with 1X lysis buffer to give a working control solution (1:10<sup>6</sup> dilution). Use freshly prepared solutions.
6. Stop solution:  
Mix thawed 5X stop solution thoroughly before diluting 1:5 with deionized water.

Fluorimeter settings:

Set fluorimeter at 365 nm excitation, 455 nm emission, 5 nm slit width. Zero the instrument using 2 ml stop solution.

### Procedure

The reaction is performed in lysis/assay buffer containing 0.85 mM 4-MU-Gal at 37  $^{\circ}\text{C}$ . The reaction is initiated by the addition of the 2X assay buffer with substrate, and is stopped by the addition of cold 1X stop solution.

#### A. Calibration Procedure

Since the activity of  $\beta$ -galactosidase in the extract is unknown, perform a preliminary assay to find the proper dilution factor so that the measured values fall within the fluorimeter's detection range.

1. Add 20  $\mu\text{l}$  of cell extract (sample volume) to a microcentrifuge tube.
2. Add 80  $\mu\text{l}$  of lysis buffer.
3. Pre-incubate 1-2 minutes at 37  $^{\circ}\text{C}$ .
4. Initiate reaction with the addition of 100  $\mu\text{l}$  of 2X assay buffer with substrate.
5. Cap tube and incubate at 37  $^{\circ}\text{C}$  for 1 hour.
6. Transfer a 20  $\mu\text{l}$  aliquot to 2 ml of cold 1X stop solution and place in an ice bath.
7. Measure the fluorescence intensity (FI) of the sample.

If the values obtained are higher than the upper limit of the instrument, dilute the sample in stop solution until measurements are on scale. Calculate the dilution factor. If values obtained are below the detection range of the instrument, repeat the reaction with a larger volume of cell extract.

### B. Multiple Reaction Kinetic Assay Procedure

Complete the Calibration Procedure in order to optimize the sample dilution or the sample volume. The procedure described below is an example for cell extract that has to be diluted by a factor determined from the calibration procedure.

REAGENT	SAMPLE ASSAY	POSITIVE ENZYME CONTROL ASSAY	NEGATIVE CONTROL ASSAY
Sample extract	20 $\mu$ l	---	---
Diluted $\beta$ -galactosidase positive control	---	20 $\mu$ l	---
Lysis buffer	80 $\mu$ l	80 $\mu$ l	100 $\mu$ l
2X assay buffer with substrate	100 $\mu$ l	100 $\mu$ l	100 $\mu$ l

1. Prepare one vial containing 2.0 ml stop solution for each time point and place in an ice bath.
2. Dilute the sample cell extract by the dilution factor determined in the calibration procedure.
3. Pipet 20  $\mu$ l of the diluted sample extract to sample assay tube and 20  $\mu$ l of diluted  $\beta$ -galactosidase positive control to positive control assay tube.
4. Add the amount of lysis buffer indicated in the above table to each assay tube. Note: In the event that a volume of cell lysate other than 20  $\mu$ l is used, adjust the amount of lysis buffer added to give a total reaction volume of 100  $\mu$ l.
5. Mix and equilibrate for 1-2 minutes at 37  $^{\circ}$ C.
6. Add 100  $\mu$ l 2X assay buffer with substrate to each assay tube to initiate the reaction.
7. Immediately remove 20  $\mu$ l aliquots (volume per test) from all three reaction mixes into 2 ml stop solution as the zero minute test points. Keep the tubes on ice.
8. Cap tubes and incubate at 37  $^{\circ}$ C.
9. Transfer additional 20  $\mu$ l aliquots at 10 minute intervals to vials containing 2 ml cold stop solution.
10. Prepare 4-MU standard dilutions of 10, 20, 40, 60, 80 and 100 nM in stop solution. Include a minimum of 5 data points in order to generate a standard curve. 10 nM equals 20 pmol in a 2 ml cuvet.
11. Measure fluorescence intensity (FI) of all samples, positive enzyme and negative controls and 4-MU standard dilutions.

### C. Calculations

1. Draw a calibration curve of 4-MU standards FI versus pmol 4-MU.
2. Calculate FI per pmol 4-MU.
3. Plot sample, negative control and positive enzyme control FI versus time.
4. Calculate FI per minute for sample, negative control and positive enzyme control.
5. Subtract the negative control FI per minute value from the sample and positive enzyme control values.
6. Determine the protein concentration of the sample extract.
7. Calculate  $\beta$ -galactosidase activity of extract in pmol 4-MU per minute per  $\mu\text{g}$  protein (units per  $\mu\text{g}$  protein), according to the following equations.
8. Positive control enzyme activity is used for confirmation that the enzymatic system is working properly.

### Sample calculations

$$\text{Activity of extract (pmol MU/min/mg protein)} = \frac{\text{FI/min}}{\text{FI/pmole MU}} \times \frac{\text{Dilution factor}}{\text{Extract conc., (mg protein/ml)}} \times 500/\text{ml}^*$$

$$500/\text{ml}^* = \frac{\text{Reaction volume (0.2 ml)}}{\text{Sample volume (0.02 ml) x Vol. per test (0.02 ml)}}$$

500/ml factor is calculated according to the following parameters: Reaction volume = 0.2 ml, Sample volume = 0.02 ml (20  $\mu\text{l}$ ), and Volume per test = 0.02 ml (20  $\mu\text{l}$ ). In the event that the sample volume is different than 20  $\mu\text{l}$  (0.02 ml), adjust the factor accordingly.

### References

1. Kain, S.R. & Ganguly, S., "Uses of Fusion Genes in Mammalian Transfection", in Current Protocols in Molecular Biology, Vol. 1, Suppl. 36, Ausubel, F.M. *et al.*, eds. (John Wiley & Sons, New York, 1996) p. 9.6.1\_
2. Young, D.C., *et al.*, *Anal. Biochem.*, **215**, 24-30 (1993)

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.

Please see reverse side of the invoice or packing slip.