

## General Information

The gene of β-galactosidase from *E. coli* is widely used as a reporter gene assay marker. Although X-gal is well known reagent to detect β-galactosidase in cell or tissue samples, the assay using these reagents require to fix cells or tissues due to the poor cell-permeability. In addition, so far developed the assay using fluorescence reagents can not clearly differentiate β-galactosidase-expressed cells or regions.

To overcome these issues, Urano, Kamiya and co-workers have successfully developed SPiDER-βGal. SPiDER-βGal ideally possesses cell-permeability and the ability to retain in intracellular region.<sup>1)</sup>

By the enzymatic reaction, SPiDER-βGal immediately forms a quinone methide that acts as electrophile when proteins containing nucleophilic functional groups nearby the molecules. By the probe undergoes the reaction with a protein, the conjugates become fluorescent compounds. Thus, SPiDER-βGal allows a single-cell analysis because it does self-immobilizing to the intracellular proteins.

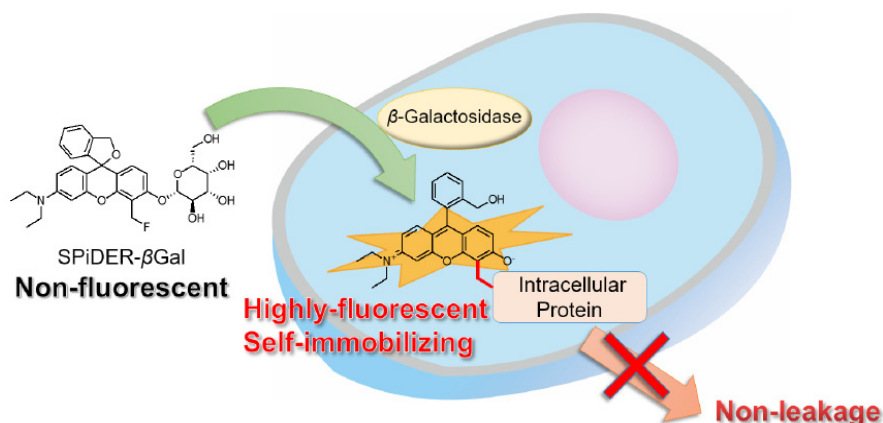


Fig. 1 Cell staining mechanism by SPiDER-βGal.

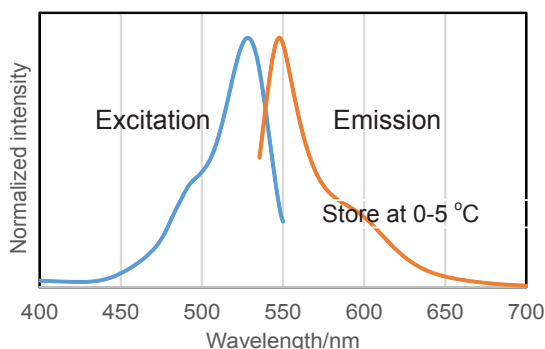


Fig. 2 Excitation and emission spectra of SPiDER-βGal after reaction with β-galactosidase

< Recommended filter >

Ex : 500 ~ 540 nm

Em : 530 ~ 570 nm

There are internal detection examples obtained using confocal microscopy and flow cytometry at an excitation wavelength of 488 nm.

## Contents

### Storage Condition

- SPiDER-βGal 20 μg x 3

Store at 0-5 °C

After a SPiDER-βGal is taken out from the seal bag, keep the unused SPiDER-βGal in the bag, seal tightly and store at 0-5°C.

### Required Equipment and Materials

- Dimethyl sulfoxide (DMSO)
- Hanks' HEPES buffer
- Micropipettes
- Microtubes

### Preparation of solutions

#### Preparation of 1 mmol/l SPiDER-βGal DMSO stock solution

Add 35 μl of DMSO to a tube of SPiDER-βGal (20 μg) and dissolve it with pipetting.

\*Store the SPiDER-βGal stock solution at -20°C.

#### Preparation of 1 μmol/l SPiDER-βGal working solution

Dilute the SPiDER-βGal DMSO stock solution with Hanks' HEPES buffer to prepare 1 μmol/l SPiDER-βGal working solution.

\*Hanks' HEPES buffer is recommended to maintain cell condition.

### General protocol

#### SPiDER-βGal staining

1. Prepare cells for the assay.
2. Discard the culture medium and wash the cells with Hanks' HEPES buffer twice.
3. Add an appropriate volume of SPiDER-βGal working solution.
4. Incubate at 37°C for 15 minutes.
5. Observe the cells under a fluorescence microscope or by a flow cytometer.

\*After staining, the cells can be observed even without washing. However, you can wash it as needed.

Fluorescence microscopic detection of  $\beta$ -galactosidase-expressed cells

1. HEK cells at  $5 \times 10^5$  cells/ml (500  $\mu$ l) and HEK/LacZ cells at  $5 \times 10^5$  cells/ml (500  $\mu$ l) were seeded in a 35 mm dish in DMEM (10% fetal bovine serum, 1% penicillin-streptomycin) and cultured overnight in a 5% CO<sub>2</sub> incubator at 37°C.
2. The cells were washed with 2 ml of Hanks' HEPES buffer twice.
3. SPiDER- $\beta$ Gal working solution (2 ml) was added to the culture dish, and the cells were incubated for 15 minutes at 37°C.
4. After the supernatant was removed, the cells were washed Hanks' HEPES buffer (2 ml) twice.
5. Hanks' HEPES buffer (2 ml) were added, and the cells were observed under a fluorescence microscope. (Fig. 3A)
6. After the supernatant was removed, 4% paraformaldehyde (PFA) /PBS solution (2 ml) was added to the culture dish, and the cells were incubated for 15 minutes at room temperature.
7. After 4% PFA/PBS solution was removed, the cells were washed Hanks' HEPES buffer (2 ml) twice.
8. Hanks' HEPES buffer (2 ml) were added, and the cells were observed under a fluorescence microscope. (Fig. 3B)

\*Filter (wavelength/band pass)

Fluorescence imaging: 550/25 nm (Ex), 605/70 nm (Em)

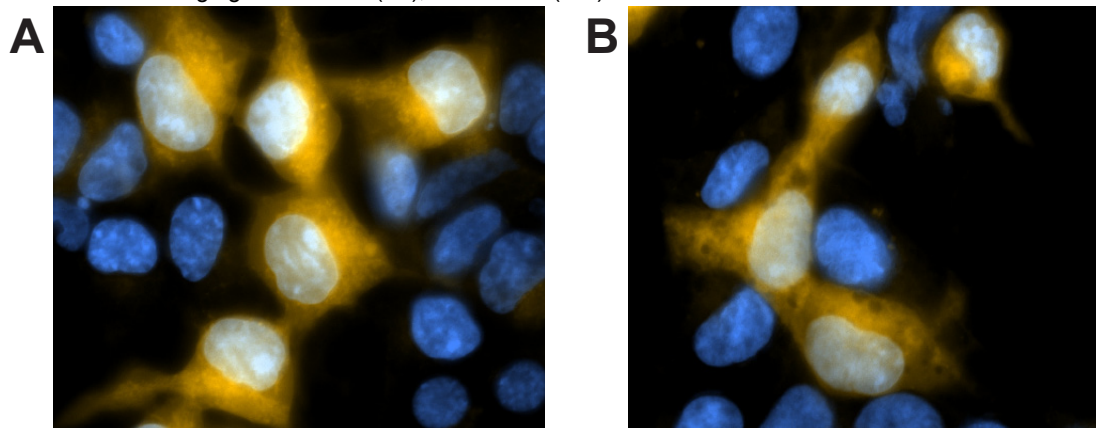


Fig. 3 Fluorescence imaging of HEK/LacZ cells and HEK cells at 1: 1 ratio.

A. living cell, B. fixing cells (4% PFA/PBS)  
(yellow: SPiDER- $\beta$ Gal, blue: Hoechst 33342)

$\beta$ -galactosidase-expressed cells (HEK/LacZ cells) were clearly observed in fluorescence imaging. In addition, the result was not changed by fixing the cells.

Flow cytometric detection of  $\beta$ -galactosidase-expressed cells

1. HEK cells at  $5 \times 10^5$  cells/ml (500  $\mu$ l) and HEK/LacZ cells at  $5 \times 10^5$  cells/ml (500  $\mu$ l) were mixed in a microtube.
2. SPiDER- $\beta$ Gal DMSO stock solution (1  $\mu$ l) was added to the tube, and the cells were incubated for 15 minutes at 37°C.
3. The cells were analyzed by a flow cytometer. (488 nm excitation, 530/30 nm bandpass filter)

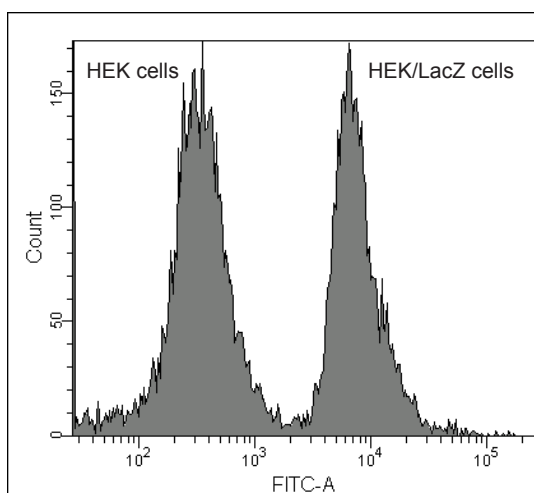


Fig. 4 Analysis of HEK/LacZ cells and HEK cells at 1: 1 ratio by flow cytometry.

$\beta$ -galactosidase-expressed cells (HEK/LacZ cells) were clearly differentiate from HEK cells in flow cytometry data analysis.

- Reference 1) T. Doura, M. Kamiya, F. Obata, Y. Yamaguchi, T. Y. Hiyama, T. Matsuda, A. Fukamizu, M. Noda, M. Miura and Y. Urano, *Angew. Chem. Int. Ed.*, **2016**, 55, 9620.

For research use only. Not for use in diagnostic procedures.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories

2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto  
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525  
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.

Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>

Dojindo EU GmbH

Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>

Dojindo China Co., Ltd

Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>