

Luminescent β -galactosidase Reporter System 3 Protocol-at-a-Glance

(PT2106-2)

Please read the *User Manual* before using this abbreviated protocol. The Protocol-at-a-Glance is provided for your convenience, but is not intended for first-time users.

A. Preparation of cell lysates for assay of β -galactosidase

Maximal levels of β -galactosidase are generally detected 48–72 hr after transfection. The volumes listed below are for adherent cells cultured on 60-mm tissue culture plates. If working with suspended cell cultures, simply collect an equivalent cell mass by centrifugation, wash three times by resuspending in 500 μ l of ice-cold PBS followed by centrifugation, and proceed from Step 6.

1. Wash cells on the plate twice with 4.0 ml of ice-cold PBS.
2. Add 1.0 ml of ice-cold PBS to each plate.
3. Scrape cells off the plate using a rubber policeman (or equivalent), and transfer cell suspension to a microcentrifuge tube on ice.
4. If residual cells are still on the plate, use another 500 μ l of PBS to collect the remaining cells and transfer to the same microcentrifuge tube.
5. Centrifuge at maximum speed for 15 sec in a microcentrifuge.
6. Aspirate supernatant, being careful not to disturb the cell pellet.
7. Gently resuspend cells in 1.0 ml of ice-cold PBS.
8. Repeat Steps 5 and 6, and carefully aspirate the last traces of PBS.
Note: The cell pellet can be stored at -20°C for 1–3 days with minimal loss of β -galactosidase activity.
9. Gently resuspend the cell pellet in 75 μ l of ice-cold lysate buffer.
10. Place the tube with cell suspension in a dry-ice/ethanol bath for 1 min (or until completely frozen).
11. Thaw the cell suspension at 37°C for 1–2 min.
12. Repeat the freeze/thaw cycle (Steps 10–11) two more times.
13. Centrifuge at 4°C for 5 min at maximum speed.
14. Transfer the supernatant to a fresh tube and keep on ice.

B. Chemiluminescent β -galactosidase assay using a tube luminometer*

1. Warm enough Reaction Buffer and Reaction Substrate to room temperature.
2. Prepare the master Reaction Buffer Mixture by adding 4 μ l of Reaction Substrate to 196 μ l of Reaction Buffer for each sample.
3. Aliquot 30–50 μ l of individual cell lysates into sample tubes.
4. Add 200 μ l of Reaction Buffer Mixture to each cell lysate and mix gently.
5. Incubate at room temperature (20 – 25°C) for 60 min.
6. Record light emission as a 5-sec integral using a tube luminometer.

Note: Since the light signal is stable for >1 hr, the emission can be recorded 0–60 min after the incubation.

*Consult the main protocol for information on detection using a plate luminometer or x-ray film.

Notice to Purchaser

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, in vitro diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without written approval of Clontech Laboratories, Inc.

Clontech, the Clontech logo and all other trademarks are the property of Clontech Laboratories, Inc. Clontech is a Takara Bio Company. ©2006.

(PR98371; published 10 September 1999)



Clontech

United States/Canada
800.662.2566

Asia Pacific
+1.650.919.7300

Europe
+33.(0)1.3904.6880

Japan
+81.(0)77.543.6116

Clontech Laboratories, Inc.
A Takara Bio Company
1290 Terra Bella Ave.
Mountain View, CA 94043
Technical Support (US)
E-mail: tech@clontech.com
www.clontech.com