

Applied Biosystems
47 Wiggins Avenue
Bedford, MA 01730
Phone: (800) 542-2369
or (781) 271-0045
FAX: (781) 275-8581
email: tropix@appliedbiosystems.com

Luciferase Assay Kit

Bioluminescent Reporter Gene Assay System for the Detection of Firefly Luciferase

Cat. Nos. BC100L, BC300L, BC2500L

Contents	Page
I. INTRODUCTION	1
II. SYSTEM COMPONENTS	1
III. LUCIFERASE DETECTION PROTOCOL	
A. PREPARATION OF CELL EXTRACTS	2
B. DIRECT LYSIS PROCEDURE FOR MICROPLATE CULTURES	2
C. BIOLUMINESCENT DETECTION PROCEDURE	3
D. PROTOCOL NOTES	3
APPENDICES	
A. PREPARATION OF CONTROLS	4
B. USE OF LUMINOMETERS	4
C. USE OF SCINTILLATION COUNTERS	4
REFERENCES	5

Applied Biosystems and Tropix are registered trademarks, and Applera, NorthStar and TR717 are trademarks of Applera Corporation or its subsidiaries in the US and certain other countries. For Research Use Only. Not for use in diagnostic procedures. Information is subject to change without notice.

I. INTRODUCTION

Reporter gene assays are widely used for studying gene regulation and function in cell biology (1,2). Luciferase is a sensitive genetic reporter due to the lack of endogenous activity in cells and tissues. Luciferase assays provide a 1000-fold increase in detection sensitivity compared to the standard assays for chloramphenicol acetyltransferase (CAT) reporter enzyme (3,4). Detection of 1 fg to 20 ng of purified luciferase is possible.

Unlike one-component luciferase assay systems, which exhibit decreased performance after long term storage, the Tropix[®] Luciferase Assay has been formulated as a two-component system in order to provide consistent results over the life of the kit. Cell lysate is mixed with Substrate A, containing ATP and inorganic salts. Light signal intensity proportional to the amount of enzyme present in an extract is measured immediately after addition of Substrate B, containing luciferin. The enhanced luciferase/luciferin reaction produces a light signal which decays with a half-life of approximately 5 min. The kinetics of light emission enable the use of luminometers with or without automatic injectors and other instrumentation in which light emission measurements can be performed, including scintillation counters (see Appendix C).

II. SYSTEM COMPONENTS

Shelf-life of all Luciferase Assay Kit components is 1 year when stored as recommended below.

	BC100L	BC300L	BC2500L
microplate assays	200	600	5,000
Lysis Solution	70 mL	210 mL	1.75 L
Substrate A	20 mL	3 x 20 mL	5 x 100 mL
Substrate B	20 mL	3 x 20 mL	5 x 100 mL

- 1. Lysis Solution:** 100 mM potassium phosphate (pH 7.8), 0.2% Triton X-100. Store at 4°C.
- 2. Substrate A:** Lyophilized ATP solution. Reconstitute in 20 mL of H₂O. Store at -20°C before reconstitution; after, store at 4°C for 1 week or store aliquots at -20°C for up to 1 year. Do not re-freeze aliquots.
- 3. Substrate B:** Lyophilized luciferin solution. Reconstitute in 20 mL of H₂O. Store at -20°C before reconstitution; after, store at 4°C for 1 week or store aliquots at -20°C for up to 1 year. Do not re-freeze aliquots.

III. LUCIFERASE DETECTION PROTOCOL

Please read the entire Protocol and Notes sections before proceeding. Perform all assays in triplicate at room temperature.

A. Preparation of Cell Extracts

See Appendix A for preparation of controls.

1. Add DTT (to 1 mM) to the required volume of Lysis Solution (see Note 1).
2. Rinse cell cultures twice with PBS.
3. Add Lysis Solution to cover the cells. Use 250 μ L per 60 mm plate.
4. Detach cells from plate with a cell scraper.
5. Transfer the cell lysate to a microfuge tube and centrifuge for 2 min to pellet debris.
6. Transfer extracts (supernatant) to a fresh tube. Use immediately or store at -70°C .

B. Direct Lysis Procedure for Microplate Cultures

This procedure is for adherent cells growing in 96-well tissue culture-treated luminometer plates.

1. Add DTT (to 1 mM) to the required volume of Lysis Solution (see Note 1).
2. Rinse cell cultures once with PBS.
3. Add 10 μ L of Lysis Solution to each well and incubate for 10 min.
4. Continue with the Bioluminescent Detection Procedure (Section C), omitting Step 2.

C. Bioluminescent Detection Procedure

Perform all assays in triplicate at room temperature.

1. Equilibrate Substrate A and B to room temperature.
2. Transfer 5-100 μL of extract to luminometer tubes or microplate wells (see Note 2).
3. Add 100 μL of Substrate A.
4. Within 10 min, inject 100 μL of Substrate B. After a 1-2 sec delay, read the signal for 0.1-1 sec (see Note 3).

D. Protocol Notes

1. The Lysis Solution provided may be substituted with alternatives. These should be compared, however, with Lysis Solution in order to ensure optimum assay performance.
2. The amount of extract required may vary depending on the luciferase expression level and the instrumentation used; the amount used should be adjusted to keep the signal within the linear range of the assay. For experiment-to-experiment consistency, the same volume of sample should be assayed every time. Lysis Solution may be added to equalize sample volumes.
3. The time between adding Substrate B and start of measurement should be as short as possible, and identical from sample to sample.

APPENDICES

A. Preparation of Controls

Positive Control

Reconstitute lyophilized luciferase (Sigma Cat. No. L-1759) to 1 mg/mL in 0.1 M sodium phosphate (pH 7.0), 0.2% BSA. The stock enzyme should be prepared fresh each time or aliquots can be stored at -70°C and used without repeated freeze/thaw. Generate a standard curve by serially diluting the stock enzyme in 10 µL of Lysis Solution. For the high-end detection limit, use 10-100 ng of enzyme.

Negative Control

Assay a volume of mock-transfected extract equivalent to that of experimental extract.

B. Use of Luminometers

We recommend using a dedicated luminometer (such as the Tropix NorthStar™ HTS workstation or TR717™ microplate luminometer) to measure the light emission from 96- or 384-well microplates. For most samples, the luminometer can be set to measure for 1-2 min/plate or 0.1-1 sec/well. The linear range of detection will vary according to cell type and on the reporter gene expression level. The number of cells or sample volume used per well should be optimized to prevent a measurement signal that is outside the linear range of the luminometer. Extremely high light signals can saturate the detector, resulting in erroneous measurements. Refer to your luminometer user's manual to determine the upper limit for your specific luminometer. Contact Tropix Technical Support for additional questions or for more information on the NorthStar™ HTS workstation or TR717™ microplate luminometer.

C. Use of Scintillation Counters

A liquid scintillation counter may be used as a substitute for a luminometer, however, sensitivity may be lower (5,6). When using a scintillation counter, it is necessary to turn off the coincident circuit in order to measure bioluminescence directly (single photon counting mode). The manufacturer of the instrument should be contacted to determine how this is done. If it is not possible to turn off the coincident circuit, a linear relationship can be established by taking the square root of the counts per minute measured minus the instrument background.

$$\text{Actual} = (\text{measured} - \text{background})^{1/2}$$

REFERENCES

1. Alam, J. and J.L. Cook. 1990. Reporter genes: Application to the study of mammalian gene transcription. *Anal. Biochem.* 188:245-254.
2. Bronstein, I., et al. 1994. Chemiluminescent and bioluminescent reporter gene assays. *Anal. Biochem.* 219:169-181.
3. Gould, S.J. and S. Subramani. 1988. Firefly luciferase as a tool in molecular and cell biology. *Anal. Biochem.* 175:5-13.
4. Brasier, A.R., et al. 1989. Optimized use of the firefly luciferase assay as a reporter gene in mammalian cell lines. *BioTechniques.* 7:1116-1122.
5. Fulton, R., and B. Van Ness. 1993. Luminescent reporter gene assays for luciferase and b-galactosidase using a liquid scintillation counter. *BioTechniques* 14:762-763.
6. Nguyen, V.T., et al. 1988. Firefly luciferase luminescence assays using scintillation counters for quantitation in transfected mammalian cells. *Anal. Biochem.* 171:404-408.

WARRANTY

Tropix, a wholly owned subsidiary of Applied Biosystems Corporation (formerly PE Corporation) ("Tropix"), warrants its products, to only the original purchaser and to no third party, against defects in materials and workmanship under normal use and application. Tropix' sole obligation and total liability under this warranty shall be to replace defective products. All products are supplied For Research Use Only as defined herein and are Not For Resale. Commercialization of products using these components requires an express license under applicable patents and intellectual property from Tropix that is not included herein. Our preparations are intended exclusively for in vitro use only. They are not for diagnostic or therapeutic use in humans or animals. Without limiting or otherwise affecting the limitations on warranty and liability stated herein: (1) those preparations with known toxicity are sent with an information sheet which describes, to our knowledge, the potential dangers in handling, (2) the absence of a toxicity warning with one of our products does not preclude a possible health hazard, (3) with all of our products, due care should be exercised to prevent human contact and ingestion, and (4) preparations should be handled by trained personnel only. This warranty is in lieu of all other warranties, express or implied and, without limitation, Tropix EXPRESSLY DISCLAIMS THE WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE. IN NO CASE OR EVENT SHALL TROPIX BE LIABLE FOR INCIDENTAL OR CONSEQUENTIAL DAMAGES, OF WHATEVER NATURE, EVEN IF TROPIX HAS BEEN ADVISED OF THE POSSIBILITY THEREOF.