

Ultrasensitive Secreted Luciferases and their use in a Dual Reporter System

Introduction to *Gaussia* and *Cypridina* Luciferases

Luciferases are enzymes that produce light as a by-product of the oxidation of substrate. This light emission can be easily detected, making luciferases ideal reporter systems. In fact, they have proven to be very useful reporters for the study of fundamental cellular processes such as the regulation of transcription and translation, mRNA and protein stability, nucleic acid and protein interactions, as well as markers of cellular physiology and responses.

The first luciferase reporters used in life science research were intracellular luciferases. These non-secreted reporters require that cells be lysed in order to measure activity. Measurement of reporter activity at a number of different time points involve complex experimental design with large numbers of wells and reagents. Additionally, the fact that each time point is taken from different cell samples can contribute to experimental variability.

The availability of secreted luciferase proteins overcomes these difficulties. New England Biolabs has recently developed reporter systems based on *Gaussia* luciferase (GLuc) and *Cypridina* luciferase (CLuc) (1,2). These luciferases produce a large number of photons per reduction cycle, making them very sensitive reporters. When expressed in mammalian cells, GLuc and CLuc are naturally secreted into the growth media (Fig. 1) (3), allowing the user to sample luciferase activity from the same source multiple times over the course of an experiment. This method of non-destructive sample collection expedites the assay protocol, simplifies experimental design and allows the cells to remain intact so they can be used in further downstream assays such as RT-PCR, Western blots, RNA expression analysis, live imaging, cell viability assays, etc.

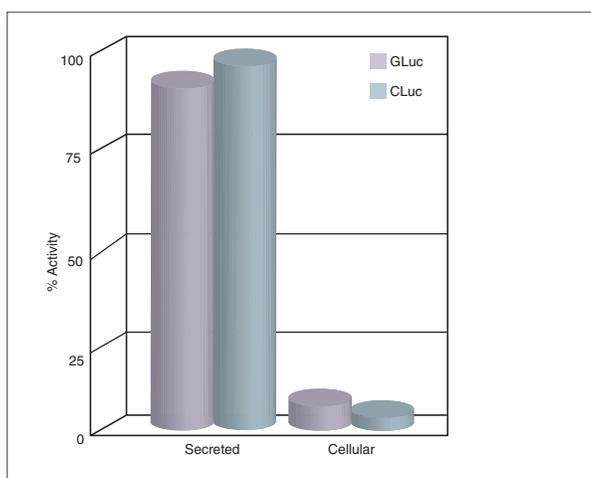


Figure 1: Both *Gaussia* and *Cypridina* luciferases are efficiently secreted from mammalian cells. GLuc and CLuc vectors were transfected into CHO-K1 cells. Supernatants were collected and cell lysates were made. Luciferase activity was measured from 20 μ l of supernatant or lysate from the same cells and the percentage of total luciferase activity in the supernatant (secreted) and in the lysate (cellular) was calculated. Activity of GLuc and CLuc was measured with its corresponding BioLux assay system.

DNA CLONING

DNA AMPLIFICATION & PCR

EPIGENETICS

RNA ANALYSIS

LIBRARY PREP FOR NEXT GEN SEQUENCING

PROTEIN EXPRESSION & ANALYSIS

CELLULAR ANALYSIS

General Protocol

Injector-equipped luminometers

1. Prepare the GLuc assay solution (e.g. 100 samples) by adding 50 μ l of BioLux[®] GLuc Substrate to 5 ml of Biolux GLuc Assay Buffer immediately before performing the assay. (Be sure to prepare enough assay solution as needed for all samples as well as for priming a particular luminometer as recommended by the manufacturer).
2. Mix well by inverting the tube several times (Do not vortex).
3. Set the luminometer with the following parameters: 50 μ l of injection, 5–15 seconds of delay & 2–10 seconds of integration.
4. Pipet samples (5–20 μ l per well) into a 96-well white (opaque) or black plate, or a luminometer tube.
5. Prime the injector with the GLuc assay solution and proceed with the measurement.

* Approximately 90% of GLuc is secreted out into the growth media after transfection and thus, the GLuc activity is typically assayed from the supernatant (i.e. growth media of GLuc-transfected cells). However, as long as the cells are alive, approximately 10% of GLuc is present inside the cells. Therefore, GLuc activity can also be assayed from the cell lysate. We recommend that the cell lysates be prepared by using Luciferase Cell Lysis Buffer (NEB #B3321), since this lysis buffer is designed to be compatible with *Cypridina*, *Gaussia*, *Renilla*, Firefly luciferase and β -gal activity assays.

(See other side)

Use of GLuc and CLuc in a Dual Reporter System

The co-transfection of an experimental reporter and a second control reporter can provide valuable information, including tracking of a reporter gene of interest, transfection efficiency and cell viability. The transfection efficiency of the secondary reporter serves as a normalizer and allows researchers to compare results within and across experiments. To take full advantage of the benefits of a secreted experimental reporter, it is best to also use a secreted control reporter, allowing the user to continuously monitor the secretion of both transfected reporters into the growth media. Additionally, it minimizes sampling error, as both reporters can be measured from a fraction of the same sample.

GLuc and CLuc utilize different substrates to carry out the light emitting oxidation reaction, and both exhibit a similar range of activity (4). Figure 2 shows that the detection of one luciferase can be monitored without any cross-reactivity from the presence of the other. Briefly, supernatants were collected from CLuc- and GLuc-expressing cells and mixed, maintaining the fraction of one supernatant constant while increasing the volume of the other supernatant. Activity of GLuc and CLuc in the mixture was measured with its corresponding BioLux assay system. These results demonstrate that GLuc and CLuc are ideal reporter partners to be used in co-expression experiments in mammalian cells.

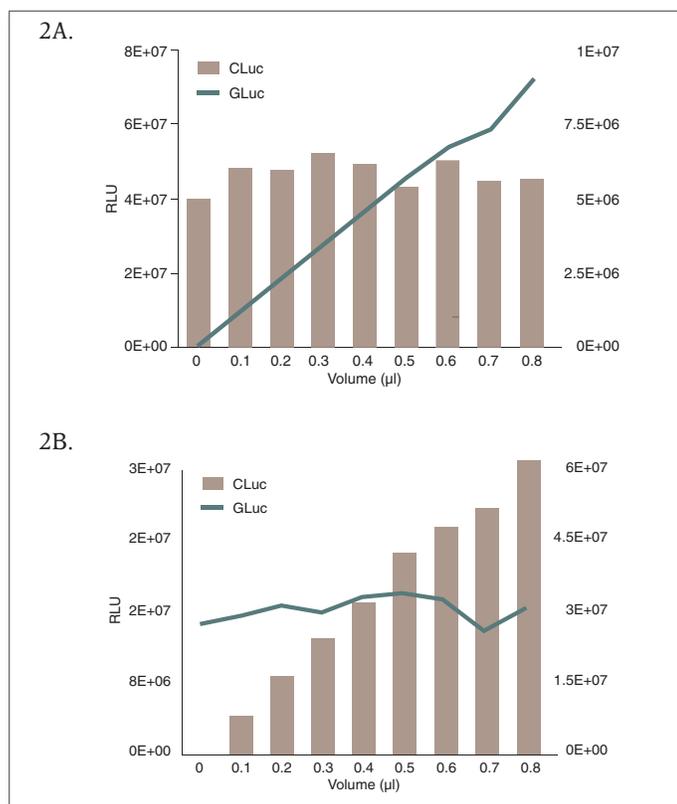


Figure 2: GLuc or CLuc activity can be measured without interference from the presence of the other luciferase. (A) supernatant from CLuc cells is held constant while volume of supernatant from GLuc cells is increased. (B) supernatant from GLuc cells is held constant while volume of supernatant from CLuc cells is increased. Activity is measured with the corresponding BioLux Assay System.

General Protocol

Luminometers without injectors

1. Prepare the GLuc assay solution (e.g. 100 samples) by adding 50 µl of BioLux GLuc Substrate to 5 ml of BioLux GLuc Assay Buffer immediately before performing the assay.
2. Mix well by inverting the tube several times (Do not vortex).
3. Set the luminometer for 2–10 seconds of integration.
4. Pipet samples (5–20 µl per well) into a 96-well white (opaque) or black plate, or a luminometer tube.
5. Add the GLuc assay solution (50 µl) to a sample (i.e. Add the assay solution to only one sample at a time) and promptly measure the luminescence.
6. Repeat Step 5 for all samples.

References

1. Verhaegen, M. and Christopoulos, T.K. (2002) *Anal. Chem.* 74, 4378–4385.
2. Nakajima, et al. (2004) *Biosci. Biotechnol. Biochem.* 68, 565–570.
3. Knappskog, S., et al. (2007) *J. Biotechnol.*, 128, 705–715.
4. Wu, C., Suzuki-Ogoh, C. and Ohmiya, Y. (2007) *Biotechniques*, 42, 290–292.

NEW ENGLAND BIOLABS® and BIOLUX® are registered trademarks of New England Biolabs, Inc.

