

Simultaneous Dual Emission detection of luciferase reporter assays

Megan Dobbs¹, Douglas Hughes¹, Janaki Narahari¹, Jae Choi¹, Georgyi Los¹, Brian Webb¹, Eric Matthews², and Carl Peters²
¹Thermo Fisher Scientific; ²BMG LABTECH

- Performing dual-luciferase assays using the microplate reader from BMG LABTECH
- Spectral resolution allows detection of two luciferase activities in one sample
- Simultaneous dual-emission detection permits one read assessment

Introduction

Luciferase-based reporter assays employ insertion of a genetic regulatory element upstream of the luciferase gene. The result is an excellent tool for monitoring gene expression in cells due to wide dynamic range and sensitive detection. The luminescence resulting from expression of the transfected luciferase reporter gene is measured to quantify the activity of the cis-acting or trans-acting components of the biological pathway. In dual luciferase assays the expression values of one luciferase is usually related to the studied subject whereas the other luciferase reporter is considered as an indicator for transfection efficiency and cell viability.

Assay Principle

The Thermo Scientific™ Pierce™ Luciferase Dual-Spectral Assays provide a simple one-step detection protocol based on the use of pairs of luciferase enzymes that are spectrally resolved. Therefore their activities can be measured with high sensitivity in the same samples (Figure 1) without the usual quenching step needed to differentiate luciferase signals. Three dual assay systems in which red firefly luciferase is paired with green *Renilla*, *Gaussia* or *Cypridina* luciferase are available.

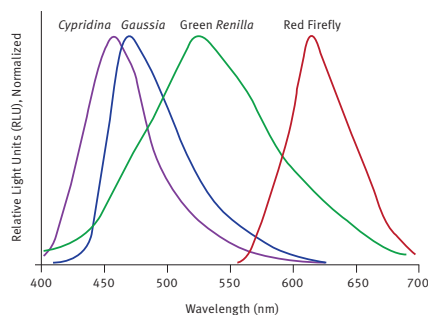


Fig. 1: Spectral emission profiles of luciferases used in Thermo Scientific™ Pierce™ Dual-Spectral Assay Kits. The emission of red firefly ($\lambda_{max} = 613\text{nm}$) allows resolution from Green *Renilla* Luc ($\lambda_{max} = 535\text{nm}$), *Gaussia* Luc ($\lambda_{max} = 470\text{nm}$) and *Cypridina* Luc ($\lambda_{max} = 463\text{nm}$).

Detecting activity from two spectrally resolved luciferases is a task for which the BMG LABTECH multidetection microplate reader is ideally suited. Light from both reporters can be measured either using the simultaneous dual-emission option which uses two photomultiplier tubes (PMT), or it can be measured

sequentially, one emission after the other. Here we demonstrate the sensitivity of the Pierce™ Luciferase Dual-Spectral Reporter Assays, over at least a 100,000-fold luciferase concentration range after normalization to a control reporter, typically red firefly luciferase.

Materials & Methods

- HEK293 stable cell lines expressing *Cypridina*, *Gaussia*, Green *Renilla* and Red Firefly luciferase under the control of CMV promoter
- Pierce™ Luciferase Cell Lysis Buffer
- Pierce™ *Cypridina*-Firefly Luciferase Dual-Spectral Assay Kit
- Pierce™ *Gaussia*-Firefly Luciferase Dual-Spectral Assay Kit
- Pierce™ *Renilla*-Firefly Luciferase Dual-Spectral Assay Kit
- White, 96-well f-bottom plates (Thermo Scientific)
- BMG LABTECH microplate reader

HEK293 cell lines stably expressing *Cypridina*, *Gaussia*, green *Renilla* and red firefly luciferases were lysed with Pierce™ Luciferase Cell Lysis Buffer. 1:10 dilutions were prepared in lysis buffer. Analysis of replicates of each serial dilution (10 $\mu\text{L}/\text{well}$) was performed. Cell lysate containing a different second reporter (10 $\mu\text{L}/\text{well}$) was added as a control. For serial dilutions of cell lysate containing *Cypridina*, *Gaussia*, or green *Renilla* luciferase, the second control reporter was red firefly. For serial dilutions of cell lysate containing red firefly, the second control reporter was *Cypridina* luciferase. The appropriate Pierce™ Luciferase Dual Assay Working Solution was prepared for each assay, and 50 $\mu\text{L}/\text{well}$ was injected into each well for a final volume of 70 $\mu\text{L}/\text{well}$. Luminescence was read through appropriate filters with a 1 second integration time. A more detailed protocol can be found online. An overview of the filter settings is given in table 1.

Table 1: Filter settings for each dual-spectral luciferase assay.

Assay	Luciferase	Filters	Gain
Cypridina-Firefly	<i>Cypridina</i>	475 +/- 30	2500
	Red Firefly	610 long pass	2000
Gaussia-Firefly	<i>Gaussia</i>	475 +/- 30	2500
	Red Firefly	610 long pass	2000
Renilla-Firefly	<i>Renilla</i>	515 +/- 30	2000
	Red Firefly	670 +/- 10	3000



Results & Discussion

Measuring the separate signals in each of the three pairs of luciferase reporter enzymes in the Pierce™ Luciferase Dual-Spectral Assay Kits was easy, using the appropriate filters (Figures 2, 3 and 4).

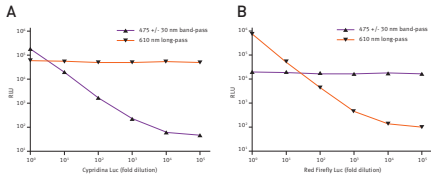


Fig. 2: Cypridina-Firefly Dual-Spectral Luciferase Assay on the POLARstar Omega using the instrument's dual-emission optics. **A.** Serial dilutions of HEK293 cell lysate expressing *Cypridina* Luc ($n = 6$). HEK293 cell lysate expressing Red Firefly Luc as a control. **B.** Serial dilutions of HEK293 cell lysate expressing red firefly luciferase ($n = 6$). HEK293 cell lysate from expressing *Cypridina* Luc as a mock "control".

By selecting filters for the luciferases in each pair that eliminate or greatly reduce interference between them, conditions for each dual-color assay system can be found that provide accurate measurement over 100,000-fold concentration range for each reporter (Figures 2 and 3).

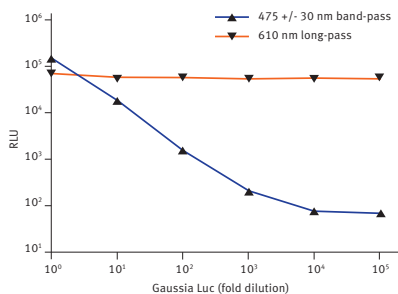


Fig. 3: Gaussia-Firefly Dual-Spectral Luciferase Assay. Serial dilutions of HEK293 cell lysate from a stable cell line expressing *Gaussia* Luc ($n = 6$). HEK293 cell lysate from a stable cell line expressing Red Firefly Luc was also added as a control.

Filter selection is critical for eliminating interference when spectral overlap is present between reporters (Figure 1). Thus for the *Renilla*-firefly assay, we used a suboptimal red-shifted 670 +/- 10nm bandpass filter for detection of red firefly, which minimizes interference from green *Renilla*. As a result, interference was only observed when

the output from green *Renilla* was significantly greater than that from red firefly (Figure 4). To separate spectral data in this region a corrective calculation is needed (see Pierce™ Dual-Spectral Calculator).

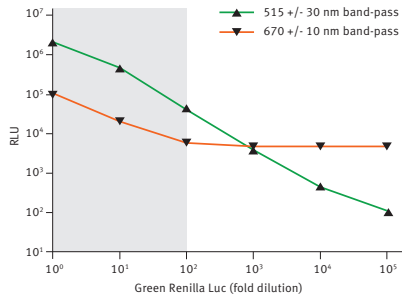


Fig. 4: Renilla-Firefly Dual-Spectral Assay. Serial dilutions of HEK293 cell lysate from a stable cell line expressing Green *Renilla* Luc ($n = 6$). HEK293 cell lysate from a stable cell line expressing red firefly luciferase was also added as a control. The shaded region indicates significant bleed-through of green *Renilla* luminescence into the red filter.

Conclusion

These experiments confirm the BMG LABTECH's microplate reader's ability to measure the activity of two luciferases that are spectrally resolved. The Thermo Scientific™ Pierce™ Luciferase Dual-Spectral Reporter Assays are sensitive over at least a 100,000-fold dilution range of each luciferase when measured using appropriate filters. Defining the limit of detection in a dual-luciferase assay is difficult and will depend upon luciferase concentration as will the dynamic range.



PHERAstar® FSX

*The PHERAstar FSX is the newest PHERAstar reader.



Omega Series