

## APPLICATION NOTE

# MONITORING CIRCADIAN GENE EXPRESSION IN MICRO-ALGAE

Luminescent imaging of luciferase reporters expressed from a circadian promoter in the algal species

Ostreococcus tauri

## Abstract

Expression of firefly luciferase from a circadian-regulated promoter is the most convenient and reliable way to image circadian gene expression in vivo over long-time series. For photosynthetic organisms, constant light conditions are required in between luminescence reads. We report how the Berthold TriStar² multimode plate reader equipped with luminescence module, in combination with an external LED illumination adaptor, can be used to image circadian rhythms in the micro algae *Ostreococcus tauri*.

## Gerben van Ooijen

Institute for Molecular Plant Sciences, School of Biological Sciences, University of Edinburgh, UK

#### **James Grand**

Berthold Technologies UK, UK



## Introduction

Most living organisms possess an endogenous timekeeper to aid efficient negotiation of daily changes to the environment. This timekeeper, the circadian clock, regulates many aspects of the metabolism and physiology of an organism to anticipate predictable changes in the external environment. In plants, this anticipation aids the responses to temporally predictable biotic and abiotic stress [1]. Conversely, disruption of circadian timekeeping severely compromises plant health and reduces yields [2-5]. It is therefore imperative that we understand the



intricate regulation of circadian rhythms in plants, including the factors that affect motion of the transcriptional clockwork itself. The extensive influence of circadian rhythms arises to a large extent from the rhythmic regulation of up to a third of the total transcriptome in the model plant Arabidopsis [6]. These rhythmic gene products are involved in a broad range of cellular processes including metabolic pathways, hormone signalling, and stress responses.

Testing circadian defects in photosynthetic organisms relies on imaging the rhythmic expression of firefly luciferase from circadian clock gene promoters, to allow assessment of circadian phenotypes in transgenic lines (cf. [7, 8]) or upon treatments such as pharmacological inhibition (cf. [9, 10]). Circadian imaging under

constant conditions allows the analysis of treatment effects on circadian period, phase, or amplitude.

Here, cells of the marine eukaryotic species *Ostreococcus tauri*, transgenically expressing the rhythmic luminescent reporter CCA1-LUC [11], are imaged under constant light conditions for 5 days in the presence of pharmacological inhibitor X or mock vehicle. In a second experiment, the effects of overexpressing transgene Y are analysed compared to the parent CCA1-LUC line. Luminescent traces directly reveal how either treatment affects circadian gene expression in these cells.

## Experimental procedures

Cells of Ostreococcus tauri transgenically expressing CCA1-LUC [11] were grown under 12h/12h light/dark cycles in a LEEC PL3 growth cabinet. Medium consists of (24 g/l NaCl, 4 g/l Na $_2$ SO $_4$ , 0.68 g/l KCl, 200mg/l NaHCO $_3$ , 100 mg/l KBr, 25 mg/l H $_3$ BO $_3$ , 3 mg/l NaF, plus hydrous salts: 50 mM MgCl $_2$ \*6H $_2$ O, 10mM CaCl $_2$ \*2H $_2$ O, 0.1  $\mu$ M SrCl\*6H $_2$ O), supplemented with Guillard's F/2 marine enrichment solution (Sigma) and 10 nM H $_2$ SeO $_3$ . Full medium was adjusted to a salinity of 30 ppt. For imaging, 40  $\mu$ l of one-week old cultures were transferred to wells of 384-well microplates (Greiner BioOne)

and 50  $\mu$ l of fresh media supplemented with 2  $\mu$ M luciferin (BioSynth AG) was added. For drugging experiments, 10  $\mu$ l of 10x inhibitor concentration or vehicle controls in media was added to the wells. Transgenic lines overexpressing transgene Y were generated as described previously [11, 12]. Plates were sealed with TopSeal-Aplus seals (Perkin-Elmer) before being transferred to the Tristar² Multimode Microplate Reader. The external light array was fitted with a single layer of Moonlight Blue filter (Lee lighting). Graphs were made using GraphPad Prism.





# TriStar<sup>2</sup> SERIES

## Fully modular microplate reading

Developed for full modularity and equipped with the proprietary ONE-4-ALL optical system, it combines the user friendliness of a multimodal optical system with the sensitivity and performance of a dedicated optical device. You can choose between the affordable Tristar<sup>2</sup> and the more advanced and flexible Tristar<sup>2</sup> S. Start with the reading technology you need to master your research today and upgrade when you need it. The TriStar<sup>2</sup> series provides you with flexibility for today, tomorrow, and beyond.

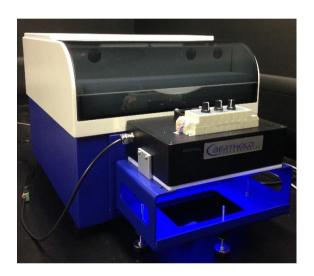
- Monochromator Technology
- High-sensitivity Luminescence
- BRET
- UV/VIS Absorbance
- UV/VIS Fluorescence
- UV/VIS FRET

- Time-Resolved Fluorescence (TRF)
- Time-Resolved FRET (TR-FRET / HTRF®)
- Fluorescence Polarization
- Top and Bottom Reading
- Incubation



## Material

- Berthold TriStar<sup>2</sup> Multimode Microplate Reader
- Luminescence module
- RGB LED illumination array for microplates



**Figure 1**: The TriStar<sup>2</sup> multimode microplate reader with external RGB LED array.



## Instrument settings

The external light was set to 2  $\mu$ mol/sec/m² as measured by a hand-held PAR radiometer (Irradian Ltd). The TriStar² was operated by the ICE software (Berthold Ltd). Microplates with 384 wells were used in this experiment, reading every well. Endpoint luminescence readings were taken for 1.5 seconds per well. An unload command was used with a delay time of 50 minutes (3000 seconds), leading to data with a time resolution of approximately 1 hour. The software was set to read 150 repeats. Screenshots of the instrument settings (left panel) and the results window (right panel) are displayed in Figure 2.

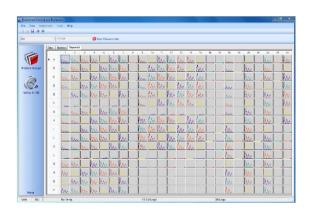


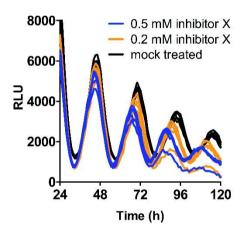
Figure 2: Screenshots of the ICE software. Left: Screenshot of instrument settings. Right: Screenshot of results window.



## Results

We have imaged CCA1-LUC circadian rhythms in constant light for 5 days in the presence of a range of concentrations of pharmacological inhibitor X or mock treatment. Stable and persistent circadian expression rhythms are visualised using the methods described here (Fig. 3, black traces). In the presence of increasing concentrations of the pharmacological inhibitor X, circadian rhythms

are increasingly fast (blue and green traces) and of decreased amplitude. Circadian gene expression was also analysed upon overexpression of a transgene, Y, in the CCA1-LUC parent background. As can be seen in Figure 4, overexpression of gene Y (blue traces) leads to clearly slower circadian rhythms compared to the parent line (black traces).



**Figure 3.** Circadian rhythms in CCA1-LUC luminescence upon pharmacological inhibition. Black lines are mock treated controls, orange and blue lines are treated with 0.2 or 0.5 mM of pharmacological inhibitor X (5 replicate wells are plotted per treatment).

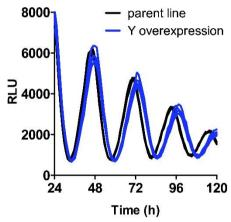


Figure 4. Circadian rhythms in parent CCA1-LUC line versus lines overexpressing transgene Y. Black lines are CCA1-LUC parent line controls, blue lines are transgenic lines overexpressing transgene Y in addition to the CCA1-LUC transgene (5 replicate wells are plotted per line).

## Conclusions

The TriStar<sup>2</sup> multimode plate reader offers a reliable and versatile platform to image luminescent rhythms from circadian reporters. The system is compatible with a range of plate formats (plates from 6 wells to 384 wells), which allows great experimental flexibility. The external light array adds further flexibility as the red, green, and blue LEDs are independently adjustable and can be equipped with an external

socket timer switch. The ICE software is intuitive and user friendly, whilst offering all the necessary flexibility to match your experimental needs. Throughout a long circadian experiment, the software interface provides a great overview of all data collected during the run. The customer service provided by Berthold has always been second to none.



## References

- 1. Harmer SL: The circadian system in higher plants. Annu Rev Plant Biol 2009, 60:357-377.
- 2. Dodd AN, Salathia N, Hall A, Kevei E, Toth R, Nagy F, Hibberd JM, Millar AJ, Webb AA: Plant circadian clocks increase photosynthesis, growth, survival, and competitive advantage. *Science* 2005, 309(5734):630-633.
- 3. Hevia MA, Canessa P, Muller-Esparza H, Larrondo LF: A circadian oscillator in the fungus Botrytis cinerea regulates virulence when infecting Arabidopsis thaliana. *Proceedings of the National Academy of Sciences of the United States of America* 2015, 112(28):8744-8749.
- 4. Wang W, Barnaby JY, Tada Y, Li H, Tor M, Caldelari D, Lee DU, Fu XD, Dong X: Timing of plant immune responses by a central circadian regulator. Nature 2011, 470(7332):110-114.
- 5. Spoel SH, van Ooijen G: Circadian redox signaling in plant immunity and abiotic stress. *Antioxid Redox Signal* 2014, 20(18):3024-3039.
- 6. Covington MF, Maloof JN, Straume M, Kay SA, Harmer SL: Global transcriptome analysis reveals circadian regulation of key pathways in plant growth and development. *Genome Biol* 2008, 9(8):R130.
- 7. van Ooijen G, Hindle M, Martin SF, Barrios-Llerena M, Sanchez F, Bouget FY, O'Neill JS, Le Bihan T, Millar AJ: Functional analysis of Casein Kinase 1 in a minimal circadian system. *PLoS One* 2013, 8(7):e70021.
- 8. Le Bihan T, Hindle M, Martin SF, Barrios-Llerena ME, Krahmer J, Kis K, Millar AJ, van Ooijen G: Label-free quantitative analysis of the casein kinase 2-responsive phosphoproteome of the marine minimal model species Ostreococcus tauri. *Proteomics* 2015.
- 9. O'Neill JS, van Ooijen G, Dixon LE, Troein C, Corellou F, Bouget FY, Reddy AB, Millar AJ: Circadian rhythms persist without transcription in a eukaryote. *Nature* 2011, 469(7331):554-558.
- 10. van Ooijen G, Dixon LE, Troein C, Millar AJ: Proteasome function is required for biological timing throughout the twenty-four hour cycle. *Curr Biol* 2011, 21(10):869-875.
- 11. Corellou F, Schwartz C, Motta JP, Djouani-Tahri el B, Sanchez F, Bouget FY: Clocks in the green lineage: comparative functional analysis of the circadian architecture of the picoeukaryote ostreococcus. *Plant Cell* 2009, 21(11):3436-3449.
- 12. van Ooijen G, Knox K, Kis K, Bouget FY, Millar AJ: Genomic Transformation of the Picoeukaryote Ostreococcus tauri. *J Vis Exp* 2012(65).

For Research Use Only. Not for use in diagnostic procedures.

© 2018 Berthold Technologies. All rights reserved. The trademarks mentioned herein are the property of Berthold Technologies or their respective owners unless otherwise specified.

Berthold Technologies GmbH & Co. KG

Calmbacher Straße 22 75323 Bad Wildbad GERMANY

Phone: +49 7081 177 0 Email: bio@berthold.com

