



MANUAL

Technology: Luminescence Assay Systems

britelite plus

Ultra-high Sensitivity Luminescence Reporter Gene Assay System

Contents	Page
1. Introduction	2
2. britelite™ plus description	3
3. Contents and storage of britelite plus	7
4. Additional requirements	8
5. Assay procedure	9
6. Stability of britelite plus reagent	10
7. Recommendations for use	10
8. Ordering information	11
9. References	12

1. Introduction

Luciferase from the North American firefly (Photinus pyralis) is one of the most frequently used enzymes for reporter gene assays. Firefly luciferase catalyzes the oxidation of the firefly-specific substrate D-luciferin to produce light. This reaction is extremely efficient and the quantum yield is the highest of any characterized bioluminescent reaction. The bright light produced from the reaction makes firefly luciferase a valuable enzyme for reporting promoter activity.^{1,2}

Light emission results from multi-step reactions. The initial reversible reaction step is the activation of D-luciferin in the presence of ATP, Mg²⁺ and luciferase enzyme which leads to enzyme bound adenyl-luciferin with the elimination of inorganic pyrophosphate (PPi).

Subsequent reaction steps involve the oxidation of adenyl- luciferin with molecular oxygen via adenyl-oxyluciferin to yield AMP, CO₂ and oxyluciferin. The oxyluciferin is generated in an electronically excited state which emits light upon transition to the ground state. The overall reaction is shown in Figure 1.

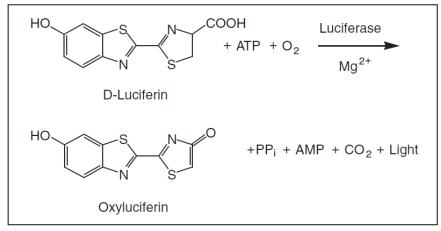


Figure 1: Reaction scheme

When light emission is initiated by the addition of luciferase to a reaction mixture containing ATP, Mg²⁺ and D-luciferin in the presence of oxygen, a fast increase in light intensity can be seen followed by a rapid decrease to a low level of sustained light (flash-type kinetics). These flash-type kinetics have been thought to be the result of the formation of intermediate product (adenyl-oxyluciferin) at the enzyme surface which inhibits the enzyme.⁴ Several substances have been described that stimulate the light production by promoting the release of the inhibitor from the enzyme^{-3,4} The enhancement in enzyme turnover yields an increase in light output.

The Revvity britelite plus assay system is a proprietary formulation that modifies the enzymatic reaction to produce a longer lasting light output at high signal intensity.

2 britelite plus description

britelite plus is a homogeneous long-lived glow type high sensitivity firefly luciferase reporter gene assay system for the quantification of firefly luciferase expression in mammalian cells. The reagent formulation contains compounds that facilitate both cell lysis as well as the luciferase enzymatic reaction. britelite plus is especially suitable for continuous process systems using microplates in ultra-high throughput environments. These in-line systems do not require a very long signal half-life since the time between addition of the reagent and reading the resulting luminescence is relatively short (minutes). In general, britelite plus has a signal half-life of approximately 30 minutes. This relative short half-life has facilitated the design of this luciferase assay system with a very high light output. britelite plus does not contain thiol compounds like dithiothreitol (DTT) and is therefore odor free.

britelite plus offers the following benefits:

- Very high assay sensitivity
- Designed for continuous process systems
- Suitable for higher density microplates such as 384-well and 1536-well plates
- Ideal for smaller-scale sample analysis when maximum sensitivity is desired
- Odor free
- Convenient storage conditions (2 8 °C)

The very high light output of britelite plus is proportional to the firefly luciferase concentration in the sample and allows detection of very low levels of luciferase in microplate formats. Figure 2 shows the assay result of a dilution series of firefly luciferase enzyme in Dulbecco's PBS/0.1% BSA (100 µL per well) using britelite plus in a white 96-well OptiPlate[™] (Revvity) measured with the Revvity TopCount[®] NXT Microplate Scintillation and Luminescence Counter. As can be seen, britelite plus allows for detection of very low levels of luciferase (low femtogram range) with excellent linearity.

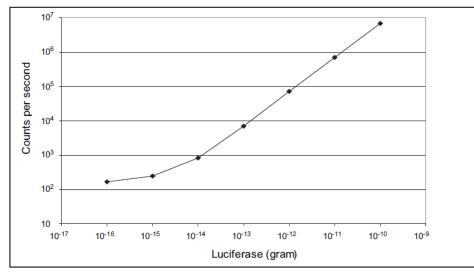


Figure 2: Sensitivity of britelite plus

Figure 3 highlights the time course of the light output of the described experiment above for 10⁻¹⁰ gram luciferase per well. After britelite plus addition, an incubation time of up to 5 minutes is needed for full signal generation.

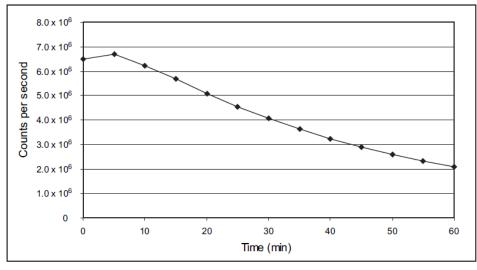


Figure 3: Time course of the luminescent signal of britelite plus

To assess the influence of different culture media on luminescence light output, the following experiment was performed using commercially available basal media in the presence and absence of Fetal Bovine Serum (FBS) and Phenol Red (PR). 100 μ L britelite plus was added to 100 μ L samples of the different media in a white 96-well OptiPlate, where each sample contained luciferase at a concentration of 5 x 10-10 gram per mL. After shaking the plate, the light output was measured using a TopCount NXT. Figure 4 illustrates the results of the

relative luminescence after 5 minutes count delay. Figure 5 shows the results of the stability of the signal as a function of the culture media used.

The results from Figure 4 show that the light output is to some extent affected by FBS but that Phenol Red has a much larger impact on the light output reflecting the concentrations of Phenol Red in the basal media. Consequently, the use of Phenol Red in culture media should be avoided to attain the highest signal. The stability of the luminescent signal is only somewhat dependent on the media as shown in Figure 5.

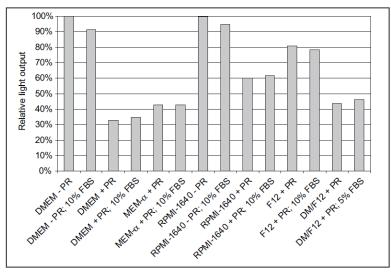


Figure 4: Relative light output in different media

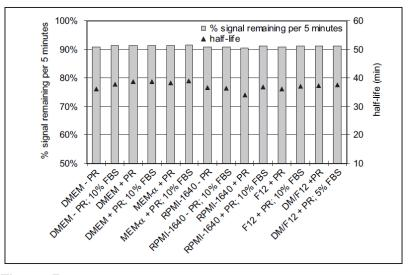


Figure 5: Signal stability in different media expressed as a percentage of signal remaining per 5 minutes and half-life (minutes)

Organic solvents are often used to introduce screening compounds, resulting in the presence of a small percentage of organic solvent in the culture medium. The effect of organic solvents on the light output of britelite plus was investigated in the following experiment. Luciferase in culture medium (DMEM without Phenol Red, supplemented with 10% FBS) was added at 100 μ L per well to a white 96-well CulturPlateTM (Revvity). The medium contained various concentrations of organic solvents (DMSO, ethanol, isopropanol and acetonitrile). Next, 100 μ L of britelite plus was added to the wells. After shaking, the plate was sealed and the luminescence was measured using TopCount NXT. The results presented in Figure 6 show that DMSO and ethanol do not affect the signal at the indicated solvent concentrations. A decrease in signal was shown for the other solvents which was dependent on the specific solvent concentration. The stability of the signal as shown in Figure 7 is not compromised when using these four solvents at the tested concentrations.

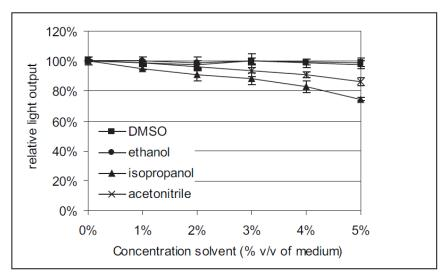


Figure 6: Relative light output in the presence of organic solvents

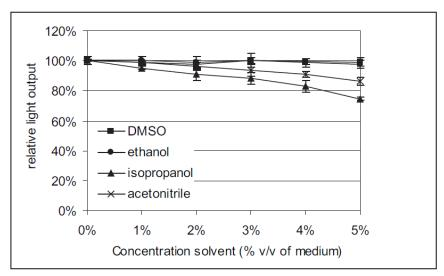


Figure 7: Signal stability in the presence of organic solvents as a percentage of signal remaining per 5 minutes

3 Contents and storage of britelite plus

6066766 – britelite plus 10 mL kit

Each 10 mL kit contains the following components:

- 1 vial britelite plus Lyophilized Substrate
- 1 bottle britelite plus Reconstitution Buffer
- Manual

6066761 – britelite plus 100 mL kit

Each 100 mL kit contains the following components:

- 10 vials britelite plus Lyophilized Substrate
- 1 bottle of britelite plus Reconstitution Buffer
- Manual

6066769 – britelite plus 1000 mL kit

Each 1000 mL kit contains the following components:

- 4 bottles britelite plus Lyophilized Substrate
- 2 bottles britelite plus Reconstitution Buffer
- Manual

britelite plus part number	Kit size	96-well plate	384-well plate	1536-well plate
6066766	10 mL	100	400	3,300
6066761	100 mL	1,000	4,000	33,000
6066769	1000 mL	10,000	40,000	330,000
* Based on recommended volumes per well:				
100 μL for 96-well, 25 μL for 384-well and 3 μL for 1536-well plate				

Number of data points per part number*

Storage conditions

britelite plus is shipped at ambient temperature and must be stored at 2 - 8 °C upon receipt.

For convenience, the britelite plus Reconstitution Buffer can be stored separately at room temperature (max. 22 °C). This allows addition of the Reconstitution Buffer to the Lyophilized Substrate without the need to equilibrate to room temperature.

If stored at the recommended conditions, the kit components are stable through the expiry date found on the kit label.

4 Additional requirements

A detection instrument such as the Revvity TopCount, MicroBeta[®], VICTOR[™] Light, VICTOR3 Multi Label Reader, EnVisionTM or EnSpire[®] is required. CCD camera systems, such as Revvity ViewLux[™] can be used for high throughput applications.

For optimum light yield, low background and minimum well-to-well crosstalk, white microplates should be used. We recommend the use of the Revvity CulturPlate, OptiPlate or ViewPlate® (when visual inspection of cells is preferred). Black plates can also be used when very high signals are expected. Black plates will reduce well-to-well crosstalk but will also quench the light output.

5 Assay procedure

1. Equilibrate the kit components to room temperature (20 - 22 °C) before reconstitution.

2. For the 10 mL and the 100 mL kit reconstitute one vial of britelite plus Lyophilized Substrate with 10 mL of britelite plus Reconstitution Buffer.

For the 1000 mL kit reconstitute one bottle of britelite plus Lyophilized Substrate with 250 mL of britelite plus Reconstitution Buffer.

Mix the contents of the vial gently by inversion and leave for 5 minutes. This should result in a clear homogeneous solution.

Keep the britelite plus reagent at room temperature (20 - 22 °C) before use.

3. Only prepare as much britelite plus reagent as needed for one day.

4. Add britelite plus reagent to each well. Equal volumes of cell culture medium and britelite plus reagent should be used.

For 96-well plates: add 100 μ L to each well containing 100 μ L of cells in medium.

For 384-well plates: add 25 μ L to each well containing 25 μ L of cells in medium.

For 1536-well plates: add 3 μ L to each well containing 3 μ L of cells in medium.

5. Mix the well contents (see section 7).

6. Seal the plate with self-adhesive TopSeal[™]-A.

For complete cell lysis and full signal generation wait at least 1 minute, but not longer than
15 minutes.

8. Measure luminescence within 15 minutes after reagent addition for maximum sensitivity.

6 Stability of britelite plus reagent

Stability of reconstituted britelite plus is approximately:

- > 85 % remaining signal after 8 hours at 20 °C
- > 95 % remaining signal after 8 hours at 4 °C
- > 90 % remaining signal after 24 hours at 4 °C
- > 80 % remaining signal after 48 hours at 4 °C

Freshly prepared reagents can be aliquoted and stored for one month at -20 °C and for three months at -80 °C.

The reagents can be subjected to at least 10 freeze - thaw cycles without significant loss of activity.

7 Recommendations for use

• Mixing of culture medium and reagent is vital to obtain low coefficients of variation between replicates. This is especially important with 384- and 1536-well plates. Optimize liquid handling procedures to attain optimal reagent/medium mixing. For detailed background information see reference 5.

• Phenol Red, as well as other colored compounds, will chemically not interfere with the luciferin/luciferase reaction, but will quench some of the emitted light, resulting in lower assay signals (see Figure 4). For optimal light output, the culture medium can be substituted with **Dulbecco's PBS containing calcium and magnesium** ions prior to the addition of britelite plus.

• The luciferase reaction requires magnesium ions. Although britelite plus does contain these ions it is strongly recommended that the sample contains sufficient amount of these ions. Normal culture media contain these ions at sufficient levels.

• When handling the plates prior to measurement, work under subdued light conditions and avoid direct sunlight or bright fluorescent light. Bright light may cause plate phosphorescence resulting in higher background levels. Plate phosphorescence has a half-life of several minutes.

• If more than one vial of substrate is reconstituted, it is advised to pool all reagents before addition to the plates.

• Optimal room and instrument temperature is 22 °C. Allow plates to adapt to room temperature after removal from the incubator and prior to the addition of the reagent. An adaptation time of 30 minutes is usually sufficient. Stacked plates will need more time to adapt to room temperature.

8 Ordering information

britelite plus	Reorder No
10 mL britelite plus assay kit	6066766
100 mL britelite plus assay kit	6066761
1,000 mL britelite plus assay kit	6066769

For further information on luminescence readers, microplates, seals and luminescence applications please contact your local Revvity representative or visit our website: http:// www.revvity.com

9 References

- 1. Alam, J. and Cook, J.L. (1990). Reporter genes: application to the study of mammalian gene transcription. Anal. Biochem. 1990 Aug 1; 188(2) 245-254
- 2. Collin Goddard (1994). Cell based screening approaches: advantages of highly automated robotics technology in HTS. Handbook for The 1994 International Forum on Advances in Screening Technologies and Data Management, p.19
- 3. Wood, K.V. Recent advantages and prospects for use of beetle luciferases as genetic reporters. In: Bioluminescence and Chemiluminescence current status. Proceedings of the VIth International Symposium on Bioluminescence and Chemiluminescence, Cambridge, September 1990. p. 543. Ed. by P. Stanley and L.J. Kricka.
- 4. Airth, R.L., Rhodes, W.C. and McElroy, W.D. (1958). The functioning of Coenzyme A in Luminescence. Biochemica and Biophysica Acta. Vol. 27, pp. 519-532
- Hancock M.K., Medina M.N., Smith B.M. and Orth A.P.. Microplate Orbital Mixing Improves High-Throughput Cell-Based Reporter Assay Readouts. J. Biomol. Screen. 2006 Nov 27. 2 2007; vol. 12: pp. 140 - 144

This product and/or the use of this product are covered by Revvity patent applications. By purchasing this product, the end user is granted a limited license to use the britelite kit and reagents for research purposes. Purchase does not include any right to use, develop or otherwise exploit this product commercially.

Limited Use License

This product is distributed and sold for life science research and commercial applications, but not for diagnostic use. Any use of this product other than for life science research and commercial applications is strictly prohibited.

GraphPad Prism is a registered trademark of GraphPad Software, Inc.

U.S. Pat. 7425422 and 7951551; European Pat. 1885874 and foreign equivalent patents



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

www.revvity.com



Revvity, Inc. 940 Winter Street Waltham, MA 02451 USA (800) 762-4000 www.revvity.com For a complete listing of our global offices, visit www.revvity.com Copyright ©2023, Revvity, Inc. All rights reserved.