



twinlite

Firefly & Renilla Luciferase
Reporter Gene Assay System



QUICK START GUIDE

A RESEARCH PRODUCT FOR RESEARCH PURPOSES ONLY

Instructions for use of product 6066709 (1000 assay kit; 96-well format)

Reagent Preparation

1. Allow all reagents to reach room temperature (20 – 22°C).
2. **Passive Lysis Solution:** PLS is a ready to use reagent.
3. **firelite plus reagent:**
 - Reconstitute one vial *firelite plus Lyophilized Substrate* with 53 mL of *firelite plus Reconstitution Buffer*.
 - Mix the contents of the vial gently by inversion and leave for 5 minutes.
 - o Unused reagent can be stored at –20°C (≤ 2 months) or –80°C (≤ 2 year).
4. **renlite plus reagent:**
 - Add 2.12 mL *renlite plus Substrate (50X)* to 106 mL *renlite plus Buffer*.
When a smaller quantity of *renlite plus reagent* is desired, add an appropriate volume of *renlite plus Substrate (50X)* to the required amount of *renlite plus Buffer* to achieve the correct concentration of 1X *renlite plus Substrate*. For example, add 300 µL *renlite plus Substrate (50X)* to 15 mL *renlite plus Buffer*.
 - Mix the contents of the bottle gently by inversion and leave for 5 minutes.
 - o Unused reagent can be stored at –20°C (≤ 3 months) or –80°C (≤ 2 year).

Cell Lysate Preparation

1. Remove cell growth medium from the cell layer.
2. Wash the cells with a sufficient amount of PBS at room temperature. Swirl briefly to remove loose cells and residual growth medium. Remove the wash solution as much as possible.
3. Add to the cell layer the recommended volume of PLS according to the table below.

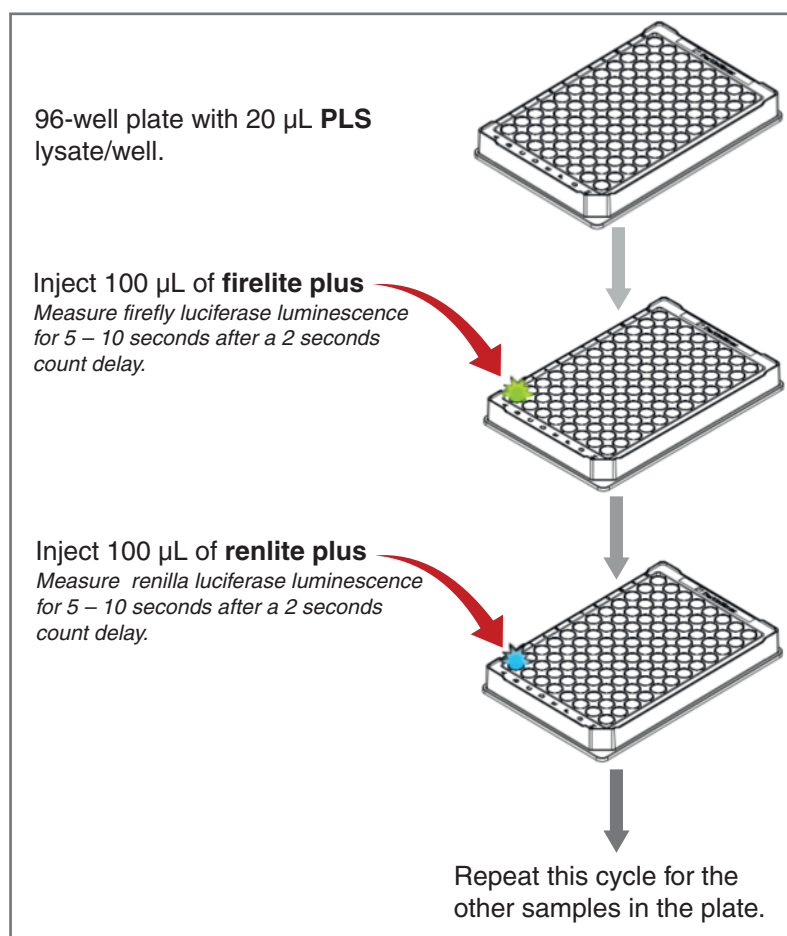
Plate type	PLS/well
6 well	500 µL
12 well	250 µL
24 well	100 µL
48 well	65 µL
96 well	20 µL

4. Place the plate on an orbital shaker or on a rocking platform so that the PLS covers the cell layer evenly for optimal lysis. Shake the plate for 15 minutes at room temperature.
5. The cell lysate can now be used in the twinlite assay. If the lysate is not needed the same day, store at –20°C (≤ 2 months) or –80°C (≤ 1 year). When the cells are cultured in an opaque 96-well plate, then the assay can be performed directly in the same plate without lysate transfer.

Measuring firefly and renilla luciferase luminescence; twinlite assay

1. Set the luminometer injectors 1 and 2 to dispense 100 μL **firelite plus** and **renlite plus** reagent respectively.
2. Set a count delay of 2 seconds between the reagent injection and measuring luminescence. Set the luminescence read time between 5 to 10 seconds.
3. Fill and rinse the designated injectors of the luminometer with the prepared reagents.
4. Load the microplate containing the samples (20 μL /well) in the luminometer, dark adapt for a few minutes to decrease plate phosphorescence (to lower plate background levels) and start the measurement.

Assay using a 96-well plate (opaque)



This product and/or its use is covered by the following patents and corresponding patent applications worldwide, owned by PerkinElmer Health Sciences B.V.: US Patent No. 8,512,968; EP Patent No. EP2222870B1; China Patent No. CN101889095B; and Australia Patent No. AU2008319571B2.



For a complete listing of our global offices visit www.perkinelmer.com

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-9254602
www.perkinelmer.com