INSTRUCTIONS

Pierce[®] Firefly Luciferase Flash Assay Kit

<u>16174 16175</u>

Number	Description	
16174	Pierce Firefly Luciferase Flash Assay Kit, sufficient reagents to perform 100 assays for Firefly luciferase activity in cultured cell lysate	
	Kit Contents:	
	Firefly Flash Assay Buffer, 5mL, store at 4°C	
	D-Luciferin, Lyophilized, 3mg, store at 4°C	
	2X Cell Lysis Buffer, 6mL, store at room temperature	
16175	Pierce Firefly Luciferase Flash Assay Kit , sufficient reagents to perform 1000 assays for Firefly luciferase activity in cultured cell lysate	
	Kit Contents:	
	Firefly Flash Assay Buffer, 50mL, store at 4°C	
	D-Luciferin, 30mg, store at 4°C	
	2X Cell Lysis Buffer, 60mL, store at room temperature	
	Storage: Upon receipt store kit at 4°C or store individual components as indicated above. Kit is shipped on dry ice.	

2367.1

Table of Contents

Introduction	1	
Procedure Summary	2	
Important Product Information	2	
Additional Materials Required	3	
Material Preparation	3	
Procedure for Firefly Luciferase Flash Assay		
A. Cell Transfection		
B. Cell Lysis	3	
C. Firefly Luciferase Flash Assay	3	
Troubleshooting		
Related Thermo Scientific Products		
General References		

Introduction

The Thermo Scientific Pierce Firefly Luciferase Flash Assay provides a highly sensitive system for detecting intracellular luciferase activity from promoter or pathway activation in mammalian cell culture experiments. The bioluminescent signal produced by firefly luciferase results from the oxidation of D-Luciferin (Figure 1). The light output correlates with the amount of firefly protein expressed, which is proportional to the activity of the promoter for firefly expression.





Figure 1. Chemical reaction of luciferin and red firefly luciferase. Light, with an emission maximum of 613nm, is produced from the oxidation of D-luciferin by red firefly luciferase in an ATP-dependent reaction.

Procedure Summary



Important Product Information

- For long-term use, store D-Luciferin (lyophilized) at 4°C protected from light. Briefly centrifuge tubes of D-Luciferin (lyophilized) before use.
- Store Firefly Luciferase Flash Assay Working Solution (Working Solution) protected from light. Working Solution must be at room temperature (20-25°C) before use and is stable for up to four hours at room temperature.
- To avoid cross-contamination, use a new disposable pipette tip for each transfer. Always use a new disposable reagent reservoir for each reagent.
- Firefly luciferase in cell lysate is subject to degradation. Store firefly luciferase-containing cell lysate on ice; perform assays immediately following cell lysis. Addition of protease inhibitors to the cell lysis buffer can increase firefly luciferase stability (see the Related Thermo Scientific Products Section).
- Avoid exposing reagents to excessive heat or light during storage and incubation.
- Do not mix reagents from different lots. Discard unused working solutions after assay completion. Do not combine leftover reagents with those reserved for additional plates.
- Individual components might contain corrosives and/or preservatives. Wear gloves while performing the assay to avoid contact with samples and reagents. Please follow proper disposal procedures.
- Dispense and equilibrate to room temperature only the reagent volumes needed for the number of plates being used.



Additional Materials Required

- Reagents and equipment for propagating mammalian cells in culture
- Reagents and materials for transfection of plasmid DNA into mammalian cells (e.g., Thermo Scientific TurboFect Transfection Reagent, Product No. R0533)
- Modified Dulbecco's Phosphate-buffered saline (DPBS) (e.g., Thermo Scientific BupH Modified Dulbecco's PBS, 8mM sodium phosphate, 2mM potassium phosphate, 140mM sodium chloride, 10mM potassium chloride; pH 7.4; Product No. 28374)
- Laboratory platform shaker
- Pipettes and/or liquid handling equipment
- Luminometer or other luminescence-monitoring instrument

Note: Use a luminometer equipped with reagent injectors to perform > 24 assays. (Optional)

• White or black opaque, 96- or 384-well microplates

Material Preparation

100X D-Luciferin Stock Solution	For 100 reactions, reconstitute lyophilized D-Luciferin pellet in 50µL of Firefly Flash Assay Buffer. Store at -20°C for up to two months.
Working Solution	For 100 reactions, add 50μ L of 100X reconstituted D-Luciferin to 5mL of Firefly Flash Assay Buffer. Use 50μ L of the Working Solution per reaction.
	Note: If using a luminometer with injectors, prepare sufficient reagent to prime the pumps in addition to reagent required for the assay.
	Note: Dilute Thermo Scientific Pierce Firefly Signal Enhancer (100X) (Product No. 16180) to 1X in Working Solution.
1X Cell Lysis Buffer	Dilute 2X Cell Lysis Buffer with an equal volume of ultrapure water.

Procedure for Firefly Luciferase Flash Assay

A. Cell Transfection

1. Plate ~10,000 cells/well in a 96-well plate. Incubate plates overnight at 37°C in 5% CO₂. If using a different plate size, adjust the cell number accordingly. Use only cells growing in log phase at a passage number \leq 15.

Note: Plate enough wells to perform the experiment in triplicate; include appropriate controls (i.e., non-transfected cell control and non-treated cell control).

- 2. Use a standard protocol to transfect mammalian cells with a Firefly luciferase plasmid.
- 3. Incubate cells for 16-72 hours at 37° C in 5% CO₂ in a cell culture incubator.
- 4. Proceed with the individual experimental protocol for cell treatment.

B. Cell Lysis

- Aspirate media from the cells, rinse once with 50-100µL/well of 1X DPBS buffer (Thermo Scientific BupH Modified Dulbecco's PBS, Product No. 28374), aspirate DPBS and add 100µL/well of 1X Cell Lysis Buffer. Do not disturb the cell monolayer during the DPBS rinse.
- 2. Shake the plate on a platform shaker at moderate speed for 15 minutes. Check for complete cell lysis using a light microscope. If lysis is incomplete, continue shaking the plate for 15 additional minutes.

C. Firefly Luciferase Flash Assay

- 1. Add 10-20 μ L/well of cell lysate to a white or black, opaque 96-well plate.
- 2. Program the luminometer; if using an injector, prime the injector with Working Solution.



- 3. Add 50µL of Working Solution to each well.
- 4. Immediately after adding the reagent, detect the light output.

Note: Follow the manufacturer's recommendations for using injector velocity to obtain a uniform coating of liquid in the well. Adjust the detector's integration time to achieve a signal within the linear range of the instrument.

Troubleshooting

Problem	Possible Cause	Solution
No signal or low signal	Low transfection efficiency	Optimize transfection conditions using a visual transfection control (e.g., a plasmid over-expressing a fluorescent protein)
		Verify plasmid DNA quality; use only transfection grade DNA
		Use actively dividing, low passage cells
		Use a different cell type
	No promoter activity	Use conditions known for promoter activation
		Incubate cells for a longer time
		Change growth conditions to improve expression
		Use a different promoter
	D-Luciferin auto-oxidized	Protect substrate from light and air and maintain 100X D-Luciferin at -20°C
		Prepare new Working Solution if used longer than 4 hours
	Low luciferase expression	Use Pierce Firefly Signal Enhancer (100X) (Product No. 16180)
		Lyse cells in a smaller volume of 1X Cell Lysis Buffer
		Use a different promoter or growth conditions to improve expression
		Increase the integration time on the instrument
		Scale-up the volume of sample and reagent per well
	Degraded luciferase protein	Store cell lysates on ice and perform assays immediately following cell lysis
High signal	High luciferase expression	Reduce incubation time before collecting samples
		Decrease the integration time on the instrument
		Dilute the sample Note: A low sample volume can increase assay variability; dilute the sample and use the recommended volume of 10-20µL per assay
High background signal	Control sample was contaminated	Incubate cells for a longer time



Related Thermo Scientific Products

See our website for a col	inplete list of related lucherase products.
16180	Pierce Firefly Signal Enhancer
16155	pMCS-Firefly Luc
16156	pCMV-Firefly Luc
16157	pTK-Firefly Luc
16189	Pierce Luciferase Cell Lysis Buffer
R0532-4	TurboFect Transfection Reagent
28374	BupH [™] Modified Dulbecco's PBS Packs, 40 packs
28344	20X Modified Dulbecco's PBS Buffer
78425	Halt Protease Inhibitor Single-use Cocktail, EDTA-free

See our website for a complete list of related luciferase products

General References

Tatsumi, H., et al. (1989). Luciferase cDNA from Japanese firefly, Luciola cruciata: cloning, structure and expression in Escherichia coli. J Biolumin Chemilumin 3(2):75-8.

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than the original purchaser of the Product ("Buyer").

No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s).

There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

Current product instructions are available at <u>www.thermoscientific.com/pierce</u>. For a faxed copy, call 800-874-3723 or contact your local distributor. © 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.