



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Steady-Luc™ Firefly HTS Assay Kit

A homogeneous firefly luciferase reporter gene assay kit with a luminescence half-life of about 3 hours, allowing batch processing in microplates.



### Product attributes

<b>Assay type/options</b>	High-throughput assay, Homogeneous assay, Luminescence (glow-type)
<b>Detection method/readout</b>	Microplate reader (luminescence)
<b>Storage Conditions</b>	Store at -60 °C or below

## Product Description

The Steady-Luc™ HTS assay system is a homogeneous, high-sensitivity firefly luciferase reporter gene assay kit. It is a glow-type assay with signal half-life of about 3 hours, allowing batch processing in microplates.

Firefly luciferase is widely used as a reporter for studying gene regulation and function, and for pharmaceutical screening. It is a very sensitive genetic reporter due to the lack of endogenous activity in mammalian cells or tissues. However, the light production resulting from the reaction leads to formation of suicidal adenylyl-luciferin at the enzyme surface. It results in very short half-life of the light emission with a flash-type kinetics. Because of this, flash-type luminescence must be detected immediately after initiating the reaction, which is not compatible with batch analysis of samples.

Biotium's Steady-Luc™ Firefly Assay Buffer contains a proprietary mixture of substances that prevent the adenylyl-luciferin reaction, producing a long lasting, glow-type signal. It produces luminescence signal with a half-life of about 3 hours, for a high-throughput microplate assay.

Glow-type luciferase assays like Steady-Luc™ have lower luminescence signal compared to flash-type assays. The sensitivity and limit of detection of the assay will depend on luciferase expression levels in your experimental system as well as luminometer sensitivity.

Biotium also offers [Steady-Luc™ Firefly HTS Assay Kit \(Lyophilized\)](#) (catalog no. 30028L), a new packaging format that includes lyophilized assay buffer for convenient room temperature shipping and storage at -20°C.

## More Luciferase Kits & Related Products

Biotium Cat. No.	Product	Description	Promega Assay	Promega Cat. No.
<a href="#">30138</a>	<a href="#">Steady-ATP™ HTS Viability Assay Kit</a>	Highly sensitive ATP-based cell luminescent viability assay with signal half-life of 5 hours for high-throughput screening in multi-well plates.	CellTiter-Glo® Luminescent Cell Viability Assay	G7570, G7571, G7572, G7573
<a href="#">30085</a>	<a href="#">Firefly Luciferase Assay Kit 2.0</a>	Flash-type firefly luciferase assay kit.	Firefly Luciferase Assay System	E1483, E1500, E1501, E4030, E4530, E4550
<a href="#">30075</a>	<a href="#">Firefly Luciferase Assay Kit (Lyophilized)</a>	Firefly luciferase assay kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20°C.		
<a href="#">30028</a>	<a href="#">Steady-Luc&amp;#153; Firefly HTS Assay Kit</a>	Glow-type firefly luciferase assay with signal half-life of about 3 hours for high-throughput screening in multi-well plates.	Steady-Glo® Assay System	E2510, E2520, E2550
<a href="#">30028L</a>	<a href="#">Steady-Luc Firefly HTS Assay Kit (Lyophilized)</a>	Steady-Luc kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20°C.		
<a href="#">30082</a>	<a href="#">Renilla Luciferase Assay Kit 2.0</a>	Flash-type Renilla luciferase assay kit.	<i>Renilla</i> Luciferase Assay System	E2810, E2820
<a href="#">30081</a>	<a href="#">Firefly and Renilla Luciferase Single Tube Assay Kit</a>	Assay kit for measuring firefly and Renilla activities sequentially in the same sample in a single tube.	Dual-Luciferase® Assay System (Firefly & <i>Renilla</i> )	E1910, E1960, E1980

This datasheet was generated on June 6, 2026 at 06:23:44 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/steady-luc-firefly-hts-assay-kit/>

