

# Luciferase Assay Systems

Our luciferase assay systems provide simple and accurate measurements for luciferase output. Our advantages include:



## Sensitivity

Highly sensitive detection of Firefly and Renilla luciferase activity that outshines competitor assays



## High-Throughput

Simple homogeneous protocol minimizes handling steps to support high throughput screening applications



## Compatibility

Works well with a variety of common media containing 0-10% serum and phenol red



## Stability

Stable luciferase signal output for over one hour, providing flexible incubation time

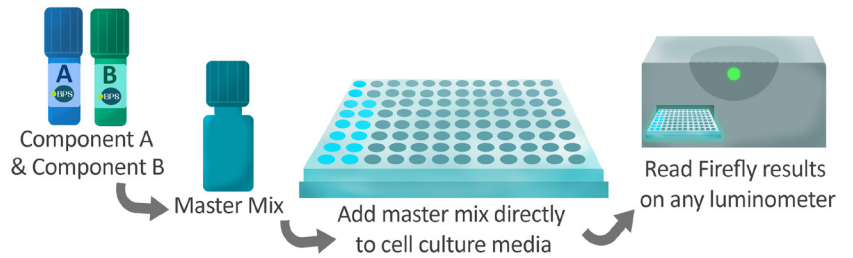
### ONE-Step™ Luciferase Assay System #60690

The ONE-Step™ product is designed to be used for high-throughput, sensitive quantitation of Firefly luciferase activity in mammalian culture.

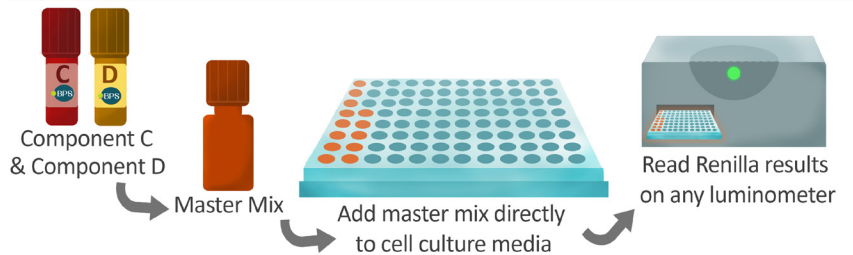
### TWO-Step (Firefly & Renilla) Luciferase Assay System #60683

The TWO-Step Luciferase product provides high-throughput, rapid quantitation of both Firefly and Renilla luciferases in order to improve signal accuracy by normalizing with the control Renilla reporter.

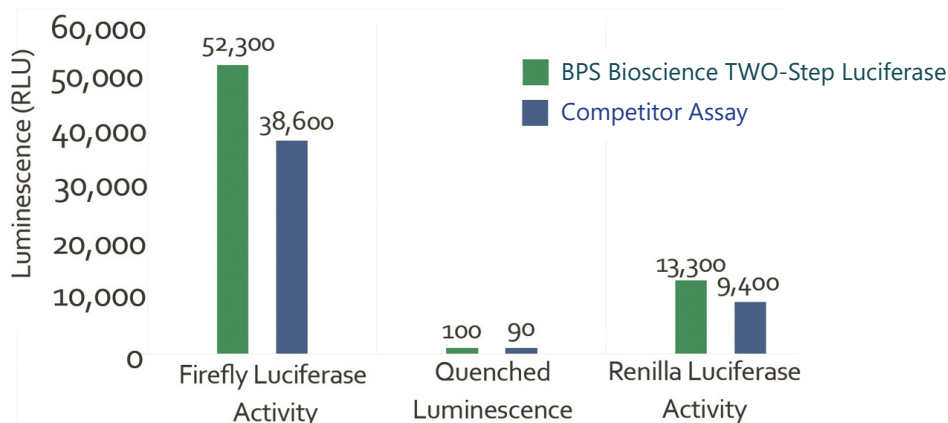
#### Firefly Luciferase Signal



#### Renilla Luciferase Signal



## Better than the Standard



Learn more at:

[bpsbioscience.com/product-types/cell-lines/media-and-reagents](https://bpsbioscience.com/product-types/cell-lines/media-and-reagents)

