

## 5X Firefly Luciferase Lysis Buffer

Additional concentrated cell lysis buffer that is designed for use with Biotium's Firefly Luciferase Assay Kits when performing large sample volume assays.



### Product attributes

**Storage Conditions** Store at -10 to -35 °C or below

## Product Description

5X Firefly Luciferase Lysis Buffer is concentrated lysis buffer designed for use with our Firefly Luciferase Assays.

- Additional 5X Firefly Luciferase Lysis Buffer for larger volume assays
- For use with Firefly Luciferase Assay Kit 2.0 ([30085](#)) and Firefly Luciferase Assay Kit (Lyophilized) ([30075](#))

Our kits are supplied with sufficient 5X Firefly Luciferase Lysis Buffer to perform the stated number of assays with cells grown in 96 – 24 well plates. 5X Firefly Luciferase Lysis Buffer is sold separately for use with the kits when more lysis buffer (e.g. >100 uL/well) is required. This Buffer is not compatible with *Renilla* Luciferase. If you require additional lysis buffer for use with *Renilla* Luciferase Assay Kit 2.0 ([30082](#)) or Firefly & *Renilla* Luciferase Single Tube Assay Kit ([30081](#)), order 1X Passive Lysis Buffer 2.0 ([99821-100mL](#)).

## More Luciferase Kits

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 <i>Renilla</i> 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 <i>Renilla</i> activities sequentially in the same sample in a single tube.	Dual-Luciferase® Assay System (Firefly & <i>Renilla</i> )	E1910, E1960, E1980