

1X Passive Lysis Buffer 2.0

Extra lysis buffer for use with Firely & Renilla Luciferase Single Tube Assay Kits.



Product attributes

Storage Conditions Store at -10 to -35 °C

Product Description

1X Passive Lysis Buffer 2.0 is a cell lysis buffer that is designed for use with Biotium's Firefly & Renilla Luciferase Single Tube Assay Kit ([30081](#)). Those kits are supplied with sufficient lysis buffer to perform the stated number of assays with cells grown in 96 – 24 well plates. 1X Passive Lysis Buffer 2.0 is sold separately for use with the kits when more lysis buffer (e.g. >100 uL/well) is required.

- Additional 1X Passive Lysis Buffer 2.0 for larger volume assays
- For use with Biotium's Firefly & Renilla Luciferase Single Tube Assay Kit ([30081](#))

Oure kits are supplied with sufficient lysis buffer to perform the stated number of assays with cells grown in 96 – 24 well plates. 1X Passive Lysis Buffer 2.0 is sold separately for use with the kits when more lysis buffer (e.g. >100 uL/well) is required.

This buffer also can also be used to dilute recombinant firefly or *Renilla* luciferase enzymes without the need to add BSA to stabilize the enzymes.

Note: If additional lysis buffer is required for the Firefly Luciferase Assay Kit 2.0 ([30085](#)) or the Firefly Luciferase Assay (Lyophilized) ([30075](#)), purchase 5X Firefly Luciferase Lysis Buffer ([99923](#)). If additional lysis buffer is required for Renilla Luciferase Assay Kit 2.0 ([30082](#)), purchase 5X Passive Lysis Buffer ([99912](#)).

More Luciferase Kits

Biotium Cat. No.	Product	Description	Promega Assay	Promega Cat. No.
30138	Steady-ATP™ HTS Viability Assay Kit	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
30085	Firefly Luciferase Assay Kit 2.0	Flash-type firefly luciferase assay kit.	Firefly Luciferase Assay System	E1483, E1500, E1501, E4030, E4530, E4550
30075	Firefly Luciferase Assay Kit (Lyophilized)	Firefly luciferase assay kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20 °C.		
30028	Steady-Luc&#153; Firefly HTS Assay Kit	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
30028L	Steady-Luc Firefly HTS Assay Kit (Lyophilized)	Steady-Luc kit with lyophilized assay buffer for economical room temperature shipping and convenient storage at -20 °C.		
30082	Renilla Luciferase Assay Kit 2.0	Flash-type Renilla luciferase assay kit.	<i>Renilla</i> Luciferase Assay System	E2810, E2820
30081	Firefly and Renilla Luciferase Single Tube Assay Kit	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