



Glowing products for science

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

## 5X Passive Lysis Buffer

Additional concentrated lysis buffer for use with our Renilla Luciferase Assay Kits in large volume assays.



### Product attributes

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

## Product Description

5X Passive Lysis Buffer is concentrated lysis buffer designed for use with our Renilla Luciferase Assays when additional volumes of lysis buffer are required.

- Additional 5X Passive Lysis Buffer for larger volume assays
- For use with Renilla Luciferase Assay Kit 2.0 ([30082](#))

Our kits are supplied with sufficient 5X Passive Lysis Buffer to complete the stated number of assays using 96-well to 24-well plates. 5X Passive Lysis Buffer is sold separately for use with the kits when more lysis buffer (e.g., >100 uL per well) is required. If you require additional lysis buffer for use with Firefly & *Renilla* Luciferase Single Tube Assay Kit ([30081](#)), use 1X Passive Lysis Buffer 2.0 ([99821-100mL](#)). If you require additional lysis buffer for use with Firefly Luciferase Assay Kit 2.0 ([30085](#)) or Firefly Luciferase Assay Kit (Lyophilized) ([30075](#)), use 5X Firefly Luciferase Lysis Buffer ([99923](#)).

## 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 Renilla luciferase assay kit.	Renilla 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 & Renilla)	E1910, E1960, E1980

## References

- Biochem J, 477, 137 (2020), [DOI: 10.1042/BCJ20190596](#)
- PLoS ONE, 14(2), e0212850 (2019), [DOI: 10.1371/journal.pone.0212850](#)
- J Tradit Complement Med, 9, 227, (2018), [DOI: 10.1016/j.itcme.2017.07.004](#)
- Int. J. Mol. Sci, 19, 774, (2018), [DOI: 10.3390/ijms19030774](#)
- Int Heart J, 59, 190, (2018), [DOI: 10.1536/ihj.16-372](#)
- Front. Immunol., 9, 1564, (2018), [DOI: 10.3389/fimmu.2018.01564](#)
- J. Orthop. Res., 35, 876, (2017), [DOI: 10.1002/jor.23342](#)
- Bulletin Chem Soc. J (2016), [DOI: 10.1246/bcsj.20150405](#)
- PLoS ONE, 8(1), e53936 (2013), [DOI: 10.1371/journal.pone.0053936](#)
- Biol Syst 1, 1, (2012), [DOI: 10.4172/bs.1000104](#)
- Int J Biochem Mol Biol, 2(4), 318, (2011), [PMCID: PMC3242428](#)

This datasheet was generated on June 17, 2026 at 06:45:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5x-passive-lysis-buffer/>