

# 1 X Passive Lysis Buffer

To be used with PromoCell's Firefly & Renilla Luciferase Single-tube Assay Kit (Cat. No. PK-CA707-30081)

## Instruction Manual

Product Name	Product Description	Size	Catalog Number
1 X Passive Lysis Buffer	To be used with PromoCell's Firefly & Renilla Luciferase Single-tube Assay Kit (Cat. No. PK-CA707-30081)	100 ml	PK-CA707-99821

### Introduction

1X Passive Lysis Buffer is a cell lysis buffer that is designed for use with PromoCell's *Firefly & Renilla Luciferase Single Tube Assay Kit* (Cat. No. PK-CA707-30081) That kit is supplied with sufficient 1X Passive Lysis Buffer to perform the stated number of assays with cells grown in 96 – 24 well plates. 1X Passive Lysis Buffer is sold separately for use with the kits when more lysis buffer (e.g. >100 µl/well) is required. The buffer also can 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 *Firefly Luciferase Assay Kit* or Firefly Luciferase Assay (lyophilized luciferase), purchase 5X Firefly Lysis Buffer (Cat. No. PK-CA707-99923). If additional lysis buffer is required for *Renilla Luciferase Assay Kit*, purchase 5X Passive Lysis Buffer (Cat. No. PK-CA707-99912).

### Storage and Stability

Store at –20°C or below. The buffer is stable for at least 12 months from date of receipt when stored as recommended. The buffer is stable to three freeze-thaw cycles.

### Cell Proliferation Assay Procedures

1. Note: 1X Passive Lysis Buffer is ready to use without dilution.

Remove the growth medium from the cultured cells and gently wash the cells once with a sufficient volume of phosphate buffered saline (PBS) to cover the surface of the culture vessel. Remove the PBS and add 1X Passive Lysis Buffer using the volume recommended below for each type of well:

Wells/plate	Lysis buffer/well
6 well	500 µl
12 well	250 µl
24 well	100 µl
48 well	65 µl
96 well	20 µl

2. Place the culture plates on a rocking platform or orbital shaker with gentle rocking/shaking to ensure complete and even coverage of the cell monolayer with 1X passive lysis buffer. Rock the culture plates at room temperature for 15 minutes.

Note: Cultures that are overgrown are often more resistant to complete lysis and typically require an increased volume of passive lysis buffer and/or an extended treatment period to ensure complete lysis and/or scraping cells off the culture plates. PromoCell offers mini cell scrapers (Cat. No. PK-CA707-22003) for harvesting lysates from 96-, 24-, and 48-well plates.

Note: 1X Passive Lysis Buffer contains protein stabilizers that may affect results of protein quantitation assays.

3. Transfer the lysate to a tube or vial. Optional: the lysate can be cleared by centrifugation for 30 seconds at top speed in a refrigerated microcentrifuge and transferred into a new tube. Place at 4°C until ready to assay. Store lysates at -20°C or -70°C if assay will not be performed on the same day.
4. To assay lysates using the PromoCell *Firefly & Renilla Luciferase Single Tube Assay*, follow the protocol provided with the kit.

## Intended Use

---

For in vitro research use only. Not for diagnostic or therapeutic procedures.

### PromoCell GmbH

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promokine.info](mailto:info@promokine.info)  
[www.promokine.info](http://www.promokine.info)

### USA/Canada

Phone: 1 – 866 – 251 – 2860 (toll free)  
Fax: 1 – 866 – 827 – 9219 (toll free)

### Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)  
Fax: 0800 – 100 83 06 (gebührenfrei)

### France

Téléphone: 0800 90 93 32 (ligne verte)  
Téléfax: 0800 90 27 36 (ligne verte)

### United Kingdom

Phone: 0800 – 96 03 33 (toll free)  
Fax: 0800 – 169 85 54 (toll free)

### Other Countries

Phone: +49 6221 – 649 34 0  
Fax: +49 6221 – 649 34 40