

# ab287865 – Luciferase Reporter Assay Kit

For the measurement of expressed luciferase in vitro.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab287865>

## Storage and Stability

The entire kit may be stored at -20°C.

## Materials Supplied

Item	Quantity	Storage Condition
Substrate A	1 vial	-20°C
Substrate B	1 vial	-20°C
Cell Lysis Buffer	100 ml	-20°C

## Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Luminescence Microplate Reader

## Reagent Preparation

- Ensure that all reagents have reached room temperature before performing assays.
- Reconstitute Substrate A and B: Add 20 ml of Cell Lysis Buffer to each vial and mix well.
- Store both substrates at -70°C after each use. Store Lysis Buffer at 4°C.

**Δ Note:** The following protocol is designed for using with adherent cultures growing in 35-mm tissue culture plates. If you are using plates of different size, adjust the volume proportionally. Read entire protocol before starting experiments.

## Assay Procedure

- Preparation of Cell Lysate: Remove media from cell culture plates and rinse once with PBS:
  - Add 1 ml of PBS and collect cells from plates by scraping and then transfer to a 1.5 ml microcentrifuge tube. Spin cells at 5000 rpm for 3 min and remove PBS.
  - Resuspend cells in 200 µl Cell Lysis Buffer and incubate on ice for 5 min. Centrifuge at 14000 rpm for 1 min. Transfer extract (supernatant) to a fresh tube and use immediately, or store at -70°C.
- Luciferase Assay:
  - Place 20 - 100 µl cell extract into an assay cuvet or microplate well. Be sure to use the same volume for each sample.
  - Add 100 µl Substrate A.
  - Within 10 min, inject 100 µl Substrate B. Read the signal immediately using a luminometer.

**Δ Note:** The amount of extract required may vary depending on the luciferase expression level and the instrumentation used; the amount used should be adjusted to keep the signal within the

linear range of the assay. The time between adding substrate B and reading signal should be as short as possible (1-2 sec) and consistent from sample to sample.

## Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)