

Product datasheet

Luciferase Reporter Assay Substrate Kit - Renilla
ab228546

[1 References](#) [1 Image](#)

Overview

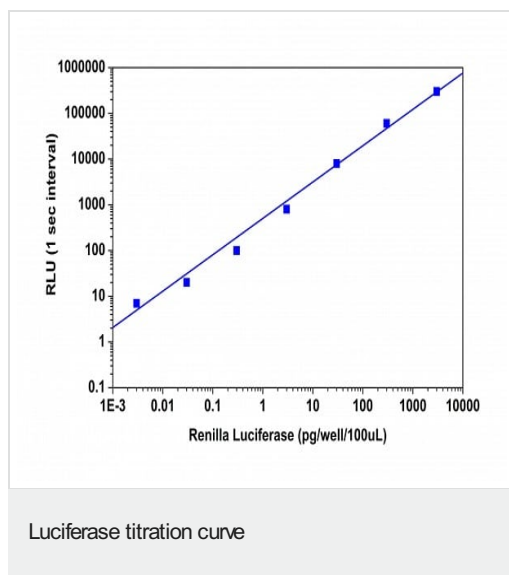
Product name	Luciferase Reporter Assay Substrate Kit - Renilla
Detection method	Luminescent
Sample type	Cell culture extracts, Adherent cells
Product overview	Luciferase Reporter Assay Substrate Kit - Renilla (ab228546) provides a fast and sensitive method to detect the luciferase from sea pansy (<i>Renilla reniformis</i>). It uses a proprietary luminogenic formulation to quantify <i>Renilla luciferase</i> activity in cell-based assays. The formulation generates a luminescent product that gives strong luminescence upon interaction with Renilla luciferase. The kit provides all the essential components. It has high sensitivity and can be performed in a convenient 96-well and 384-well microtiter-plate format. The “glow-type” signal with a half-life of one hour provides a consistent signal across large number of assay plates. The assay is compatible with standard cell growth media. This kit enables the measurement of primary expression or gene expression with wild type and the synthetic hRluc genes.
Notes	Common reporter genes include beta-galactosidase, beta-glucuronidase and luciferase. The most versatile reporter gene is the firefly luciferase. Recently there is steadily increasing use of other luciferases, such as Renilla luciferase since these reporters are smaller and do not require the presence of ATP.
Platform	Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 units
100X Luciferase Substrate	1 x 100µl
Assay Buffer	1 x 10ml

Images



Renilla Luciferase was measured with Luciferase Reporter Assay Substrate Kit – Renilla (ab228546) in a white 96-well plate.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors