

# Calcium Assay Kit (Luminometric) ab112114

### 画像数 1

#### 製品の概要

<b>製品名</b>	Calcium Assay Kit (Luminometric)
<b>検出方法</b>	Luminescent
<b>サンプルの種類</b>	Adherent cells, Suspension cells
<b>アッセイタイプ</b>	Cell-based
<b>製品の概要</b>	<p>Calcium flux assays are preferred methods in drug discovery for screening G protein coupled receptors (GPCR). This kit uses a highly calcium-sensitive and membrane-permeable coelenterazine analog as a calcium indicator for the cells that are transfected with apoaequorin gene. Aequorin is a calcium-sensitive bioluminescent protein from the jellyfish <i>Aequorea victoria</i> that has been used extensively as a calcium indicator in cells. The aequorin complex emits blue light when bound to calcium ions. The luminescence intensity is directly proportional to the <math>Ca^{2+}</math> concentration.</p> <p>ab112114 is much more sensitive than the fluorescence-based calcium assays. This kit provides an optimized assay method for monitoring the G-protein-coupled receptors and calcium channels. The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation. It might be useful for monitoring the intracellular calcium mobilization in a specified compartment given that recombinant apoaequorin proteins can now be targeted to specific organelles, cells and tissues. ab112114 is more sensitive than the fluorescent calcium assays.</p> <p>Visit our <a href="#">FAQs page</a> for tips and troubleshooting.</p>
<b>特記事項</b>	ab112114 should be stored desiccated.
<b>試験プラットフォーム</b>	Microplate reader

#### 製品の特性

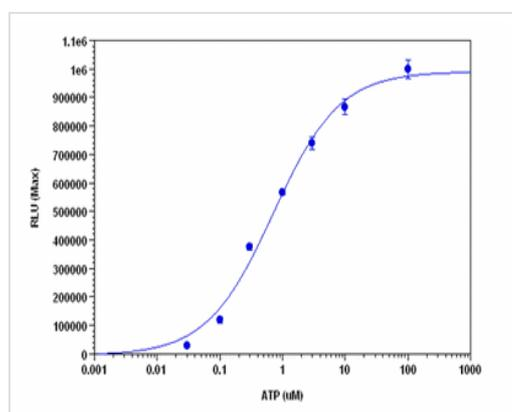
**保存方法** Store at -20°C. Please refer to protocols.

内容	10 x 96 tests	100 x 96 tests
100% ETOH	1 x 500µl	1 x 5000µl
Assay Buffer	1 x 100ml	1 x 100ml

内容	10 x 96 tests	100 x 96 tests
Coelenterazine Analog	1 vial	10 vials

**関連性** Calcium is essential for all living organisms, where  $\text{Ca}^{2+}$  sequestration and release into and out of the cytoplasm functions as a signal for many cellular processes. 99% of calcium is found in bones and teeth with the remaining 1% found in the blood and soft tissue. Serum calcium levels are tightly controlled (8.4-11.4 mg/dL) and any variation outside this range can have serious effects. Calcium plays a role in mediating the constriction and relaxation of blood vessels, nerve impulse transmission, muscle contraction, and hormone secretion. Calcium ion channels control the migration of calcium ions across cell membranes, permitting the activation and inhibition of a wide variety of enzymes. Causes of low calcium levels include chronic kidney failure, vitamin D deficiency, and low blood magnesium levels that can occur in severe alcoholism.

## 画像



ATP Dose Response on CHO-aeq cells. CHO cells stably transfected with apoaequorin were seeded overnight at 50,000 cells/100  $\mu\text{L}$ /well in a white wall/clear bottom 96-well plate. The growth medium was removed and the cells were incubated with 100  $\mu\text{L}$  of dye-loading solution using ab112114 for 3 hours at room temperature and protected from light. ATP (25  $\mu\text{L}$ /well) was added to achieve the final indicated concentrations. The  $\text{EC}_{50}$  of ATP is about 0.8  $\mu\text{M}$ .

Functional Studies - Calcium Assay Kit  
(Luminometric) (ab112114)

**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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

---

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