

# Fluo-4 Assay Kit (Calcium) ab228555

[2 References](#) [画像数 1](#)

### 製品の概要

製品名	Fluo-4 Assay Kit (Calcium)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
製品の概要	Fluo-4 Assay Kit (Calcium) (ab228555) provides a homogeneous fluorescence-based assay for detecting the intracellular calcium mobilization. The kit offers an optimized assay method for monitoring G-protein-coupled receptors (GPCRs) and calcium channels.

The assay can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation.

**特記事項** This product is intended to be used for monitoring calcium fluctuations in vivo in live cells using the following HTS imaging plate readers: FLIPR™, FDSS, BMG NOVOstar™, FLEXStation, ViewLux, IN Cell Analyzer or Arrayscan.

Calcium flux assays are preferred methods in drug discovery for screening G protein coupled receptors (GPCR). Cells expressing a GPCR of interest that signals through calcium are pre-loaded with Fluo-4 AM which can cross cell membrane. Once inside the cell, the lipophilic blocking groups of Fluo-4 AM are cleaved by non-specific cell esterase, resulting in a negatively charged Fluo-4 dye that stays inside cells, and its fluorescence is greatly enhanced upon binding to calcium. When cells are stimulated with screening compounds, the receptor signals release of intracellular calcium, which greatly increase the fluorescence of Fluo-4.

**試験プラットフォーム** Microplate reader

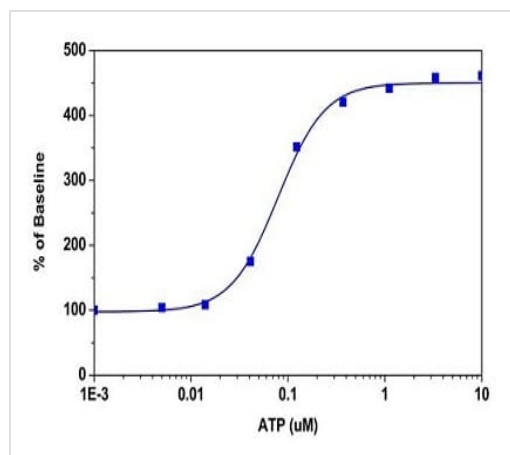
### 製品の特性

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

内容	10 units
10X F127 Plus	1 x 10ml
Fluo-4 AM	1 vial

内容	10 units
HHBS	1 x 100ml

## 画像



ATP dose response was measured in CHO-K1 cells with Fluo-4 Assay Kit (Calcium)

CHO-K1 cells were seeded overnight at 60,000 cells/100  $\mu$ L/well in a Costar black wall/clear bottom 96-well plate. The cells were incubated with 100  $\mu$ L of dye-loading solution using the Fluo-4 Assay Kit (Calcium) for 1 hour at 37°C, and then at room temperature for another 15 minutes. ATP (50 $\mu$ L/well) was added by Flexstation 3 to achieve the final indicated concentrations.

**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