

Calcium Assay kit (Fluorometric) ab112115

[6 References](#) [画像数 2](#)

製品の概要

製品名	Calcium Assay kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Serum, Plasma, Cell Lysate
アッセイタイプ	Quantitative
検出感度	0.03 mM
検出範囲	0.04 mM - 1.5 mM
全工程の試験時間	0h 30m
製品の概要	<p>Calcium Assay Kit (Fluorometric) ab112115 provides a simple method for detecting calcium in physiological solutions by using a red fluorescence probe. The fluorescence signal can be easily read by a fluorescence microplate reader at Ex/Em = 540/590 nm.</p> <p>The kit can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step. The calcium assay can be completed within 30 minutes.</p> <p>The calcium assay kit can be used to detect as little as 0.03 mM calcium. The kit has a broad dynamic range (30 µM to 1 mM).</p> <p>Calcium assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix - incubate for 5-30 min whilst measuring the fluorescence intensity with a microplate reader
特記事項	<p>This product is intended to be used <i>in vitro</i> with solution-based samples such as cell extracts. If you would like to detect calcium <i>in vivo</i> using live cells, please check our Fluo-8, Rhod-4 or Fura-2 based assays.</p> <p>Previously called Calcium Quantification Kit.</p>
試験プラットフォーム	Microplate reader

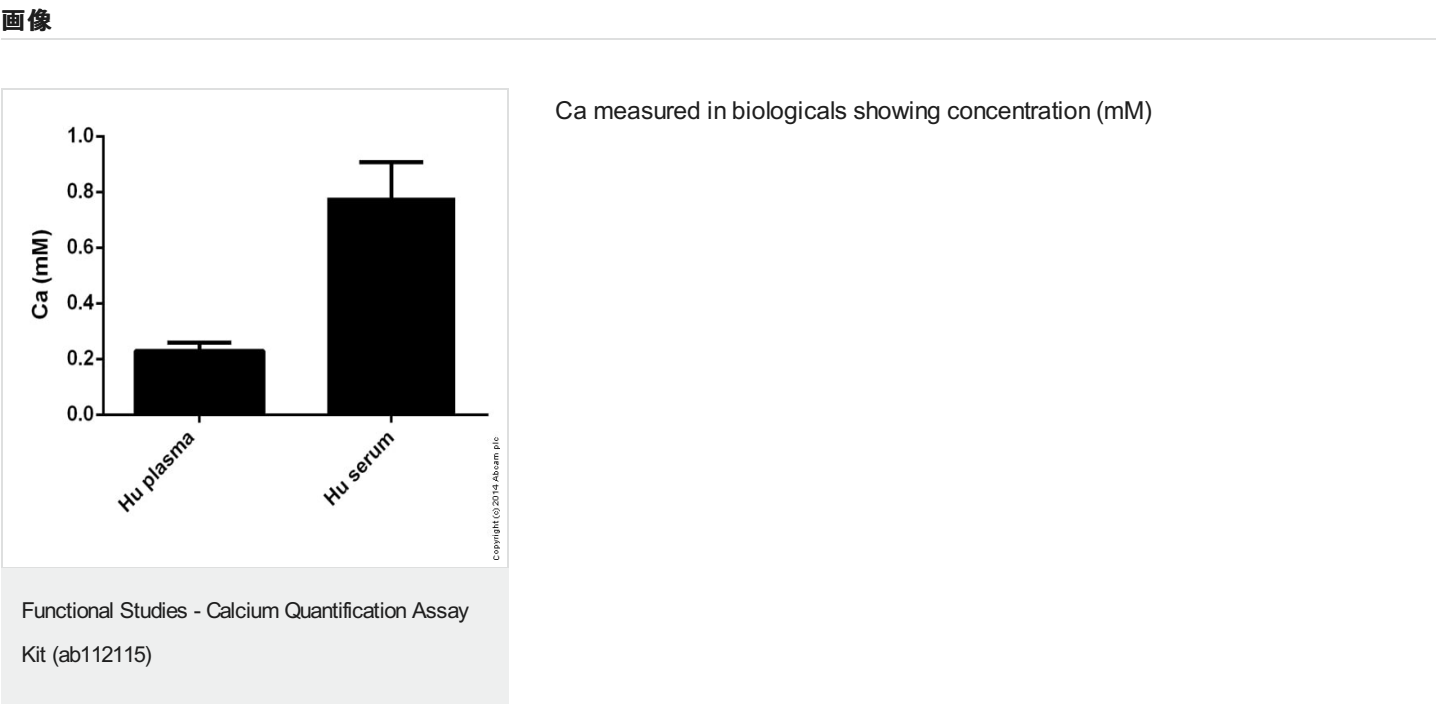
製品の特性

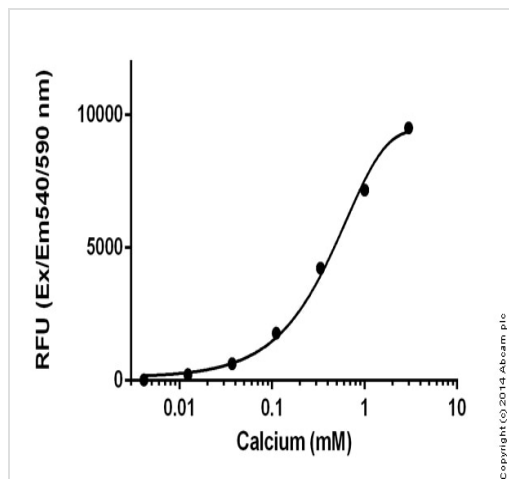
保存方法 Store at -20°C. Please refer to protocols.

内容	200 tests
300 mM Calcium Standard	1 x 0.5ml
Assay Buffer	1 x 10ml
Rhod Red Indicator	2 vials

関連性

Calcium is essential for all living organisms, where 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.





Standard curve (fluorimetric): mean of duplicates (+/-SD) with background readings subtracted

Functional Studies - Calcium Quantification Assay
Kit (ab112115)

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