

Phosphate Assay Kit (Fluorometric) ab102508

3 References [画像数 1](#)

製品の概要

製品名	Phosphate Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate
アッセイタイプ	Quantitative
検出感度	> 0.1 nmol/well
検出範囲	2 μM - 10 μM
全工程の試験時間	2h 00m
製品の概要	<p>Abcam's Phosphate Assay Kit (Fluorometric) provides a sensitive, easy, quick means of assessing phosphate over a wide range of concentrations. In the assay, inorganic phosphate reacted with maltose to produce glucose in the present of a special enzyme. The glucose is specifically oxidized to generate a product that reacts with the OxiRed probe to generate fluorescence (Ex/Em=535/587nm). The kit can be used to detect Pi produced through reactions involving ATPases, GTPases, 5'-nucleotidase, protein phosphatases, acid and alkaline phosphatases, and phosphorylase, etc. from a variety of samples. Glucose background can be eliminated by doing a glucose control.</p>

Visit our [FAQs page](#) for tips and troubleshooting.

特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K420 Phosphate Assay Kit (Fluorometric). K420-100 is the same size as the 100 test size of ab102508.</p> <p>Inorganic phosphate (Pi) is one of the most important ions in biological systems. It functions in a variety of roles. One of the most important roles is as a molecular switch, turning enzyme activity on and off through the mediation of the various protein kinases and phosphatases in biological systems. A highly sensitive assay is desired to monitor Pi in variety samples or monitor Pi changes during kinase and phosphatase reactions.</p>
------	--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

試験プラットフォーム	Microplate reader
------------	-------------------

製品の特性

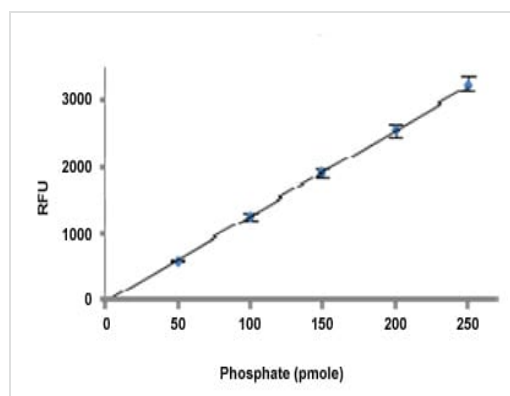
保存方法	Store at -20°C. Please refer to protocols.
------	--------------------------------------------

内容	100 tests
Converter	1 vial
Developer	1 vial
Phosphate Assay Buffer	1 x 25ml
Phosphate Standard	1 x 50µl
Phosphate Substrate	1 x 0.2ml
PicoProbe (in DMSO)	1 x 200µl

関連性

Phosphate is one of the most important of the inorganic ions in biological systems. It functions in a variety of roles. One of the most important roles is as a molecular switch, turning enzyme activity on and off through the mediation of the various protein kinases and phosphatases in biological systems. Phosphate is also of great importance in mineralization processes and is a primary stimulus of algal blooms frequently found in bodies of fresh water, due to run-off from areas of high fertilizer use.

画像



Phosphate Standard Curve - 2 hours

Functional Studies - Phosphate Fluorescent Assay
Kit (ab102508)

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