

Fructose-6-Phosphate Assay Kit (Fluorometric) ab204720

8 References 画像数 2

製品の概要

製品名	Fructose-6-Phosphate Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Cell Lysate, Tissue Lysate
検出感度	1 µM
製品の概要	Fructose-6-Phosphate Assay Kit (ab204720) is a fluorescence based simple, highly sensitive and rapid means of quantifying fructose-6-phosphate (F6P) in a variety of samples. In the assay, F6P is converted to glucose-6-phosphate (G6P) which is subsequently oxidized with the generation of a fluorescent product. This assay kit can detect F6P in the range of 0.01 to 0.5 nmoles with detection sensitivity ~ 1 µM of F6P.
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K689 PicoProbe™ Fructose-6-Phosphate Fluorometric Assay Kit. K689-100 is the same size as the 100 test size of ab204720.</p> <p>Fructose-6-phosphate (F6P) is an important intermediate in the glycolytic pathway which leads from glucose to pyruvate. It is formed from glucose-6-phosphate and is further phosphorylated to fructose-1,6-diphosphate which is subsequently cleaved to glyceraldehyde phosphate and dihydroxyacetone phosphate. The transformation of glucose-6-phosphate is controlled by phosphoglucose isomerase, a very interesting enzyme in that it possesses multiple functions, as an isomerase, a neuroleukin, an autocrine motility factor and a differentiation and maturation mediator.</p>
試験プラットフォーム	Microplate reader

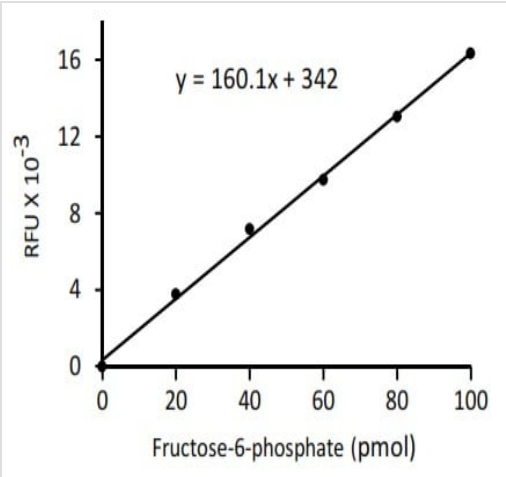
製品の特性

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

内容	100 tests
Assay Buffer V	1 x 25ml
Converter Enzyme X	1 vial
Developer IX	1 vial

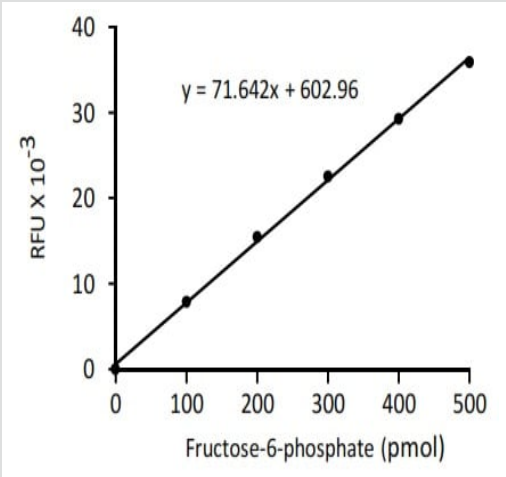
内容	100 tests
Development Enzyme Mix IX	1 vial
F6P Standard	1 vial
PicoProbe I	1 x 400µl

画像



Typical F6P Standard calibration curve

Fructose-6-Phosphate Assay Kit (ab204720)



Typical F6P Standard calibration curve

Fructose-6-Phosphate Assay Kit (ab204720)

**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