

Starch Assay Kit ab83393

★★★★★ 1 Abreviews 5 References 画像数 1

製品の概要

製品名	Starch Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell Lysate, Tissue Lysate, Food samples
アッセイタイプ	Quantitative
検出感度	> 4 µg/ml
検出範囲	4 µg/ml - 2000 µg/ml
全工程の試験時間	0h 40m
製品の概要	Abcam's Starch Assay Kit provides an easy, accurate assay to measure starch levels in a variety of samples. In the assay, starch is hydrolyzed to glucose which is oxidized to generate color (570 nm) and fluorescence (Ex/Em = 535/587 nm). The assay can detect starch at 0.0004 to 2 mg/ml.

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 K647 Starch Colorimetric/Fluorometric Assay Kit. K647-100 is the same size as the 100 test size of ab83393.</p> <p>Starch is a complex carbohydrate consisting of a large number of glucose units. All plants contain starch, present as amylose, (linear α-1,4 linked polymer) and amylopectin, (highly α-1,6 branched α-1,4 polymer). Starch generally contains 0-25% amylose and 75-100% amylopectin.</p>
------	---

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

製品の特性

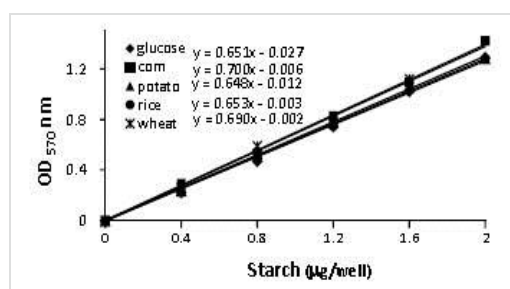
保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	100 tests
Development Buffer	1 x 25ml
Development Enzyme Mix (Lyophilised)	1 vial
Hydrolysis Buffer	1 x 25ml

内容	100 tests
Hydrolysis Enzyme Mix (Lyophilised)	1 vial
OxiRed Probe in DMSO	1 vial
Starch Standard (2.0 mg/ml)	1 x 100µl

**関連性** Starch is a complex carbohydrate consisting of a large number of glucose units. All plants contain starch, present as amylose, (linear alpha-1,4 linked polymer) and amylopectin, (highly alpha-1,6 branched alpha-1,4 polymer). Starch generally contains 0-25% amylose and 75-100% amylopectin.

## 画像



Starch Standard Curve: Different types of pure starch were extracted with 10N KOH/H<sub>3</sub>PO<sub>4</sub> as described. Assays were performed following the kit protocol.

Functional Studies - Starch Assay Kit (ab83393)

**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