

Maltose Assay Kit ab83388

1 References 画像数 2

製品の概要

製品名	Maltose Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
全工程の試験時間	1h 00m
製品の概要	Abcam's Maltose Assay Kit provides a fast, easy and sensitive method for quantifying maltose in various biological samples (e.g. serum, plasma, body fluids, food, growth medium, etc.).

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 K628 Maltose Colorimetric/Fluorometric Assay Kit. K628-100 is the same size as the 100 test size of ab83388.</p> <p>Maltose (C<sub>12</sub>H<sub>22</sub>O<sub>11</sub>; FW: 342.3), one of the main fuel sources to generate the universal energy molecule ATP, is the major disaccharide that is generated from hydrolysis of starch in food. Maltose contains two glucose units joined by a α-1,4-glycosidic linkage, which can be easily converted to two glucoses by α-D-glucosidase. The generated glucose can be specifically oxidized to produce a product that interacts with the probe to generate color and fluorescence. Thus, maltose can be determined by either colorimetric (spectrophotometry at λ = 570 nm) or fluorometric (Ex/Em = 535/587 nm) methods.</p>
------	--

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

製品の特性

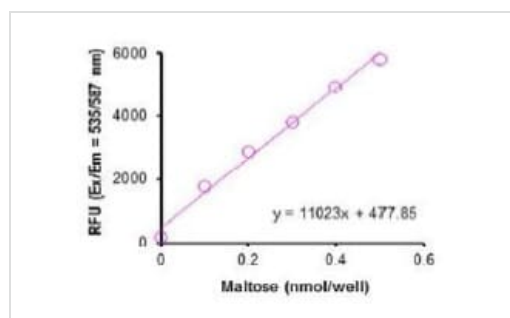
保存方法	Store at -20°C. Please refer to protocols.
------	--

内容	100 tests
a-D-Glucosidase (Lyophilized)	1 vial
Enzyme Mix (Lyophilized)	1 vial
Maltose Assay Buffer	1 x 25ml

内容	100 tests
Maltose Probe in DMSO	1 x 200μl
Maltose Standard (100 nmol/μl)	1 x 100μl

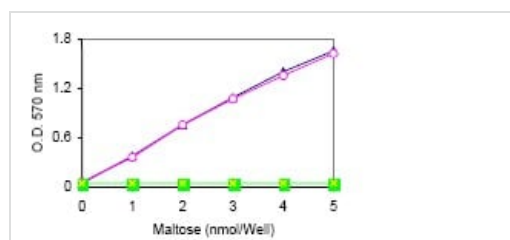
**関連性** Maltose ( $C_{12}H_{22}O_{11}$ ; FW: 342.3), one of the main fuel sources to generate the universal energy molecule ATP, is the major disaccharide that is generated from hydrolysis of starch in food.

## 画像



Example of Fluorometric Standard Curve

Maltose with  $\alpha$ -D-Glucosidase Standard Curve. Assays were performed following kit instructions



Functional Studies - Maltose Assay Kit (ab83388)

Maltose Standard Curve. Assays were performed following kit instructions. Open circle is maltose with  $\alpha$ -D-Glucosidase. Solid square is maltose without  $\alpha$ -D-Glucosidase. Triangle is free glucose (divided by 2).

**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