

Maltose and Glucose Assay Kit ab65335

2 References [画像数 1](#)

製品の概要

| | |
|----------|--|
| 製品名 | Maltose and Glucose 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 |
| 検出感度 | > 0.01 nmol/well |
| 検出範囲 | 0.01 nmol/well - 10 nmol/well |
| 全工程の試験時間 | 1h 00m |
| 製品の概要 | <p>Maltose and Glucose Assay Kit (ab65335) can measure glucose or maltose levels in a variety of biological samples (e.g. serum, plasma, body fluids, food, growth medium, etc.). The assay is based on the ability of glucose oxidase to specifically oxidize free glucose, producing a product that interacts with the glucose probe. The amount of product can be either colorimetric (spectrophotometry at OD= 570 nm) or fluorometric (Ex/Em = 535/587 nm) method.</p> <p>This kit can detect 10 pmol to 10 nmol glucose per assay. It can also be used for isomaltose detection.</p> |

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

特記事項 This product is manufactured by BioVision, an Abcam company and was previously called K618 Maltose and Glucose Colorimetric/Fluorometric Assay Kit. K618-100 is the same size as the 100 test size of ab65335.

Glucose (C₆H₁₂O₆; FW: 180.16) and Maltose (C₁₂H₂₂O₁₁; FW: 342.3) are the main fuel sources to generate the universal energy molecule ATP. Maltose is the major disaccharide that 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.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

試験プラットフォーム Microplate reader

製品の特性

保存方法

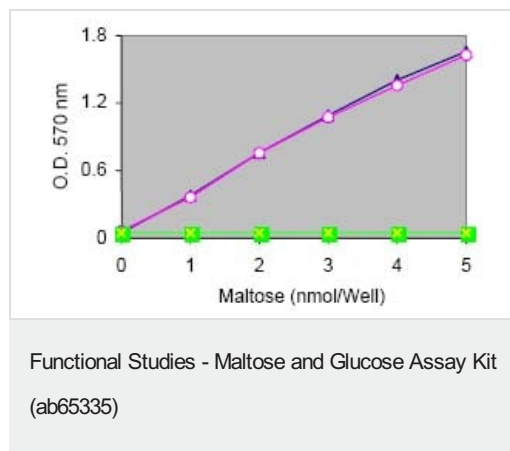
Store at -20°C. Please refer to protocols.

| 内容 | 100 tests |
|-----------------------------------|-----------|
| alpha-D-Glucosidase (Lyophilized) | 1 vial |
| Assay Buffer | 1 x 25ml |
| Glucose Enzyme Mix (lyophilized) | 1 vial |
| Glucose Probe | 1 x 0.2ml |
| Maltose Standard (100 nmol/μl) | 1 x 100μl |

関連性

Maltose ($C_{12}H_{22}O_{11}$; FW:342.30) is a disaccharide formed from two units of glucose ($C_6H_{12}O_6$; FW: 180.16) joined via a α -1,4-glycosidic linkage. Maltose is the major disaccharide generated from hydrolysis of starch and in living organisms, maltose is hydrolyzed by maltase. Once inside the cells, glucose will be processed and enter the glycolysis pathway to ultimately produce energy in form of ATP.

画像



Maltose Standard Curve. Assays were performed using various amounts of Maltose standard according to kit instructions. Circle is maltose with α -D-Glucosidase. Square is maltose without α -D-Glucosidase. Triangle is glucose with 2 times nmol of Maltose.

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