

Glucose and Sucrose Assay Kit ab65334

3 References [画像数 2](#)

製品の概要

製品名	Glucose and Sucrose 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	0h 40m
種交差性	交差種: Mammals
製品の概要	<p>Abcam's Glucose and Sucrose Assay Kit provides a convenient means for measuring glucose and sucrose levels from various biological samples (e.g. serum, plasma, body fluids, food, growth medium, etc.). To measure glucose level, glucose oxidase specifically oxidizes free-glucose generating a compound that reacts with the glucose probe to produce resorufin, which can be detected colorimetrically (OD570nm) or fluorometrically (Ex/Em 535/587). To measure sucrose, invertase can be added to the reaction to convert sucrose to free glucose and fructose, so total glucose level can be measured. Then the sucrose level = Total Glucose – Free Glucose. Visit our FAQs page for tips and troubleshooting.</p>
特記事項	<p>Glucose (C₆H₁₂O₆; FW: 180.16) and sucrose (C₁₂H₂₂O₁₁; FW:342.3) are the important fuel sources to generate universal energy molecule ATP. Measurement of glucose or sucrose level can be very important in both research and development process. Sucrose is a disaccharide which can be converted into one glucose and one fructose when adding Invertase.</p>
試験プラットフォーム	Microplate reader

製品の特性

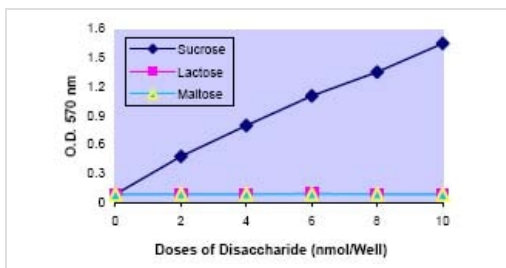
保存方法 Store at -20°C. Please refer to protocols.

内容	100 tests
Assay Buffer II	1 x 25ml
Development Enzyme Mix II	1 vial
Invertase	1 vial

内容	100 tests
OxiRed Probe	1 x 200µl
Sucrose Standard	1 x 100µl

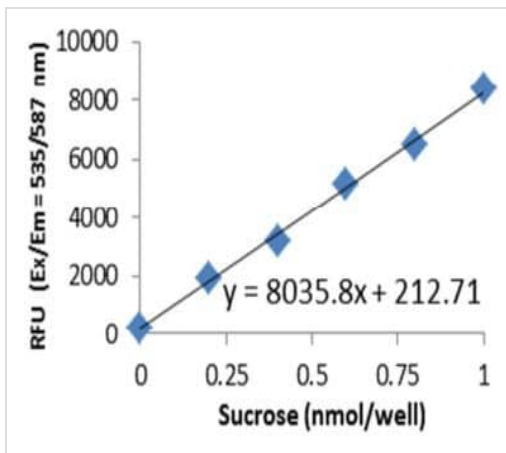
関連性 Sucrose ($C_{12}H_{22}O_{11}$; FW:342.3) is a disaccharide composed of glucose ($C_6H_{12}O_6$; FW: 180.16) and fructose ($C_6H_{12}O_6$; FW:180.16). Sucrose can be broken down by different In the human body, sucrose can be broken down by sucrase and/or invertase into its components, which are rapidly absorbed into the bloodstream. Once inside the cells, both monosaccharides will be processed and enter the glycolysis pathway to ultimately produce energy in form of ATP.

画像



Colorimetric standard calibration Curve of Sucrose.

Functional Studies - Glucose and Sucrose Assay Kit (ab65334)



Fluorometric Standard Curves of Sucrose

Glucose and Sucrose Assay Kit Fluorometric Standard Curve

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