

Product datasheet

Glucose Assay Kit ab65333

15 References 4 Images

Overview

Product name	Glucose Assay Kit
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids
Assay type	Quantitative
Sensitivity	1 μ M
Range	1 μ M - 10000 μ M
Assay time	0h 40m
Species reactivity	Reacts with: Mammal
Product overview	<p>Glucose Assay Kit (ab65333) provides direct measurement of glucose in various biological samples (e.g., serum, plasma, body fluid, food, growth medium, etc.). Glucose Enzyme Mix specifically oxidizes glucose to generate a product which reacts with a dye to generate color ($\lambda = 570$ nm) and fluorescence (Ex/Em = 535/587 nm). The generated color and fluorescence is proportionally to the glucose amount. The method is rapid, simple, sensitive, and suitable for high throughput. The assay is also suitable for monitoring glucose level during fermentation and glucose feeding in protein expression processes. The kit detects 1-10000 μM glucose samples.</p> <p>If you have reducing substances in your samples, we recommend using Glucose Detection Kit II (ab102517).</p>

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

Notes	<p>Glucose ($C_6H_{12}O_6$; FW: 180.16) is a very important fuel source to generate universal energy molecules ATP. Glucose level is a key diagnostic parameter for many metabolic disorders. Measurement of glucose can be very important in both research and drug discovery processes.</p>
--------------	--

Tested applications	Suitable for: Functional Studies
----------------------------	---

Properties

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

Components	Identifier	100 tests
Assay Buffer	WM	1 x 25ml
Glucose Enzyme Mix (lyophilized)	Green	1 vial

Components	Identifier	100 tests
Glucose Probe (in DMSO)	Red	1 x 200µl
Glucose Standard (100 nmol/µl)	Yellow	1 x 100µl

Relevance

Glucose (C₆H₁₂O₆; FW: 180.16) is a ubiquitous energy source in most organisms, from bacteria to humans. The breakdown of carbohydrates produces mono- and disaccharides, most of which is glucose. Through glycolysis and TCA (citric acid cycle), glucose is oxidized to eventually form CO₂ and water, generating the universal energy molecule ATP. Glucose is a primary source of energy for the brain and a critical component in the production of proteins and in lipid metabolism and therefore measurement of glucose level is a key diagnostic parameter for many metabolic disorders.

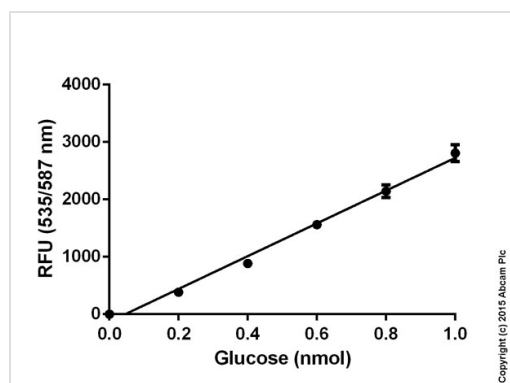
Applications

Our [Abpromise guarantee](#) covers the use of **ab65333** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

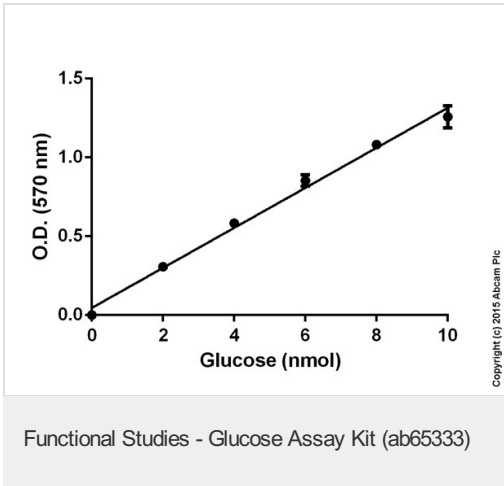
Application	Abreviews	Notes
Functional Studies		Use at an assay dependent dilution.

Images

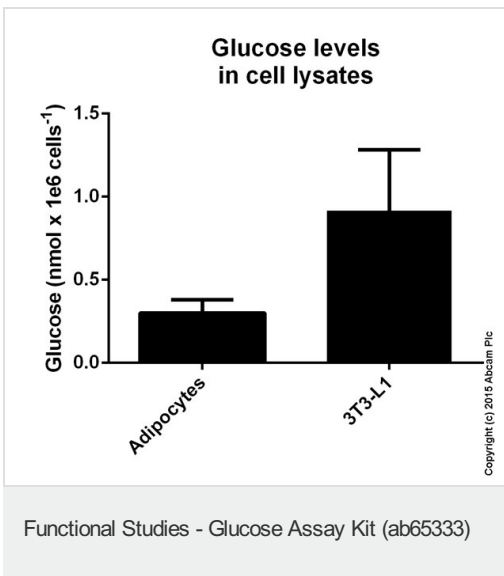


Standard curve: mean of duplicates (+/- SD)
with background reads subtracted

Functional Studies - Glucose Assay Kit (ab65333)

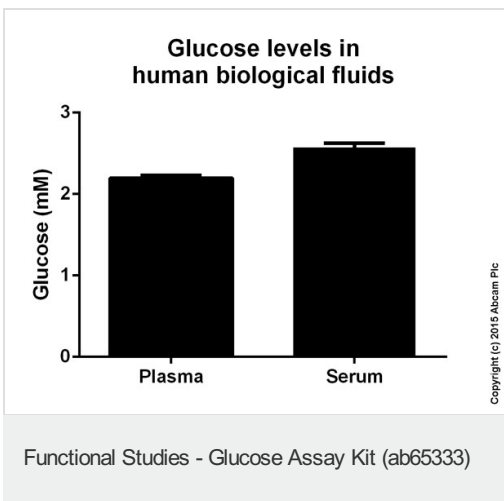


Standard curve: mean of duplicates (+/- SD) with background reads subtracted



Glucose measured in cell lysates showing quantity (nmol) per million cells.

Samples with the concentration of 2×10^7 cells/mL were used. Samples were diluted 1.5-13.5 fold and measured colorimetrically.



Glucose measured in human biological fluids showing quantity (μmol) per mL of tested sample. Samples were diluted 13.5 fold and measured colorimetrically.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors