

# Glucose Assay Kit (Fluorometric, High Sensitivity)

## ab169559

[9 References](#) [2 Images](#)

Overview

Product name	Glucose Assay Kit (Fluorometric, High Sensitivity)
Detection method	Fluorescent
Sample type	Serum, Plasma, Other biological fluids, Tissue, Adherent cells, Suspension cells, Tissue Culture Media
Assay type	Quantitative
Sensitivity	< 0.5 µM
Species reactivity	<b>Reacts with:</b> Mammals, Other species
Product overview	Picoprobe Glucose Assay Kit (ab169559) is a simple, rapid, ultra-sensitive assay to measure glucose. It is suitable for high-throughput use. In this assay, D-glucose is enzymatically oxidized to form a product which reacts with a colorless probe to generate the fluorescence (Ex/Em = 535/587 nm). The fluorescence generated is directly proportional to the amount of glucose. This assay kit can detect less than 0.5 µM glucose in various biological samples.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K688 PicoProbe™ Glucose Fluorometric Assay Kit. K688-100 is the same size as the 100 test size of ab169559.</p> <p>Glucose is the main energy source for virtually all living organisms. 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>
Platform	Microplate reader

Properties

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

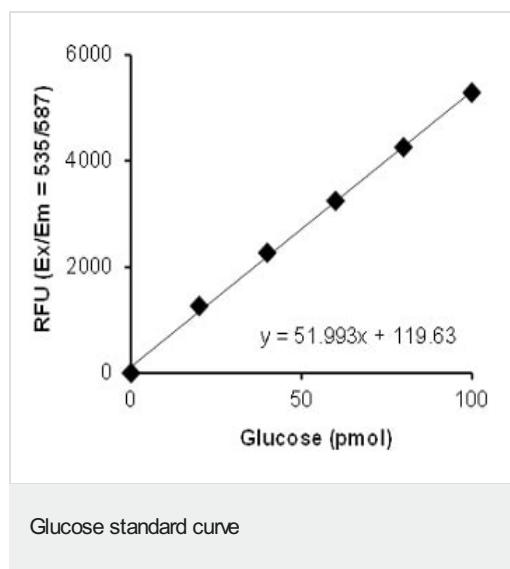
Components	100 tests
Assay Buffer II	1 x 25ml
Glucose Enzyme Mix	1 vial

Components	100 tests
Glucose Standard	1 x 100µl
PicoProbe I	1 x 0.4ml
Substrate Mix I	1 vial

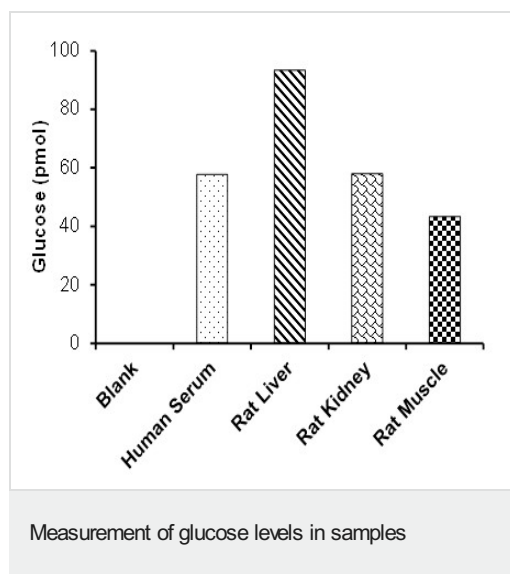
## Relevance

Glucose ( $C_6H_{12}O_6$ ; 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  $CO_2$  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.

## Images



Representative standard curve.



Measurement of glucose levels in human serum (1 µL of 1:10 diluted) and rat tissue lysates from liver, kidney and muscle (0.14 µg, 0.19 µg and 0.93 µg respectively).

**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.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