

Product datasheet

Glucose-6-Phosphate Assay Kit - High Sensitivity (Fluorometric) ab107923

[3 Images](#)

Overview

Product name	Glucose-6-Phosphate Assay Kit - High Sensitivity (Fluorometric)
Detection method	Fluorescent
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Quantitative
Sensitivity	> 1 μM
Range	1 μM - 500 μM
Assay time	0h 20m
Product overview	<p>Abcam's Glucose-6-Phosphate Assay Kit - High Sensitivity (Fluorometric) (ab107923) is a simple, sensitive and rapid means of quantifying glucose-6-Phosphate (G6P) in a variety of samples. In the assay, glucose-6-phosphate is oxidized with the generation of a product that converts a nearly colorless probe to an intensely fluorescent product (Ex/Em 535/587 nm). This assay can detect G6P in the range of 10 to 500 pmoles which is equivalent to the range of 1-500 μM in the original sample assuming a dilution of 5X during processing. This Glucose-6-Phosphate Assay Kit - High Sensitivity is more sensitive than our colorimetric assay, Glucose-6-Phosphate Assay Kit (ab83426).</p> <p>Visit our FAQs page for tips and troubleshooting.</p>
Notes	<p>Glucose-6-phosphate (G6P) is a key intermediate for glucose transport into cells which then enters either metabolic pathways or storage. G6P can enter the glycolytic pathway, the pentose phosphate shunt or be stored as glycogen or starch. G6P is utilized by its dehydrogenase to generate reducing equivalents in the form of NAD(P)H. This is particularly important in red blood cells where a G6PDH deficiency leads to hemolytic anemia.</p>
Platform	Microplate

Properties

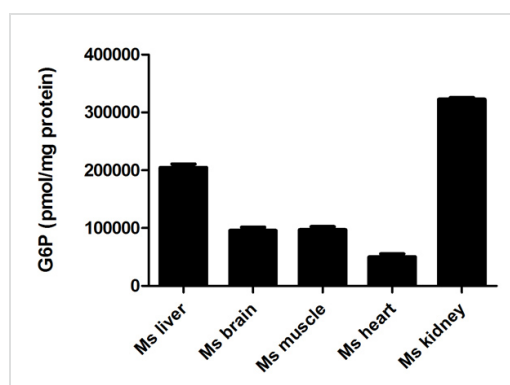
Storage instructions	Store at -20°C. Please refer to protocols.
Storage buffer	Preservative: None Constituents: DMSO

Components	Identifier	100 tests
Glucose 6 phosphate assay buffer	WM	1 x 25ml
Glucose 6 phosphate Enzyme Mix (lyophilized)	Green	1 vial
Glucose 6 phosphate Standard (10 μ mol; lyophilized)	Yellow	1 vial
Glucose 6 phosphate Substrate Mix (lyophilized)	Red	1 vial
Probe (in DMSO)	Blue	1 x 0.4ml

Relevance

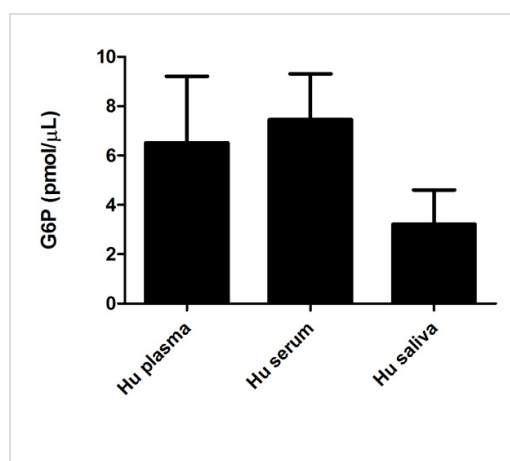
Glucose 6 phosphate (G6P) is a key sugar intermediate for glucose to get into cells, and then enter either metabolic pathways or storage. G6P can enter the glycolytic pathway, the pentose phosphate shunt or be stored as glycogen or starch. G6P is utilized by its dehydrogenase to generate reducing equivalents in the form of NADPH. This is particularly important in red blood cells where a G6PDH deficiency leads to hemolytic anemia.

Images



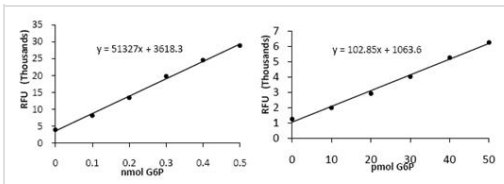
Glucose 6 phosphate measured in mouse tissue lysates showing quantity (pmol) per mg protein of tested sample

Functional Studies - Glucose 6 phosphate Assay Kit(ab107923)



Glucose 6 phosphate measured in biological fluids showing quantity (pmol) per microliter of tested sample

Functional Studies - Glucose 6 phosphate Assay Kit (ab107923)



Glucose 6 phosphate standard curves (0 – 500 and 0 – 50 pmol range) generated using ab107923.

Functional Studies - Glucose 6 phosphate Assay Kit
(ab107923)

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 <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