

# Amplite<sup>™</sup> Fluorimetric Glucose-6-Phosphate Assay Kit

Catalog number: 13804 Unit size: 200 Tests

Component	Storage	Amount
Component A: Enzyme Probe	Freeze (<-15 °C), Minimize light exposure	2 bottles (lyophilized powder)
Component B: Assay Buffer	Freeze (<-15 °C)	1 bottle (10 mL)
Component C: NADP	Freeze (<-15 °C), Minimize light exposure	1 vial
Component D: Glucose-6-Phosphate Standard	Freeze (<-15 °C), Minimize light exposure	3.04 mg/vial

# OVERVIEW

Glucose-6-phosphate (G6P) is a key intermediate for glucose transport into cells. G6P may also be converted to glycogen or starch for storage in the liver and muscles. G6P is utilized by glucose-6-phosphate dehydrogenase (G6PD) to generate the reducing equivalents in the form of NADPH. This is particularly important in red blood cells where G6PD deficiency leads to hemolytic anemia. AAT Bioquest's Amplite™ Fluorimetric Glucose-6-Phosphate Assay Kit provides a simple, sensitive and rapid fluorescence-based method for detecting G6P in biological samples such as serum, plasma, urine, as well as in cell culture samples. In the coupled enzyme assay, the G6P concentration is proportionally related to NADPH that is specifically monitored by a fluorescence microplate reader. With the Amplite™ G6P Assay Kit, we were able to detect as little as 0.3 µM G6P in a 100 µL reaction volume.

# AT A GLANCE

### **Protocol summary**

- 1. Prepare G6P working solution (50 μL)
- 2. Add G6P standards or test samples (50 µL)
- 3. Incubate at room temperature for 30 minutes 2 hours
- 4. Monitor fluorescence increase at Ex/Em = 540/590 nm (Cutoff = 570 nm)

**Important** Thaw kit components at room temperature before starting the experiment.

## **KEY PARAMETERS**

Instrument:	Fluorescence microplate reader
Excitation:	540 nm
Emission:	590 nm
Cutoff:	570 nm
Recommended plate:	Solid black
Instrument:	Absorbance microplate reader
Absorbance:	575/605 nm
Recommended plate:	Clear bottom

## PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

#### 1. NADP stock solution (100X):

Add 100  $\mu L$  of  $H_2O$  into the vial of NADP (Component C) to make 100X NADP stock solution.

#### 2. G6P standard solution (100 mM):

Add 100  $\mu$ L of H<sub>2</sub>O or 1x PBS buffer into the vial of G6P Standard (Component D) to make 100 mM G6P standard solution.

# PREPARATION OF STANDARD SOLUTION

#### G6P standard

For convenience, use the Serial Dilution Planner: https://www.aatbio.com/tools/serial-dilution/13804

Add 10  $\mu$ L of 100 mM G6P standard solution into 990  $\mu$ L 1x PBS buffer to generate 1 mM G6P standard solution. Then, add 100  $\mu$ L of 1 mM G6P standard solution into 900  $\mu$ L 1 x PBS buffer to make 100  $\mu$ M G6P standard solution (G6P7). Take 100  $\mu$ M G6P standard solution (G6P7) and perform 1:3 serial dilutions to get serially diluted G6P standards (G6P6 - G6P1) with 1x PBS buffer.

**Note** Diluted G6P standard solution is unstable, and should be used within 4 hours.

## PREPARATION OF WORKING SOLUTION

1. Add 5 mL of Assay Buffer (Component B) into one bottle of Enzyme Probe (Component A), and mix well.

2. Add 50  $\mu L$  of 100X NADP stock solution into the bottle of Component A+B, and mix well to make G6P working solution.

Note This G6P working solution is enough for one 96-well plate.

## SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of G6P standards and test samples in a solid black 96-well microplate. G6P=G6P Standards (G6P1 - G6P7, 0.14 to 100  $\mu$ M), BL=Blank Control, TS=Test Samples.

BL	BL	TS	TS
G6P1	G6P1		
G6P2	G6P2		
G6P3	G6P3		
G6P4	G6P4		
G6P5	G6P5		
G6P6	G6P6		
G6P7	G6P7		

Table 2. Reagent composition for each well.

Well	Volume	Reagent
G6P1 - G6P7	50 µL	Serial Dilution (0.14 to 100 µM)
BL	50 μL	Dilution Buffer
TS	50 µL	test sample

Tel: 408-733-1055 | Fax: 408-733-1304 | Email: support@aatbio.com

©2012 AAT Bioquest, Inc. Last revised October 2018. For more information and tools, please visit https://www.aatbio.com

- 1. Prepare G6P standards (G6P), blank controls (BL), and test samples (TS) according to the layout provided in Tables 1 and 2. For a 384-well plate, use 25  $\mu$ L of reagent per well instead of 50  $\mu$ L.
- 2. Add 50  $\mu$ L of GGP working solution to each well of GGP standard, blank control, and test samples to make the total GGP assay volume of 100  $\mu$ L/well. For a 384-well plate, add 25  $\mu$ L of GGP working solution into each well instead, for a total volume of 50  $\mu$ L/well.
- 3. Incubate the reaction at room temperature for 30 minutes to 2 hours, protected from light.
- Monitor the fluorescence increase with a fluorescence plate reader at Excitation = 530 - 570 nm, Emission = 590 - 600 nm (optimal Ex/Em = 540/590 nm, Cutoff = 570nm).

**Note** The contents of the plate can also be transferred to a white clear bottom plate and read by absorbance microplate reader at the ratio of  $A_{575nm}/A_{605nm}$ . The absorption detection has lower sensitivity compared to fluorescence reading.

# EXAMPLE DATA ANALYSIS AND FIGURES

The reading (RFU) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the baseline corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate Glucose-6-Phosphate samples. We recommend using the Online Linear Regression Calculator which can be found at:

https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-onlinecalculator

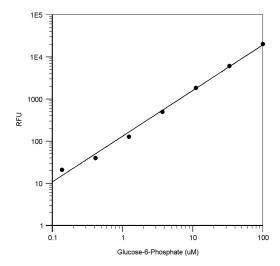


Figure 1. G6P dose response was measured with Amplite<sup>™</sup> Fluorimetric G6P Assay Kit in a 96-well solid black plate using a Gemini (Molecular Devices) microplate reader.

# DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.