# Amplite<sup>TM</sup> Fluorimetric Glucose Quantitation Kit

\*Red Fluorescence\*

Ordering InformationStorage ConditionsInstrument PlatformProduct Number: 40005 (500 assays)Keep at -20 °C and protect from lightFluorescence microplate readers

### Introduction

Glucose, a monosaccharide, is the most important carbohydrate in biology. It is a source of energy and metabolic intermediate for cell growth. As one of the main products of photosynthesis, glucose starts cellular respiration in both prokaryotes and eukaryotes. Glucose level is a key diagnostic parameter for many metabolic disorders, e.g., diabetes.

This Amplite<sup>TM</sup> Fluorimetric Glucose Quantification Kit provides a quick and sensitive method for the measurement of glucose. It uses glucose oxidase-based enzyme coupled reactions to detect glucose through the production of hydrogen peroxide, which is monitored by our Amplite<sup>TM</sup> Red peroxidase substrate. Amplite<sup>TM</sup> Red peroxidase substrate can be read by a fluorescence microplate reader at Ex/Em = 540/590 nm. The assay is robust, and can be readily adapted for a wide variety of applications that require the measurement of glucose. With the Amplite<sup>TM</sup> Fluorimetric Glucose Quantitation Kit, we can detect as little as  $0.1~\mu M$  D-glucose.

## **Kit Key Features**

Sensitive: Detect as low as 0.1 μM D-glucose in solution.

**Continuous:** Easily adapted to automation without a separation step.

**Convenient:** Formulated to have minimal hands-on time. No wash is required.

**Non-Radioactive:** No special requirements for waste treatment.

## Kit Components

Components	Amount		
Component A: Amplite <sup>TM</sup> Red (light-sensitive)	1 vial		
Component B: Assay Buffer	1 bottle (50 mL)		
Component C: Horseradish Peroxidase (HRP)	1 vial (10 units)		
Component D: Glucose Oxidase	1 vial (100 units)		
Component E: DMSO	1 vial (200 μL)		
Component F: Glucose	1 vial (144 mg)		

## **Assay Protocol for One 96-Well Plate**

## **Brief Summary**

Prepare assay reaction mixture (50  $\mu$ L)  $\rightarrow$  Add Glucose standards or test samples (50  $\mu$ L)  $\rightarrow$  Incubate at 37 °C for 10-30 minutes  $\rightarrow$  Monitor fluorescence intensity at Ex/Em = 540/590 nm

Note: Thaw all the kit components to room temperature before starting the experiment.

#### 1. Prepare stock solutions:

1.1 <u>250X Amplite<sup>™</sup> Red stock solution:</u> Add 100 μL of DMSO (Component E) into the vial of Amplite<sup>™</sup> Red substrate (Component A). The stock solution should be used promptly. Any remaining solution should be aliquoted and refrozen at -20 °C.

Note 1: Avoid repeated freeze-thaw cycles.

Note 2: The Amplite<sup>TM</sup> Red substrate is unstable in the presence of thiols such as dithiothreitol (DTT) and 2-mercaptoethanol. The final concentration of DTT or 2-mercaptoethanol in the reaction should be no higher than 10  $\mu$ M. The Amplite<sup>TM</sup> Red substrate is also unstable at high pH (> 8.5). Therefore, the reaction should be performed at pH 7–8. The provided assay buffer (pH 7.4) is recommended.

- 1.2 <u>10 U/mL HRP stock solution:</u> Add 1 mL of assay buffer (Component B) into the vial of horseradish peroxidase (Component C).
  - Note: The unused HRP solution should be divided into single use aliquots and stored at -20°C.
- 1.3 100 U/mL glucose oxidase solution: Add 1 mL of assay buffer (Component B) into the vial of glucose oxidase (Component D).
  - Note: The unused glucose oxidase solution should be divided into single use aliquots and stored at -20°C.
- 1.4 <u>800 mM glucose stock solution:</u> Add 1 mL of assay buffer (Component B) into the vial of glucose (Component F).

*Note: The unused glucose solution should be stored at -20 °C.* 

## 2. Prepare assay reaction mixture:

Prepare Assay reaction mixture according to the following tables, protected from light.

**Table 1** Assay reaction mixture for one 96-well plate (2X)

Components	Volume
250X Amplite™ Red Stock Solution (from Step 1.1)	20 μL
10 U/mL HRP Stock Solution (from Step 1.2)	100 μL
100 U/mL Glucose Oxidase Solution (from Step 1.3)	100 μL
Assay Buffer (Component B)	4.78 mL
Total volume	5 mL

**Table 2** Layout of glucose standards and test samples in a solid black 96-well microplate

BL	BL	TS	TS	 			
GS1	GS1			 			
GS2	GS2						
GS3	GS3						
GS4	GS4						
GS5	GS5						
GS5 GS6	GS6						
GS7	GS7						

*Note: GS= Glucose standards. BL=Blank control. TS=test samples.* 

**Table 3**. Reagent composition for each well

Glucose Standard	Blank Control	Test Sample	
Serial Dilutions*: 50 μL	Assay Buffer (Component B): 50 µL	50 μL	

\*Note 1: Add the serially diluted glucose standards from approximately 0.03  $\mu$ M to 30  $\mu$ M into each well from GS1 to GS7 in duplicate.

Note 2: High concentration of glucose (e.g.,  $100 \mu M$  in test sample or standard) may cause reduced fluorescence signal due to the overoxidation of Amplite<sup>TM</sup> red substrate (to a non-fluorescent product).

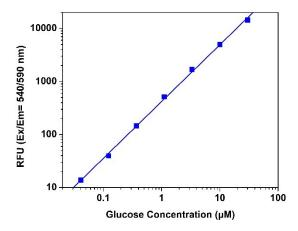
## 3. Run Glucose assay:

- 3.1 Prepare a glucose standard by diluting the appropriate amount of the 800 mM glucose stock solution (from Step 1.4) into assay buffer (Component B) to produce glucose concentrations of 30  $\mu$ M. Then perform 1:3 serial dilutions in assay buffer (Component B) to get approximately 10, 3, 1, 0.3, 0.1 and 0.03  $\mu$ M serially diluted glucose standards. A non-glucose buffer control is included as blank control.
- 3.2 Add 50  $\mu$ L of assay reaction mixture (from Step 2) into each well of glucose standard, blank control, and test samples (see Step 2, Table 3) to make the total glucose assay volume of 100  $\mu$ L/well *Note: For a 384-well plate, add 25 \muL of sample and 25 \muL of assay reaction mixture into each well.*
- 3.3 Incubate the reaction for 10 to 30 minutes at 37 °C, protected from light.
- 3.4 Monitor the fluorescence intensity with a fluorescence plate reader at Ex/Em = 530-570 nm/590-600 nm (optimal Ex/Em = 540/590 nm).

## **Data Analysis**

The fluorescence in blank wells (with the assay buffer only) is used as a control, and is subtracted from the values for those wells with glucose reactions. A glucose standard curve is shown in Figure 1.

Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.



**Figure 1**. Glucose dose response was measured with Amplite<sup>TM</sup> Fluorimetric Glucose Quantitation Kit (Cat #40005) on a 96-well black plate using a Gemini microplate reader (Molecular Devices). As low as 0.1 μM glucose was detected with 30 minutes incubation (n=3).

## **References**

- 1. Delva P, Degan M, Trettene M, Lechi A. (2006) Insulin and glucose mediate opposite intracellular ionized magnesium variations in human lymphocytes. J Endocrinol, 190, 711.
- 2. Delva P, Degan M, Pastori C, Faccini G, Lechi A. (2002) Glucose-induced alterations of intracellular ionized magnesium in human lymphocytes. Life Sci, 71, 2119.
- 3. Wang XT, Au SW, Lam VM, Engel PC. (2002) Recombinant human glucose-6-phosphate dehydrogenase. Evidence for a rapid-equilibrium random-order mechanism. Eur J Biochem, 269, 3417.
- 4. Leira F, Louzao MC, Vieites JM, Botana LM, Vieytes MR. (2002) Fluorescent microplate cell assay to measure uptake and metabolism of glucose in normal human lung fibroblasts. Toxicol In Vitro, 16, 267.

Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at info@aatbio.com if you have any questions.