

PhosphoWorks™ Colorimetric MESG Phosphate Assay Kit *UV Absorption*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 21659 (200 assays)	Keep in freezer and protect from light	Spectrophotometer Absorbance microplate reader

Introduction

Phosphate is involved in many biological processes. For example, phosphatases, ATPases and several other enzymes catalyze biochemical reactions in which inorganic phosphate (Pi) is released from a phosphoester substrate. The detection of many phosphoester-metabolizing enzymes is difficult because suitable substrates are not available. It is usually necessary to determine inorganic phosphate release using tedious colorimetric assays or radioisotope-based methods.

This PhosphoWorks™ Colorimetric MESG Phosphate Assay Kit has been developed for measuring the activity of any Pi-generating enzyme using MESG reagent. The measurement of Pi is based on absorbance change of MESG by phosphate. In the presence of inorganic phosphate, MESG is converted to 2-amino-6-mercapto-7-methylpurine by purine nucleoside phosphorylase with absorption wavelength shift to red. This feature has been used to develop our convenient MESG phosphate assay kit, an alternative to hazardous radioactive methods. The MESG substrate gives an absorbance increase at 360 nm on phosphorylation at pH 6.5-8.5. The assay is shown to quantitate phosphate at the final concentration as low as 0.2 μM. The kit has been used for monitoring ATPase activities. It can also be used for monitoring phosphatase activities.

Kit Components

Components	Amount
Component A: Assay Buffer	1 Bottle (10 mL)
Component B: MESG Substrate	1 vial (lyophilized powder)
Component C: Purine Nucleoside Phosphorylase (PNP)	1 vial (lyophilized powder)
Component D: 1 mM KH ₂ PO ₄	1 vial (1 mL)

Assay Protocol for One 96-Well Plate

Brief Summary

Prepare 50 μL of test samples and/or phosphate standards → Add 50 μL of assay solution → Incubate at room temperature for 30 minutes → Monitor absorbance at 360 nm

1. Prepare assay reagents:

- 1.1 Thaw all the four components at room temperature before use.
- 1.2 Prepare MESG Substrate (Component B) Solution: Add 500 μL of ddH₂O to the vial of MESG Substrate (Component B). Mix well by vortexing to get MESG Substrate Solution.
Note: 250 ul is enough for one plate make single used aliquos and store it at -20°C Immediately.,
- 1.3 Prepare Purine Nucleoside Phosphorylase (Component C) Solution: Add 100 μL of ddH₂O to the vial of Purine Nucleoside Phosphorylase (PNP; Component C). Mix well by vortexing to get Purine Nucleoside Phosphorylase Solution.
- 1.4 Prepare Assay Solution: Add the whole volume of MESG Substrate Solution (from Step 1.2) and Purine Nucleoside Phosphorylase Solution (from Step 1.3) into the bottle of Assay Buffer (Component A), mix well to get the assay solution. Place the assay solution on ice.
Note 1: This Assay Solution is stable for at least 4 hours on ice. It is not recommended to freeze the assay solution for another assay.
Note 2: To achieve the desirable results, UV-transparent plates or cuvettes are required.
Note 3: Due to the high sensitivity of this assay to Pi, it is extremely important to use Pi-free laboratory ware and

2. Prepare serially diluted phosphate standards and/or test samples:

- 2.1 Prepare Phosphate Standard: Add 50 μL of 1 mM KH₂PO₄ (Component D) into 950 μL of deionized water or enzyme reaction buffer to give 50 μM phosphate standard solution.
- 2.2 Take 200 μL of 50 μM phosphate standard solution to perform 1:2 serial dilutions to give 25, 12.5, 6.25, 3.125, 1.56, and 0.78 μM serially diluted phosphate standards.
- 2.3 Add phosphate-containing test samples and/or phosphate standards into a clear UV-transparent 96-well microplate according to Tables 1 and 2.

