

PhosphoWorks™ Fluorimetric Pyrophosphate Assay Kit *Blue Fluorescence*

Catalog number: 21611
Unit size: 200 Tests

Component	Storage	Amount (Cat No. 21611)
Component A: Assay Buffer	Freeze (< -15 °C)	1 bottle (25 mL)
Component B: PPI Sensor	Freeze (< -15 °C), Minimize light exposure	1 vial (lyophilized powder)
Component C: Pyrophosphate Standard	Freeze (< -15 °C), Minimize light exposure	1 vial (1 mL, 50 mM)
Component D: DMSO	Freeze (< -15 °C)	1 vial (200 µL)

OVERVIEW

Pyrophosphate (PPI) is produced by a number of biochemical reactions, such as ATP hydrolysis, DNA and RNA polymerizations, cyclic AMP formation by the enzyme adenylate cyclase and the enzymatic activation of fatty acids to form their coenzyme A esters. Our PhosphoWorks™ Pyrophosphate Assay Kit provides the most robust spectrophotometric method for measuring pyrophosphate. This kit uses our proprietary fluorogenic pyrophosphate sensor that has its fluorescence intensity proportionally dependent upon the concentration of pyrophosphate. Our assay is much easier and more robust than the enzyme-coupling pyrophosphate methods that require at least two enzymes for their pyrophosphate detections. The kit provides all the essential components for assaying pyrophosphate. This kit has been successfully used in high throughput screening (HTS). Please inquire special HTS bulk package discount for the screening of >10,000 assays.

AT A GLANCE

Protocol Summary

1. Prepare Pyrophosphate standards and/or test samples (50 µL)
2. Add Pyrophosphate working solution (50 µL)
3. Incubate at room temperature for 10 to 30 minutes
4. Monitor fluorescence intensity at Ex/Em = 316/456 nm (Cutoff = 420 nm)

Important Note

Thaw all the four components at room temperature before use.

KEY PARAMETERS

Fluorescence microplate reader

Cutoff	420 nm
Emission	456 nm
Excitation	316 nm
Recommended plate	Solid black

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

PPI Sensor stock solution (200X)

Add 50 µL of DMSO (Component D) into the vial of PPI Sensor (Component B) to make 200X PPI Sensor stock solution. Protect from light.

Note 25 µL of the PPI Sensor Stock Solution is enough for one 96-well plate.

Pyrophosphate standard solution (1 mM)

Add 10 µL of 50 mM Pyrophosphate Standard (Component C) into 490 µL of ddH₂O or 50 mM Hepes buffer (pH 7) to make 1 mM Pyrophosphate standard solution.

PREPARATION OF STANDARD SOLUTIONS

For convenience, use the Serial Dilution Planner:
<https://www.aatbio.com/tools/serial-dilution/21611>

Pyrophosphate standard

Add 50 µL of 1 mM Pyrophosphate standard solution into 450 µL of ddH₂O or 50 mM Hepes buffer to get 100 µM Pyrophosphate standard solution (PS7). Take 100 µM Pyrophosphate standard solution and perform 1:3 serial dilutions in ddH₂O or 50 mM Hepes buffer to get serially diluted Pyrophosphate standards (PS6 - PS1).

PREPARATION OF WORKING SOLUTION

Add 25 µL of 200X PPI Sensor stock solution to 5 mL of Assay Buffer (Component A) and mix well to make PPI working solution.

Note Due to the high sensitivity of this assay to PPI, it is important to use PPI-free labware and reagents. DTT ≥ 1 mM will increase the background, MgCl₂ ≥ 2 mM will decrease the response.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of Pyrophosphate standards and test samples in a solid black 96-well microplate. PS = Pyrophosphate Standard (PS1 - PS7, 0.14 to 100 µM), BL = Blank Control, TS = Test Sample.

BL	BL	TS	TS
PS1	PS1
PS2	PS2
PS3	PS3		
PS4	PS4		
PS5	PS5		
PS6	PS6		
PS7	PS7		

Table 2. Reagent composition for each well.

Well	Volume	Reagent
PS1 - PS7	50 µL	Serial Dilutions (0.14 to 100 µM)
BL	50 µL	Assay Buffer

TS	50 μ L	test sample
----	------------	-------------

1. Prepare Pyrophosphate standards (PS), blank controls (BL), and test samples (TS) according to the layout provided in Tables 1 and 2. For a 384-well plate, use 25 μ L of reagent per well instead of 50 μ L.
2. Add 50 μ L of PPi working solution to each well of Pyrophosphate standard, blank control, and test samples to make the total Pyrophosphate assay volume of 100 μ L/well. For a 384-well plate, add 25 μ L of PPi working solution into each well instead, for a total volume of 50 μ L/well. Mix the reagents thoroughly.
3. Incubate at room temperature for 10 to 30 minutes.
4. Monitor the fluorescence increase with a fluorescence plate reader at Ex/Em = 316/456 nm (Cutoff = 420 nm).

EXAMPLE DATA ANALYSIS AND FIGURES

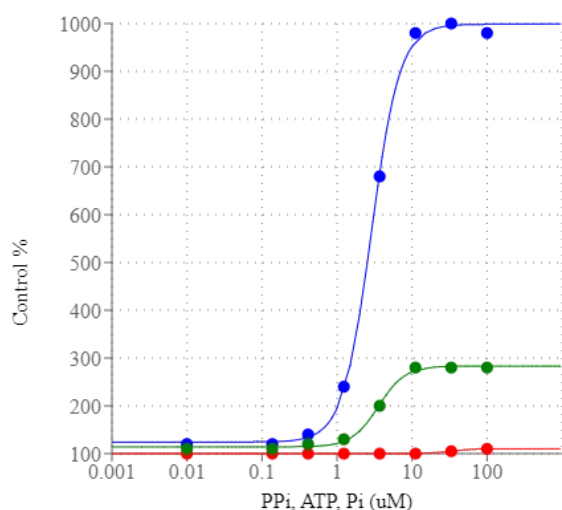


Figure 1. Pyrophosphate, ATP and phosphate dose responses were measured with PhosphoWorks™ Fluoremetric Pyrophosphate Assay Kit in a solid black 96-well plate using a fluorescence microplate reader.

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.