
Product Manual

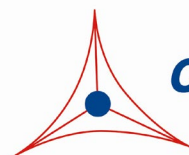
Adenosine Monophosphate Assay Kit

Catalog Number

MET-5160

100 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Adenosine monophosphate (AMP) is a nucleotide made up of a phosphate group, the sugar ribose, and the nucleobase adenine. AMP is an important part of various metabolic reactions since AMP is often converted to adenosine diphosphate (ADP) or adenosine triphosphate (ATP). AMP is a constituent of RNA strand synthesis. AMP regulates AMP-activated protein kinase (AMPK), an enzyme that controls cell migration, calcium signaling, and the secretion of cytokines, and t-cell priming capacity in immune cells such as dendritic cells.

Cell Biolabs' Adenosine Monophosphate Assay Kit is a simple fluorometric assay that measures the amount of adenosine monophosphate present in biological samples in a 96-well microtiter plate format. Each kit provides sufficient reagents to perform up to 100 assays*, including blanks, adenosine monophosphate standards, and unknown samples. Sample adenosine monophosphate concentrations are determined by comparison with a known adenosine monophosphate standard. The kit has a detection sensitivity limit of 1.6 μ M adenosine monophosphate.

****Note: Each sample replicate requires 2 assays, one treated with pyruvate phosphate dikinase (+PPDK) and one without (-PPDK). Adenosine monophosphate is calculated from the difference in RFU readings from the 2 wells.***

Assay Principle

Cell Biolabs' Adenosine Monophosphate Assay Kit measures total adenosine monophosphate within biological samples. Adenosine monophosphate (with pyrophosphate and phosphoenolpyruvate) is converted by PPDK to pyruvate + adenosine triphosphate (ATP) and phosphate. Pyruvate is converted by pyruvate oxidase in the presence of phosphate and oxygen into acetyl phosphate, carbon dioxide, and hydrogen peroxide. The resulting hydrogen peroxide is then detected with a highly specific fluorometric probe. Horseradish peroxidase catalyzes the reaction between the probe and hydrogen peroxide, which bind in a 1:1 ratio. Samples are compared to a known concentration of adenosine monophosphate standard within the 96-well microtiter plate format. Samples and standards are incubated for 30 minutes and then read with a standard 96-well fluorometric plate reader (Figure 1).

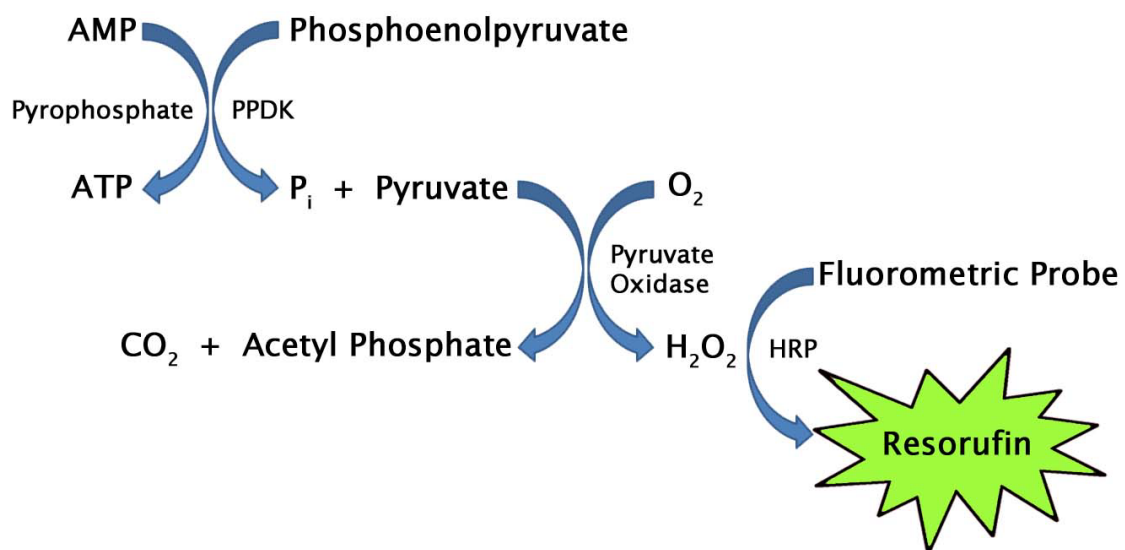


Figure 1. Adenosine monophosphate Assay Principle

Related Products

1. MET-5159: Pyrophosphate Assay Kit
2. MET-5029: Pyruvate Assay Kit (Fluorometric)
3. MET-5125: Pyruvate Assay Kit (Colorimetric)
4. MET-5158: Methionine Assay Kit
5. MET-5152: s-Adenosylmethionine ELISA Kit
6. MET-5151: s-Adenosylhomocysteine ELISA Kit
7. MET-5001: Lactose Assay Kit
8. MET-5013: Lactate Assay Kit (Fluorometric)
9. MET-5023: Glycogen Assay Kit (Fluorometric)
10. STA-399: Free Glycerol Assay Kit (Fluorometric)
11. STA-674: Glutamate Assay Kit
12. STA-675: Hydroxyproline Assay Kit
13. STA-681: Glucose Assay Kit (Fluorometric)
14. STA-682: Total Carbohydrate Assay Kit

Kit Components

Box 1 (shipped on blue ice packs)

1. AMP Standard (Part No. 51601C): One 50 μ L tube of adenosine monophosphate at 10 mM.
2. Pyrophosphate (Part No. 51602C): One 50 μ L tube of pyrophosphate at 20 mM.
3. PEP (Part No. 51585C): One 50 μ L tube of phosphoenolpyruvate (PEP) at 100 mM.
4. Na₂HPO₄ (Part No. 51586A): One 50 μ L tube of sodium phosphate dibasic (Na₂HPO₄) at 100 mM.
5. Fluorometric Probe (Part No. 50231C): One 50 μ L tube in DMSO.
6. HRP (Part No. 234402-T): One 10 μ L tube of a 100 U/mL solution in glycerol.
7. FAD (Part No. 50293C): One 50 μ L tube of 2 mM Flavin Adenine Dinucleotide (FAD).
8. TPP (Part No. 50294C): One 50 μ L tube of 2 mM Thiamine Pyrophosphate (TPP).
9. Pyruvate Oxidase (Part No. 50295C): One 300 μ L tube.
10. PPDK (Part No. 51588D): One 200 μ L tube of phosphate pyruvate dikinase (PPDK).

Box 2 (shipped on blue ice packs)

1. 10X Assay Buffer (Part No. 51582A): One 25 mL bottle.

Materials Not Supplied

1. Distilled or deionized water
2. 10 kDa molecular weight cutoff (MWCO) centrifuge spin filter (e.g., Amicon Ultra 0.5mL)
3. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
4. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
5. Standard 96-well fluorescence black microtiter plate and/or black cell culture microplate
6. Multichannel micropipette reservoir
7. Fluorescence microplate reader capable of reading excitation in the 530-570 nm range and emission in the 590-600 nm range.

Storage

Upon receipt, store the 10X Assay Buffer at room temperature. Store the PPDK at -80°C. Store all other components at -20°C. The Fluorometric Probe is light sensitive and must be stored accordingly. Avoid multiple freeze/thaw cycles.

Preparation of Reagents

- 1X Assay Buffer: Dilute the 10X Assay Buffer 1:10 with deionized water for a 1X solution. Stir or vortex to homogeneity. Store at 4°C.

- Reaction Mix and Negative Control Mix: Prepare two separate mixtures according to the table below. The PPDK is omitted from the Negative Control Mix.

Component	Reaction Mix (20 assays)	Negative Control Mix (20 assays)
PPDK	40 μ L	-----
Pyrophosphate	10 μ L	10 μ L
PEP	10 μ L	10 μ L
HRP	2 μ L	2 μ L
Pyruvate Oxidase	60 μ L	60 μ L
FAD	10 μ L	10 μ L
TPP	10 μ L	10 μ L
Na ₂ HPO ₄	10 μ L	10 μ L
Fluorometric Probe	10 μ L	10 μ L
1X Assay Buffer	838 μ L	878 μ L
Total	1000 μL	1000 μL

Note: Prepare only enough for immediate use and scale proportionally as needed.

Preparation of Samples

- Cell culture supernatants: Cell culture media formulated with pyruvate should be avoided. To remove insoluble particles, centrifuge at 10,000 x g for 5 min. Collect the supernatant and filter the solution with a 10kDa spin filter to deproteinate the sample. Collect flow through. The flow through may be assayed directly or diluted as necessary into PBS.

Note: Maintain pH between 7 and 8 for optimal working conditions as the Fluorometric Probe is unstable at high pH (>8.5).

- Tissue lysates: Sonicate or homogenize tissue sample in PBS and centrifuge at 10,000 x g for 10 minutes at 4°C. Collect the supernatant and filter the solution with a 10kDa spin filter to deproteinate the sample. Collect flow through. The flow through may be assayed directly or diluted as necessary into PBS.
- Cell lysates: Resuspend cells in PBS. Homogenize or sonicate the cells on ice. Centrifuge to remove debris. Collect the supernatant and filter the solution with a 10kDa spin filter to deproteinate the sample. Collect flow through. The flow through may be assayed undiluted or diluted as necessary into PBS.
- Serum, plasma, saliva, or urine: To remove insoluble particles, centrifuge at 10,000 x g for 5 min. Collect the supernatant and filter the solution with a 10kDa spin filter to deproteinate the sample. Collect flow through. The flow through may be assayed directly or diluted as necessary into PBS.

Notes:

- All samples should be assayed immediately or stored at -80°C for up to 1-2 months. Run proper controls as necessary. Optimal experimental conditions for samples must be determined by the investigator. Always run a standard curve with samples.
- Samples with NADH concentrations above $10\ \mu\text{M}$ and glutathione concentrations above $50\ \mu\text{M}$ will oxidize the Fluorometric Probe and could result in erroneous readings. To minimize this interference, it is recommended that superoxide dismutase (SOD) be added to the reaction at a final concentration of $40\ \text{U/mL}$ (Votyakova and Reynolds, Ref. 2).
- Avoid samples containing DTT or β -mercaptoethanol since the Fluorometric Probe is not stable in the presence of thiols (above $10\ \mu\text{M}$).

Preparation of Standard Curve

Prepare fresh Adenosine Monophosphate standards by diluting into 1X Assay Buffer according to Table 1 below.

Standard Tubes	10 mM Adenosine monophosphate Solution (μL)	1X Assay Buffer (μL)	Adenosine monophosphate (μM)
1	5	495	100
2	250 of Tube #1	250	50
3	250 of Tube #2	250	25
4	250 of Tube #3	250	12.5
5	250 of Tube #4	250	6.25
6	250 of Tube #5	250	3.125
7	250 of Tube #5	250	1.56
8	0	250	0

Table 1. Preparation of Adenosine monophosphate Standards.

Assay Protocol

1. Prepare and mix all reagents thoroughly before use. Each sample, including unknowns and standards, should be assayed in duplicate or triplicate.
Note: Each unknown sample replicate requires two paired wells, one to be treated with PPDK (+PPDK) and one without the enzyme (-PPDK) to measure endogenous.
2. Add $50\ \mu\text{L}$ of each adenosine monophosphate standard or unknown sample into wells of a 96-well microtiter plate.
3. Add $50\ \mu\text{L}$ of Reaction Mix to the standards and to one half of the paired sample wells, and mix the well contents thoroughly.
4. Add $50\ \mu\text{L}$ of Negative Control Mix to the other half of the paired sample wells.
5. Mix the well contents thoroughly and incubate for 30 minutes at 37°C protected from light.
Note: This assay is continuous (not terminated) and therefore may be measured at multiple time points to follow the reaction kinetics.

6. Read the plate with a fluorescence microplate reader equipped for excitation in the 530-570 nm range and for emission in the 590-600 nm range.

Calculation of Results

1. Determine the average Relative Fluorescence Unit (RFU) values for each sample, control, and standard.
2. Subtract the average zero standard value from itself and all standard values.
3. Graph the standard curve (see Figure 2).
4. Subtract the sample well values without PPDK (-PPDK) from the sample well values containing PPDK (+PPDK) to obtain the difference. The fluorescence difference is due to the PPDK activity.

$$\text{Net RFU} = (\text{RFU}_{+\text{PPDK}}) - (\text{RFU}_{-\text{PPDK}})$$

5. Compare the net RFU of each sample to the standard curve to determine and extrapolate the quantity of AMP present in the sample. Only use values within the range of the standard curve.

Example of Results

The following figures demonstrate typical Adenosine Monophosphate Assay Kit results. One should use the data below for reference only. This data should not be used to interpret or calculate actual sample results.

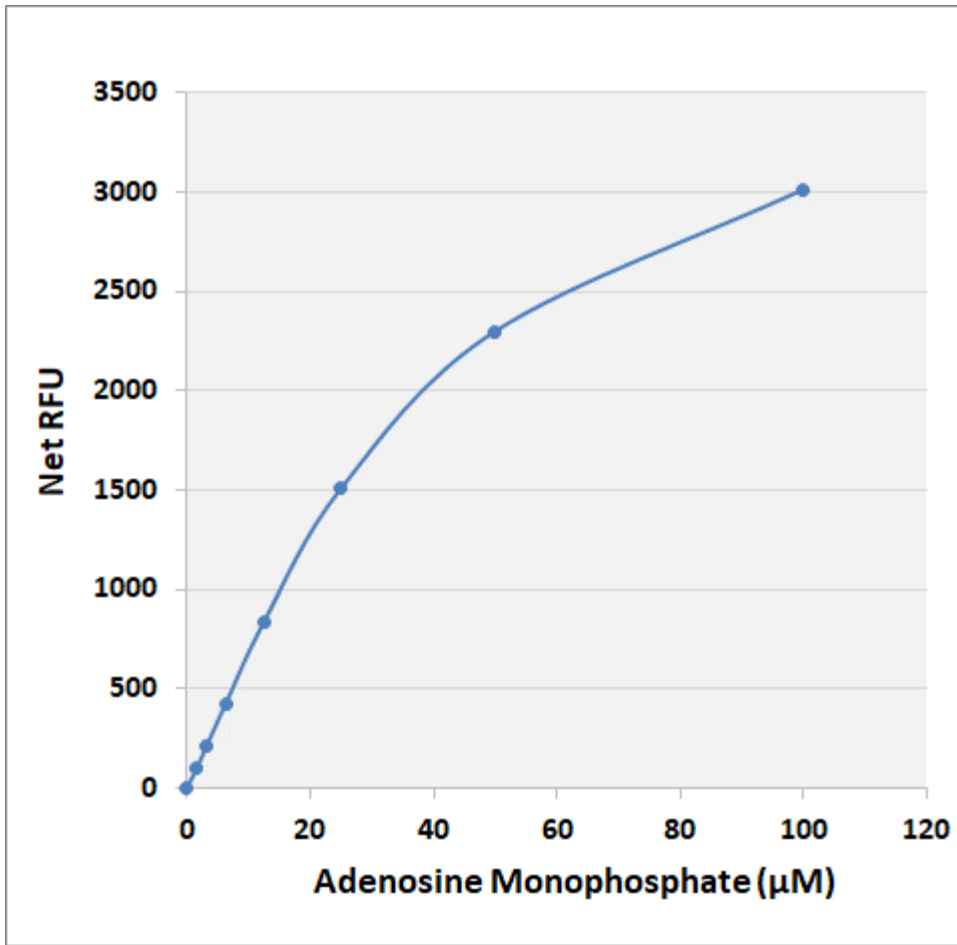


Figure 2: Adenosine monophosphate Standard Curve.

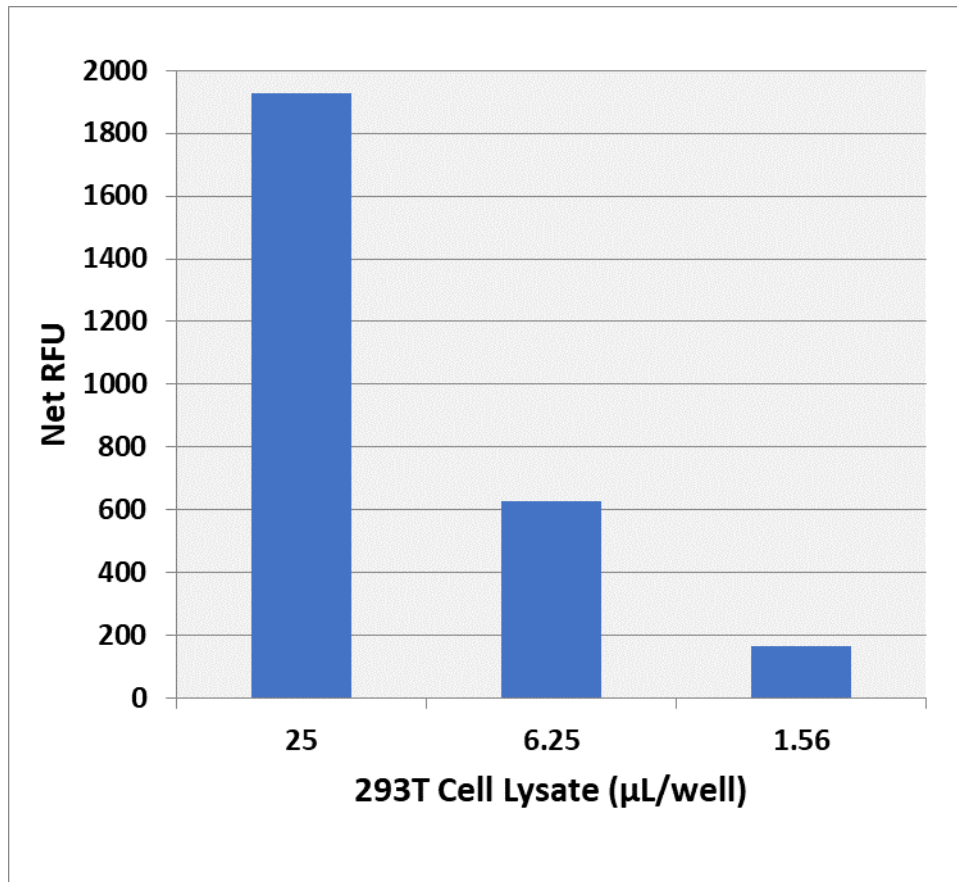


Figure 3: Adenosine monophosphate Detection in 293T cells using the Adenosine monophosphate Assay Kit. 293T cells were homogenized and deproteinated according to the preparation of samples section above.

References

1. Jauker M, Griesser H, and Richert C (2015). *Angewandte Chemie*. **54**: 14564–14569
2. Votyakova TV, and Reynolds IJ (2001) *Neurochem*. **79**:266.
3. Bressen C and Saghatelian A (2020) *Autophagy*. **17**: 828-829.
4. Long YC and Zierath JR (2006) *J. Clin Invest*. **116**:1776-1783.
5. Steinberg GR, Watt MJ, and Febbraio MA (2009) *Front Biosci (Landmark Ed)* **14**:1902-1916.
6. Wculek SK, Khouili SC, Priego E, Heras-Murillo I, and Sancho D (2019) *Front Immunol* **10**:775.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS’s sole obligation and purchaser’s exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In

no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858 271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2022: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.