

---

Product Manual

# Total Phosphatidic Acid Assay Kit

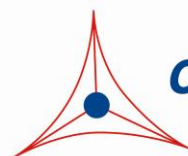
Catalog Number

MET-5019

100 assays

**FOR RESEARCH USE ONLY**  
**Not for use in diagnostic procedures**

---



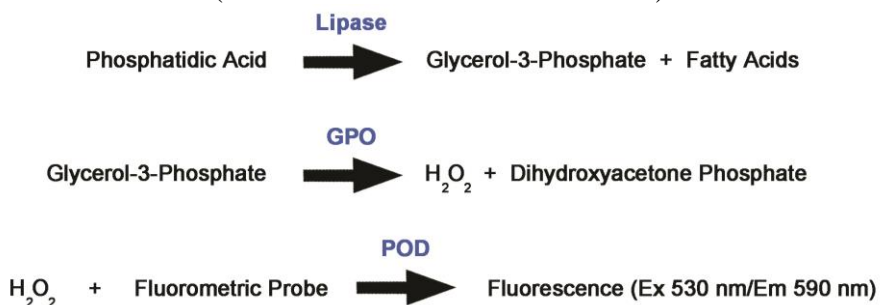
**CELL BIOLABS, INC.**

*Creating Solutions for Life Science Research*

## **Introduction**

Phosphatidic Acid (PA) is a critical precursor for the biosynthesis of many lipids in the cell. PA plays a critical role in membrane structure and acts as a signaling lipid, recruiting cytosolic proteins to the cell membrane. Within the cell, PA concentrations are maintained at extremely low levels by the activity of potent phospholipid phosphatases, converting phosphatidic acid to diacylglycerol (DAG). Because DAG is another important lipid precursor, it too is rapidly metabolized into other membrane lipid components (e.g., phosphatidylcholine, phosphatidylserine, phosphatidylethanolamine). Ultimately, measurement of cellular PA is useful for monitoring lipid synthesis and metabolism.

Cell Biolabs' Total Phosphatidic Acid Assay Kit measures total phosphatidic acid content, including lysophosphatidic acid (LPA), in samples by a coupled enzymatic reaction system. First, lipase is used to hydrolyze phosphatidic acid in samples to glycerol-3-phosphate. Next, the glycerol-3-phosphate product is oxidized by glycerol-3-phosphate oxidase (GPO), producing hydrogen peroxide which reacts with the kit's Fluorometric Probe (Ex. 530-560 nm/Em. 585-595 nm).



The Total Phosphatidic Acid Assay Kit is a simple, fluorometric assay that quantitatively measures total PA (PA and LPA) in samples using a 96-well microtiter plate format. Each kit provides sufficient reagents to perform up to 100 assays, including blanks, standards and unknown samples. The kit contains an L- $\alpha$ -Phosphatidic Acid Standard and has a detection sensitivity limit of ~5  $\mu\text{M}$ .

*Note: This kit is not intended for urine, plasma, or serum samples.*

## **Related Products**

1. STA-397: Serum Triglyceride Quantification Kit (Fluorometric)
2. STA-399: Free Glycerol Assay Kit (Fluorometric)
3. STA-600: Phosphatidylcholine Assay Kit
4. STA-619: Free Fatty Acid Assay Kit (Fluorometric)
5. MET-5129- Lysine Assay Kit (Fluorometric)

## **Kit Components**

1. Phosphatidic Acid Standard (Part No. 50191D): One 200  $\mu\text{L}$  vial of 1 mM L- $\alpha$ -Phosphatidic Acid.
2. 10X Assay Buffer (Part No. 50192D): One 1.5 mL vial.
3. Lipase Solution (Part No. 50193D): Three 1.4 mL vials.
4. Enzyme Mixture (Part No. 50194D): Three 1.75 mL vials.
5. Fluorometric Probe (Part No. 261901): One 110  $\mu\text{L}$  amber vial.

## **Materials Not Supplied**

1. Standard 96-well fluorescence black microtiter plate
2. PBS (containing Magnesium and Calcium)
3. Extraction reagents for cellular lipids (methanol, chloroform, 1M NaCl)

## **Storage**

Store the entire kit at -80°C. Avoid multiple freeze/thaws by aliquoting. The Fluorometric Probe is light sensitive and should be maintained in amber tubes.

## **Preparation of Reagents**

- Phosphatidic Acid Standard: Thaw at room temperature. Once homogeneous and mixed well, maintain the standard at room temperature during assay preparation. The solution is stable for 1 week at room temperature. For longer term storage, the solution should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- 1X Assay Buffer: 10X Assay Buffer should be thawed/maintained at 4°C during assay preparation. Dilute the 10X Assay Buffer with deionized water. Stir to homogeneity. The 1X solution is stable for 1 month at 4°C. For longer term storage, any unused 10X stock material should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- Lipase Solution and Enzyme Mixture: Thaw at 4°C. Once homogeneous and mixed well, maintain the solution at 4°C during assay preparation. The solution is stable for 1 week at 4°C. For longer term storage, the solution should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.  
*Note: These components are provided in multiple tubes to minimize multiple freeze/thaws.*
- Fluorometric Probe: Thaw and maintain at room temperature during assay preparation. Any unused material should be aliquoted and frozen at -80°C to avoid multiple freeze/thaws.
- PEU (pre-equilibrated upper phase) Solution: Mix 50 mL of chloroform, 50 mL of methanol, and 45 mL of 1M NaCl in a glass container. Shake or mix the solution well, then allow it to separate into 2 phases. Use the upper phase for washing during the extraction.

## **Preparation of Phosphatidic Acid Standard**

Thaw the Phosphatidic Acid Standard at room temperature. Mix well by vortexing to ensure the solution is homogeneous. Freshly prepare a dilution series of standard in the concentration range of 0  $\mu$ M – 250  $\mu$ M by diluting the standard stock solution (provided at 1 mM) in 1X Assay Buffer (see Table 1). Standards should be prepared fresh, vortexed well and used immediately.

<b>Standard Tubes</b>	<b>1 mM Phosphatidic Acid Standard (<math>\mu</math>L)</b>	<b>1X Assay Buffer (<math>\mu</math>L)</b>	<b>Final Phosphatidic Acid Standard (<math>\mu</math>M)</b>
1	25	75	250
2	50 of Tube #1	50	125
3	50 of Tube #2	50	62.5
4	50 of Tube #3	50	31.2
5	50 of Tube #4	50	15.6
6	50 of Tube #5	50	7.8
7	50 of Tube #6	50	3.9
8	0	50	0

**Table 1. Preparation of Phosphatidic Acid Standards**

## **Preparation of Samples**

- Urine, plasma and serum: This kit is not recommended for these samples.
- Cell Lysates:  
For adherent cells, remove media and wash cells twice with cold PBS. Harvest  $\sim 1 \times 10^7$  cells by using a rubber policeman. Do not use proteolytic enzymes. Centrifuge at 1500 x g for 10 minutes. Carefully remove the supernatant and resuspend in 1 mL of cold PBS. Proceed to step 1 of the extraction procedure below.  
For suspension cells, collect  $\sim 1 \times 10^7$  cells by centrifugation at 1500 x g for 10 minutes. Carefully remove the supernatant and wash the cell pellet with cold PBS. Repeat PBS wash/centrifugation once more. Carefully discard the supernatant and resuspend in 1 mL of cold PBS. Proceed to step 1 of the extraction procedure below:

### **Extraction Procedure**

1. Sonicate the 1 mL of cell suspension on ice.
2. Add 1.5 mL of methanol to the sonicated sample.
3. Add 2.25 mL of 1 M NaCl and 2.5 mL of chloroform to the sample. Vortex well.
4. Centrifuge at 1500 x g for 10 minutes at 4°C to separate the phases.
5. Carefully remove the upper aqueous phase and discard.
6. Wash the lower chloroform phase 2 times with 2 mL of pre-equilibrated upper phase (PEU) (see Preparation of Reagents Section). Separate the phases each time by centrifuging at 1500 x g for 10 minutes at 4°C. Carefully remove the upper phase and discard each time.
7. After the final wash, carefully transfer the lower organic phase to a glass vial or tube (a syringe works well). Avoid transferring any remaining upper, aqueous phase.
8. Dry the lower phase in a speedvac or under a gentle stream of nitrogen.
9. Resuspend the dried sample with 50  $\mu$ L of 1X Assay Buffer (1:20 of the original volume). Samples may be stored at -80°C for up to a month.

## **Assay Protocol**

*Note: Each PA standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed. Maintain the Lipase Solution and Enzyme Mixture at 4°C during assay preparation.*

1. Add 10  $\mu$ L of the PA standards, samples or blanks to the 96-well microtiter plate.
2. Add 40  $\mu$ L of Lipase Solution to each well.
3. Incubate at 37°C for 30 minutes.
4. During the step 3 incubation, separately prepare the desired volume of Detection Enzyme Mixture according to Table 2, based on the number of tests to be performed. Maintaining all components and mixtures at 4°C throughout this step, add components in the following sequence:
  - a. In a tube, add the appropriate volume of Enzyme Mixture.
  - b. To the Enzyme Mixture, add the corresponding volume of Fluorometric Probe. Mix well and immediately use.

Note: Detection Enzyme Mixture may appear slightly pink in color. This is normal background and should be subtracted from all absorbance values (see step 9 for calculation).

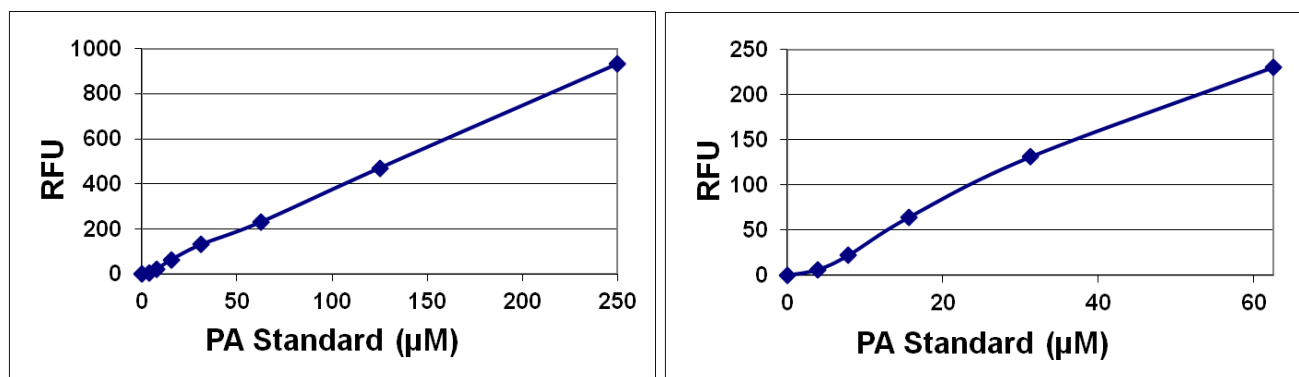
Enzyme Mixture (mL)	Fluorometric Probe ( $\mu$ L)	Total Volume of Detection Enzyme Mixture (mL)	# of Tests in 96-well Plate (100 $\mu$ L/test)
5	50	5.05	100
2.5	25	2.525	50
1.25	13	1.263	25

**Table 2. Preparation of Detection Enzyme Mixture**

5. Transfer 50  $\mu$ L of the above Detection Enzyme Mixture (from step 4) to each well.
6. Cover the plate wells to protect the reaction from light.
7. Incubate at room temperature for 10 minutes.
8. Read the plate with a fluorescence microplate reader equipped for excitation in the 530-560 nm range and for emission in the 585-595 nm range.
9. Calculate the concentration of phosphatidic acid within samples by comparing the sample fluorescence to the standard curve. Negative controls (without PA) should be subtracted from each PA sample and standard.

### Example of Results

The following figures demonstrate typical Total Phosphatidic Acid Assay Kit results. One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1: Total Phosphatidic Acid Assay Standard Curve.** PA standard curve was performed according to the Assay Protocol. Background has been subtracted.

### References

1. Wang, X. et al. (2006) *Prog. Lipid Res.* **45**, 250-278.
2. Merida, I. et al. (2008) *Biochem. J.* **409**, 1-18.
3. Jenkins, G.M. and Frohman, M.A. (2005) *Cell Mol. Life Sci.* **62**, 2305-2316.
4. Foster, D.A. and Xu, L. (2003). *Mol. Cancer Res.* **1**, 789-800.
5. Fang, Y. et al. (2001) *Science* **294**, 1942-1945.
6. Avila-Flores, A. et al. (2005). *J. Biol. Chem.* **280**, 10091-10099.
7. Schmelzle, T. and Hall, M.N. (2000) *Cell* **103**, 253-262.
8. Lim, H.K. et al. (2003) *J. Biol. Chem.* **278**, 45117-45127.

## **Recent Product Citations**

1. Liang, B. et al. (2021). TBX3 functions as a tumor suppressor downstream of activated CTNNB1 mutants during hepatocarcinogenesis. *J Hepatol*. doi: 10.1016/j.jhep.2021.01.044.
2. Uji, T. et al. (2022). Phospholipase D activation is required for 1-aminocyclopropane 1-carboxylic acid signaling during sexual reproduction in the marine red alga *Neopyropia yezoensis* (Rhodophyta). *BMC Plant Biol*. **22**(1):181. doi: 10.1186/s12870-022-03575-z.
3. Nath, A.S. et al. (2022). Modulation of the cell membrane lipid milieu by peroxisomal  $\beta$ -oxidation induces Rho1 signaling to trigger inflammatory responses. *Cell Rep*. **38**(9):110433. doi: 10.1016/j.celrep.2022.110433.
4. Li, R.H. et al. (2021). A phosphatidic acid-binding lncRNA SNHG9 facilitates LATS1 liquid-liquid phase separation to promote oncogenic YAP signaling. *Cell Res*. doi: 10.1038/s41422-021-00530-9.
5. Nadeem, A. et al. (2021). Phosphatidic acid-mediated binding and mammalian cell internalization of the *Vibrio cholerae* cytotoxin MakA. *PLoS Pathog*. **17**(3):e1009414. doi: 10.1371/journal.ppat.1009414.
6. Loh, J. et al. (2019). An acute decrease in plasma membrane tension induces macropinocytosis via PLD2 activation. *J Cell Sci*. **132**(17). pii: jcs232579. doi: 10.1242/jcs.232579.
7. Kumar Sah, R. et al. (2019). Phosphatidic acid homeostasis regulated by a type-2 phosphatidic acid phosphatase represents a novel druggable target in malaria intervention. *Cell Death Discov*. **5**:107. doi: 10.1038/s41420-019-0187-1.
8. Zhang, P. et al. (2019). Lipin 2/3 phosphatidic acid phosphatases maintain phospholipid homeostasis to regulate chylomicron synthesis. *J Clin Invest*. **129**(1):281-295. doi: 10.1172/JCI122595.
9. Han, H. et al. (2018). Regulation of the Hippo Pathway by Phosphatidic Acid-Mediated Lipid-Protein Interaction. *Mol Cell*. **72**(2):328-340.e8. doi: 10.1016/j.molcel.2018.08.038.
10. Lutkewitte, A.J. et al. (2018). Lipin deactivation after acetaminophen overdose causes phosphatidic acid accumulation in liver and plasma in mice and humans and enhances liver regeneration. *Food Chem Toxicol*. **115**:273-283. doi: 10.1016/j.fct.2018.03.014.

## **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.  
5628 Copley Drive  
San Diego, CA 92111  
Worldwide: +1 858 271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2016-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.