



# Phospholipid Assay Kit

Catalog Number KA1635

100 assays

Version: 06

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	4
<b>Assay Protocol .....</b>	<b>5</b>
Assay Procedure .....	5
<b>Data Analysis.....</b>	<b>7</b>
Calculation of Results.....	7

## Introduction

### Intended Use

- ✓ Applications:
  - Assays: phospholipid in biological samples such as serum and non-EDTA plasma.
  - Drug Discovery/Pharmacology: effects of drugs on choline-containing phospholipid metabolism.
  
- ✓ Features
  - Sensitive: Use 20  $\mu$ L samples. Linear detection range: colorimetric assay 3 - 200  $\mu$ M, fluorimetric assay 0.6 - 20  $\mu$ M phospholipid.

### Background

Phospholipids are a class of lipids which constitute a major component of cell membranes and play important roles in signal transduction. Most phospholipids contain one diglyceride, a phosphate group, and one choline. Phospholipid Assay Kit provides a simple, direct and high-throughput assay for measuring choline-containing phospholipids in biological samples. In this assay, phospholipids (such as lecithin, lysolecithin and sphingomyelin) are enzymatically hydrolyzed to choline which is determined using choline oxidase and a H<sub>2</sub>O<sub>2</sub> specific dye. The optical density of the pink colored product at 570 nm or fluorescence intensity (530/585 nm) is directly proportional to the phospholipid concentration in the sample.

## General Information

### Materials Supplied

List of component

Component	Amount
Assay Buffer	10 mL
PLD Enzyme	120 $\mu$ L
Enzyme Mix (Dried)	1 vial
Dye Reagent	120 $\mu$ L
Standard: 2 mM phosphatidylcholine	400 $\mu$ L

### Storage Instruction

Store all components at -20°C. Shelf life of six months after receipt.

### Materials Required but Not Supplied

- ✓ Pipetting devices
- ✓ Centrifuge tubes
- ✓ clear flat-bottom uncoated 96-well plates
- ✓ Optical density plate reader
- ✓ black flat-bottom uncoated 96-well plates
- ✓ Fluorescence plate reader.

### Precautions for Use

- ✓ Precautions  
Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

## Assay Protocol

### Assay Procedure

✓ Colorimetric assay

Liquid samples such as serum and plasma can be assayed directly. Solid samples can be homogenized in the assay buffer.

*Note: SH-containing reagents (e.g.  $\beta$ -mercaptoethanol, dithiothreitol,  $> 5 \mu\text{M}$ ), sodium azide, EDTA, and sodium dodecyl sulfate are known to interfere in this assay and should be avoided in sample preparation.*

1. Equilibrate all components to room temperature. Briefly centrifuge the tubes before opening. Keep thawed tubes on ice during assay. Reconstitute Enzyme Mix with 120  $\mu\text{L}$  Assay Buffer. Reconstituted Enzyme Mix is stable for 1 month when stored at  $-20^\circ\text{C}$ . *Note: a yellow precipitate may form after thawing reconstituted Enzyme Mix. If a precipitate forms, pellet it by centrifuging for 2 min at 14000 rpm and use the clear supernatant.*
2. Standards: mix 24  $\mu\text{L}$  2 mM Standard with 216  $\mu\text{L}$   $\text{dH}_2\text{O}$  (final 200  $\mu\text{M}$ ). Dilute standard in  $\text{dH}_2\text{O}$  as follows.

No	200 $\mu\text{M}$ STD + $\text{H}_2\text{O}$	Vol ( $\mu\text{L}$ )	Standard ( $\mu\text{M}$ )
1	100 $\mu\text{L}$ + 0 $\mu\text{L}$	100	200
2	60 $\mu\text{L}$ + 40 $\mu\text{L}$	100	120
3	30 $\mu\text{L}$ + 70 $\mu\text{L}$	100	60
4	0 $\mu\text{L}$ + 100 $\mu\text{L}$	100	0

Transfer 20  $\mu\text{L}$  diluted standards into separate wells of a clear flat-bottom 96-well plate.

Samples: transfer 20  $\mu\text{L}$  of each sample into separate wells of the plate.

*Note: if a sample is known to contain choline, prepare an extra sample blank well with 20  $\mu\text{L}$  of the sample.*

3. Color reaction. Prepare enough Working Reagent by mixing, for each well, 85  $\mu\text{L}$  Assay Buffer, 1  $\mu\text{L}$  PLD Enzyme, 1  $\mu\text{L}$  Enzyme Mix and 1  $\mu\text{L}$  Dye Reagent. Add 80  $\mu\text{L}$  Working Reagent to each well.

For samples that contain choline, prepare a blank control reagent with no PLD Enzyme (i.e., 85  $\mu\text{L}$  Assay Buffer, 1  $\mu\text{L}$  Enzyme Mix and 1  $\mu\text{L}$  Dye Reagent). Add 80  $\mu\text{L}$  of the Control Reagent to the Sample Blank well.

Tap plate to mix. Incubate 30 min at room temperature.

*Note: if precipitation occurs with certain samples, carry out the reaction in centrifuge tubes. After the 30 min incubation, centrifuge 5 min at 14,000 rpm. Transfer the supernatant into the wells for OD reading.*

4. Read optical density at 570 nm (550-585 nm).

✓ Fluorimetric assay

The fluorimetric assay procedure is similar to the colorimetric procedure except that (1) 0, 6, 12 and 20  $\mu\text{M}$  phospholipid standards and (2) a black 96-well plate are used. Read fluorescence intensity at  $\lambda_{\text{ex}} = 530 \text{ nm}$  and  $\lambda_{\text{em}} = 585 \text{ nm}$ .

*Note: if the calculated phospholipid concentration of a sample is higher than 200  $\mu\text{M}$  in the Colorimetric Assay or 20  $\mu\text{M}$  in the Fluorimetric Assay, dilute sample in 0.5% Triton X-100 and repeat the assay. Multiply result by the dilution factor  $n$ .*

## Data Analysis

### Calculation of Results

Subtract blank value (#4) from the standard values and plot the  $\Delta OD$  or  $\Delta F$  against standard concentrations. Determine the slope and calculate the phospholipid concentration of Sample,

$$[\text{Phospholipid}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{\text{Slope}(\mu\text{M}^{-1})} \times n \quad (\mu\text{M})$$

$R_{\text{SAMPLE}}$  and  $R_{\text{BLANK}}$  are optical density or fluorescence intensity readings of the Sample and H<sub>2</sub>O Blank (or Sample Blank if sample contains choline), respectively.  $n$  is the sample dilution factor.

