

## EnzyChrom™ Phospholipid Assay Kit (EPLP-100)

### Quantitative Colorimetric/Fluorimetric Phospholipid Determination

#### DESCRIPTION

**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. BioAssay Systems' method 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.

#### KEY FEATURES

**Sensitive.** Use 20  $\mu$ L samples. Linear detection range: colorimetric assay 3 - 200  $\mu$ M, fluorimetric assay 0.6 - 20  $\mu$ M phospholipid.

#### APPLICATIONS

**Assays:** phospholipid in biological samples such as serum and non-EDTA plasma.

**Drug Discovery/Pharmacology:** effects of drugs on choline-containing phospholipid metabolism.

#### KIT CONTENTS

**Assay Buffer:** 10 mL  
**Enzyme Mix:** Dried  
**Standard:** 400  $\mu$ L 2 mM phosphatidylcholine

**PLD Enzyme:** 120  $\mu$ L  
**Dye Reagent:** 120  $\mu$ L

**Storage conditions.** The kit is shipped on ice. Store all components at -20°C. Shelf life of six months after receipt.

**Precautions:** reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### 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  $\mu$ 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  $\mu$ L Assay Buffer. Reconstituted Enzyme Mix is stable for 1 month when stored at -20°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  $\mu$ L 2 mM Standard with 216  $\mu$ L dH<sub>2</sub>O (final 200  $\mu$ M). Dilute standard in dH<sub>2</sub>O as follows.

No	200 $\mu$ M STD + H <sub>2</sub> O	Vol ( $\mu$ L)	Standard ( $\mu$ M)
1	100 $\mu$ L + 0 $\mu$ L	100	200
2	60 $\mu$ L + 40 $\mu$ L	100	120
3	30 $\mu$ L + 70 $\mu$ L	100	60
4	0 $\mu$ L + 100 $\mu$ L	100	0

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

**Samples:** transfer 20  $\mu$ 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  $\mu$ L of the sample.*

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

*For samples that contain choline, prepare a blank control reagent with no PLD Enzyme (i.e., 85  $\mu$ L Assay Buffer, 1  $\mu$ L Enzyme Mix and 1  $\mu$ L Dye Reagent). Add 80  $\mu$ 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  $\mu$ M phospholipid standards and (2) a black 96-well plate are used. Read fluorescence intensity at  $\lambda_{ex}$  = 530 nm and  $\lambda_{em}$  = 585 nm.

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

#### CALCULATION

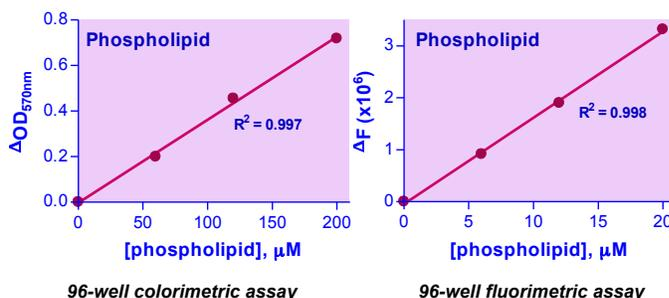
Subtract blank value (#4) from the standard values and plot the  $\Delta$ OD or  $\Delta$ F against standard concentrations. Determine the slope and calculate the phospholipid concentration of Sample,

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

$R_{\text{SAMPLE}}$  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.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

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.



#### PUBLICATIONS

1. Kim S et al (2012) Aberrant upregulation of astroglial ceramide potentiates oligodendrocyte injury. *Brain Pathol* 22(1):41-47.
2. Mukherjee A et al (2012) Lysophosphatidic acid activates lipogenic pathways and de novo lipid synthesis in ovarian cancer cells. *J Biol Chem* 287(3):24990-50000.
3. Stoll C et al (2011). Liposomes alter thermal phase behavior and composition of red blood cell membranes. *Biochim Biophys Acta* 1808(1):474-481.