

Product datasheet

Phospholipid Assay Kit (Colorimetric/Fluorometric)
ab234050

4 References 5 Images

Overview

Product name	Phospholipid Assay Kit (Colorimetric/Fluorometric)
Detection method	Colorimetric/Fluorometric
Sample type	Serum, Plasma
Assay type	Quantitative
Product overview	Phospholipid Assay Kit (Colorimetric/Fluorometric) (ab234050) is a simple, accurate and reproducible way to measure total choline-containing phospholipids in biological samples such as serum and plasma.

This kit provides a high-throughput assay of inhibitors/inducers affecting phospholipid metabolism. The signal is directly proportional to the phospholipid concentration of phospholipids in the sample and can detect phospholipids in the range of 0.1 to 10 nmol per sample.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K351 Phospholipid Assay Kit (Colorimetric/Fluorometric). K351-100 is the same size as the 100 test size of ab234050.

Phospholipids are a major component of lipid bilayer membranes. In addition, phospholipids are biologically important molecules as they are closely involved in signal transduction. Most phospholipids contain one diglyceride, a phosphate group, and one choline moiety. In human serum and plasma, almost 95% of the phospholipids are choline-containing phospholipids such as phosphatidylcholine and sphingomyelin. The remaining 5% is mainly choline-free phospholipids (i.e. phosphatidylethanolamine and phosphatidylserine). Phospholipid concentration is considered a disease biomarker as it can be altered in many lung, liver and heart diseases.

Platform Microplate reader

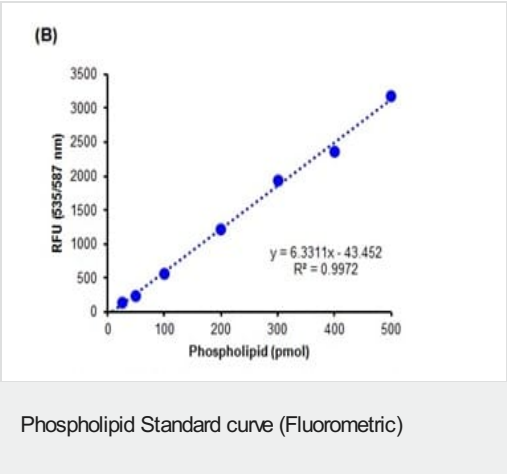
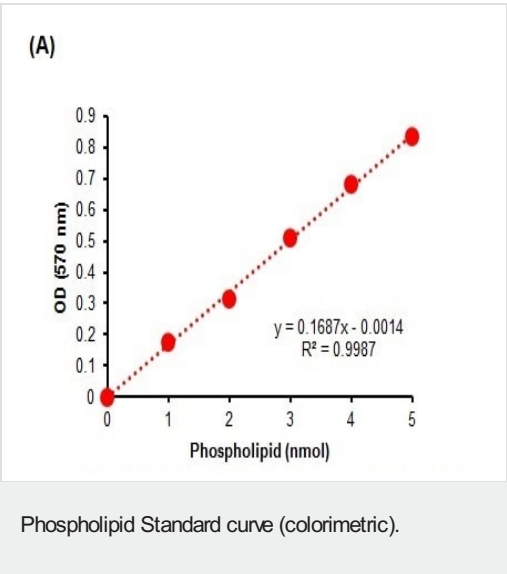
Properties

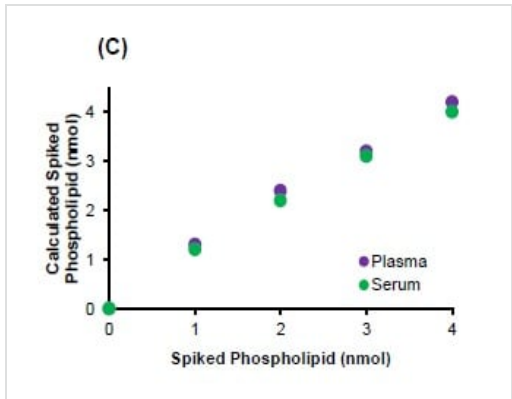
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Assay Buffer I	1 x 25ml

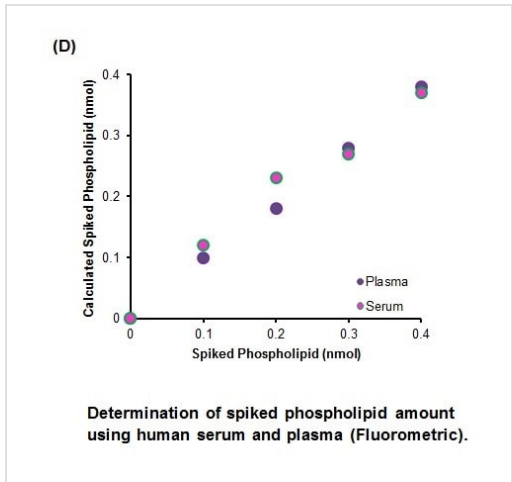
Components	100 tests
Enzyme Mix IV	1 vial
OxiRed Probe	1 x 0.2ml
PC Hydrolysis Enzyme	1 vial
Phospholipid Standard	1 vial

Images



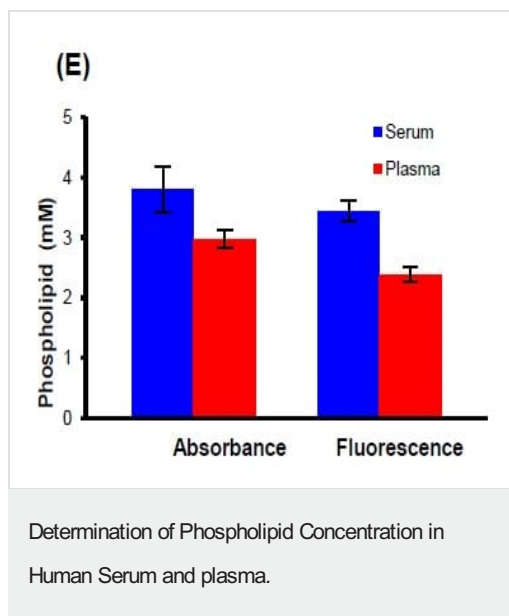


Determination of spiked phospholipid amount using human serum and plasma (colorimetric)



Determination of spiked phospholipid amount using human serum and plasma (Fluorometric).

Determination of spiked phospholipid amount using human serum and plasma (Fluorometric)



Normal concentrations in human ranges between 1 and 4 mM.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors