

Phosphatidylcholine Assay Kit (Colorimetric/Fluorometric) ab83377

23 References 4 图像

概述

产品名称	Phosphatidylcholine Assay试剂盒(Colorimetric/Fluorometric)
检测方法	Colorimetric/Fluorometric
样品类型	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
检测类型	Quantitative
灵敏度	> 0.1 nmol/well
范围	0.1 nmol/well - 10 nmol/well
检测时间	0h 40m
产品概述	Phosphatidylcholine Assay Kit (Colorimetric/Fluorometric) (ab83377) is a simple convenient means of measuring phosphatidylcholine in a variety of biological samples. This assay uses an enzyme-coupled reaction to hydrolyze phosphatidylcholine to release choline, which subsequently oxidizes the OxiRed probe in order to generate fluorescence (Ex/Em 535 nm 587 nm) and absorbance (570 nm). This assay measures phosphatidylcholine in the range of 0.1 to 10 nmol per sample.

Visit our [FAQs page](#) for tips and troubleshooting.

说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K576 Phosphatidylcholine Colorimetric/Fluorometric Assay Kit. K576-100 is the same size as the 100 test size of ab83377.</p> <p>Phosphatidylcholine (PC) is a phospholipid which incorporates choline as the headgroup of the lipid. PC is a major constituent of biological membranes and is involved in cell signaling through release of choline by phospholipase D leaving the second messenger phosphatidic acid.</p> <p>PC is present in serum at ~0.2-2.5 mM (~50-200 mg/dL).</p>
----	---

平台	Microplate reader
----	-------------------

性能

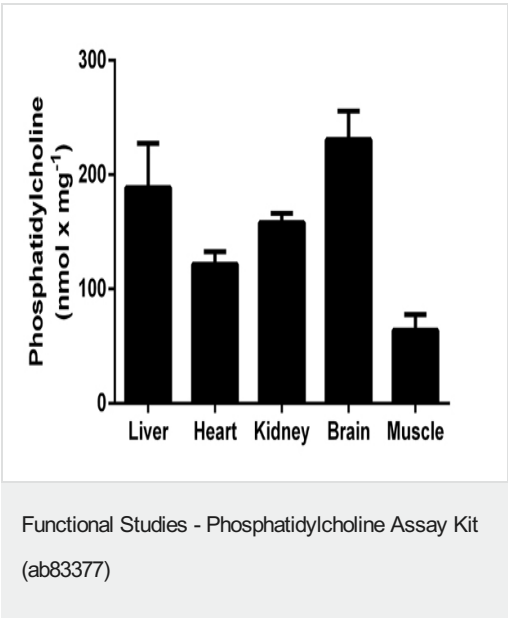
存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	100 tests
Assay Buffer I	1 x 25ml
Enzyme Mix IV	1 vial
OxiRed Probe	1 x 200µl
PC Hydrolysis Enzyme	1 vial
Phospholipid Standard	1 vial

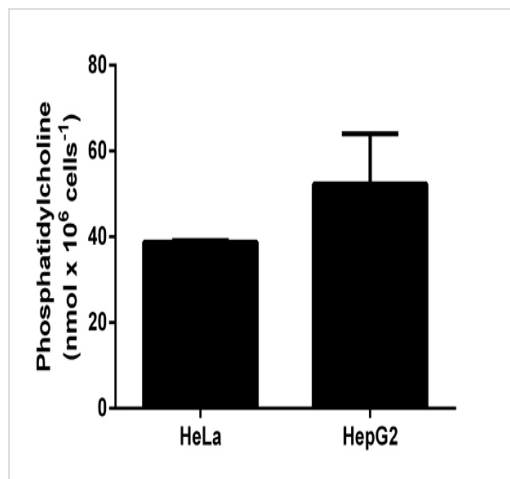
相关性

Phosphatidylcholine (PC) is a phospholipid which incorporates choline as the headgroup of the lipid. PC is a major constituent of biological membranes and is involved in cell signaling through release of choline by phospholipase D leaving the second messenger phosphatidic acid.

图片

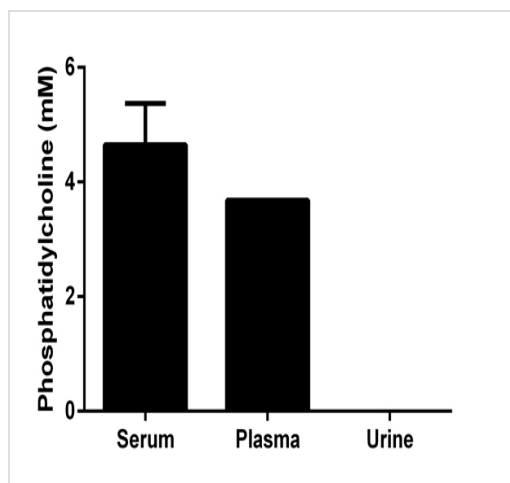


Phosphatidylcholine levels measured fluorometrically in mouse tissue lysates (mg of extracted protein; background signal subtracted, mean of duplicates; +/- SD).



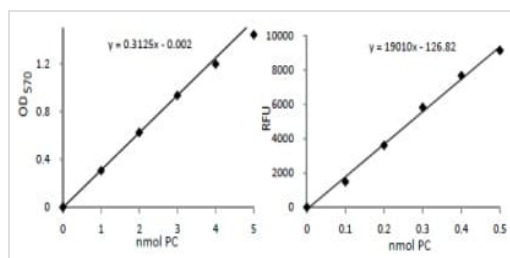
Phosphatidylcholine levels colorimetrically measured in cell lysates (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Phosphatidylcholine Assay Kit (ab83377)



Phosphatidylcholine levels colorimetrically measured in rat biological fluids (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Phosphatidylcholine Assay Kit (ab83377)



Examples of Colorimetric and Fluorometric Phosphatidylcholine standard curves using ab83377.

Functional Studies - Phosphatidylcholine Assay Kit (ab83377)

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.cn/abpromise> or contact our technical team.

Terms and conditions

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