

## Product datasheet

# Phosphatidylethanolamine Assay Kit ab241005

1 References 2 Images

### Overview

**Product name** Phosphatidylethanolamine Assay Kit

**Detection method** Fluorescent

**Product overview** The Phosphatidylethanolamine assay kit (ab241005) is a microplate based enzymatic assay for the quantitation of PE in cells and tissues. PE Converter hydrolyses PE to an intermediate, which converts a colorless probe to a fluorescent product via enzymatic reaction (Ex/Em: 535/587).

The intermediate formed through PE converter hydrolysis is specific to phosphatidylethanolamine. Thus no other phospholipids (i.e. phosphatidylcholine, phosphatidylinositol or phosphatidic acid) will be detected, making the kit highly specific. This assay kit can detect as low as 0.2 nmol per well.

**Platform** Microplate reader

### Properties

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

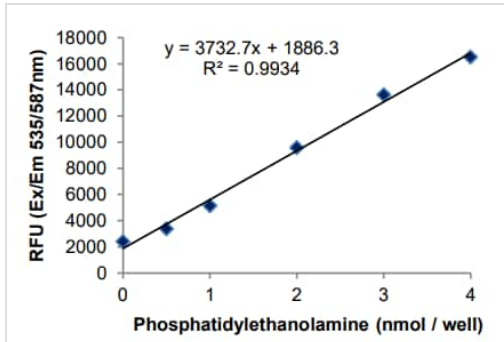
Components	Identifier	100 tests
PE Assay Buffer		1 x 25ml
PE Converter	Cap colour is white.	1 vial
PE Developer		1 x 600µl
PE Enzyme Mix		1 vial
PE Probe		1 x 200µl
PE Standard (1 mM)		1 x 100µl

### Relevance

Phosphatidylethanolamine (PE), also known as cephalin is the second most abundant phospholipid in animal and plant tissues, and is present on the cytoplasmic side of the plasma membrane. Phosphatidylethanolamine is a neutral phospholipid consisting of a phosphatidyl group ester linked to an ethanolamine molecule. Its functions include membrane fission/fusion, maintenance of membrane curvature and stabilization of membrane proteins, since it can form

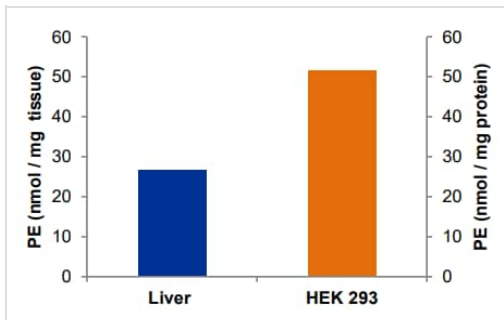
hydrogen bonds with proteins through an ionizable amine group. It acts as a chaperone during assembly of membrane proteins and aids in their translocation from the cytoplasm to the membrane. It is also involved in secretion of very low density lipoproteins in the liver.

## Images



Phosphatidylethanolamine standard curve

Phosphatidylethanolamine standard curve



Phosphatidylethanolamine content in rat liver (100 µg wet tissue) and HEK 293 cells (25 µg protein). Sample preparation and assay was carried out according to kit protocol.

Phosphatidylethanolamine content in rat liver (100 µg wet tissue) and HEK 293 cells (25 µg protein).

**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