

Phosphatidylglycerol Assay Kit (Fluorometric) ab273288

画像数 4

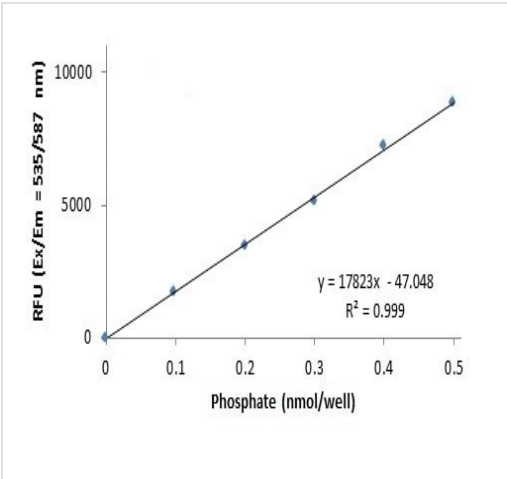
製品の概要

| | |
|------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 製品名 | Phosphatidylglycerol Assay Kit (Fluorometric) |
| 検出方法 | Fluorescent |
| サンプルの種類 | Cell culture extracts, Tissue Extracts |
| アッセイタイプ | Quantitative |
| ステップ | Multiple steps standard assay |
| 製品の概要 | Phosphatidylglycerol Assay Kit (Fluorometric) (ab273288) utilizes specific enzymes to generate an intermediate that reacts with a probe. This yields a signal that can be quantified fluorometrically and is proportional to the amount of Phosphatidylglycerol (PG) present in the sample, unaffected by the presence of phosphatidic acid, cardiolipin, or other lipids. When used as described, the assay is highly selective against other phospholipids and capable of detecting as little as 20 pmoles of phosphatidylglycerol. |
| 特記事項 | This product is manufactured by BioVision, an Abcam company and was previously called K488 Phosphatidylglycerol Assay Kit (Fluorometric). K488-100 is the same size as the 100 test size of ab273288. |
| 試験プラットフォーム | Microplate |

製品の特性

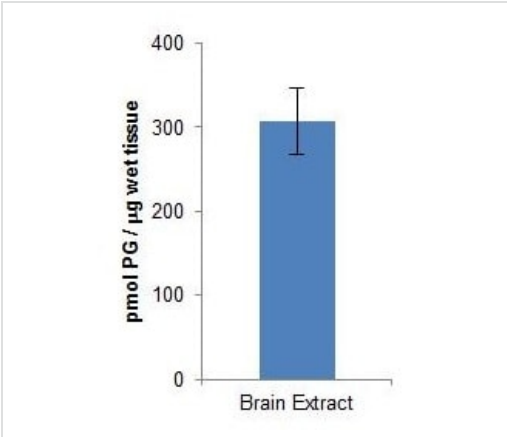
保存方法 Store at -20°C. Please refer to protocols.

| 内容 | 100 tests |
|-------------------------------|-----------|
| Assay Buffer I | 1 x 25ml |
| Enzyme Mix VI | 1 vial |
| OxiRed Probe | 1 x 200µl |
| PG Converter Mix | 1 vial |
| Phosphatidylglycerol Standard | 1 vial |

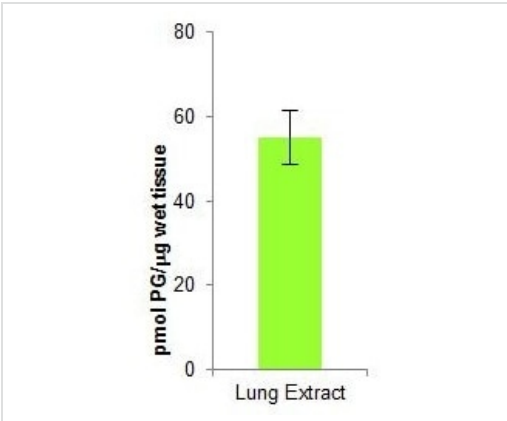


PG standard curve.

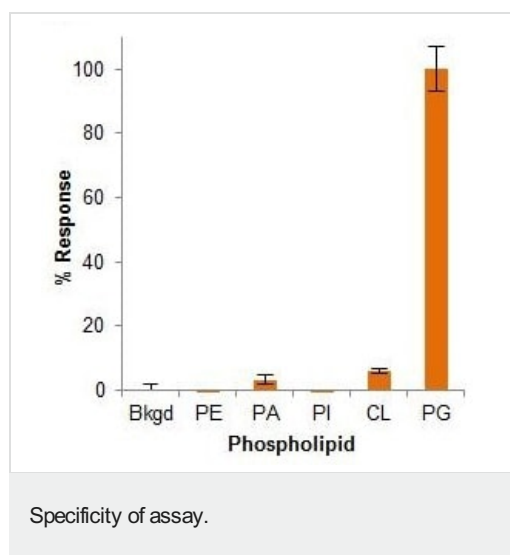
Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.



Determination of total PG concentration in lipid extracted from brain lysate.



Determination of Phosphatidylglycerol in lipid extracted from lung tissue.



Specificity of assay, when used as described, for Phosphatidylglycerol (PG) versus Phosphatidylethanolamine (PE), Phosphatidic Acid (PA), Phosphatidylinositol (PI), and cardiolipin (CL). The amount of phospholipid for each sample was 1 nmole.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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