

# Phosphate Assay Kit (Colorimetric) ab65622

**81 References**   **5 图像**

### 概述

产品名称	Phosphate Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Serum, Plasma, Other biological fluids, Tissue Extracts, Cell Lysate, Cell culture media
检测类型	Enzyme activity
灵敏度	> 1 $\mu$ M
范围	0.001 mM - 1 mM
检测时间	2h 00m
产品概述	<p>Phosphate Assay Kit (Colorimetric) (ab65622) provides an easy, quick and simple method for measuring phosphate levels.</p> <p>The phosphate assay protocol uses a proprietary formulation of malachite green and ammonium molybdate which forms a chromogenic complex with phosphate ion giving an intense absorption band around OD = 650nm.</p> <p>Phosphate assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add reaction mix to sample and standard wells</li> <li>- incubate for 30 min</li> <li>- analyze with a microplate reader</li> </ul>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K410 Phosphate Colorimetric Assay Kit. K410-500 is the same size as the 500 test size of ab65622.</p> <p>This assay can be used with biological fluids but also inorganic samples such as algal blooms and water from run-off areas of high fertilizer use.</p>
平台	Microplate reader

### 性能

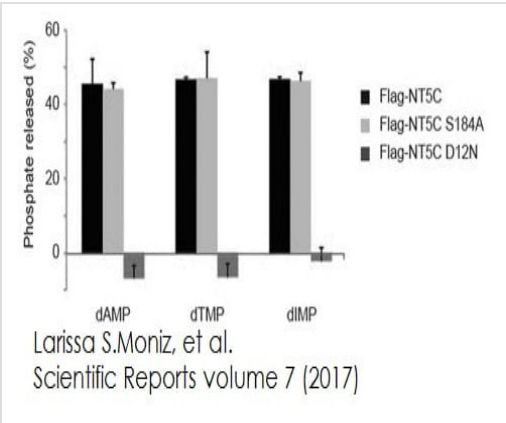
存放说明	Store at room temperature. Please refer to protocols.
------	---

组件	500 tests
Phosphate Reagent	1 x 15ml
Phosphate Standard	1 x 500µl

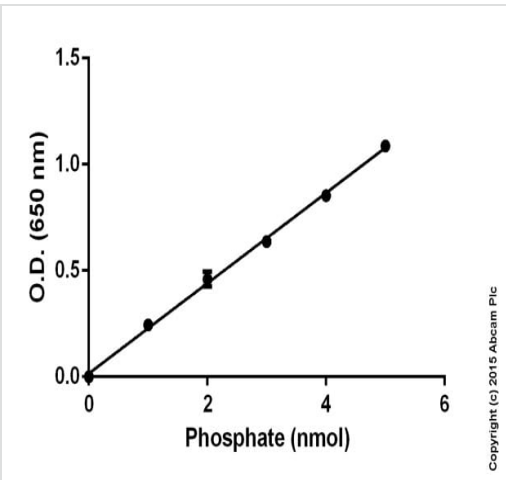
相关性

Phosphate is one of the most important of the inorganic ions in biological systems. It functions in a variety of roles. One of the most important roles is as a molecular switch, turning enzyme activity on and off through the mediation of the various protein kinases and phosphatases in biological systems. Phosphate is also of great importance in mineralization processes and is a primary stimulus of algal blooms frequently found in bodies of fresh water, due to run-off from areas of high fertilizer use.

图片



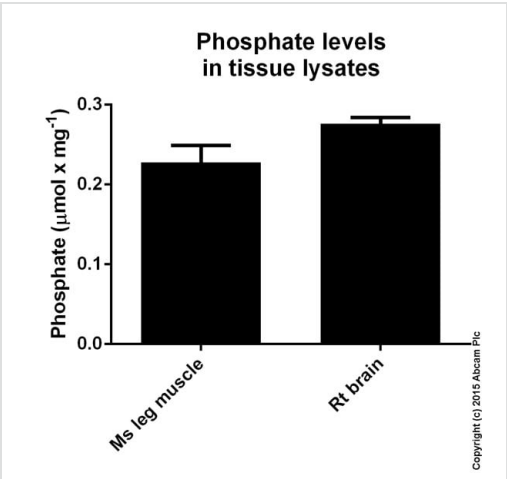
Functional Studies - Phosphate Assay Kit  
 (Colorimetric) (ab65622)  
 Larissa S.Moniz, et al., Scientific Reports volume 7, Fig 5, Article number: 39985 (2017)



Functional Studies - Phosphate Colorimetric Assay  
 Kit (ab65622)

Immunoprecipitates of Flag-NT5C ectopically expressed in HEK293 cells were incubated with 5 mM of the indicated nucleotides. Phosphate release was measured using a malachite green colorimetric assay (ab65622) and expressed as a percent of total nucleotide. The experiment was performed in duplicate and repeated 3 times independently. Error bars are sem.

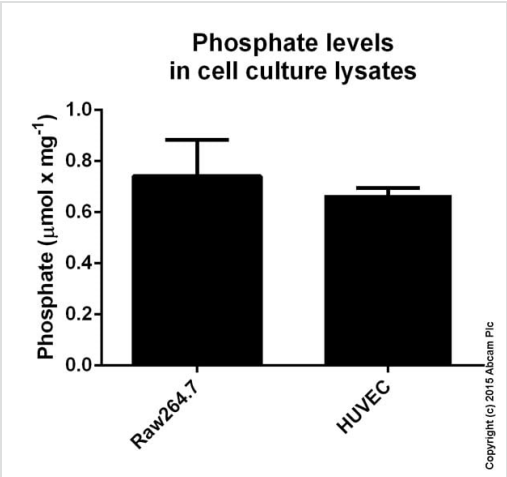
Standard curve: mean of duplicates (+/- SD) with background reads subtracted



Functional Studies - Phosphate Colorimetric Assay  
Kit (ab65622)

Phosphate measured in tissue lysates showing quantity (μmol) per mg of extracted protein.

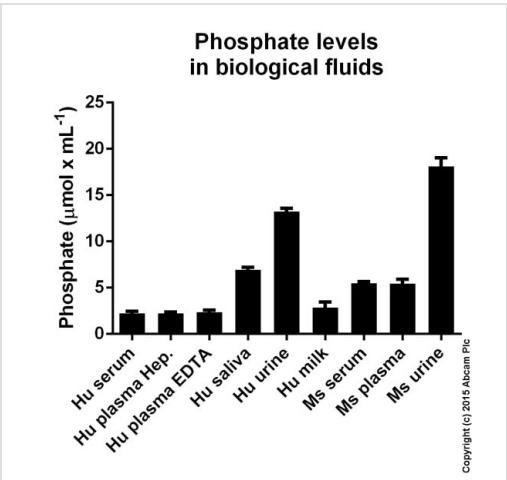
Protein concentration for samples varied from 9 mg/mL to 14 mg/mL. Samples were diluted 400-4000 fold.



Functional Studies - Phosphate Colorimetric Assay  
Kit (ab65622)

Phosphate measured in cell culture lysates showing quantity (μmol) per mg of extracted protein.

Samples with the concentration of 1.3e7 cells/mL were used. Samples were diluted 400-4000 fold.



Functional Studies - Phosphate Colorimetric Assay  
Kit (ab65622)

Phosphate measured in biological fluids showing quantity (μmol) per mL of tested sample. Samples were diluted 100-1000 fold.

**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