

Pyruvate Kinase Assay Kit ab83432

27 References **画像数 8**

製品の概要

製品名	Pyruvate Kinase Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids, Tissue
アッセイタイプ	Enzyme activity (quantitative)
検出感度	= 0.1 mU/ml
全工程の試験時間	0h 40m
製品の概要	<p>Pyruvate Kinase Assay Kit ab83432 provides a simple, direct and automation-ready assay for measuring pyruvate kinase activity in cell culture supernatant, plasma, serum, tissue, urine and other biological fluids (not validated for use with whole blood).</p> <p>In the pyruvate kinase assay protocol, PEP and ADP are catalyzed by pyruvate kinase to generate pyruvate and ATP. The generated pyruvate is oxidized by pyruvate oxidase to produce color (absorbance 570 nm) and fluorescence (Ex/Em 535/587 nm). Since the increase in color or fluorescence intensity is proportional to the increase in pyruvate amount, the PK activity can be accurately measured.</p> <p>The kit detects 0.1 mU/ml pyruvate kinase.</p> <p>Pyruvate kinase assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix - analyze with microplate reader immediately and after 10-20 min
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K709 Pyruvate Kinase Activity Colorimetric/Fluorometric Assay Kit. K709-100 is the same size as the 100 test size of ab83432.</p> <p>Pyruvate kinase (PK, EC 2.7.1.40) is an enzyme involved in glycolysis. It catalyzes the transfer of a phosphate group from phosphoenolpyruvate (PEP) to ADP, yielding one molecule of pyruvate and one molecule of ATP. Lack of pyruvate kinase will slow down the process of glycolysis which causes the disease known as pyruvate kinase deficiency.</p>
試験プラットフォーム	Microplate reader

製品の特性

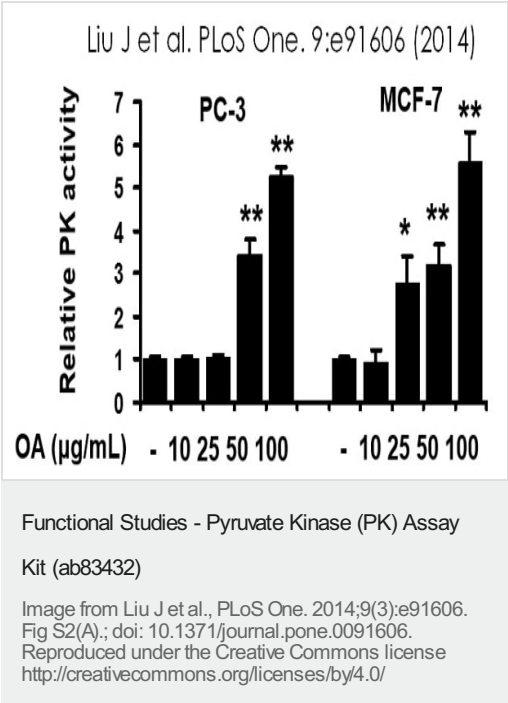
保存方法Store at -20°C. Please refer to protocols.

内容	100 tests	2000 tests
Development Enzyme Mix I	1 vial	20 vials
OxiRed Probe	1 x 200µl	20 x 200µl
PK Assay Buffer	1 x 25ml	20 x 25ml
PK Positive Control	1 vial	20 vials
PK Substrate Mix	1 vial	20 vials
Pyruvate Standard	1 x 100µl	20 x 100µl

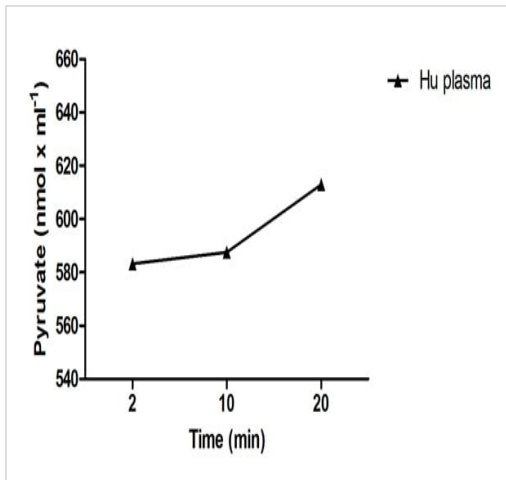
関連性Pyruvate kinase (PK, EC 2.7.1.40) is an enzyme involved in glycolysis. It catalyzes the transfer of a phosphate group from phosphoenolpyruvate (PEP) to ADP, yielding one molecule of pyruvate and one molecule of ATP. Lack of pyruvate kinase will slow down the process of glycolysis which causes the disease known as pyruvate kinase deficiency.

細胞内局在Cytoplasmic and Nuclear

画像

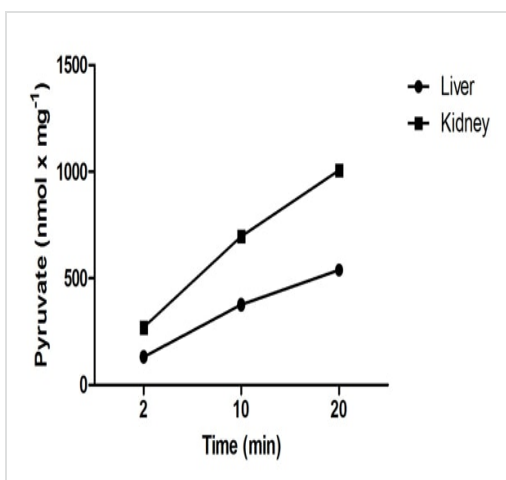


PC-3 and MCF-7 cells were treated with OA at the indicated doses and PK activity was determined 12 hr after OA treatment using ab83432. The average values of three independent experiments were shown as Mean ± SD. *, P<0.05, **, P<0.01.



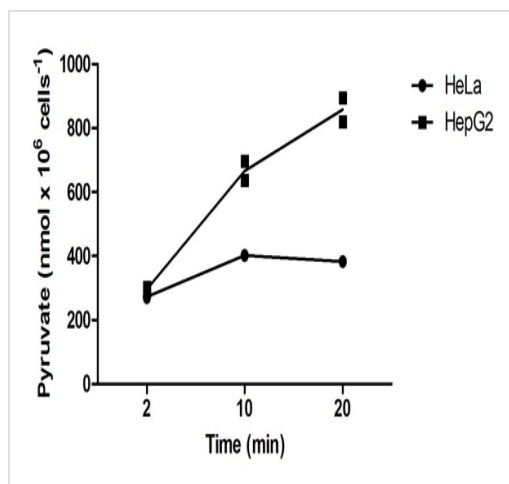
Pyruvate measured in biological fluid after 2 min, 10 min and 20 min incubation time showing quantity (nmol) per ml of tested sample.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



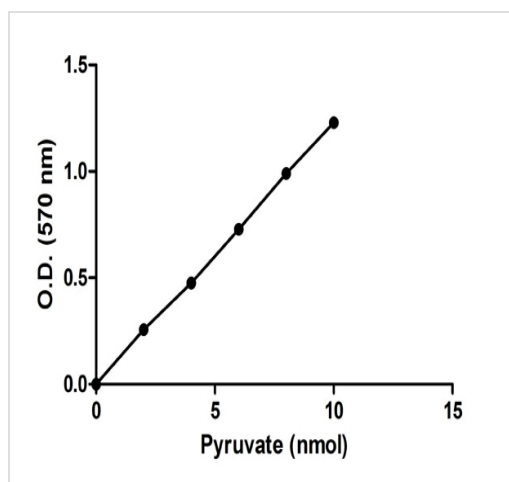
Pyruvate measured in mouse tissue lysates after 2 min, 10 min and 20 min incubation time showing quantity (nmol) per mg of tested sample.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



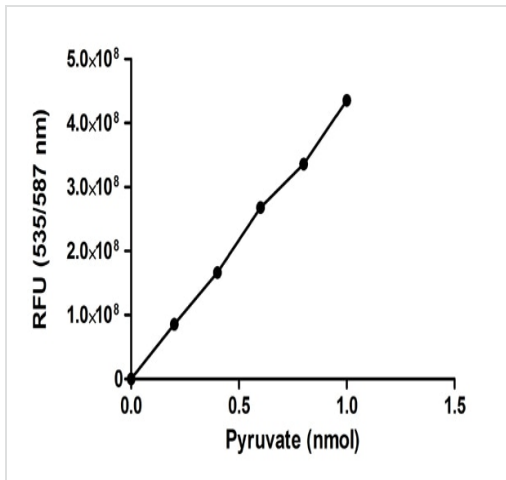
Pyruvate measured in cell lysates after 2 min, 10 min and 20 min incubation time showing quantity (nmol) per 1 mln of tested cells.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



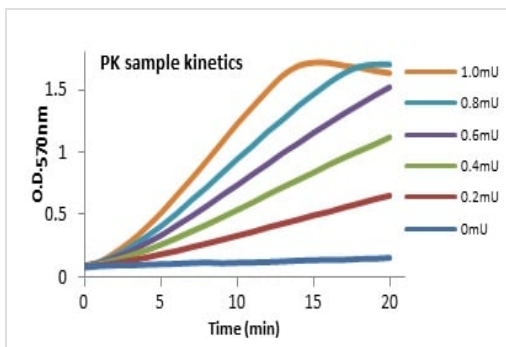
Colorimetric standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



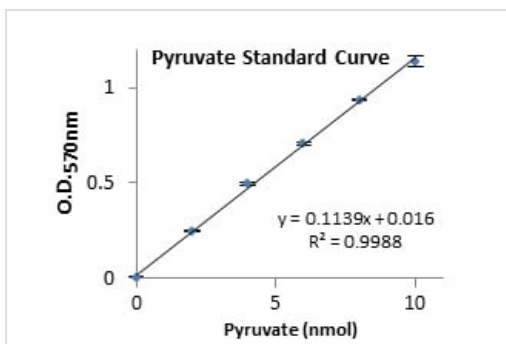
Fluorometric standard curve: mean of duplicates (+/- SD) with background reads subtracted.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



Pyruvate kinase samples time line using ab83432.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)



Pyruvate standard curve using ab83432.

Functional Studies - Pyruvate Kinase (PK) Assay
Kit (ab83432)

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