

Pyruvate Assay Kit ab65342

★★★★★ [2 Abreviews](#) [81 References](#) [画像数 6](#)

製品の概要

製品名	Pyruvate Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Urine, Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	> 1 μ M
検出範囲	1 μ M - 200 μ M
全工程の試験時間	0h 40m
製品の概要	<p>Pyruvate Assay Kit (Colorimetric/Fluorometric) ab65342 provides a simple, direct and automation-ready procedure for measuring pyruvate concentration in various biological samples such as blood, cells, culture and fermentation media, etc.</p> <p>In the pyruvate assay protocol, pyruvate is oxidized by pyruvate oxidase via enzyme reactions to generate color (λ= 570 nm) and fluorescence (at Ex/Em = 535/587 nm). Since the color or fluorescence intensity is proportional to pyruvate content, the pyruvate concentration can be accurately measured.</p> <p>The kit detects 1-200 μM pyruvate.</p> <p>Pyruvate assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix and incubate for 30 min at room temp - analyze with microplate reader
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K609 Pyruvate Colorimetric/Fluorometric Assay Kit. K609-100 is the same size as the 100 test size of ab65342.</p>
試験プラットフォーム	Microplate reader

製品の特性

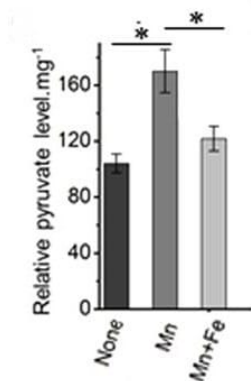
保存方法 Store at -20°C. Please refer to protocols.

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

関連性

Pyruvate is a key intersection molecule in the network of metabolic pathways. Pyruvate can be made from glucose through glycolysis converted back to carbohydrates via gluconeogenesis, or to fatty acids or energy via acetyl CoA. It can also be used to construct the amino acid alanine and be converted to ethanol. Glucose is converted to pyruvate which is then converted into acetyl CoA, the main input for the citric acid cycle (TCA). If insufficient oxygen is available, pyruvate breaks down anaerobically, creating lactate in animals and ethanol in plants and microorganisms.

画像

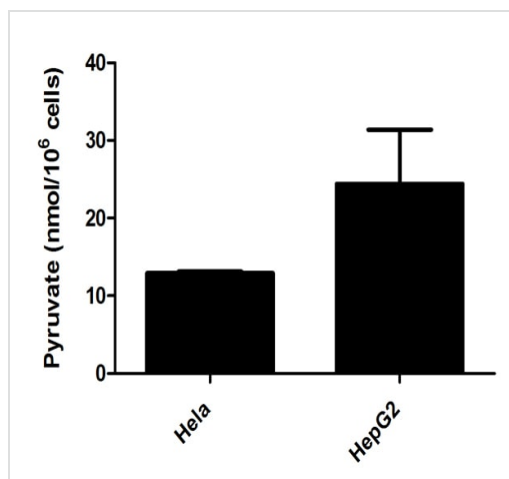


Kaur, Gursharan et al.
Scientific reports vol. 7,1
11645. (2017)

Pyruvate levels under manganese stress and iron supplementation were measured with ab65342 and the results were plotted as a function of normalized absorbance at 450 nm per milligram of cellular proteins.

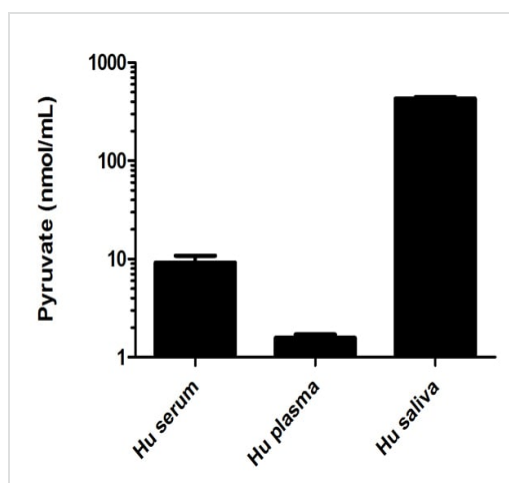
Functional Studies - Pyruvate Assay Kit (ab65342)

Kaur, Gursharan et al., Scientific reports vol. 7,1 11645.,
Fig 2, doi:10.1038/s41598-017-12004-3



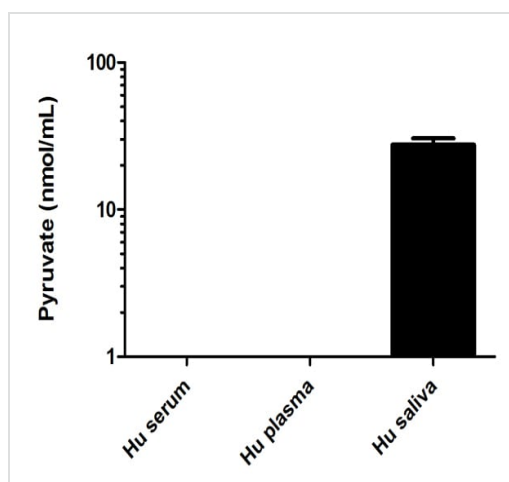
Pyruvate measured fluorometrically in biological fluids showing quantity (nmol) per mln of tested cells

Functional Studies - Pyruvate Assay Kit (ab65342) - Fluorometric



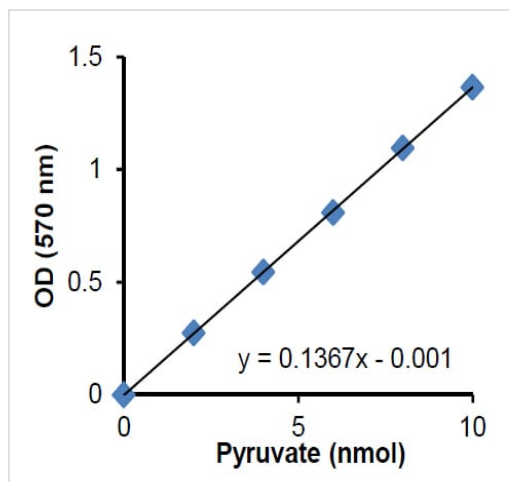
Pyruvate measured fluorometrically in biological fluids showing quantity (nmol) per mL of tested sample

Functional Studies - Pyruvate Assay Kit (ab65342) - Fluorometric



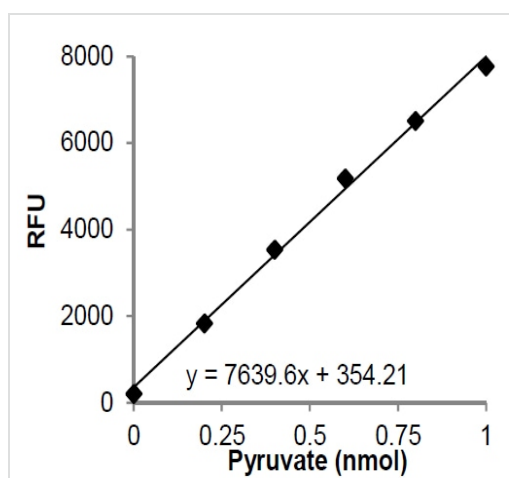
Pyruvate measured colourimetrically in biological fluids showing quantity (nmol) per mL of tested sample

Functional Studies - Pyruvate Assay Kit (ab65342) - Colourimetric



Pyruvate colorimetric standard curve.

Functional Studies - Pyruvate Assay Kit (ab65342)



Pyruvate fluorometric standard curve.

Functional Studies - Pyruvate Assay Kit (ab65342)

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