

Hexokinase Activity Assay Kit (Colorimetric) ab136957

44 References **画像数 5**

製品の概要

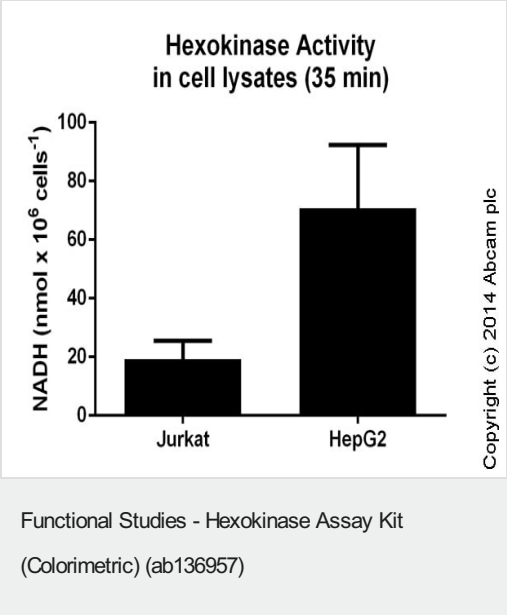
製品名	Hexokinase Activity Assay Kit (Colorimetric)
検出方法	Colorimetric
サンプルの種類	Serum, Tissue, Adherent cells, Suspension cells, Tissue Extracts
アッセイタイプ	Enzyme activity (quantitative)
検出感度	< 0.1 mU/well
全工程の試験時間	1h 00m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Hexokinase Activity Assay Kit (ab136957) uses a simple, sensitive and rapid procedure which can detect hexokinase activity to below 0.1 mU/well.</p> <p>In the hexokinase assay protocol, glucose is converted to glucose-6-phosphate by hexokinase. The glucose-6-phosphate is then oxidized by glucose-6-phosphate dehydrogenase to form NADH. The NADH reduces a colorless probe to a colored product with strong absorbance at 450 nm.</p> <p>Hexokinase activity assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix and incubate for 20-60 min - analyze repeatedly with a microplate reader during the incubation
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K789 Hexokinase Colorimetric Assay Kit. K789-100 is the same size as the 100 test size of ab136957.</p> <p>Hexokinases (HK) are found in many organisms including bacteria, plants and mammals and play an important role in glucose metabolism. Hexokinases phosphorylate glucose and generate glucose-6-phosphate for glycolysis. Hexokinases have four isoforms (HK-I, II, III and IV). HK-I, HK-II and HK-III have low Km, while HK-IV has 100 fold high Km.</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
Assay Buffer LX	1 x 25ml
Developer Solution III	1 vial
Development Enzyme Mix IX	1 vial
Hexokinase Positive Control	1 vial
HK Coenzyme	1 vial
HK Substrate	1 x 1ml
NADH Standard I	1 vial

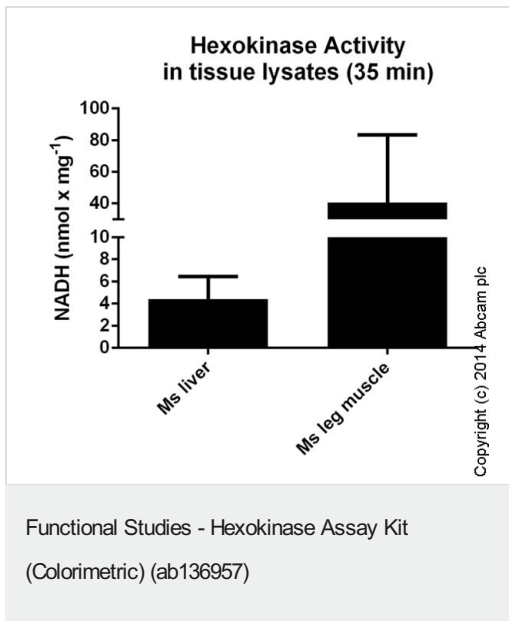
画像



Hexokinase Activity measured in cell lysates showing quantity (nmol) per 1 mln cells after 35 min of incubation.

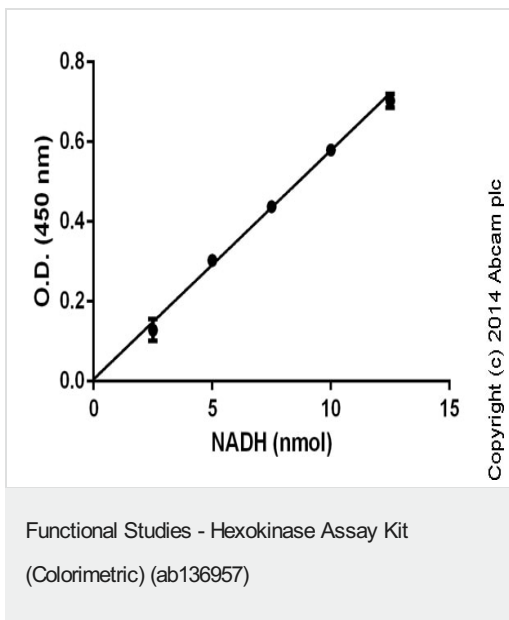
Samples with the concentration of 5e6 cells/mL were used.

Samples were diluted 1-27 fold.

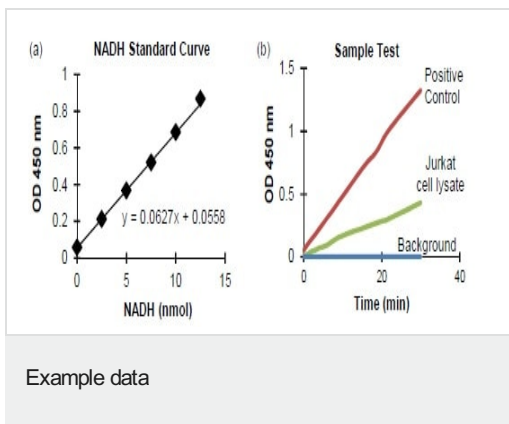


Hexokinase Activity measured in tissue lysates showing quantity (nmol) per mg of extracted protein after 35 min of incubation.

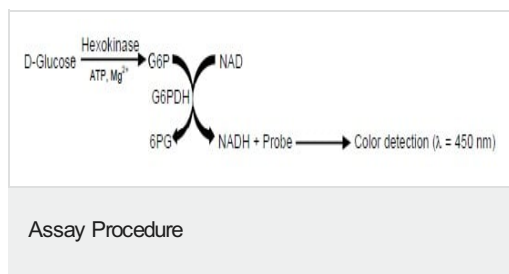
Protein concentration for samples varied from 13 mg/mL to 43 mg/mL. Samples were diluted 1-9 fold.



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



NADH standard curve (a). Hexokinase activity in Positive Control and Jurkat cell lysate (40 µg) (b).



In this kit glucose is converted to glucose-6-phosphate by hexokinase; the glucose-6-phosphate is oxidized by glucose-6-phosphate dehydrogenase to form NADH, which reduces a colorless probe to a colored product with strong absorbance at 450 nm.

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