

Hexokinase Activity Assay Kit (Fluorometric) ab211103

1 References 3 图像

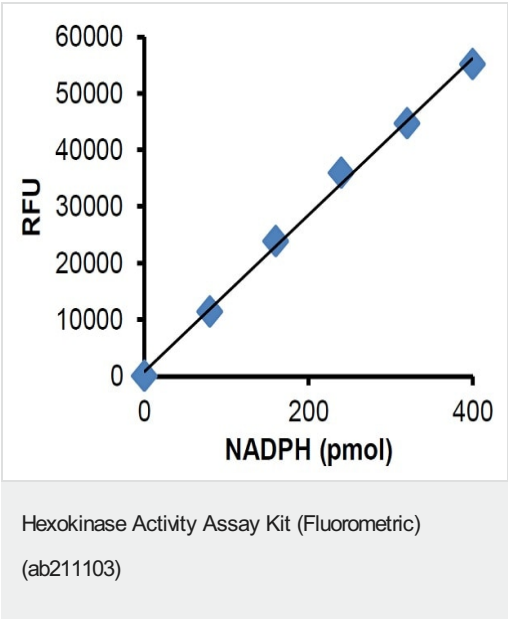
概述	
产品名称	Hexokinase Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell Lysate, Tissue Lysate
检测类型	Enzyme activity (quantitative)
灵敏度	2 μU
种属反应性	与反应: Mammals, Other species
产品概述	<p>The Hexokinase Activity Assay Kit (Fluorometric) (ab211103) provides a simple, sensitive and quick method for monitoring hexokinase (HK) activity in cells, serum, and animal or plant tissues. In this assay, HK converts glucose into glucose-6-phosphate, which in turn undergoes a series of reactions and reduces the sensitive probe to generate an intense fluorescent product that can be easily detected at Ex/Em = 535/587 nm.</p> <p>This assay can detect as low as 2 μU of HK activity.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K769 PicoProbe™ Hexokinase Activity Assay Kit (Fluorometric). K769-100 is the same size as the 100 test size of ab211103.</p> <p>Hexokinase (HK, 6-Phosphate glucose kinase, ATP:D-Hexose 6- Phosphotransferase, ATP-dependent hexokinase, EC 1.1.1.49) is responsible for phosphorylating hexoses (six-carbon sugars) to form hexose phosphate. Hexokinases play an important role in glucose metabolism, as glucose is the most important substrate of hexokinases. Hexokinases are found in many organisms including bacteria, plants and mammals. In mammals, there are four isoforms (HK-I, II, III and IV). HK-I, HK-II, and HK-III are referred as “low Km” because of their high affinity for glucose (Km < 1mM), while HK-IV (also known as Glucokinase) has a Km for glucose 100-fold higher and can only phosphorylate glucose when the substrate concentration is high enough.</p> <p>Hexokinase deficiency leads to diseases such as X-linked muscular dystrophy and rare autosomal recessive hemolytic anemia. On the other hand, increased hexokinase activity is detected in various human tumors and is associated with metastasis.</p>
平台	Microplate reader
性能	

存放说明

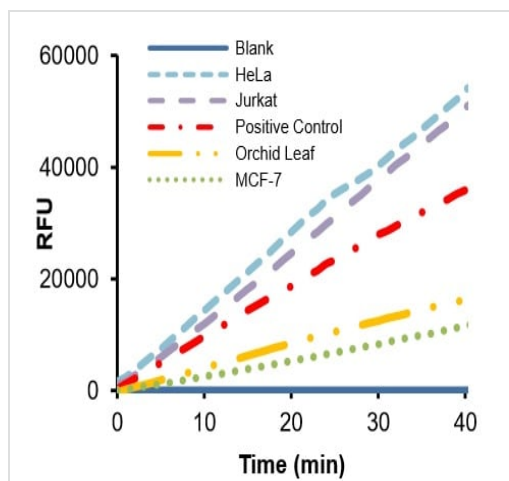
Store at -20°C. Please refer to protocols.

组件	100 tests
Assay Buffer LX	1 x 25ml
ATP II	1 vial
Developer IX	1 vial
Development Enzyme Mix IX	1 vial
HK Positive Control	1 vial
HK Substrate	1 x 1ml
NADPH Standard	1 vial
PicoProbe I	1 x 400µl

图片

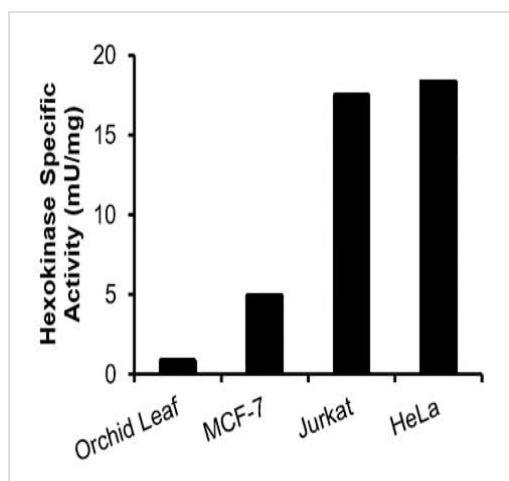


Typical NADPH standard calibration curve.



Kinetic curves showing Hexokinase activity in positive control (included in kit), lysates from HeLa (0.62 μ g), Jurkat (0.65 μ g) and MCF-7 cells (0.56 μ g), and lysates from orchid leaf (4 μ g).

Hexokinase Activity Assay Kit (Fluorometric)
(ab211103)



Hexokinase specific activity in lysates from orchid leaf (4 μ g), MCF-7 (0.56 μ g), Jurkat (0.65 μ g) and HeLa cells (0.62 μ g).

Hexokinase Activity Assay Kit (Fluorometric)
(ab211103)

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