

Hexokinase Inhibitor Assay Kit (Colorimetric) ab211107

1 References 1 图像

概述

产品名称	Hexokinase抑制剂Assay 试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Inhibitor compounds
检测类型	Semi-quantitative
检测时间	0h 30m
产品概述	Hexokinase Inhibitor Screening Kit (Colorimetric) (ab211107) provides a rapid, simple and adaptable test for high-throughput screening of Hexokinase inhibitors within 30 minutes. 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 OD = 450 nm. In the presence of a Hexokinase inhibitor, the reaction is impeded/abolished resulting in decrease or total loss of absorbance.

This simple and high-throughput adaptable assay kit can be used to screen, study or characterize potential inhibitors of Hexokinase.

**说明** This product is manufactured by BioVision, an Abcam company and was previously called K828 Hexokinase (HK) Inhibitor Screening Kit (Colorimetric). K828-100 is the same size as the 100 test size of ab211107.

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.

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.

平台 Microplate reader

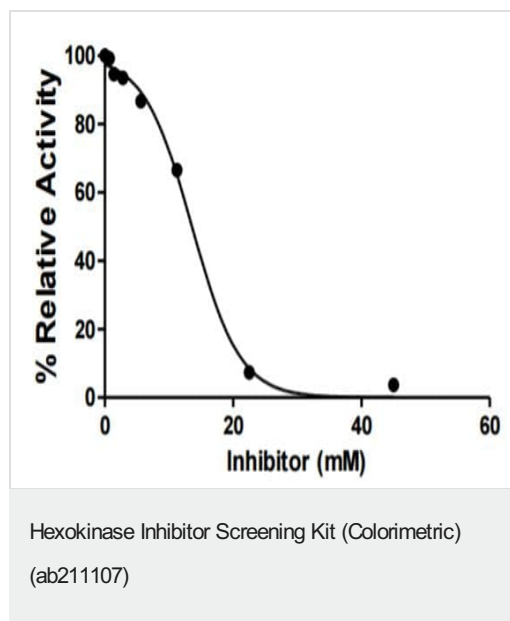
性能

## 存放说明

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

组件	100 tests
HK Assay Buffer	1 x 25ml
Hexokinase (0.5 U)	1 vial
HK Coenzyme (25 mg)	1 vial
HK Converter (10 U)	1 vial
HK Developer (30 mg)	1 vial
HK Inhibitor Control (40 $\mu$ mole)	1 vial
HK Substrate	1 x 1ml

## 图片



Typical inhibition curve of Hexokinase activity by Hexokinase Inhibitor (Bromopyruvic Acid). Assay was performed following the kit protocol.

**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