

# Hexokinase II Inhibitor Assay Kit (Colorimetric) ab211114

[3 References](#) [1 Image](#)

### Overview

---

<b>Product name</b>	Hexokinase II Inhibitor Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Assay time</b>	0h 30m
<b>Product overview</b>	Hexokinase II Inhibitor Screening Kit (Colorimetric) (ab211114) provides a sensitive, quick, and easy method for screening potential inhibitors of Hexokinase II (HK-II) in less than 30 minutes. The assay is based in the ability of HK-II to convert glucose into glucose-6-phosphate. Glucose-6-phosphate is oxidized by glucose-6-phosphate dehydrogenase to form NADH, which reduces a probe that shows strong absorbance at OD 450 nm. In the presence of a HK-II inhibitor such as Bromopyruvic Acid, the reaction is impeded, decreasing the rate or extent of generation of HK-II-dependent absorbance at OD 450 nm.

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

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K713 Human Hexokinase (HK) Inhibitor Screening Kit (Colorimetric). K713-100 is the same size as the 100 test size of ab211114.

Hexokinases (HK, EC: 2.7.1.1) are found in many organisms including bacteria, plants and mammals and play an important role in glucose metabolism. The Hexokinase family phosphorylates glucose and generates glucose-6-phosphate for glycolysis. Four Hexokinase isoforms (HK-I, II, III and IV) are found in numerous species. HK-I, HK-II and HK-III have low affinity for glucose, one of their natural substrates, while HK-IV show high affinity for this monosaccharide. Recent studies have found increased hexokinase activity in various human metastatic tumors. Moreover, Hexokinase II (HKII) is the main isoform of the Hexokinases and is responsible for malignant phenotypes. HKII binds to the outer mitochondrial membrane via the Voltage-Dependent Anion Channel (VDAC), a Porin-like protein. HKII has become an attractive therapeutic target for its role in cancer metastasis.

**Platform** Microplate reader

### Properties

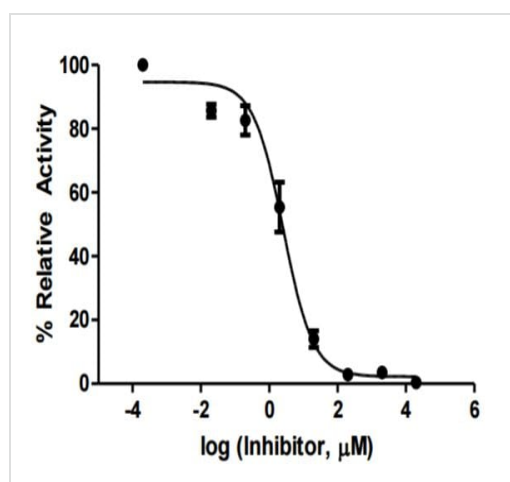
---

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	100 tests
Hexokinase II (Human)	1 x 10 $\mu$ l
HK Assay Buffer	1 x 25ml
HK Coenzyme (25 mg)	1 vial
HK Converter (10 U)	1 vial
HK Developer (30 mg)	1 vial
HK Inhibitor Control (Bromopyruvic Acid) (50 $\mu$ mol)	1 vial
HK Substrate	1 x 1ml

<b>Tissue specificity</b>	Predominant hexokinase isozyme expressed in insulin-responsive tissues such as skeletal muscle.
<b>Pathway</b>	Carbohydrate metabolism; hexose metabolism.
<b>Sequence similarities</b>	Belongs to the hexokinase family. Contains 2 hexokinase domains.
<b>Domain</b>	The N- and C-terminal halves of this hexokinase show extensive sequence similarity to each other. The catalytic activity is associated with the C-terminus while regulatory function is associated with the N-terminus. Each domain can bind a single glucose and Gluc-6-P molecule.
<b>Cellular localization</b>	Mitochondrion outer membrane. Its hydrophobic N-terminal sequence may be involved in membrane binding.

## Images



Typical inhibition curve of human Hexokinase II activity by the hexokinase inhibitor Bromopyruvic Acid.  $IC_{50}$  was determined to be 3  $\mu$ M. Assay was performed following the kit protocol.

Hexokinase II Inhibitor Screening Kit (Colorimetric)  
(ab211114)

**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.com/abpromise> or contact our technical team.

## **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors