

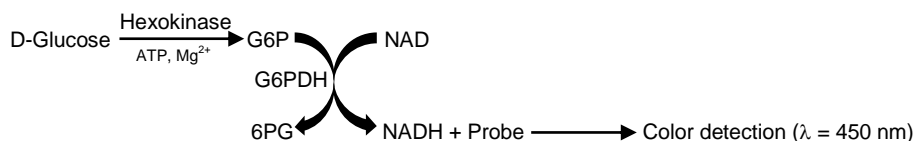
Hexokinase Colorimetric Assay Kit

rev. 06/18

(Catalog # K789-100; 100 assays; Store at -20°C)

I. Introduction:

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 K_m , while HK-IV has 100 fold high K_m . Hexokinase deficiency leads to severe human diseases such as X-linked muscular dystrophy and a rare autosomal recessive hemolytic anemia. On the other hand, increased hexokinase activity is detected in various human tumors and is associated with metastasis. Early detection of abnormal hexokinase activity is crucial for diagnosis, prediction and treatment of the disease. In BioVision's Hexokinase Assay 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. The assay is simple, sensitive and rapid and can detect hexokinase activity even less than 0.1 mU/well.



II. Application:

- Measurement of Hexokinase activity in various tissues/cells.
- Analysis of glucose metabolism and cell signaling in various cell types.
- Screening anti-diabetic drugs.

III. Sample Type:

- Serum
- Animal tissues: Liver, Heart, Kidney etc.
- Cell culture: Adherent or suspension cells.

IV. Kit Contents:

Components	K789-100	Cap Code	Part Number
HK Assay Buffer	25 ml	WM	K789-100-1
HK Substrate	1 ml	White	K789-100-2
HK Coenzyme (Lyophilized)	1 vial	Purple	K789-100-3
HK Enzyme Mix (Lyophilized)	1 vial	Green	K789-100-4
HK Developer (Lyophilized)	1 vial	Red	K789-100-5
NADH Standard (Lyophilized)	1 vial	Yellow	K789-100-6
HK Positive Control (Lyophilized)	1 vial	Orange	K789-100-7

V. User Supplied Reagents and Equipment:

- 96-well plate with flat clear bottom
- Multi-well spectrophotometer (ELISA reader)

VI. Storage and Handling:

Store kit at -20°C, protected from light. Warm all buffers to room temperature before use. Briefly centrifuge all small vials prior to opening.

VII. Reagent Preparation and Storage Conditions:

- **HK Coenzyme:** Reconstitute with 220 μ l HK Assay Buffer to generate 0.2 M solution. Store at -20°C. Use within two months. Keep on ice while in use.
- **HK Enzyme Mix:** Reconstitute with 220 μ l HK Assay Buffer. Pipette up and down to dissolve completely. Aliquot and store at -20°C. Avoid repeated freeze/thaw cycles. Use within two months. Keep on ice while in use.
- **HK Developer:** Reconstitute with 220 μ l dH₂O. Pipette up and down to dissolve completely. Store at -20°C. Use within two months.
- **NADH Standard:** Reconstitute with 400 μ l dH₂O to generate 1.25 mM (1.25 nmol/ μ l) NADH Standard solution. Store at -20°C. Use within two months. Keep on ice while in use.
- **HK Positive Control:** Reconstitute with 100 μ l HK Assay Buffer and mix thoroughly. Aliquot and store at -20°C.

VIII. Hexokinase Assay Protocol:

1. **NADH Standard Curve:** Add 0, 2, 4, 6, 8 and 10 μ l of 1.25 mM NADH Standard into a series of wells in duplicate in 96 well plate to generate 0, 2.5, 5.0, 7.5, 10 and 12.5 nmol/well of NADH Standard. Adjust volume to 50 μ l/well with HK Assay Buffer.
2. **Sample Preparation:** Rapidly homogenize tissue (10 mg) or cells (1×10^6) with 200 μ l ice cold HK Assay Buffer for 10 minutes on ice. Centrifuge at 12000 rpm for 5 min. Collect the supernatant. Add 1-50 μ l sample (40 μ g) per well, adjust final volume to 50 μ l with HK Assay Buffer. Prepare a parallel sample well as the background control to avoid interference from the NADH in the sample. **Note:** For unknown samples, we suggest testing several doses to ensure the readings are within the standard curve range.

FOR RESEARCH USE ONLY!

3. HK Positive Control: Dilute Positive control solution 1:99 in HK Assay Buffer. Use 1-10 μl of diluted Positive Control into the desired well(s) and adjust the final volume to 50 μl with HK Assay Buffer.

4. Reaction Mix: Mix enough reagents for the number of assays to be performed. For each well, prepare 50 μl Mix containing:

	Reaction Mix	Background Control Mix
HK Assay Buffer	34 μl	44 μl
HK Enzyme Mix	2 μl	2 μl
HK Developer	2 μl	2 μl
HK Coenzyme	2 μl	2 μl
HK Substrate	10 μl	----

Add 50 μl of the reaction mix to each well containing the Standard, Positive Control and test samples and 50 μl of background control mix to each well containing the background control sample. Mix well.

5. Measurement: Incubate for 20-60 min at room temperature and measure $\text{OD}_{450\text{nm}}$. **Note:** Incubation time depends on the Hexokinase activity in the samples. We recommend measuring the OD in a kinetic mode, and choose two time points (T_1 & T_2) in the linear range to calculate the hexokinase activity of the samples. The NADH standard curve can read in endpoint mode (i.e., at the end of incubation time).

6. Calculation: Subtract the 0 standard reading from all standard readings. Plot the NADH standard curve. Correct sample background by subtracting the value derived from the background control from all sample readings. Calculate the hexokinase activity of the test sample: $\Delta\text{OD} = A_2 - A_1$. Apply the ΔOD to the NADH standard curve to get B nmol of NADH generated by hexokinase during the reaction time ($\Delta T = T_2 - T_1$).

$$\text{Sample Hexokinase activity} = \frac{B}{(\Delta T \times V)} \times \text{Dilution Factor} = \text{nmol/min/ml} = \text{mU/ml}$$

Where: **B** is the NADH amount from standard curve (nmol).

ΔT is the reaction time (min).

V is the sample volume added into the reaction well (ml).

Unit Definition: One unit of hexokinase is the amount of enzyme that will generate 1.0 μmol of NADH per min at pH 8 at room temperature.

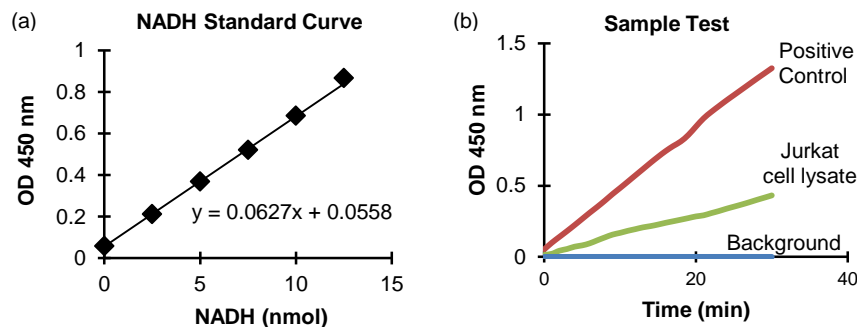


Figure 1: NADH standard curve (a). Hexokinase activity in Positive Control and Jurkat cell lysate (40 μg) (b). Assays were performed following kit protocol.

IX. RELATED PRODUCTS:

Glucose and Sucrose Assay Kit
 Glucose Uptake Colorimetric Assay Kit
 Glucose Uptake Fluorometric Assay Kit
 Maltose and Glucose Assay Kit
 PicoProbe™ Glucose-6-Phosphate Assay Kit

NAD/NADH Quantification Kit
 Glucose-6-Phosphate Dehydrogenase Assay Kit
 Glucose Assay kit
 NADP/NADPH Quantification Kit
 Glucose Dehydrogenase Activity Assay Kit

FOR RESEARCH USE ONLY! Not to be used on humans.