



## **ELISA PRODUCT INFORMATION & MANUAL**

Hexokinase 1  
***NBP2-60571***

Enzyme-linked Immunosorbent Assay for quantitative detection of Human Hexokinase 1. For research use only.  
Not for diagnostic or therapeutic procedures.

## Assay Summary

**Step 1.** Add 50  $\mu\text{l}$  of Standard or Sample per well.  
Incubate 2 hours.

**Step 2.** Wash, then add 50  $\mu\text{l}$  of Biotinylated Antibody per well.  
Incubate 2 hours.

**Step 3.** Wash, then add 50  $\mu\text{l}$  of SP Conjugate per well.  
Incubate 30 minutes.

**Step 4.** Wash, then add 50  $\mu\text{l}$  of Chromogen Substrate per well.  
Incubate 30 minutes.

**Step 5.** Add 50  $\mu\text{l}$  of Stop Solution per well.  
Read at 450 nm immediately.

## Symbol Key



Consult instructions for use.



# Human Hexokinase-1 ELISA Kit

Catalog No. NBP2-60571

*Sample insert for reference use only*

## Introduction

Hexokinases (HK) catalyse the phosphorylation of hexose to produce hexose 6-phosphate. The four mammalian hexokinases are designated HK1, HK2, HK3, and HK4. Although highly conserved in amino acid sequence, these enzymes differ in molecular mass, tissue distribution, regulation, and catalytic properties. HK1 – HK3 have molecular masses of approximately 100 kDa and show a broad but distinct tissue distribution. They have a relatively high affinity for glucose and are subject to feedback regulation by physiological levels of glucose 6-phosphate (G6P). The 100-kDa HK originates from a 50-kDa precursor via gene duplication and tandem ligation (1). Hexokinase-1 (HK1) is the most ubiquitously expressed isoform out of the four hexokinases. It is constitutively expressed in the brain, kidney, and red blood cells. It is located at the mitochondria outer membrane. HK1 gene encodes a 102-kDa and 917-amino acid protein (2). It is comprised of two similar domains fused into a single polypeptide chain. The C-terminal half of HK1 is catalytically active, whereas the N-terminal half is necessary for the relief of product inhibition by phosphate (3). It plays important roles in glucose catabolism and cell survival.

## Principle of the Assay

The Human Hexokinase-1 ELISA (Enzyme-Linked Immunosorbent Assay) Kit is designed for detection of hexokinase-1 in **cell lysate samples**. This assay employs a quantitative **sandwich enzyme immunoassay** technique that measures human hexokinase-1 in approximately 5 hours. A polyclonal antibody specific for human hexokinase-1 has been pre-coated onto a 96-well microplate with removable strips. Hexokinase-1 in standards and samples is sandwiched by the immobilized antibody and a biotinylated polyclonal antibody specific for human hexokinase-1, which is recognized by a streptavidin-peroxidase (SP) conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

## Caution and Warning

- This product is for **Research Use Only** and is not intended for use in diagnostic procedures.

- Prepare all reagents (diluent buffer, wash buffer, standard, biotinylated antibody, and SP conjugate) as instructed, prior to running the assay.
- Prepare all samples prior to running the assay. The dilution factors for the samples are suggested in this insert. However, the user should determine the optimal dilution factor.
- Spin down the SP conjugate vial, the biotinylated antibody vial, and the standard diluent vial before opening and using contents.
- The Stop Solution is an acidic solution.
- The kit should not be used beyond the expiration date.

## Reagents

- **Human Hexokinase-1 Microplate:** A 96-well polystyrene microplate (12 strips of 8 wells) coated with a polyclonal antibody against human hexokinase-1.
- **Sealing Tapes:** Each kit contains 3 precut, pressure sensitive sealing tapes that can be cut to fit the format of the individual assay.
- **Human Hexokinase-1 Standard:** Human hexokinase-1 in a buffered protein base (1 µg, lyophilized, 2 vials).
- **Biotinylated Human Hexokinase-1 Antibody (50x):** A 50-fold concentrated biotinylated polyclonal antibody against human hexokinase-1 (120 µl).
- **EIA Diluent Concentrate (10x):** A 10-fold concentrated buffered protein base (20 ml).
- **Standard Diluent (1x):** A buffered protein base with stabilizer (2 ml).
- **Wash Buffer Concentrate (20x):** A 20-fold concentrated buffered surfactant (30 ml, 2 bottles).
- **SP Conjugate (100x):** A 100-fold concentrate (80 µl).
- **Chromogen Substrate (1x):** A stabilized peroxidase chromogen substrate tetramethylbenzidine (8 ml).
- **Stop Solution (1x):** A 0.5 N hydrochloric acid solution to stop the chromogen substrate reaction (12 ml).

## Storage Condition

- Upon arrival, immediately store components of the kit at recommended temperatures up to the expiration date.
- Store Standard, SP Conjugate, and Biotinylated Antibody at -20°C.
- Store Microplate, Diluent Concentrate (10x), Standard Diluent (1x), Wash Buffer, Stop Solution, and Chromogen Substrate at 2-8°C.
- Unused microplate wells may be returned to the foil pouch with the desiccant packs and resealed. May be stored for up to 30 days in a vacuum desiccator.

## Other Supplies Required

- Microplate reader capable of measuring absorbance at 450 nm
- Pipettes (1-20  $\mu\text{l}$ , 20-200  $\mu\text{l}$ , 200-1000  $\mu\text{l}$ , and multiple channel)
- Deionized or distilled reagent grade water

## Sample Collection, Preparation, and Storage

- **Cell Lysate:** Rinse cell with cold PBS and then scrape the cell into a tube with 5 ml of cold PBS and 0.5 M EDTA. Centrifuge suspension at 1500 rpm for 10 minutes at 4°C and aspirate supernatant. Re-suspend pellet in ice-cold Lysis Buffer (10 mM Tris, pH 8.0, 130 mM NaCl, 1% Triton X-100, protease inhibitor cocktail). For every  $1 \times 10^6$  cells, add approximately 100  $\mu\text{L}$  of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at 4°C and collect supernatant.

*Applicable samples may also include biofluids, cell culture, and tissue homogenates. If necessary, user should determine optimal dilution factor depending on application needs.*

**Refer to Dilution Guidelines for further instruction.**

<b>Guidelines for Dilutions of 100-fold or Greater</b> <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
<b>100x</b>	<b>10000x</b>
A) 4 $\mu\text{l}$ sample: 396 $\mu\text{l}$ buffer (100x) = 100-fold dilution  <i>Assuming the needed volume is less than or equal to 400 <math>\mu\text{l}</math>.</i>	A) 4 $\mu\text{l}$ sample : 396 $\mu\text{l}$ buffer (100x) B) 4 $\mu\text{l}$ of A : 396 $\mu\text{l}$ buffer (100x) = 10000-fold dilution  <i>Assuming the needed volume is less than or equal to 400 <math>\mu\text{l}</math>.</i>
<b>1000x</b>	<b>100000x</b>
A) 4 $\mu\text{l}$ sample : 396 $\mu\text{l}$ buffer (100x) B) 24 $\mu\text{l}$ of A : 216 $\mu\text{l}$ buffer (10x) = 1000-fold dilution  <i>Assuming the needed volume is less than or equal to 240 <math>\mu\text{l}</math>.</i>	A) 4 $\mu\text{l}$ sample : 396 $\mu\text{l}$ buffer (100x) B) 4 $\mu\text{l}$ of A : 396 $\mu\text{l}$ buffer (100x) C) 24 $\mu\text{l}$ of B : 216 $\mu\text{l}$ buffer (10x) = 100000-fold dilution  <i>Assuming the needed volume is less than or equal to 240 <math>\mu\text{l}</math>.</i>

## Reagent Preparation

- Freshly dilute all reagents and bring all reagents to room temperature before use.
- EIA Diluent Concentrate (10x):** If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the EIA Diluent Concentrate 10-fold with reagent grade water to produce a 1x solution. Store for up to 30 days at 2-8°C.
- Human Hexokinase-1 Standard:** Reconstitute the Human Hexokinase-1 Standard (1 µg) with 0.5 ml of **Standard Diluent** to generate a 2 µg/ml standard stock solution. Allow the vial to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare duplicate or triplicate standard points by serially diluting from the standard stock solution (2 µg/ml) 2-fold with equal volume of **EIA Diluent** to produce 1000, 500, 250, 125, 62.5, 31.25, and 15.625 ng/ml solutions. EIA Diluent serves as the zero standard (0 ng/ml). Aliquot remaining stock solution to limit repeated freeze-thaw cycles. This solution should be stored at -20°C and used within 48 hours.

Standard Point	Dilution	[Hexokinase-1] (ng/ml)
P1	1 part Standard (2 µg/ml) + 1 part EIA Diluent	1000
P2	1 part P1 + 1 part EIA Diluent	500
P3	1 part P2 + 1 part EIA Diluent	250
P4	1 part P3 + 1 part EIA Diluent	125
P5	1 part P4 + 1 part EIA Diluent	62.5
P6	1 part P5 + 1 part EIA Diluent	31.25
P7	1 part P6 + 1 part EIA Diluent	15.625
P8	EIA Diluent	0.0

- Biotinylated Human Hexokinase-1 Antibody (50x):** Spin down the antibody briefly and dilute the desired amount of the antibody 50-fold with EIA Diluent to produce a 1x solution. The undiluted antibody should be stored at -20°C.
- Wash Buffer Concentrate (20x):** If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the Wash Buffer Concentrate 20-fold with reagent grade water to produce a 1x solution.
- SP Conjugate (100x):** Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 100-fold with EIA Diluent to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

## Assay Procedure

- Prepare all reagents, standard solutions, and samples as instructed. Bring all reagents to room temperature before use. The assay is performed at room temperature (20-25°C).
- Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
- Add 50  $\mu$ l of Human Hexokinase-1 Standard or sample to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
- Wash five times with 200  $\mu$ l of Wash Buffer manually. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a machine, wash six times with 300  $\mu$ l of Wash Buffer and then invert the plate, decanting the contents; hit 4-5 times on absorbent material to completely remove the liquid.
- Add 50  $\mu$ l of Biotinylated Human Hexokinase-1 Antibody to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours.
- Wash the microplate as described above.
- Add 50  $\mu$ l of SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
- Wash the microplate as described above.
- Add 50  $\mu$ l of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate for 30 minutes or until the optimal blue color density develops.
- Add 50  $\mu$ l of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
- Read the absorbance on a microplate reader at a wavelength of 450 nm **immediately**. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.



## Data Analysis

- Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
- To generate a standard curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best fit line can be determined by regression analysis using log-log or four-parameter logistic curve fit.
- Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.

## Typical Data

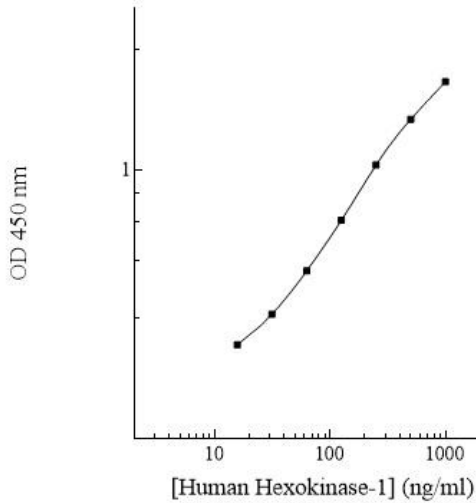
- The typical data is provided for reference only. Individual laboratory means may vary from the values listed. Variations between laboratories may be caused by technique differences.

Standard Point	ng/ml	OD	Average OD
P1	1000	2.140 2.114	2.127
P2	500	1.547 1.533	1.540
P3	250	1.048 1.036	1.042
P4	125	0.652 0.645	0.649
P5	62.5	0.425 0.421	0.423
P6	31.25	0.291 0.289	0.290
P7	15.625	0.225 0.221	0.223
P8	0.0	0.129 0.123	0.126

## Standard Curve

- The curve is provided for illustration only. A standard curve should be generated each time the assay is performed.

### Human Hexokinase-1 Standard Curve



### Performance Characteristics

- The minimum detectable dose of human hexokinase-1 as calculated by 2SD from the mean of a zero standard was established to be 9.5 ng/ml.

### Recovery

Standard Added Value	62.5 – 500 ng/ml
Recovery %	89 – 113%
<b>Average Recovery %</b>	<b>96%</b>

### Cross-Reactivity

Protein	Cross-Reactivity (%)
Hexokinase-2	45%
Hexokinase-3	65%
Hexokinase-4	30%
Ketohexokinase	None

## Troubleshooting

Issue	Causes	Course of Action
Low Precision	Use of expired components	<ul style="list-style-type: none"> <li>Check the expiration date listed before use.</li> <li>Do not interchange components from different lots.</li> </ul>
	Improper wash step	<ul style="list-style-type: none"> <li>Check that the correct wash buffer is being used.</li> <li>Check that all wells are empty after aspiration.</li> <li>Check that the microplate washer is dispensing properly.</li> <li>If washing by pipette, check for proper pipetting technique.</li> </ul>
	Splashing of reagents while loading wells	<ul style="list-style-type: none"> <li>Pipette properly in a controlled and careful manner.</li> </ul>
	Inconsistent volumes loaded into wells	<ul style="list-style-type: none"> <li>Pipette properly in a controlled and careful manner.</li> <li>Check pipette calibration.</li> <li>Check pipette for proper performance.</li> </ul>
	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none"> <li>Thoroughly agitate the lyophilized components after reconstitution.</li> <li>Thoroughly mix dilutions.</li> </ul>
	Improperly sealed microplate	<ul style="list-style-type: none"> <li>Check the microplate pouch for proper sealing.</li> <li>Check that the microplate pouch has no punctures.</li> <li>Check that three desiccants are inside the microplate pouch prior to sealing.</li> </ul>
Unexpectedly Low or High Signal Intensity	Microplate was left unattended between steps	<ul style="list-style-type: none"> <li>Each step of the procedure should be performed uninterrupted.</li> </ul>
	Omission of step	<ul style="list-style-type: none"> <li>Consult the provided procedure for complete list of steps.</li> </ul>
	Steps performed in incorrect order	<ul style="list-style-type: none"> <li>Consult the provided procedure for the correct order.</li> </ul>
	Insufficient amount of reagents added to wells	<ul style="list-style-type: none"> <li>Check pipette calibration.</li> <li>Check pipette for proper performance.</li> </ul>
	Wash step was skipped	<ul style="list-style-type: none"> <li>Consult the provided procedure for all wash steps.</li> </ul>
	Improper wash buffer	<ul style="list-style-type: none"> <li>Check that the correct wash buffer is being used.</li> </ul>
	Improper reagent preparation	<ul style="list-style-type: none"> <li>Consult reagent preparation section for the correct dilutions of all reagents.</li> </ul>
	Insufficient or prolonged incubation periods	<ul style="list-style-type: none"> <li>Consult the provided procedure for correct incubation time.</li> </ul>
Deficient Standard Curve Fit	Non-optimal sample dilution	<ul style="list-style-type: none"> <li>Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay.</li> <li>Competitive ELISA: If samples generate OD values lower than the highest standard point (P1), dilute samples further and repeat the assay.</li> <li>User should determine the optimal dilution factor for samples.</li> </ul>
	Contamination of reagents	<ul style="list-style-type: none"> <li>A new tip must be used for each addition of different samples or reagents during the assay procedure.</li> </ul>
	Contents of wells evaporate	<ul style="list-style-type: none"> <li>Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.</li> </ul>
	Improper pipetting	<ul style="list-style-type: none"> <li>Pipette properly in a controlled and careful manner.</li> <li>Check pipette calibration.</li> <li>Check pipette for proper performance.</li> </ul>

	Insufficient mixing of reagent dilutions	<ul style="list-style-type: none"><li>• Thoroughly agitate the lyophilized components after reconstitution.</li><li>• Thoroughly mix dilutions.</li></ul>
--	--	---

## References

- (1) Printz RL *et al.* (1997) *Biochemical Society Transactions*. 25(1):107-112.
- (2) Nishi S *et al.* (1988) *Biochem Biophys Res Commun*. 157(3):937-943.
- (3) Aleshin AE *et al.* (1998) *J Mol Biol*. 282(2):345-357.

Version 1.0