

# PRODUCT INFORMATION & MANUAL

# Phosphofructokinase/PFK Activity Assay Kit (Colorimetric) NBP3-25908

For research use only.

Not for diagnostic or therapeutic procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

# Phosphofructokinase/PFK Activity Assay Kit (Colorimetric)

Catalog No: NBP3-25908

Method: Colorimetric method

Specification: 96T (Can detect 96 samples without duplication)

Instrument: Microplate reader

Sensitivity: 0.27 U/L

Detection range: 0.27-32.29 U/L

Average intra-assay CV (%): 3.0

Average inter-assay CV (%): 8.0

Average recovery rate (%): 100

- ▲ This kit is for research use only.
- ▲ Instructions should be followed strictly, changes of operation may result in unreliable results.
- ▲ Please kindly provide us the lot number (on the outside of the box) of the kit for more efficient service.

# General information

#### ▲ Intended use

This kit can be used to measure phosphofructokinase (PFK) activity in serum, plasma, tissue and cell samples.

# **▲ Detection principle**

Phosphofructokinase (PFK), also known as 6-phosphofructokinase, is a class of kinases that can act on fructose-6-phosphate. Phosphofructokinase catalyzes fructose-6-phosphate and ATP to produce fructose-1, 6-diphosphate and ADP, then pyruvate kinase and lactate dehydrogenase further catalyze the oxidation of NADH to NAD<sup>+</sup>. The activity of PFK can be calculated by measuring the change of absorbance value at 340 nm.

# ▲ Kit components & storage

| Item      | Component       | Specification    | Storage                          |  |
|-----------|-----------------|------------------|----------------------------------|--|
| Reagent 1 | Buffer Solution | 20 mL × 1 vial   | -20°C , 12 months, shading light |  |
| Reagent 2 | Substrate A     | Powder × 2 vials | -20°C , 12 months shading light  |  |
| Reagent 3 | Substrate B     | Powder × 2 vials | -20°C , 12 months, shading light |  |
| Reagent 4 | Enzyme Reagent  | Powder × 2 vials | -20°C , 12 months, shading light |  |
|           | UV Microplate   | 96 wells         | No requirement                   |  |
|           | Plate Sealer    | 2 pieces         |                                  |  |

Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other.

# ▲ Materials prepared by users



# **1** Instruments

Microplate reader (330-350 nm, optimum wavelength: 340nm)



#### Reagents:

Normal saline (0.9% NaCl)

#### **▲** Safety data

Some of the reagents in the kit contain dangerous substances. It should be avoided to touch the skin and clothing. Wash immediately with plenty of water if touching it carelessly. All the samples and waste material should be treated according to the relevant rules of laboratory's biosafety.

#### **▲ Precautions**

Before the experiment, please read the instructions carefully, and wear gloves and work clothes.

#### ▲ The key points of the assay

- 1.The working solution should be stored at -20°C with shading light and avoid repeated freeze-thaw.
- 2. All reagents should be stored with shading light strictly.

# **Pre-assay preparation**

#### ▲ Reagent preparation

- 1. Bring all reagents to room temperature before use.
- 2. Preparation of working solution:

Dissolve a vial of reagent 2 powder and reagent 3 powder with 10 mL reagent 1. The prepared solution can be divided into smaller packages at -20°C with shading light for 3 days.

3. Preparation of reagent 4 working solution:

Dissolve a vial of reagent 4 with 1.2 mL of double distilled water. Preserve it on the ice box with shading light for use and the prepared solution can be stored at -20°C with shading light for 7 days.

#### **▲** Sample preparation

#### 1. Serum and plasma samples:

Detect directly (If the sample is turbid, centrifuge at 12000 g for 10 min before detection).

#### 2. Tissue sample:

Weigh the tissue accurately and add normal saline (0.9% NaCl) at a ratio of weight (g): volume (mL) =1: 9, homogenize the tissue in ice bath, centrifuge at 10000 g for 10 min at 4°C, then take the supernatant for measurement. The supernatant after centrifugation must be clarified, and if there is turbidity, it must be centrifuged again. Meanwhile, determine the protein concentration of supernatant.

#### 3. Cell sample:

Collect the  $1\times10^6$  cells, add 200 µL normal saline (0.9% NaCl). Homogenize the cells sample with homogenizer on ice. Centrifuge the homogenized cells at 10000 g for 10 min, then take the supernatant and preserve it on ice for detection. Meanwhile, determine the protein concentration of supernatant.

# **▲ Dilution of sample**

It is recommended to take 2~3 samples with expected large difference to do pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (0.27-32.29 U/L).

The recommended dilution factor for different samples is as follows (for reference only)

| Sample type                        | Dilution factor |  |  |  |
|------------------------------------|-----------------|--|--|--|
| 10% Rat liver tissue homogenate    | 1               |  |  |  |
| 10% Rat kidney tissue homogenate   | 1               |  |  |  |
| 10% Rat brain tissue homogenate    | 1               |  |  |  |
| 10% Mouse liver tissue homogenate  | 1               |  |  |  |
| 10% Mouse spleen tissue homogenate | 1               |  |  |  |
| 10% Mouse kidney tissue homogenate | 1               |  |  |  |
| 10% Mouse heart tissue homogenate  | 1               |  |  |  |
| Rat serum                          | 1               |  |  |  |
| Rat plasma                         | 1               |  |  |  |
| 1×10 <sup>6</sup> Jurkat cell      | 1               |  |  |  |

Note: The diluent is normal saline (0.9% NaCl).

# **Assay protocol**

# ▲ Plate set up

|   | 1  | 2   | 3   | 4   | 5   | 6   | 7   | 8   | 9   | 10  | 11  | 12  |
|---|----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| Α | S1 | S9  | S17 | S25 | S33 | S41 | S49 | S57 | S65 | S73 | S81 | S89 |
| В | S2 | S10 | S18 | S26 | S34 | S42 | S50 | S58 | S66 | S74 | S82 | S90 |
| С | S3 | S11 | S19 | S27 | S35 | S43 | S51 | S59 | S67 | S75 | S83 | S91 |
| D | S4 | S12 | S20 | S28 | S36 | S44 | S52 | S60 | S68 | S76 | S84 | S92 |
| E | S5 | S13 | S21 | S29 | S37 | S45 | S53 | S61 | S69 | S77 | S85 | S93 |
| F | S6 | S14 | S22 | S30 | S38 | S46 | S54 | S62 | S70 | S78 | S86 | S94 |
| G | S7 | S15 | S23 | S31 | S39 | S47 | S55 | S63 | S71 | S79 | S87 | S95 |
| Н | S8 | S16 | S24 | S32 | S40 | S48 | S56 | S64 | S72 | S80 | S88 | S96 |

Note:S1-S96, sample wells.

# ▲ Detailed operation steps

- (1) Sample well: Add 10 µL of sample to the wells.
- (2) Add 20  $\mu$ L of reagent 4 working solution and 170  $\mu$ L of working solution into each well.
- (3) Measure the OD value of each well at 20 s and 5 min 20 s respectively at 340 nm with microplate reader, recorded as  $A_1$ ,  $A_2$ ,  $\Delta A = A_1 A_2$ .

# **▲** Summary operation table

|  | Sample well |  |  |  |
|--|-------------|--|--|--|
| Sample (µL)  | 10          |  |  |  |
| Reagent 4 working solution (µL)  | 20          |  |  |  |
| Working solution (µL)  | 170         |  |  |  |
| Measure the OD value of each well at 20 s and 5 min 20 s respectively, |             |  |  |  |
| recorded as $A_1$ , $A_2$ , $\Delta A = A_1 - A_2$ .                   |             |  |  |  |

#### **▲** Calculation

#### 1. Serum/plasma samples:

Definition: The enzyme amount of 1 µmol of NADH consumed by 1 L of liquid sample per minute at room temperature is defined as 1 unit.

PFK activity (U/L) = 
$$\Delta A_{340} \div (6220 \times d) \div T \times f \times 10^6$$

#### 2. Tissue and cell samples:

Definition: The enzyme amount of 1 µmol of NADH consumed by 1 g sample protein per minute at room temperature is defined as 1 unit.

PFK activity (U/gprot) = 
$$\Delta A_{340} \div (6220 \times d) \div C_{pr} \div T \times f \times 10^6$$

#### Note:

 $\Delta A_{340}$ :  $A_1 - A_2$ .

6220: The molar extinction coefficient of NADH, L/mol•cm

d: Optical path, 0.6 cm

C<sub>pr</sub>: Concentration of protein in sample, gprot/L.

T: The time of reaction, 5 min.

f: Dilution factor of sample before test.

10^6: 1 moL/L=1000000 μmol/L

# **Appendix I Data**

#### **▲ Example analysis**

For rat liver tissue, take 10  $\mu$ L of 10% rat liver tissue homogenate, and carry the assay according to the operation table.

#### The results are as follows:

the  $A_1$  of the sample is 1.245, after 5 minutes of reaction, the  $A_2$  of the sample is 0.849, the concentration of protein in sample is 14.45 gprot/L, and the calculation result is:

PFK activity (U/gprot) = 
$$(1.245 - 0.849) \div (6220 \times 0.6) \div 14.45 \div 5 \times 10^6$$
  
= 1.47 U/gprot