

β-N-Acetylglucosaminidase (NAG) Activity Assay Kit

Instruction Manual

Catalog Number	PK-CA577-K733		
Description	<p>β-N-Acetylglucosaminidase (NAG, EC 3.2.1.52) is a lysosomal enzyme that is expressed in various tissues, including kidney, liver and lungs. NAG can cleave N-acetyl-glucosamine, a monosaccharide derivative of glucose. Its concentration in urine is minimal due to its inability to cross the glomerular basal membrane. Increased concentration of NAG in urine indicates renal tubular cell breakdown. Acute Kidney Injury (AKI) is the sudden loss of kidney functions, causing electrolyte imbalance, and retention of urea and other nitrogenous products. NAG has become one of the most studied and used biomarkers for the detection and diagnosis of AKI.</p> <p>PromoKines's NAG Activity Assay Kit provides a simple and sensitive method for monitoring NAG enzymatic activity. In this assay, NAG uses a synthetic p-nitrophenol derivative (R-pNP) as a NAG substrate and releases pNP which can be measured at absorbance (OD 400 nm). The assay can detect as low as 50 μU of NAG activity in a variety of samples.</p> <p style="text-align: center;"> $\text{R-pNP} \xrightarrow{\text{NAG; [H+]}} \text{R + pNP} \xrightarrow{\text{Stop Solution}} \text{pNP (OD: 400 nm)}$ </p>		
Quantity	100 assays		
Applications	<ul style="list-style-type: none"> • Measurement of NAG in biological samples from different mammalian species 		
Sample Type	<ul style="list-style-type: none"> • Biological fluids such as urine, serum etc. • Tissues such as kidney, liver, brain etc. • Cells such as HeLa, HepG2, CHO etc. 		
	<p>Components</p> <p>NAG Assay Buffer</p> <p>NAG Substrate</p> <p>p-Nitrophenol (pNP) (20 mM)</p> <p>NAG Stop Solution</p> <p>NAG Positive Control</p>	<p>Quantity</p> <p>35 ml</p> <p>6 ml</p> <p>0.1 ml</p> <p>3 ml</p> <p>1 vial</p>	<p>Color Code</p> <p>NM</p> <p>Amber</p> <p>Yellow</p> <p>Clear</p> <p>Blue</p>
User Supplied Reagents & Equipment	<ul style="list-style-type: none"> • Multiwell absorbance reader (ELISA reader) • 96-well clear plate with flat bottom 		
Storage and Reagents Preparation	<p>Store kit at -20°C, protected from light. Briefly centrifuge all small vials prior to opening. Read entire protocol before performing the assay procedure.</p> <ul style="list-style-type: none"> • NAG Assay Buffer: Bring to room temperature (RT) before use. Store at 4°C or -20°C. • NAG Substrate: Light sensitive. Store at -20°C. Once opened, use within two months. Bring to RT before use. Mix well. If precipitation is observed, sonicate the contents in a water bath sonicator (interval: 2 minutes). Repeat if necessary. • p-Nitrophenol (pNP): Light sensitive. Bring to RT before use. Store at -20°C. Once opened, use within two months. • NAG Stop Solution: Bring to RT before use. Store at -20 °C. Once opened, use within two months. • NAG Positive Control: Reconstitute with 40 μl NAG Assay Buffer. Mix well. Store at -20°C. Use within two months. 		
Assay Protocol	<p>1. Sample Preparation: Serum samples can be used directly. Centrifuge urine samples at 10,000 x g, 4°C for 3 minutes, if precipitation is observed. Collect supernatant. Homogenize tissues (~10-20 mg) or cells (~1x10⁶) on ice using 100-200 μl NAG Assay Buffer. Centrifuge homogenate at 10,000 x g, 4°C for 3 minutes. Collect supernatant. Add 1-70 μl sample into desired well(s) in a 96-well plate. Prepare a parallel sample well as sample background control. For Positive Control, add 1-5 μl of Positive Control into desired well(s). Adjust the volume of Positive Control, background control and sample wells to 70 μl/well with NAG Assay Buffer.</p> <p>Note:</p> <p>NAG enzymatic activity varies over a wide range depending on the sample. For unknown samples, we recommend doing a pilot experiment and testing several doses to ensure the</p>		

readings are within the Standard Curve linear range.

2. Standard Curve Preparation: Prepare 2 mM pNP Standard by adding 10 µl of 20 mM pNP Standard into 90 µl of NAG Assay Buffer. Add 0, 2, 4, 6, 8, and 10 µl of 2 mM pNP Standard into a series of wells in a 96-well plate to generate 0, 4, 8, 12, 16 and 20 nmol of pNP/well respectively. Adjust the volume to 125 µl/well with NAG Assay Buffer.

3. Substrate Hydrolysis: Add 55 µl NAG Substrate into each well containing samples, and positive controls. To the sample background control well, add 55 µl NAG Assay Buffer. Mix well. Incubate for 5-30 minutes at 37°C. After incubation, add 25 µl NAG Stop Solution into each well containing samples, Positive Control, background control and Standards. Mix well.

Note:

Incubation time depends on the NAG enzymatic activity in samples. Longer incubation time may be required for samples having low NAG activity.

4. Measurement: To stabilize the reaction, incubate plate at 37°C for 10 minutes, protected from light. Measure absorbance (OD 400 nm) with end point setting.

5. Calculation: Subtract 0 Standard reading from all readings. Plot the pNP Standard Curve. If sample background control reading is significant, then subtract sample background control reading from sample reading. Apply sample corrected OD to pNP Standard Curve to get B nmol of pNP generated by NAG activity in the sample well.

Sample NAG Activity = (BXD)/(TXV) nmol/min/ml or mU/ml or U/L

Where: B is amount of pNP in the sample well from Standard Curve (nmol)

T is incubation time for substrate hydrolysis (5-30 minutes)

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

D is sample dilution factor

Sample NAG activity can also be expressed as mU/mg of protein or mU/mg creatinine (urine).

Unit Definition: One unit of NAG activity is the amount of enzyme that generates 1 µmol of pNP per minute at pH 4.2 at 37°C.

Intended Use

For in vitro research use only. Not for diagnostic or therapeutic procedures.

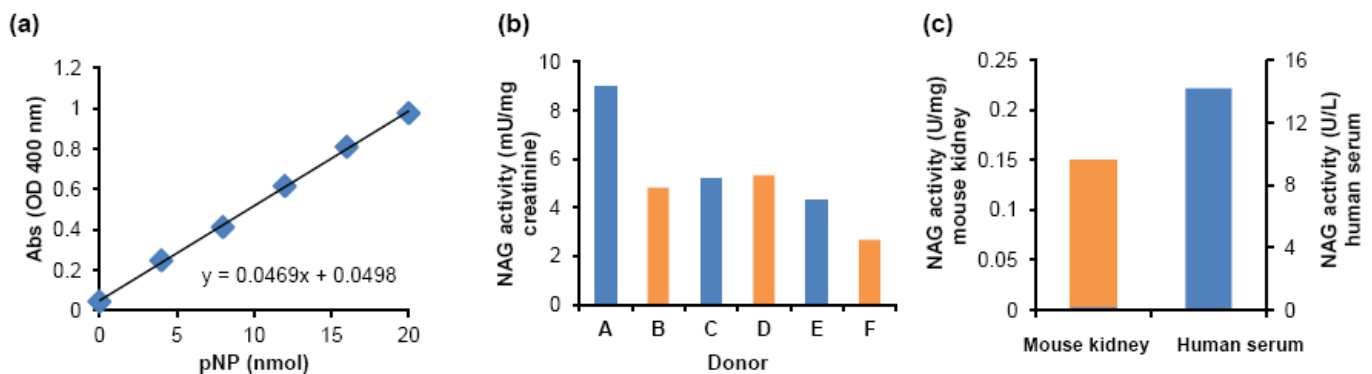


Figure: (a) pNP Standard Curve. (b) Measurement of NAG activity in human urine from different donors. Undiluted samples (70 µl) were incubated for 30 minutes with NAG substrate. (c) Measurement of NAG activity in mouse kidney (10 µg) and human serum (20 µl). Samples were incubated for 30 minutes with NAG substrate. All assays were performed following the kit protocol.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

USA/Canada

Phone: 1 – 866 – 251 – 2860 (toll free)
Fax: 1 – 866 – 827 – 9219 (toll free)

Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)
Fax: 0800 – 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 – 96 03 33 (toll free)
Fax: 0800 – 169 85 54 (toll free)

Other Countries

Phone: +49 6221 – 649 34 0
Fax: +49 6221 – 649 34 40