

Product Information

β -N-Acetylglucosaminidase Assay Kit

Product Code **CS0780**

Storage Temperature 2–8 °C

TECHNICAL BULLETIN

Product Description

β -N-Acetylglucosaminidase (NAG, N-acetyl- β -D-glucosaminidase) is a lysosomal enzyme in mammalian cells, which takes part in intracellular degradation of glycolipids and glycoproteins.¹ High activities of this enzyme have been detected in human kidney, lung, and liver lysosomes. In humans, different genetically determined deficiencies of the enzyme result in lipid storage disorders.^{2,3} Moreover, elevated levels of serum NAG are associated with certain forms of cancer,⁴ and the increase in total NAG activity in the urine is associated with renal disorder.⁵

The assay is based on the hydrolysis of the NAG substrate, 4-Nitrophenyl N-acetyl- β -D-glucosaminide (NP-GlcNAc), by the enzyme. This enzymatic hydrolysis of the substrate releases *p*-nitrophenol, which upon ionization in basic pH, can be measured colorimetrically at 405 nm.

The β -N-Acetylglucosaminidase Assay Kit provides all the reagents required for efficient detection of β -N-acetylglucosaminidase activity in cell lysates, tissue homogenates, lysosomal isolated fractions, or purified enzyme preparations.

The kit was tested on CHO, HeLa, HEK 239T, A431, HepG2, U937, and Jurkat cells, and on rat brain, spleen, kidney, and liver tissues.

Components

The kit is sufficient for fifty 1 ml reactions or 500 multiwell plate reactions of 100 μ l.

Dilution Buffer Product Code D2318	8 ml
4-Nitrophenyl N-acetyl- β -D-glucosaminide Product Code N9376	50 mg
Citrate Buffer Solution, 0.09 M Product Code C2488	100 ml

<i>p</i> -Nitrophenol Standard Solution, 10 mM Product Code N7660	1 ml
--	------

β -N-Acetylglucosaminidase from Jack beans Product Code A2264	1 vial
---	--------

Sodium Carbonate Product Code S2127	5 g
--	-----

Equipment required but not provided

- Spectrophotometer or ELISA reader (405 nm)
- Cuvettes (3 ml, Product Code C5291), or 96 well plates (flat bottom, Product Code P7366)
- 37 °C water bath

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Use ultrapure water (17 M Ω .cm or equivalent) for preparation of reagents.

Substrate Solution (1 mg/ml) - Dissolve 10 mg of the 4-Nitrophenyl N-acetyl- β -D-glucosaminide (NP-GlcNAc, Product Code N9376) in 10 ml of the 0.09 M Citrate Buffer Solution (Product code C2488). Mix the Substrate Solution with rocking/shaking on a horizontal shaker at room temperature.

Note: The substrate does not easily dissolve in the buffer. It may take ~1 hour of shaking to completely dissolve the substrate. Use of a larger container (50 ml) may aid dissolution.

The Substrate Solution should be stored on ice during the experiment. For longer storage of at least one month, store at –20 °C.

NAG Control Enzyme - Just before use, dilute a portion of the β -N-Acetylglucosaminidase (NAG, Product Code A2264) 1:100 in the Dilution Buffer (Product Code D2318). The enzyme is supplied as an ammonium sulfate suspension, therefore, vortex the tube vigorously and pipette up and down a few times to ensure a uniform suspension before diluting the enzyme. Store the NAG Control Enzyme on ice. Do not freeze.

Note: If the signal obtained with the NAG Control Enzyme is too high/low, adjust the dilution accordingly.

Stop Solution (sodium carbonate solution) - Add 118 ml of ultrapure water to the sodium carbonate bottle (Product Code S2127) and mix well with a magnetic stirrer until completely dissolved. Store the Stop Solution at room temperature.

Standard Solution - Before performing the assay in 96 well plates, dilute 5 μ l of the 10 mM *p*-Nitrophenol Standard Solution (Product Code N7660) with 995 μ l of Stop Solution. Vortex briefly and store on ice.

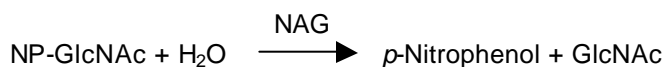
NAG Sample – The use of the CellLytic™ M (Product Code C2978) or CellLytic MT (Product Code C3228) Cell Lysis reagents is recommended for cell or tissue sample preparation, respectively.

Storage/Stability

The kit is shipped on wet ice and storage at 2–8 °C is recommended. Upon arrival it is recommended to store the NP-GlcNAc (Product Code N9376) at –20 °C and the Sodium Carbonate (Product Code S2127) at room temperature.

Procedure

The NAG enzyme hydrolyzes the NP-GlcNAc substrate under acidic conditions (pH ~4.7) at 37 °C. The enzymatic hydrolysis of the substrate releases *p*-nitrophenol.



Addition of the basic Stop Solution causes ionization of the *p*-nitrophenol to form the yellow *p*-nitrophenylate ion. The absorbance of the *p*-nitrophenylate ion is measured at 405 nm.^{6,7}

I. Assay in Tubes (3 ml final volume)

1. Equilibrate the Substrate Solution to 37 °C by incubating for several minutes.
2. Set the spectrophotometer at 405 nm.
3. Add the reaction components to the tubes according to Table 1 and vortex the tubes briefly.

Table 1.

Reaction Scheme for Tube Assays

	Substrate Solution	Sample
Blank*	1 ml	----
Positive control	995 μ l	5 μ l of NAG Control Enzyme
Test	900-990 μ l	10-100 μ l of sample

* A blank reaction (Substrate Solution without enzyme) should be run, since a portion of the NP-GlcNAc may hydrolyze spontaneously during the incubation time.

Note: If required (the absorbance is too low), the volume of the sample/NAG Control Enzyme may be as large as 10% of the reaction volume.

4. Incubate the tubes for 5-30 minutes at 37 °C. The incubation time required for the test sample(s) depends on the enzyme concentration. The incubation time required for the positive control is usually 5-10 minutes.
5. Stop the reaction by adding 2 ml of Stop Solution. Vortex briefly and transfer the reaction mixture to a cuvette. The reaction mixture should be yellow in color.
6. Use the blank reaction mixture to zero the spectrophotometer and then read the absorbance of the samples.

Note: Read the samples immediately after the incubation, within 30 minutes.

II. 96 Well Plate Assay

Note: It is recommended to perform the assays in triplicates.

1. Equilibrate the Substrate Solution and the Standard Solution to 37 °C by incubating for several minutes.
2. Set the plate reader at 405 nm.
3. Add the reaction components to the 96 well plates according to Table 2 and mix using a horizontal shaker or by pipetting.

Table 2.

Reaction Scheme for 96 Well Plate Assays

	Substrate Solution	Sample	Standard Solution
Blank*	100 µl	-	-
Standard	-	-	300 µl
Positive control**	98 µl	2 µl of NAG Control Enzyme	-
Test	90-99 µl	1-10 µl of sample	-

* A blank reaction (Substrate Solution without enzyme) should be run, since a portion of the NP -GlcNAc may hydrolyze spontaneously during the incubation time.

** If required (the absorbance is too low), the volume of the enzyme can be increased up to 10% (10 µl) of the total reaction volume.

4. Incubate the plate for 5-10 minutes at 37 °C. If you suspect that the NAG activity of the test sample may be low, the incubation time can be extended up to 30 minutes.
5. Stop the reactions by adding 200 µl of Stop Solution to each well, except to wells containing the Standard Solution.
6. Measure the absorption at 405 nm.

Calculations

Unit definition: 1 unit will hydrolyze 1.0 µmole of 4-Nitrophenyl N-acetyl-β-D-glucosaminide to *p*-nitrophenol and N-acetyl-β-D-glucosaminide per 1 minute at pH 4.7 at 37 °C.

A. Activity calculation for cuvette assay:

$$\text{Units/ml} = \frac{A_{405\text{sample}} \times 3 \times \text{DF}}{18.3 \times \text{time} \times V_{\text{enz}}}$$

$A_{405\text{sample}}$ – absorbance of the sample at 405 nm

3 – final volume in milliliters in the reaction cuvette after addition of stop solution

DF – enzyme dilution factor

18.3 – millimolar extinction coefficient for *p*-nitrophenolate ion

time – minutes

V_{enz} – volume of the sample in milliliter

B. Activity calculation for 96 well plate assay:

$$\text{Units/ml} = \frac{(A_{405\text{sample}} - A_{405\text{blank}}) \times 0.05 \times 0.3 \times \text{DF}}{A_{405\text{standard}} \times \text{time} \times V_{\text{enz}}}$$

$A_{405\text{sample}}$ – absorbance of the sample at 405 nm

$A_{405\text{blank}}$ – absorbance of the blank at 405 nm

0.05 – µmole/ml of 4-nitrophenol in the Standard Solution

0.3 – final volume in milliliters of the 96 well plate reaction after addition of the stop solution

DF – enzyme dilution factor

$A_{405\text{standard}}$ – absorbance of the Standard Solution at 405 nm

time – minutes

V_{enz} – volume of the sample in milliliter

References

1. Legler, G., *et al.*, Bovine N-acetyl-beta-D-glucosaminidase: affinity purification and characterization of its active site with nitrogen containing analogs of N-acetylglucosamine. *Biochim. Biophys. Acta*, **1080**, 89-95 (1991).
2. Myerowitz, R., *et al.*, Molecular pathophysiology in Tay-Sachs and Sandhoff diseases as revealed by gene expression profiling. *Hum. Mol. Genet.*, **11**, 1343-1350 (2002).
3. Olah, A.V., *et al.*, Age dependence of serum β -N-acetylhexosaminidase (NAG) activity. *Clin. Chem. Lab. Med.*, **42**, 305-306 (2004).
4. Gil-Martin, E., *et al.*, N-acetyl-beta-hexosaminidase activity and isoenzymes in human gastric adenocarcinoma. *Oncology*, **56**, 142-154 (1999).
5. Price, R.G., Urinary N-acetyl-beta-D-glucosaminidase (NAG) as an indicator of renal disease. *Curr. Probl. Clin. Biochem.*, **9**, 150-163 (1979).
6. Horak, E., *et al.*, Spectrophotometric assay for urinary N-acetyl-beta-D-glucosaminidase activity. *Clin. Chem.*, **27**, 1180-1185 (1981).
7. Shibata, H., and Yagi, T., Rate assay of N-acetyl-beta-D-hexosaminidase with 4-nitrophenyl N-acetyl-beta-D-glucosaminide as an artificial substrate. *Clin. Chim. Acta*, **251**, 53-64 (1996).

EB,MAM 10/05-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.