

Product Information

Chitinase Assay Kit

Catalog Number **CS0980**

Storage Temperature 2–8 °C

TECHNICAL BULLETIN

Product Description

Chitinase catalyzes the hydrolytic cleavage of the $\beta(1\rightarrow4)$ -glycoside bonds present in biopolymers of N-acetylglucosamine, primarily in chitin. Chitinases are widely distributed in living organisms and are found in fungi, bacteria, parasites, plants, and animals.¹ They are classified in families based on amino acid sequence similarities.²

The chitinolytic enzymes are also categorized based on their enzymatic action on chitin substrates. Endochitinases are defined as the enzymes catalyzing the random cleavage at internal points in the chitin chain. Exochitinases catalyze the progressive release of acetylchitobiose or N-acetylglucosamine from the non-reducing end of chitin, and thus, are referred to as chitobiosidase and β -N-acetylglucosaminidase, respectively.^{3,4}

Chitinases perform different functions in different organisms. In bacteria they are mainly involved in nutritional processes, while in yeast and various fungi, these enzymes participate in morphogenesis. In animals and plants, chitinases primarily play a role in the defense of the organism against pathogen attack.

Human chitotriosidase (Chit), a chitinolytic enzyme, is a member of the chitinase family. In human plasma, Chit activity has been proposed as a biological marker of macrophage activity in several lysosomal diseases, and was found at higher levels in patients with *Plasmodium falciparum* malaria infection. This suggests that Chit induction may reflect an immunological response.⁵

Another member of the chitinase family is the acid mammalian chitinase (AMCase), thought to play a role in inflammatory disorders. Elevated levels of ACMase were observed in lung tissue of asthmatic patients, suggesting a role for this enzyme in asthma pathophysiology.⁶

The kit assay is based on the enzymatic hydrolysis of chitinase substrates. This hydrolysis releases *p*-nitrophenol (4-nitrophenol), which upon ionization in basic pH, can be measured colorimetrically at 405 nm.⁷

The Chitinase Assay Kit provides all the reagents required for efficient detection of chitinase activity in fungal and bacterial growth media, macrophage lysates, and purified enzyme preparations. In addition, the kit provides three different substrates for the detection of the various types of the chitinolytic activity:³

4-Nitrophenyl N,N'-diacetyl- β -D-chitobioside – a substrate suitable for exochitinase activity detection (chitobiosidase activity)

4-Nitrophenyl N-acetyl- β -D-glucosaminide – a substrate suitable for exochitinase activity detection (β -N-acetylglucosaminidase activity)

4-Nitrophenyl β -D-N,N',N''-triacetylchitotriose – a substrate suitable for endochitinase activity detection

The kit was tested on *Trichoderma viride*, *Streptomyces griseus*, and human macrophages.

Components

Assay Buffer	20 ml
Catalog Number A4855	
4-Nitrophenyl N-acetyl- β -D-glucosaminide	10 mg
Catalog Number N9376	
4-Nitrophenyl N,N'-diacetyl- β -D-chitobioside	5 mg
Catalog Number N6133	

4-Nitrophenyl β -D-N,N',N''-triacetylchitotriose Catalog Number N8638	1 mg
Chitinase from <i>Trichoderma viride</i> Catalog Number C6242	1 mg
<i>p</i> -Nitrophenol Solution, 10 mM Catalog Number N7660	1 ml
Sodium Carbonate Catalog Number S2127	1 g

The kit is sufficient for the number of multiwell plate reactions indicated for each substrate:

Catalog Number N9376 at a substrate concentration of 1 mg/ml is sufficient for 100 reactions.

Catalog Number N6133 at a substrate concentration of 1 mg/ml is sufficient for 50 reactions.

Catalog Number N8638 at a substrate concentration of 1 mg/ml is sufficient for 10 reactions.

Note: For other substrate concentrations see Substrate Solution(s) Notes under Preparation Instructions.

Reagents and Equipment Required but Not Provided

- Dulbecco's Phosphate Buffered Saline (PBS, Catalog Number D8537)
- Ultrapure water
- Spectrophotometer or ELISA reader (405 nm)
- 96 well plates (flat bottom, Catalog Number P7366)
- 37 °C water bath
- For macrophages lysis: CellLytic™ M Cell Lysis Reagent (Catalog Number C2978).

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.

Stop Solution (sodium carbonate solution) – Add 24 ml of ultrapure water to the contents of the sodium carbonate bottle (Catalog Number S2127) and mix well with a magnetic stirrer until completely dissolved. Store the Stop Solution at room temperature.

Substrate Solution(s) (1 mg/ml) – Dissolve 1 mg of the appropriate substrate in 1 ml of Assay Buffer (Catalog Number A4855):

- 4-Nitrophenyl N-acetyl- β -D-glucosaminide, Catalog Number N9376, or
- 4-Nitrophenyl β -D-N,N',N''-triacetylchitotriose, Catalog Number N8638 or
- 4-Nitrophenyl N,N'-diacetyl- β -D-chitobioside, Catalog Number N6133

1 ml of Substrate Solution is sufficient for ~10 reactions. Mix the Substrate Solution(s) with rocking/shaking on a horizontal shaker at room temperature.

Notes: A 1 mg/ml substrate solution is recommended. If required, the 4-Nitrophenyl β -D-N,N',N''-triacetylchitotriose and the 4-Nitrophenyl N,N'-diacetyl- β -D-chitobioside Substrate Solutions can be used at lower concentrations in the reaction (0.5 or 0.2 mg/ml, respectively).

The substrates do not dissolve easily in the buffer. It may take ~1 hour of shaking to completely dissolve the substrates. Use of a larger container (15 ml) may aid dissolution. The Substrate Solution(s) should be stored on ice during the experiment. For long term storage, up to one month, store at –20 °C.

Chitinase Control Enzyme – Add 5 ml of PBS to the contents of the chitinase bottle (Catalog Number C6242) to give a final chitinase concentration of 0.2 mg/ml. Vortex until dissolved. The chitinase dissolves immediately to give a slightly hazy solution. For long term storage, store in working aliquots at –20 °C (stable for at least 3 months at –20 °C). Just before use, dilute the chitinase 20-fold with PBS and keep on ice.

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

Sample preparation – *Trichoderma viride* and *Streptomyces griseus* growth medium can be sampled directly from the growing culture (since the chitinase is secreted into the growth medium) and can be used in the assay after a brief centrifugation to remove particles from the medium. Human macrophage proteins can be extracted with CellLytic M Cell Lysis Reagent (Catalog Number C2978).

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 4-Nitrophenyl N-acetyl- β -D-glucosaminide (Catalog Number N9376) and the Chitinase from *Trichoderma viride* (Catalog Number C6242) at –20 °C, and the Sodium Carbonate (Catalog Number S2127) at room temperature.

Procedure

The chitinase hydrolysis is performed in an acidic environment (pH ~4.8) at 37 °C.⁶ The enzymatic hydrolysis liberates *p*-nitrophenol.⁸ 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.³

In order to quantitate the total chitinolytic activity, separate reactions should be run with the three substrates supplied in the kit. Profiling of the chitinolytic enzymes can be determined after separation of the chitinolytic enzymes by SDS-PAGE, using an agarose overlay containing fluorescent substrates.^{3,9} Note that in crude preparations there may be additive/synergist activity of different chitinases.

It is recommended to perform the assays in duplicates. For each substrate, perform a separate activity assay according to the following instructions.

1. Equilibrate the Substrate Solution(s) and the Standard Solution to 37 °C by incubating for several minutes in a 37 °C water bath.
2. Set the plate reader at 405 nm.
3. Add the reaction components to the 96 well plate according to Table 1 and mix using a horizontal shaker or by pipetting. The substrate should be added first and the enzyme should be added last.

Table 1.

Reaction Scheme for 96 Well Plate Assays

	Substrate Solution	Sample	Standard Solution
Blank*	100 μ l	–	–
Standard**	–	–	300 μ l
Positive control***	90–99 μ l	1–10 μ l of Chitinase Control Enzyme	–
Test	90–99 μ l	1–10 μ l of sample	–

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

** A standard should be run when activity calculations are required.

*** The volume of the enzyme can range between 1–10 μ l, depending on the reaction duration (i.e., for a shorter time a higher volume of the enzyme is required). The positive control enzyme is the 20-fold diluted enzyme. If required, the concentrated, non-diluted enzyme may be used.

4. Incubate the plate for 30 minutes at 37 °C. If required, the incubation time for highly active samples can be reduced to as low as 5 minutes.
5. Stop the reactions by adding 200 μ l of Stop Solution to each well, except for the wells containing the Standard Solution. After the addition of the Stop Solution the reaction mixture will develop a yellow tint.
6. Measure the absorption at 405 nm no later than 30 minutes after ending the reaction.

Calculations

Unit definition: One unit will release 1.0 μ mole of *p*-nitrophenol from the appropriate substrate per minute at pH 4.8 at 37 °C.

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

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

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

0.05 – μ mole/ml of *p*-nitrophenol in the Standard Solution

0.3 – final volume of the 96 well plate reaction after addition of the Stop Solution (ml)

DF - Dilution Factor - fold dilution of the original chitinase enzyme or biological solution to prepare sample for the test

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

time – minutes

V_{enz} – volume of the sample (ml)

References

1. Kasprzewska, A., Plant Chitinases - Regulation and Function. *Cell. Mol. Biol. Lett.*, **8**, 809-824 (2003).
2. Henrrisat, B., New families in the classification of glycosyl hydrolases based on amino acid sequence similarities. *Biochem. J.*, **280**, 309-316 (1991).
3. Tronsmo, A., and Harman, G.E., Detection and quantification of N-acetyl-beta-D-glucosaminidase, chitobiosidase, and endochitinase in solutions and on gels. *Anal. Biochem.*, **208**, 74-79 (1993).
4. Sahai, A.S., and Manocha, M.S., Chitinases of fungi and plants: their involvement in morphogenesis and host-parasite interaction. *FEMS Microbiol. Rev.*, **11**, 317-338 (1993).
5. Malaguamera, L., *et al.*, Interferon- γ , tumor necrosis factor- α , and lipopolysaccharide promote chitotriosidase gene expression in human macrophages. *J. Clin. Lab. Anal.*, **19**, 128-132 (2005).
6. Donnelly, L.E., and Barnes, P.J., Acidic mammalian chitinase - a potential target for asthma therapy. *Trends Pharmacol. Sci.*, **25**, 509-511 (2004).
7. Duo-Chuan, L.I., *et al.*, Purification and partial characterization of two chitinases from the mycoparasitic fungus *Talaromyces flavus*. *Mycopathologia*, **159**, 223-229 (2005).
8. Frandberg, E., and Schnurer, J., Evaluation of chromogenic chito-oligosaccharide analogue, *p*-nitrophenyl- β -D-N,N'-diacetylchitobiose, for the measurements of the chitinolytic activity of bacteria. *J. Appl. Bacteriol.*, **76**, 259-263 (1994).
9. Chernin, L.S., *et al.*, Chitinolytic Activity in *Chromobacterium violaceum*: Substrate analysis and regulation by quorum sensing. *J. Bacteriol.*, **180**, 4435-4441 (1998).

CellLytic is a trademark of Sigma-Aldrich®
Biotechnology LP and Sigma-Aldrich Co.

KAA,ESS,EB,MAM 09/08-1

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.