

## Technical Bulletin

# Lysozyme Inhibitor Screening Kit

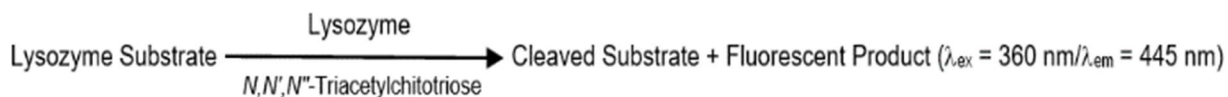
**Catalog Number MAK393****Product Description**

Lysozyme, also known as muramidase or N-acetylmuramide glycanhydrolase, is a glycoside hydrolase. It is ubiquitously found in a wide range of biological fluids such as tears, saliva, and serum, where it serves as a key effector of the innate immune system. It is also synthesized by certain carcinomas. The excessive production of lysozymes by cancer cells (especially myelomonocytic leukemia) results in higher levels of lysozyme which leads to renal failure. It has been demonstrated that the expression level of lysozyme is an independent prognostic factor which can be used to predict both relapse-free survival and overall survival in

patients with breast carcinomas. Therefore, lysozyme inhibitors may constitute attractive, potential targets for developing anti-inflammatory and/or anti-tumor drugs.

The Lysozyme Inhibitor Screening Kit uses *N,N',N''*-Triacetylchitotriose, a competitive inhibitor that binds to the active site of lysozyme. The inhibitor decreases the catalytic ability of lysozyme to hydrolyze a fluorogenic substrate. The kit provides a simple, rapid, sensitive and reliable test suitable for screening of lysozyme inhibitors.

The kit is suitable for the screening, studying and characterizing of lysozyme inhibitors.

**Components**

The kit is sufficient for 100 fluorometric assays in 96 well plates.

- Lysozyme Assay Buffer 25 mL  
Catalog Number MAK393A
- Lysozyme Substrate (in DMSO) 65  $\mu\text{L}$   
Catalog Number MAK393B
- Lysozyme 1 vial  
Catalog Number MAK393C
- *N,N',N''*-Triacetylchitotriose 1 vial  
Catalog Number MAK393D

**Reagents and Equipment  
Required but Not Provided**

- Pipetting devices and accessories (including multichannel pipettor)
- 96-well opaque flat-bottom plate. Cell culture or tissue culture treated plates are **not** recommended.
- Fluorescence multiwell plate reader

## Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

## Storage/Stability

The kit is shipped on wet ice. Store components at -20 °C, protected from light. Briefly centrifuge small vials at low speed prior to opening.

## Preparation Instructions

Lysozyme Assay Buffer: Warm to 37 °C before use. Store at either 2-8 °C or -20 °C.

Lysozyme Substrate (in DMSO): Aliquot and store at -20 °C. Bring to room temperature prior to use.

Lysozyme (lyophilized): Reconstitute vial with 1060 µL of Lysozyme Assay Buffer and pipette up and down to mix thoroughly. Aliquot and store at -20 °C. Avoid repeated freeze/thaw cycles. Keep on ice while in use. Stable for two months after reconstitution.

*N,N',N''*-Triacetylchitotriose: Reconstitute vial with 60 µL of purified water and store at -20 °C. Use within two months of reconstitution.

## Procedure

**Note:** To achieve better kinetic progress curves, pre-incubate the 96-well plate and assay buffer at 37 °C prior to using.

### Test Compound Preparation

1. Dissolve test inhibitors to 100× in appropriate solvent.
2. Further dilute 10-fold (10×) with Lysozyme Assay Buffer.
3. Add 10 µL of the diluted test inhibitor into designated wells (S).
4. Additional wells with serial dilutions of the test inhibitors may be prepared at this time, if desired, containing 10 µL in

each candidate well if IC<sub>50</sub> values need to be estimated.

### Lysozyme Enzyme Control

Add 10 µL of purified water to a well designated as Enzyme Control (EC).

### Inhibitor Control

Add 10 µL of *N,N',N''*-Triacetylchitotriose to a well designated as Inhibitor Control (IC).

### Solvent Control

High solvent concentration might affect lysozyme activity. Prepare parallel well(s) as Solvent Control (SC) to test the effect of the solvent on lysozyme activity where water is substituted with the final solvent concentration in the samples.

### Lysozyme Enzyme Solution Preparation

1. Mix enough reagents for the number of assays to be performed. For each well prepare 40 µL of Lysozyme Enzyme Solution according to Table 1. Mix well.

**Table 1.**

Preparation of Lysozyme Enzyme Solution

Reagent	Volume
Lysozyme Assay Buffer	30 µL
Lysozyme	10 µL

2. Add 40 µL of the Lysozyme Enzyme Solution into all wells, mix well.

### Lysozyme Substrate Solution Preparation

1. Prepare an 80-fold dilution of Lysozyme Substrate by diluting 4 µL of Lysozyme Substrate with 316 µL of Lysozyme Assay Buffer. Vortex briefly and keep on ice.
2. Add 50 µL of the freshly diluted substrate to each well. Mix well. Always use freshly prepared substrate solutions. Do **not** store the diluted substrate solution.



### Measurement

Measure fluorescence at  $\lambda_{Ex} = 360 \text{ nm}$ /  
 $\lambda_{Em} = 445 \text{ nm}$  in kinetic mode at  $37^\circ\text{C}$  for  
60 minutes. Choose two time points  
( $T_1$  and  $T_2$ ) in the linear range of the plot and  
obtain the corresponding fluorescence values  
( $RFU_1$  and  $RFU_2$ ).

### Results

Calculate the slope for all samples, including  
Enzyme Control (EC), by dividing the net  
 $\Delta RFU$  ( $RFU_2 - RFU_1$ ) values by the time  
 $\Delta T$  ( $T_2 - T_1$ ).

For Solvent Controls that differ substantially  
from the EC, use their values in the  
equations below instead of EC. Calculate %  
Relative Inhibition as follows:

% Relative Inhibition =

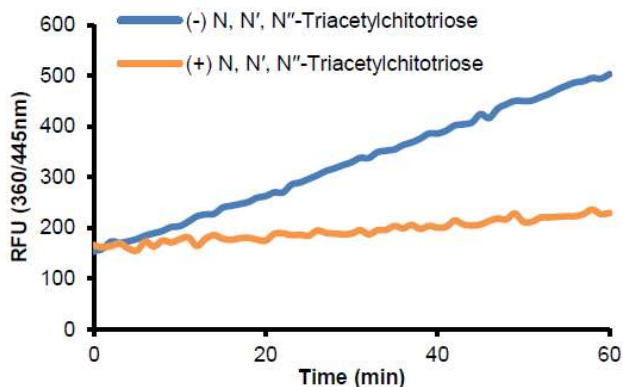
$$\frac{\text{Slope of [EC]} - \text{Slope of [S]}}{\text{Slope of [EC]}} \times 100\%$$

% Relative Activity =

$$\frac{\text{Slope of [S]}}{\text{Slope of [EC]}} \times 100\%$$

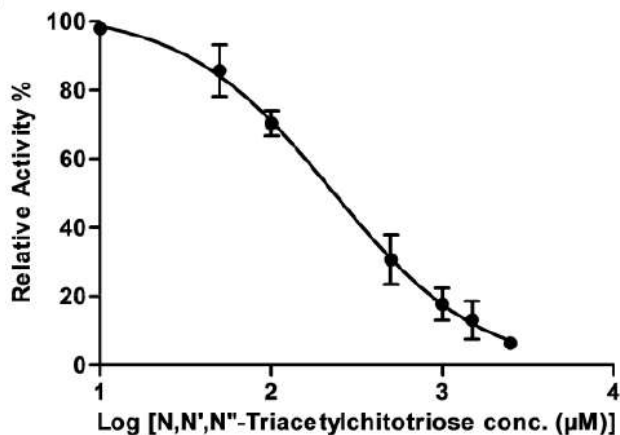
**Figure 1.**

Progress curve of lysozyme activity in the  
presence or absence of the inhibitor  
 $N,N',N''$ -Triacetylchitotriose.



**Figure 2.**

$IC_{50}$  of  $N,N',N''$ -Triacetylchitotriose was  
calculated to be  $227.5 \pm 1.2 \mu\text{M}$ . Assay was  
performed following the kit protocol.



**MERCK**

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

## Technical Service

Visit the tech service page on our web site at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

## Standard Warranty

The applicable warranty for the products listed in this publication may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).