

## Product Information

### Chymotrypsin Activity Assay Kit

Catalog Number **MAK345**Storage Temperature  $-20^{\circ}\text{C}$ 

## TECHNICAL BULLETIN

### Product Description

Chymotrypsin is a key serine protease involved in dietary protein digestion in mammals. It is primarily produced by the pancreas, but may be expressed in other tissues, including the spleen and liver. In the pancreas, chymotrypsin is initially expressed as the inactive proenzyme chymotrypsinogen, which is cleaved by other proteases to chymotrypsin, its active form. Chymotrypsin specifically cleaves peptide bonds at the C-terminal end of bulky hydrophobic or aromatic amino acids (such as tyrosine, tryptophan, or phenylalanine).

The Chymotrypsin Activity Assay kit uses a synthetic fluorogenic substrate, enabling kinetic measurement of chymotrypsin activity in cell and tissue lysates. A chymotrypsin activator cleaves chymotrypsinogen to form active chymotrypsin, which then hydrolyzes the non-fluorescent substrate to release a stable fluorophore ( $\lambda_{\text{ex}} = 380 \text{ nm}/\lambda_{\text{em}} = 460 \text{ nm}$ ). The kit includes a selective chymotrypsin inhibitor that can be used to measure specific chymotrypsin activity in samples containing non-specific proteases and endopeptidases that may also metabolize the substrate. The assay can detect as low as 0.01 mU of Chymotrypsin.

### Components

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

Chymotrypsin Assay Buffer Catalog Number MAK345A	25 mL
Chymotrypsin Substrate Catalog Number MAK345B	200 $\mu\text{L}$
Chymotrypsin Activator Catalog Number MAK345C	1 vial
Chymotrypsin Inhibitor Catalog Number MAK345D	80 $\mu\text{L}$

Coumarin Standard (1 mM) Catalog Number MAK345E	100 $\mu\text{L}$
--	-------------------

Chymotrypsin Positive Control Catalog Number MAK345F	1 vial
---	--------

### Reagents and Equipment Required but Not Provided.

- Pipetting devices and accessories (e.g., multichannel pipettor)
- 96-well clear or white plate with flat bottom
- Fluorescence multiwell plate reader
- Anhydrous Dimethylsulfoxide (DMSO) (Catalog Number 276855)

### 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  $-20^{\circ}\text{C}$ , protected from light upon receiving. Briefly centrifuge small vials prior to opening.

### Preparation Instructions

Allow Chymotrypsin Assay Buffer, Chymotrypsin Substrate, Chymotrypsin Inhibitor, and Coumarin Standard to thaw completely before use. Mix well.

### Reagent Preparation

Chymotrypsin Activator – Reconstitute with 220  $\mu\text{L}$  of Chymotrypsin Assay Buffer immediately before use. Aliquot remainder and store at  $-80^{\circ}\text{C}$ . Once reconstituted, use within 2 months.

Chymotrypsin Positive Control – Reconstitute with 22  $\mu\text{L}$  of Chymotrypsin Assay Buffer immediately before use. Remaining positive control should be aliquoted and stored at  $-80^{\circ}\text{C}$ . Once reconstituted, use within 2 months.

## Procedure

### Sample Preparation

1. Homogenize cells ( $1 \times 10^6$ ) or tissue (20 mg) with 100  $\mu\text{L}$  of ice-cold Chymotrypsin Assay Buffer, and keep on ice for 10 minutes.
2. Centrifuge at  $10,000 \times g$  for 10 minutes at  $4^\circ\text{C}$  and transfer the supernatant to a fresh tube.
3. Determine protein concentration. Protein concentration should range between 5–20 mg/mL. Samples with a protein concentration  $>20$  mg/mL may be diluted with Chymotrypsin Assay Buffer to obtain the desired range.
4. Aliquot and store lysates at  $-80^\circ\text{C}$  unless being used immediately.
5. Use 5–20  $\mu\text{L}$  sample per well using a clear 96 well plate.
6. Prepare two identical wells for each sample labelled “Sample Background Control” (SBC), and “Sample” (S).
7. An additional well called “Sample + Inhibitor” (SI) may be prepared for samples in which nonspecific chymotrypsin-like protease activity is likely to be present. For SI, add 2  $\mu\text{L}$  of chymotrypsin inhibitor in addition to sample.
8. Adjust volume in each well to 50  $\mu\text{L}$  with Chymotrypsin Assay Buffer.
9. For positive control (PC), add 1–4  $\mu\text{L}$  of Chymotrypsin Positive Control into desired well(s) and adjust the final volume to 50  $\mu\text{L}$  with Chymotrypsin Assay Buffer.
10. For reagent background control (BC), add 50  $\mu\text{L}$  of Chymotrypsin Assay Buffer to a well.
11. For unknown samples, it is suggested to test several concentrations to ensure the readings are within the Standard Curve range.

### Coumarin Standards

1. Add 0, 2, 4, 6, 8, and 10  $\mu\text{L}$  from the provided 1 mM Coumarin Standard stock solution into a series of wells in a clear 96 well plate.
2. Bring the total volume up to 100  $\mu\text{L}$  per well with Chymotrypsin Assay Buffer to generate 0, 2, 4, 6, 8, and 10 nmol/well of Coumarin Standard.
3. Mix by pipetting, making sure that no bubbles are introduced in the wells.
4. If sample activity is low (outside standard curve RFU values), another standard curve ranging from 0.1 to 1 nmol/well may be generated. For this, dilute the provided Coumarin Standard 10-fold with DMSO to obtain a 100  $\mu\text{M}$  Coumarin Standard solution. Add 0, 2, 4, 6, 8, and 10  $\mu\text{L}$  of the 100  $\mu\text{M}$  solution into a series of wells in a 96 well plate and bring the total volume up to 100  $\mu\text{L}$  with Chymotrypsin Assay Buffer to generate 0, 0.2, 0.4, 0.6, 0.8, and 1 nmol/well of Coumarin Standard.

### Reaction Mix

Prepare reaction mixes for test sample and corresponding sample background control wells according to Table 1. Mix enough reagents for the number of assays to be performed (50  $\mu\text{L}$ /well).

**Table 1.**

Reagent	SBC Mix	Reaction Mix
Chymotrypsin Assay Buffer	48 $\mu\text{L}$	46 $\mu\text{L}$
Chymotrypsin Substrate	–	2 $\mu\text{L}$
Chymotrypsin Activator	2 $\mu\text{L}$	2 $\mu\text{L}$

### Assay Reaction

1. Add 50  $\mu\text{L}$  of the SBC Mix to each of the “Sample Background Control” wells.
2. Add 50  $\mu\text{L}$  of the Reaction Mix to wells containing samples (S), sample + inhibitor (SI), positive control (PC), and reagent background control (BC) for a final volume of 100  $\mu\text{L}$  per well. Turbidity upon addition of Chymotrypsin Substrate to Chymotrypsin Assay buffer is normal and will disappear following vortexing.
3. Immediately start recording the fluorescence ( $\lambda_{\text{ex}} = 380 \text{ nm}/\lambda_{\text{em}} = 460 \text{ nm}$ ) in kinetic mode (i.e., at 30 second intervals) for 30–60 minutes at  $25^\circ\text{C}$ . Ideal measurement time depends on the chymotrypsin activity in samples. It is recommended to run the assay in kinetic mode to ensure that the linear reaction phase is recorded. The Coumarin Standard curve can be read in endpoint mode.

## Results

1. Subtract the 0 pmol/well reading from all other Coumarin Standard readings and plot the standard curve.
2. For sample reaction wells (including paired inhibitor control wells), choose two time points ( $t_1$  and  $t_2$ ) in the linear phase of the reaction progress curve.
3. Obtain the corresponding fluorescence values at those points ( $RFU_1$  and  $RFU_2$ ) and determine the change in fluorescence over the time interval:  
 $\Delta F = RFU_2 - RFU_1$ .
4. Subtract the reagent background control (BC)  $\Delta F$  value from the respective sample (S)  $\Delta F$  values. If sample background control (SBC)  $\Delta F$  values are higher than BC, subtract the SBC from the corresponding sample (rather than subtracting the reagent BC).
5. Chymotrypsin specific activity is obtained by applying the background-corrected  $\Delta F$  values to the Coumarin Standard curve to get B pmol of substrate metabolized during the reaction time.

Chymotrypsin Specific Activity ( $\mu U/mg$ ) =

$$\Delta B / (\Delta t \times p) = \text{pmol/min/mg}$$

where:

$\Delta B$  = change in coumarin concentration during reaction (in pmol)

$\Delta t = t_2 - t_1$  (in minutes)

$p$  = sample protein content added to well (in mg)

**Note:** If chymotrypsin inhibitor is being used, calculate chymotrypsin activity as follows:

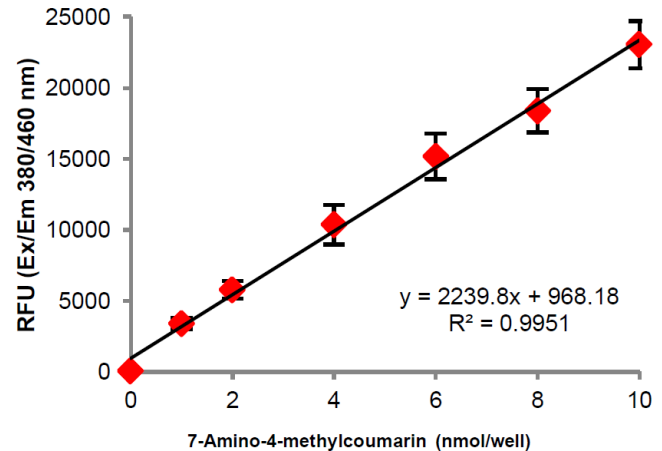
Chymotrypsin activity =

Total activity in sample – Activity in presence of Chymotrypsin Inhibitor

Unit Definition: One unit of Chymotrypsin is the amount of enzyme that generates 1.0  $\mu\text{mol}$  of coumarin per minute at pH 8 at 25 °C.

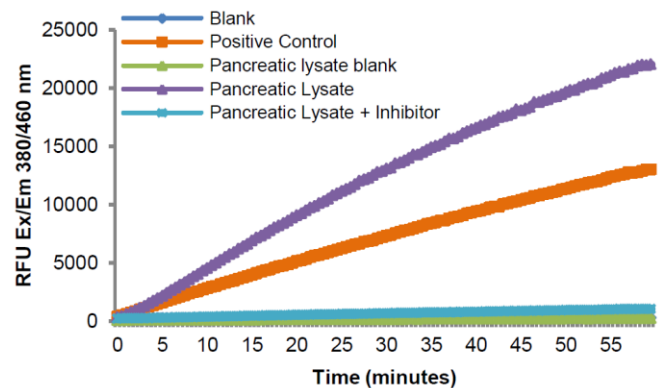
**Figure 1.**

Typical Coumarin Standard Curve Showing a Range of 0–10 nmol/well.

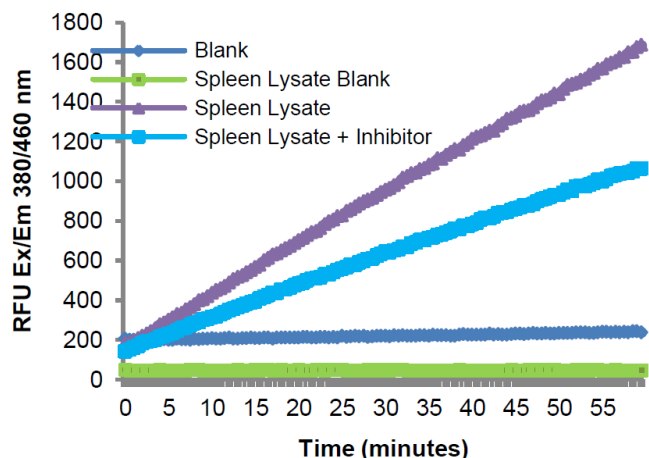


**Figure 2.**

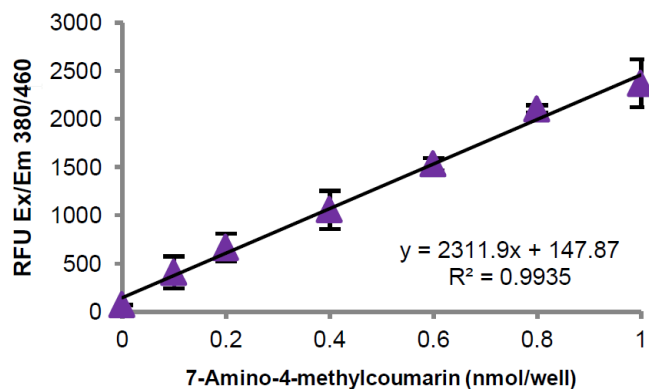
Reaction Kinetics for Positive Control and Rat Pancreatic Lysate (9  $\mu\text{g}$ ).



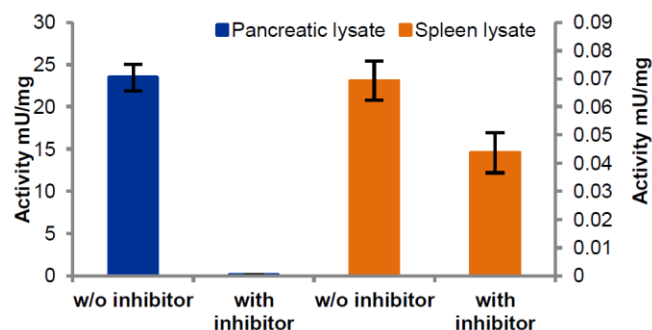
**Figure 3.**  
Reaction Kinetics for Rat Spleen Lysate (160 µg)



**Figure 4.**  
Typical Coumarin Standard Curve Showing a Range of 0-1 nmol/well.

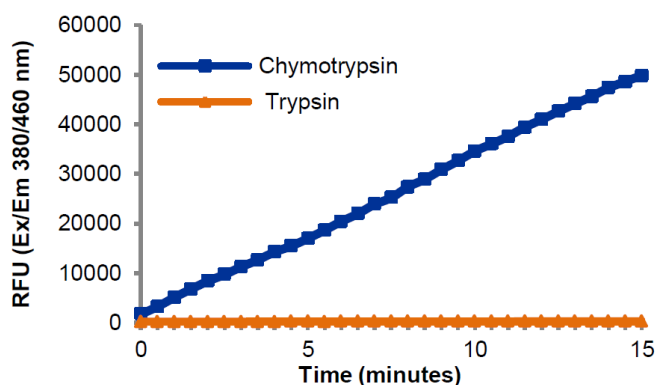


**Figure 5.**  
Chymotrypsin Activity With and Without Inhibitor in Rat Pancreas and Rat Spleen.



The presence of non-specific chymotrypsin-like proteases in spleen leads to some activity in presence of the selective chymotrypsin inhibitor.

**Figure 6.**  
Reaction Kinetics Using Substrate in The Presence of Chymotrypsin or Trypsin.



The substrate is cleaved by chymotrypsin, but not trypsin, indicating the assay is free from trypsin interference.

SS,HM,MAM 02/19-1