

## Product Information

### Cathepsin B Inhibitor Screening Kit

Catalog Number **MAK200**

Storage Temperature  $-20^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Cathepsin B (CTSB; EC 3.4.22.1) is a lysosomal cysteine proteinase and is involved in intracellular degradation and turnover of proteins.<sup>1</sup> It is involved in the proteolytic processing of amyloid precursor protein (APP). Incomplete proteolytic processing of APP is a causative factor in Alzheimer's disease.<sup>2</sup> CTSB has also been implicated in the invasion and metastasis of certain cancers.<sup>3</sup>

The Cathepsin B Inhibitor Screening Kit is a simple assay for screening potential inhibitors of CTSB. CTSB activity is measured by cleaving a synthetic 7-amino-4-trifluoromethylcoumarin (AFC)-based substrate to yield AFC, a fluorescent product ( $\lambda_{\text{ex}} = 400/$   $\lambda_{\text{em}} = 505 \text{ nm}$ ), proportional to the enzymatic activity present.

### Components

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

|                                                            |                   |
|------------------------------------------------------------|-------------------|
| CTSB Reaction Buffer<br>Catalog Number MAK200A             | 15 mL             |
| CTSB Reagent<br>Catalog Number MAK200B                     | 100 $\mu\text{L}$ |
| Cathepsin B, human<br>Catalog Number MAK200C               | 1 vial            |
| CTSB Substrate, Ac-RR-AFC, 10 mM<br>Catalog Number MAK200D | 0.2 mL            |
| CTSB Inhibitor, F-F-FMK, 1 mM<br>Catalog Number MAK200E    | 20 $\mu\text{L}$  |

### Reagents and Equipment Required but Not Provided

- 96 well flat-bottom plate – It is recommended to use black plates with clear bottoms for fluorescence assays.
- Fluorescence multiwell plate reader

### Precautions and Disclaimer

This product is 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.

### Preparation Instructions

Briefly centrifuge vials before opening. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

CTSB Reaction Buffer – Warm to room temperature before use.

CTSB Reagent – Aliquot and store at  $-20^{\circ}\text{C}$ .

Cathepsin B, human – Reconstitute with 110  $\mu\text{L}$  of CTSB Reaction Buffer. Mix well by pipetting. Aliquot and store at  $-80^{\circ}\text{C}$ .

CTSB Substrate and CTSB Inhibitor – Ready to use. Store at  $-20^{\circ}\text{C}$ .

### Storage/Stability

The kit is shipped on wet ice and storage at  $-20^{\circ}\text{C}$ , protected from light, is recommended. Briefly centrifuge the vials at low speed before opening.

### Procedure

#### Sample Preparation

Prepare a 10 $\times$  Sample Inhibitor Solution by diluting sample inhibitors with CTSB Reaction Buffer to 10 $\times$  the final testing concentration. An initial concentrated inhibitor solution may be in a different solvent if the inhibitor is minimally soluble in aqueous CTSB Reaction Buffer.

To correct for background in samples, include a Sample Blank by omitting the Cathepsin B. The Sample Blank readings can then be subtracted from the sample readings.

For unknown samples, it is suggested to test several sample dilutions.

Prepare an Enzyme Control (uninhibited enzyme) by using CTSB Reaction Buffer in place of sample inhibitor.

An Inhibitor Control may be prepared by diluting 1  $\mu\text{L}$  of CTSB Inhibitor with 9  $\mu\text{L}$  of CTSB Reaction Buffer.

Add 10  $\mu\text{L}$  of sample inhibitor (10 $\times$  Sample Inhibitor Solution), Sample Blank (10 $\times$  Sample Inhibitor Solution), Enzyme Control (CTSB Reaction Buffer), or Inhibitor Control into duplicate wells of a 96 well plate.

#### Assay Reaction

1. Set up Inhibition Reaction Mixes according to the scheme in Table 1. 50  $\mu\text{L}$  of the appropriate Inhibition Reaction Mix is required for each reaction (well).

**Table 1.**  
Inhibition Reaction Mixes

| Reagent              | Samples and Controls | Sample Blank     |
|----------------------|----------------------|------------------|
| CTSB Reaction Buffer | 48 $\mu\text{L}$     | 49 $\mu\text{L}$ |
| CTSB Reagent         | 1 $\mu\text{L}$      | 1 $\mu\text{L}$  |
| Cathepsin B          | 1 $\mu\text{L}$      | –                |

2. Add 50  $\mu\text{L}$  of the appropriate Inhibition Reaction Mix to each of the wells. Mix well using a horizontal shaker or by pipetting. Incubate the plate at room temperature for 10–15 minutes. Protect the plate from light during the incubation.
3. Set up an Enzymatic Reaction Mix according to the scheme in Table 2. 40  $\mu\text{L}$  of the Enzymatic Reaction Mix is required for each reaction (well).

**Table 2.**  
Enzymatic Reaction Mix

| Reagent              | Samples, Controls, and Sample Blank |
|----------------------|-------------------------------------|
| CTSB Reaction Buffer | 38 $\mu\text{L}$                    |
| CTSB Substrate       | 2 $\mu\text{L}$                     |

4. Add 40  $\mu\text{L}$  of the Enzymatic Reaction Mix to each reaction well. Mix well using a horizontal shaker or by pipetting.
5. Measure the fluorescence (FLU,  $\lambda_{\text{ex}} = 400/\lambda_{\text{em}} = 505 \text{ nm}$ ) in a microplate reader in kinetic mode for 30–60 minutes at 37 °C. It is recommended to take fluorescent readings every minute.

## Results

### Calculations

Plot the fluorescence (FLU) for each well versus time.

Choose two time points (T1 and T2) in the linear range of the plot and obtain the slope for each well between T1 and T2. Determine the FLU at each time (FLU1 and FLU2) and use them to determine the slope of the plot ( $\Delta\text{FLU}/\text{minute}$ ).

**Note:** The Enzymatic Control must be set up each time the assay is run.

Subtract the slope of the Sample Blank from the slope of the samples to obtain the corrected measurement. Use the corrected measurement to determine the % Relative Inhibition.

### % Relative Inhibition

$$\text{Slope} = (\text{FLU}_2 - \text{FLU}_1)/(\text{T}_2 - \text{T}_1) = \Delta\text{FLU}/\text{minute}$$

$$\% \text{ Relative Inhibition} = \frac{(\text{Slope}_{\text{EC}} - \text{Slope}_{\text{SM}})}{\text{Slope}_{\text{EC}}} \times 100\%$$

where:

$\text{Slope}_{\text{SM}}$  = the slope of the Sample Inhibitor

$\text{Slope}_{\text{EC}}$  = the slope of the Enzyme Control

**Note:** Irreversible inhibitors that completely inhibit Cathepsin B activity will have  $\Delta\text{FLU} = 0$ . The % Relative Inhibition will be 100%.

### Sample Calculation

$$\text{Slope}_{\text{SM}} = 0.435 \text{ FLU}/\text{min}$$

$$\text{Slope}_{\text{EC}} = 0.755 \text{ FLU}/\text{min}$$

$$\% \text{ Relative Inhibition} = \frac{(0.755 - 0.435)}{0.755} \times 100\% = 42.4\%$$

## References

1. Reiser, J. et al., Specialized roles for cysteine cathepsins in health and disease. *J. Clin. Invest.*, **120**, 3421–3431 (2010).
2. Cho, K. et al., CA-074Me, a cathepsin B inhibitor, decreases APP accumulation and protects primary rat cortical neurons treated with okadaic acid. *Neurosci. Lett.*, **548**, 222–227 (2013).
3. Bao, W., et al., Silencing of Cathepsin B suppresses the proliferation and invasion of endometrial cancer. *Oncol. Rep.*, **30**, 723–730 (2013).

**Troubleshooting Guide**

| <b>Problem</b>                                 | <b>Possible Cause</b>                                     | <b>Suggested Solution</b>                                                                  |
|------------------------------------------------|-----------------------------------------------------------|--------------------------------------------------------------------------------------------|
| Assay not working                              | Cold assay buffer                                         | Assay Buffer must be at room temperature                                                   |
|                                                | Omission of step in procedure                             | Refer and follow Technical Bulletin precisely                                              |
|                                                | Plate reader at incorrect wavelength                      | Check filter settings of instrument                                                        |
|                                                | Type of 96 well plate used                                | For Fluorometric assays, use black plates with clear bottoms                               |
| Samples with erratic readings                  | Samples prepared in different buffer                      | Use the Assay Buffer provided or refer to Technical Bulletin for instructions              |
|                                                | Cell/Tissue culture samples were incompletely homogenized | Repeat the sample homogenization, increasing the length and extent of homogenization step. |
|                                                | Samples used after multiple freeze-thaw cycles            | Aliquot and freeze samples if needed to use multiple times                                 |
|                                                | Presence of interfering substance in the sample           | If possible, dilute sample further                                                         |
|                                                | Use of old or inappropriately stored samples              | Use fresh samples and store correctly until use                                            |
| Lower/higher readings in samples and standards | Improperly thawed components                              | Thaw all components completely and mix gently before use                                   |
|                                                | Use of expired kit or improperly stored reagents          | Check the expiration date and store the components appropriately                           |
|                                                | Allowing the reagents to sit for extended times on ice    | Prepare fresh Reaction Mixes before each use                                               |
|                                                | Incorrect incubation times or temperatures                | Refer to Technical Bulletin and verify correct incubation times and temperatures           |
|                                                | Incorrect volumes used                                    | Use calibrated pipettes and aliquot correctly                                              |
| Non-linear standard curve                      | Use of partially thawed components                        | Thaw and resuspend all components before preparing the Reaction Mixes                      |
|                                                | Pipetting errors in preparation of standards              | Avoid pipetting small volumes                                                              |
|                                                | Pipetting errors in the Reaction Mix                      | Prepare Reaction Mixes whenever possible                                                   |
|                                                | Air bubbles formed in well                                | Pipette gently against the wall of the plate well                                          |
|                                                | Standard stock is at incorrect concentration              | Refer to the standard dilution instructions in the Technical Bulletin                      |
|                                                | Calculation errors                                        | Recheck calculations after referring to Technical Bulletin                                 |
|                                                | Substituting reagents from older kits/lots                | Use fresh components from the same kit                                                     |
| Unanticipated results                          | Samples measured at incorrect wavelength                  | Check the equipment and filter settings                                                    |
|                                                | Samples contain interfering substances                    | If possible, dilute sample further                                                         |
|                                                | Sample readings above/below the linear range              | Concentrate or dilute samples so readings are in the linear range                          |

GD,KVG,LS,MAM 10/14-1