

SensoLyte® 390 Cathepsin D Assay Kit *Fluorimetric*

Revision Number: 1.1	Last updated: October 2014	
Catalog #	AS-72098	
Kit Size	100 Assays (96-well plate)	

- Optimized Performance: This kit is optimized to detect Cathepsin D activity.
- Enhanced Value: It provides enough reagents to perform 100 assays in a 96-well format.
- *High Speed:* The entire process can be completed in one hour
- Assured Reliability: Detailed protocol and references are provided.

Kit Components, Storage and Handling

Component	Description	Quantity
Component A	Mca/Dnp, Cathepsin D substrate, Ex/Em=330 nm / 390 nm upon cleavage	1.6 mM, 50 μL
Component B	Mca fluorescence reference standard, Ex/Em=330 nm/390 nm	2 mM, 10 μL
Component C	Cathepsin D, bovine spleen	$0.1 \text{ mg/mL}, 20 \mu\text{L}$
Component D	Assay Buffer	20 mL
Component E	Pepstatin A	4 mM, 100 μL
Component F	DTT	1 M, 100 μL

Other Materials Required (but not provided)

- <u>96-well microplate</u>: Black, flat-bottom, non-binding 96-well plate.
- <u>Fluorescence microplate reader</u>: Capable of detecting emission at 390 nm with excitation at 330 nm.

Storage and Handling

- Store all kit components at -20°C, except for Component C
- Store Component C at -80°C
- Protect Components A and B from light and moisture.
- Component D can be stored at room temperature for convenience.

Introduction

Cathepsin D is the lysosomal aspartic proteinase implicated in intracellular protein degradation. It is involved in several pathological processes, such as inflammatory states, atherosclerosis, thrombosis, apoptosis, neoplastic proliferation or Alzheimer disease. ¹⁻⁵

The SensoLyte[®] 390 Cathepsin D Assay Kit provides a convenient assay for screening of enzyme inhibitors and for continuous assay of Cathepsin D activity using a Mca/Dnp FRET peptide. The sequence of this FRET peptide is derived from the cleavage site of Cathepsin D.^{6,7} In the FRET peptide, the fluorescence of Mca is quenched by Dnp. Upon cleavage into two separate fragments by Cathepsin D, the fluorescence of Mca is recovered, and can be monitored at excitation/emission = 330 nm/390 nm.

Protocol

Note 1: For standard curve, please refer to Appendix II (optional).

Note 2: Please use protocol A or B based on your needs.

Protocol A. Screening Cathepsin D inhibitors using purified enzyme.

1. Prepare working solutions.

Note: Warm all kit components until thawed to room temperature before starting the experiments.

1.1 Prepare assay buffer: Prepare fresh assay buffer for each experiment according Table 1. Use this DTT-containing assay buffer in all the following steps.

Table 1. Assay buffer for one 96-well plate (100 assays)

	1 \ /
Components	Volume
Assay buffer (Component D)	9.95 mL
1 M DTT (Component F)	50 μL
Total volume	10 mL

1.2 Cathepsin D substrate solution: Dilute cathepsin D substrate (Component A) 1:100 in assay buffer. For each experiment, prepare fresh substrate solution.

Table 2. Cathepsin D substrate solution for one 96-well plate (100 assays)

Components	Volume
Cathepsin D substrate (100X, Component A)	50 μL
Assay buffer	4.95 mL
Total volume	5 mL

1.3 <u>Cathepsin D diluent</u>: Dilute the enzyme (Component C) 1:800 in assay buffer. This amount of enzyme is enough for a full 96-well plate. If not using the entire plate, adjust the amount of enzyme to be diluted accordingly.

<u>Note:</u> Prepare enzyme diluents immediately before use. Do not vortex the enzyme solutions. Prolonged storage or vigorous agitation of the diluted enzyme will cause denaturation. Store the enzyme solution on ice.

1.4 Pepstatin A (Cathepsin D inhibitor): Dilute the 4 mM inhibitor solution (Component E) 1:10 in assay buffer. The diluted pepstatin solution has a concentration of 400 μM. Add 10 μl of the diluted pepstatin into each of the inhibitor control well.

2. Set up the enzymatic reaction.

 $\underline{2.1}$ Add test compounds and diluted enzyme solution to the microplate wells. The suggested volume of enzyme solution for one well of a 96-well plate is 40 μ L and test compound is 10 μ L.

- <u>2.2</u> Simultaneously establish the following control wells, as deemed necessary:
 - ➤ <u>Positive control</u> contains the enzyme without test compound.
 - ➤ <u>Inhibitor control</u> contains Cathepsin D enzyme and pepstatin.
 - ➤ <u>Vehicle control</u> contains Cathepsin D enzyme and vehicle used in delivering test compound (e.g. DMSO, concentration not to exceed 1%).
 - ➤ <u>Test compound control</u> contains assay buffer and test compound. Some test compounds have strong autofluorescence and may give false results.
 - Substrate control contains assay buffer.
- 2.3 Using the assay buffer, bring the total volume of all controls to 50 µL.
- 2.4 Optional: Pre-incubate the plate for 10 min. at assay temperature. Any temperature (the *assay temperature*) from room temperature to 37°C may be used, as long as the subsequent incubations are performed at the same temperature.

3. Run the enzymatic reaction.

- 3.1 Add 50 μL of Cathepsin D substrate solution into each well. For best accuracy, it is advisable to have the substrate solution equilibrated to the assay temperature. Mix the reagents completely by shaking the plate gently for 30 sec.
- 3.2 Measure fluorescence signal:
 - <u>For kinetic reading</u>: Immediately start measuring fluorescence intensity at Ex/Em=330 nm/390 nm continuously and record data every 5 min. for 30 to 60 min.
 - <u>For end-point reading</u>: Incubate the reaction for 30 to 60 min. Keep plate from direct light. Then measure fluorescence intensity at Ex/Em=330 nm/390 nm.
- 3.3 For methods of data analysis: Refer to Appendix I.

<u>Protocol B.</u> Measuring Cathepsin D activity in biological samples.

1. Prepare cathepsin containing biological samples.

1.1 Prepare cell extracts:

- Suspension cells are collected by centrifugation at 500 X g for 5 min. For adherent cells, simply aspirate the growth medium.
- Add an appropriate amount of assay buffer (see step 2.1) to cells or cell pellet. Collect the cell suspension in a microcentrifuge tube.
- Incubate the cell suspension at 4°C for 10 minutes.
- Centrifuge the cell suspension for 10 minutes at 2,500X g, 4°C. Collect the supernatant and store at -70°C until use.

 $\underline{\text{Note}}$: If activity of cell extracts is low, prepare more concentrated samples with centrifugation at 20,000 x g. Use precipitated pellets.

2. Prepare working solutions.

Note: Warm all kit components until thawed to room temperature before starting the experiments.

<u>2.1</u> <u>Prepare assay buffer:</u> Prepare fresh assay buffer for each experiment according Table 1. Use this DTT-containing assay buffer in all the following steps.

Table 1. Assay buffer for one 96-well plate (100 assays)

Components	Volume
Assay buffer (Component D)	9.95 mL
1 M DTT (Component F)	50 μL
Total volume	10 mL

2.2 Cathepsin D substrate solution: Dilute cathepsin D substrate (Component A) 1:100 in assay buffer. For each experiment, prepare fresh substrate solution.

Table 2. Cathepsin D substrate solution for one 96-well plate (100 assays)

Components	Volume
Cathepsin D substrate (100X, Component A)	50 μL
Assay buffer	4.95 mL
Total volume	5 mL

2.3 Cathepsin D diluent: If using purified cathepsin D as a positive control, then dilute the enzyme 1:250 in assay buffer. Add 50 μl of the diluted enzyme into each of the positive control well.

3. Set up enzymatic reaction.

- 3.1 Add 5-50 μL of cathepsin D containing biological sample.
- 3.2 Set up the following control wells at the same time, as deemed necessary:
 - Positive control contains purified active cathepsin D.
 - > Substrate control contains assay buffer.
- 3.3 Using the assay buffer, bring the total volume of all controls to 50 µL.
- 3.3 Optional: Pre-incubate the plate for 10 min. at assay temperature. Any temperature (the *assay temperature*) from room temperature to 37°C may be used, as long as the subsequent incubations are performed at the same temperature.

4. Run the enzymatic reaction.

- 4.1 Add 50 μL of cathepsin D substrate solution into each well. For best accuracy, it is advisable to have the substrate solution equilibrated to the assay temperature. Mix the reagents completely by shaking the plate gently for 30 sec.
- 4.2 Measure fluorescence signal:
 - <u>For kinetic reading</u>: Immediately start measuring fluorescence intensity at Ex/Em=330 nm/390 nm continuously and record data every 5 min. for 30 to 60 min.
 - <u>For end-point reading</u>: Incubate the reaction for 30 to 60 min. Keep plate from direct light. Then measure fluorescence intensity at Ex/Em=330 nm/390 nm.
- 4.3 For methods of data analysis: Refer to Appendix I.

Appendix I. Data Analysis

- The fluorescence reading from the substrate control well is used as the background fluorescence. This background reading should be subtracted from the readings of the other wells containing substrate. All fluorescence readings are expressed in relative fluorescence units (RFU).
- For kinetics analysis:
 - ➤ Plot data as RFU versus time for each sample. If you want to convert RFUs to the concentration of the product of the enzymatic reaction, please refer to <u>Appendix II</u> for establishing a fluorescence reference standard.
 - ➤ Determine the range of initial time points during which the reaction is linear. Typically, the first 10-15% of the reaction will be the optimal range.
 - ➤ Obtain the initial reaction velocity (V_o) in RFU/min by determining the slope of the linear portion of the data plot.
 - ➤ A variety of data analyses can be done, e.g., determining inhibition %, EC₅₀, IC₅₀, K_m, K_i, etc.
- For endpoint analysis:
 - ➤ Plot data as RFU versus concentration of test compounds.
 - ➤ A variety of data analyses can be done, e.g., determining inhibition %, EC₅₀, IC₅₀, etc.

Appendix II. Instrument Calibration

- Mca fluorescence reference standard: Dilute 2 mM Mca (Component B) to 20 μM in assay buffer. Do 2-fold serial dilutions to get concentrations of 10, 5, 2.5, 1.2, 0.6 and 0.3 μM, include an assay buffer blank. Add 50 μL/well of these serially diluted Mca reference solutions.
- Add 50 μL/well of the diluted cathepsin substrate solution (refer to Protocol A, step 1.1 for preparation).

<u>Note</u>: The cathepsin substrate solution is added to the Mca reference standard to correct for the absorptive quenching by the FRET peptide. If multiple concentrations of substrate are used, this step must be performed for each concentration.

- Measure the fluorescence intensity of the reference standard and substrate control wells at Ex/Em=330 nm/390 nm. Use the same setting of sensitivity as used in the enzyme reaction.
- Plot the Mca fluorescent reference standard curve as RFU (relative fluorescent units) versus concentration as shown in Figure 1.
- The final concentrations of Mca reference standard are 10, 5, 2.5, 1.2, 0.6, 0.3, 0.15 and 0 μM. This reference standard is used to calibrate the variation of different instruments and different experiments. It is also an indicator of the amount of final product of the enzymatic reaction.

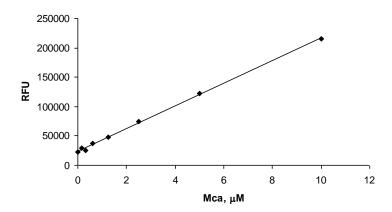


Figure 1. Mca reference standard, Mca was serially diluted in assay buffer, containing cathepsin D substrate, and the fluorescence recorded at Ex/Em=330 nm/390 nm. (Flexstation 384II, Molecular Devices)

References

- 1. Berchem, G. et al, Oncogene 21, 5951(2002).
- 2. Liaudet-Coopman, E. et al, Cancer Lett. 237, 167 (2006).
- 3. Davidson, Y. et al, J. Neurol. Neurosurg Psychiat. 77, 515 (2006).
- 4. Laurent-Matha, V. et al, J. Cell Biol. 168, 489 (2005).
- 5. Simon, DI. et al, *Biochemistry* **33**, 6555 (1994).
- 6. Baechle, D. et al, *J. Peptide Sci.* **11**, 166 (2005).
- 7. Yasuda, Y. et al, *J. Biochem.* **125**, 1137 (1999).