



## SensoLyte<sup>®</sup> 440 Deubiquitination Assay Kit \*Fluorimetric\*

|                             |  |                              |  |
|-----------------------------|--|------------------------------|--|
| <i>Revision Number: 1.1</i> |  | <i>Last Updated: 3/26/15</i> |  |
| <b>Catalog #</b>            |  | <b>AS-72204</b>              |  |
| <b>Kit Size</b>             |  | 100 Assays (96-well plate)   |  |

- **Optimized Performance:** This kit is optimized to detect deubiquitin protease activity
- **Enhanced Value:** It provides ample reagents to perform 100 assays in a 96-well plate 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 | 440 DUB fluorogenic substrate, Ex/Em = 354/442 nm upon cleavage | 20 $\mu$ M, 50 $\mu$ L |
| Component B | Fluorescence reference standard, Ex/Em=354/442 nm               | 20 $\mu$ M, 20 $\mu$ L |
| Component C | Deubiquitin protease, Recombinant Human UCH-L3                  | 20 $\mu$ L             |
| Component D | 2X Assay Buffer   | 25 mL                  |
| Component E | DTT   | 1 M, 500 $\mu$ L       |
| Component F | UCH-L3 Inhibitor  | 20 mM, 10 $\mu$ L      |

#### Other Materials Required (but not provided)

- 96-well microplate: Black, flat-bottom, non-binding 96-well plate.
- Fluorescence microplate reader: Capable of detecting emission at 354 nm with excitation at 442 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.

[Type text]

---

## Introduction

Modification of proteins with ubiquitin is a key step in protein degradation controlling many intracellular processes such as transcription, cell cycle progression, receptor internalization, and DNA repair.<sup>1</sup> Recent studies have demonstrated that regulation also occurs at the level of deubiquitination.<sup>2,3</sup> Deubiquitinating enzymes (DUBs) are proteases that reverse ubiquitin modifications.

Over 100 DUBs have been identified in humans. They are grouped into five distinct families based on their sequence similarities and mechanisms of action. Four of the families are cysteine proteases, while the fifth is a novel type of zinc-dependent metalloprotease.<sup>4,5</sup> The majority of human DUBs belong to two cysteine proteases families, the ubiquitin specific proteases (USPs) and the ubiquitin carboxyterminal hydrolases (UCHs). Both protease families hydrolyze the peptide bond downstream of the C terminus of ubiquitin, which is either a classic peptide bond in the proforms of ubiquitin or an isopeptide bond to an  $\epsilon$ -amino group of a lysine residue within an ubiquitin modified protein. DUBs have been implicated in the pathogenesis of many human diseases, such as neurodegenerative disorders and cancer.<sup>6,7</sup> Consequently they have become actively studied molecular targets for drug discovery.<sup>8,9</sup>

The Sensolyte<sup>®</sup> 440 Deubiquitination Assay Kit provides a convenient assay for inhibitors/activators screening in deubiquitinylation or for continuous assay of enzyme activity. A natural substrate, ubiquitin, coupled with the AMC (7-amino-4-methylcoumarin) fluorophore is used in the kit. Upon the cleavage by a deubiquitin protease, this substrate generates the blue fluorophore that can be detected at excitation/emission=354/442 nm. The substrate supplied in the kit is exquisitely sensitive for ubiquitin hydrolases (UCHs).

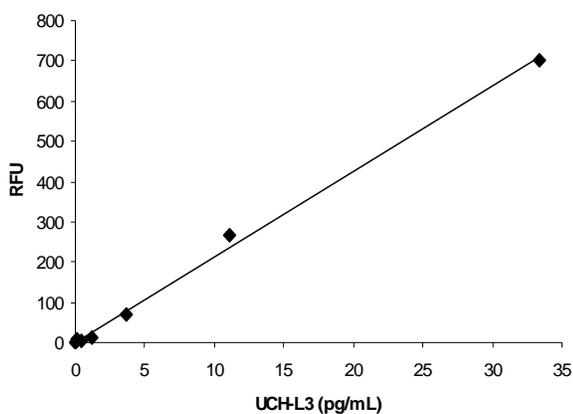


Figure 1. Detection of UCH-L3 activity with the Sensolyte<sup>®</sup> 440 Deubiquitination Assay Kit. The detection limit can reach as low as 3.7 pg/mL.

---

## Protocol

Note 1: For instrument calibration, refer to Appendix II (optional).

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

### Protocol A. Screening compounds using purified enzyme.

#### 1. Prepare working solutions.

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

- 1.1 Prepare 1X assay buffer: Prepare fresh DTT-containing 1X assay buffer for each experiment. Refer to Table 1. **Use this DTT-containing assay buffer in all the following steps.** If not using the entire plate, adjust the amount of assay buffer to be diluted accordingly.

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

| Components                    | Volume      |
|-------------------------------|-------------|
| 2X assay buffer (Component D) | 5 mL        |
| 1 M DTT (Component E)         | 100 $\mu$ L |
| Deionized water               | 4.9 mL      |
| Total volume                  | 10 mL       |

- 1.2 The 440 DUB substrate solution: Dilute deubiquitin substrate (Component A) 1:100 in 1X assay buffer. Refer to Table 2. For each experiment, prepare fresh substrate solution. If not using the entire plate, adjust the amount of substrate to be diluted accordingly.

**Table 2. The 440 DUB substrate solutions for one 96-well plate (100 assays)**

| Components                      | Volume     |
|---------------------------------|------------|
| 440 DUB substrate (Component A) | 50 $\mu$ L |
| 1X assay buffer                 | 4.95 mL    |
| Total volume                    | 5 mL       |

- 1.3 Protease diluents: Dilute the enzyme, UCH-L3 (Component C) 1:200 in 1X 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.

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

- 1.4 UCH-L3 inhibitor: Dilute the 20 mM inhibitor solution (Component F) 1:10 in 1X assay buffer to get a concentration of 2 mM. Add 10  $\mu$ L of the diluted inhibitor into each of the inhibitor control well.

## 2. Set up the enzymatic reaction.

- 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  $\mu$ L and test compound is 10  $\mu$ L.

- 2.2 Simultaneously set up the following control wells, as deemed necessary:

- Positive control contains the diluted UCH-L3 without test compound.
- Inhibitor control contains the diluted UCH-L3 and inhibitor.
- Vehicle control contains enzyme and vehicle used in delivering test compound (e.g. DMSO, concentration not to exceed 1%).
- Test compound control contains 1X assay buffer and test compound. Some test compounds have strong autofluorescence and may give false results.
- Substrate control contains 1X assay buffer.

- 2.3 Using 1X assay buffer, bring the total volume of all controls to 50  $\mu$ L.

- 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  $\mu\text{L}$  of the 4400 DUB substrate solutions 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:

- For kinetic reading: Immediately start measuring fluorescence at  $\text{Ex/Em}=354/442$  nm continuously and record data every 5 min. for 30 to 60 min.
- For end-point reading: Incubate the reaction for 30 to 60 min. Keep plate from direct light. Mix the reagents and measure fluorescence intensity at  $\text{Ex/Em}=354/442$  nm.

3.3 For methods of data analysis: Refer to Appendix I.

## **Protocol B. Measuring deubiquitination activity in biological samples.**

### 1. Prepare working solutions.

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

1.1 The 440 DUB substrate solution: Dilute deubiquitin substrate (Component A) in 2X assay buffer (Component D). Refer to Table 1. If not using the entire plate, adjust the amount of substrate to be diluted accordingly.

**Table 1. The 440 DUB substrate solution for one 96-well plate (100 assays)**

| <b>Components</b>               | <b>Volume</b>     |
|---------------------------------|-------------------|
| 440 DUB substrate (Component A) | 50 $\mu\text{L}$  |
| 1 M DTT (Component E)           | 100 $\mu\text{L}$ |
| 2X assay buffer (Component D)   | 4.85 mL           |
| Total volume                    | 5 mL              |

### 2. Set up the enzymatic reaction.

2.1 Add 50  $\mu\text{L}$  of DUBs containing sample.

2.2 Set up the following control wells at the same time, as deemed necessary:

- Positive control contains purified active UCH-L3.
- Substrate control contains deionized water.

2.3 Bring the total volume of all controls to 50  $\mu\text{L}$ .

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  $\mu\text{L}$  of the 440 DUB substrate solutions 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.

3.4 Measure fluorescence signal:

- For kinetic reading: Immediately start measuring fluorescence at Ex/Em=354 nm/442 nm continuously and record data every 5 min. for 30 to 60 min.
- For end-point reading: Incubate the reaction for 30 to 60 min. Keep plate from direct light. Mix the reagents and measure fluorescence intensity at Ex/Em=354 nm/442 nm.

3.5 For methods of data analysis: Refer to Appendix I.

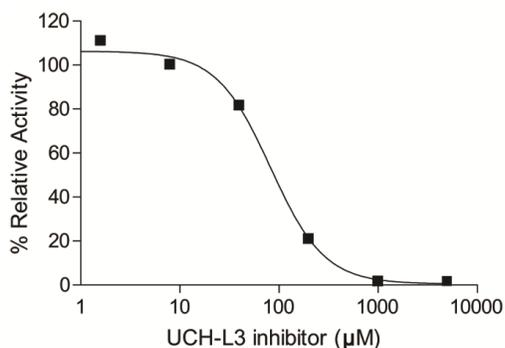


Figure 2. Inhibition of UCH-L3 activity as measured with SensoLyte® 440 Deubiquitination Assay Kit.

---

## 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. To convert RFUs to the concentration of the product of the enzymatic reaction, please refer to [Appendix II](#) 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 %,  $EC_{50}$ ,  $IC_{50}$ ,  $K_m$ ,  $K_i$ , 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 %,  $EC_{50}$ ,  $IC_{50}$ , etc.

---

## Appendix II. Instrument Calibration

- Fluorescence reference standard: Dilute the 20 µM fluorescence standard (Component B) to 200 nM in 1X DTT-containing assay buffer. Do 2-fold serial dilutions to get concentrations

of 100, 50, 25, 12.5, 6.25, and 3.13 nM, include an assay buffer blank. Add 50  $\mu$ L/well of these serially diluted reference solutions.

- Add 50  $\mu$ L/well of the diluted 440 DUB substrate solutions (refer to Protocol A, Step 1.2 for preparation).

Note: The 440 DUB substrate solution is added to the reference standard to normalize for the intrinsic substrate fluorescence. If multiple concentrations of substrate are used, this step must be performed for each concentration.

- Measure the fluorescence of the reference standard and substrate control wells at Ex/Em=354/442 nm. Use the same setting of sensitivity as used in the enzyme reaction.
- Plot the reference standard curve as RFU (relative fluorescent units) versus concentration.
- The final concentrations of fluorescence reference standard are 100, 50, 25, 12.5, 6.25, 3.13, 1.56 and 0 nM. 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 in the enzymatic reaction.

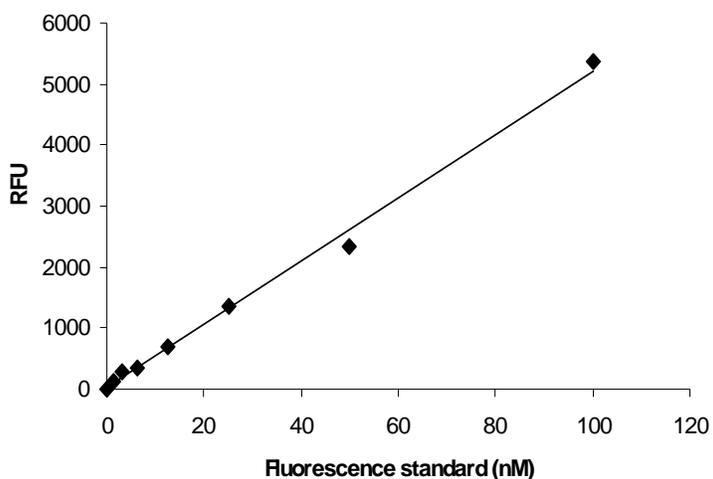


Figure 3. AMC reference standard. AMC was serially diluted in assay buffer containing substrate, and the fluorescence was recorded at Ex/Em=354/442 nm. (FLx800, Bio-Tek Instruments).

---

## References :

1. Hershko, A. and Ciechanover, A. *Annu Rev Biochem* **67**, 425 (1998).
2. Amerik, AY. and M. Hochstrasser, *Biochim Biophys Acta* **1695**, 189 (2004)
3. Kim, J.H. et al. *J Biochem.* **134**, 9 (2003).
4. Soboleva, T.A. *Curr Protein Pepy Sci* **5**, 191 (2004).
5. Karen, H. et al. *Biochem J* **414**, 161 (2008).
6. Setsuie, R. et al. *Neurochem Int.* **51**, 105 (2007).
7. Hussain, S. et al. *Cell Cycle* **8**, 1688 (2009).
8. Nicholson, B. et al. *Future Oncol* **3**, 191 (2007).
9. Colland, F. *Biochem Soc Trans* **38**, 137 (2010).

Revised: July 26, 2011