



## SensoLyte<sup>®</sup> FDP Alkaline Phosphatase ELISA Assay Kit \**Fluorimetric*\*

|                      |                            |
|----------------------|----------------------------|
| Revision Number: 1.1 | Last Updated: October 2014 |
| Catalog #            | AS-71101-R                 |
| Kit Size             | 500 Assays (96-well plate) |

- **Convenient Format:** Complete kit includes all the assay components.
- **Optimized Performance:** Optimal conditions for AP-labeled secondary antibody detection.
- **Enhanced Value:** Less expensive than the sum of individual components.
- **High Speed:** Minimal hands-on time.
- **Assured Reliability:** Detailed protocol and references are provided.

### Kit Components, Storage and Handling

| Component   | Description  | Quantity |
|-------------|--|----------|
| Component A | FDP, fluorogenic alkaline phosphatase substrate      | 1 vial   |
| Component B | Assay buffer   | 60 mL    |
| Component C | Stop solution  | 30 mL    |
| Component D | 10X Wash buffer                                      | 60 mL    |
| Component E | DMSO   | 0.3 mL   |
| Component F | Alkaline phosphatase-conjugated goat anti-rabbit IgG | 50 µL    |

### Other Materials Required (but not provided)

- Microplate: Black, high-binding 96-well plate for ELISA.
- Fluorescence microplate reader: Capable of detecting emission at 520 nm with excitation at 420 nm.

### Storage and Handling

- Except for Component A, store all components at 4°C.
- Store Component A at -20°C.

---

## Introduction

The SensoLyte<sup>®</sup> FDP Alkaline Phosphatase Assay Kit is used in ELISA with alkaline phosphatase (AP) conjugated antibody or streptavidin. FDP (3,6-fluorescein diphosphate) is a sensitive fluorogenic substrate for alkaline phosphatase. The final hydrolytic product of FDP is fluorescein, a fluorophore that has a very high emission quantum yield. This characteristics of fluorescein makes FDP a superior substrate for alkaline phosphatase and increases the assay sensitivity by 100 times compared to the colorimetric substrate, *p*NPP<sup>1</sup>. The signal can easily be read by a fluorescence plate reader at Ex/Em=485±20/528±20 nm.

---

## Protocol

Note 1: Prepare the ELISA assay plate according to standard ELISA procedures (refer to [Appendix](#)). Alkaline phosphatase-conjugated goat anti-rabbit IgG (Component E) is provided in this kit.

Note 2: Warm up all kit components to room temperature when the ELISA plate is ready for detection.

### 1. Prepare FDP stock solution.

- FDP stock solution (200X): Reconstitute the substrate by adding 250 µL of DMSO (Component E) into the FDP vial (Component A). The stock solution is good for 3-4 months if stored at -20°C.

### 2. Prepare the FDP reaction mixture.

- Prepare the FDP reaction mixture according to the following table. Keep the reaction mixture away from light.

Table 1. FDP reaction mixture for one 96-well plate (100 assays).

| Components                 | Volume |
|----------------------------|--------|
| FDP stock solution (200X)  | 50µL   |
| Assay Buffer (Component B) | 10 mL  |
| Total volume               | 10mL   |

### 3. Optional: if phosphate-buffered saline was used in ELISA procedures, the microplate must be washed with 1X wash buffer.

- Add 10 mL of 10 X wash buffer (Component D) to 90 mL deionized water to get 1X wash buffer.
- Wash microplate with 200 µL/well of 1X wash buffer three times, then pad dry on paper towels. For better sensitivity, we recommend using the buffers described in [Appendix](#).

### 4. Start the alkaline phosphatase reaction.

- Add 100µL/well of FDP reaction mixture prepared in Step 2.
- Incubate the reaction for 10 to 30 minutes, away from light.

Note: The reaction can be stopped by adding 50µL/well of stop solution (Component C). The signal is stable for at least 45 minutes.

- Read the plate using a fluorescence microplate reader with a filter set of excitation/emission at  $485\pm 20/528\pm 20$  nm.
- The results can be plotted as shown in Figure 1.

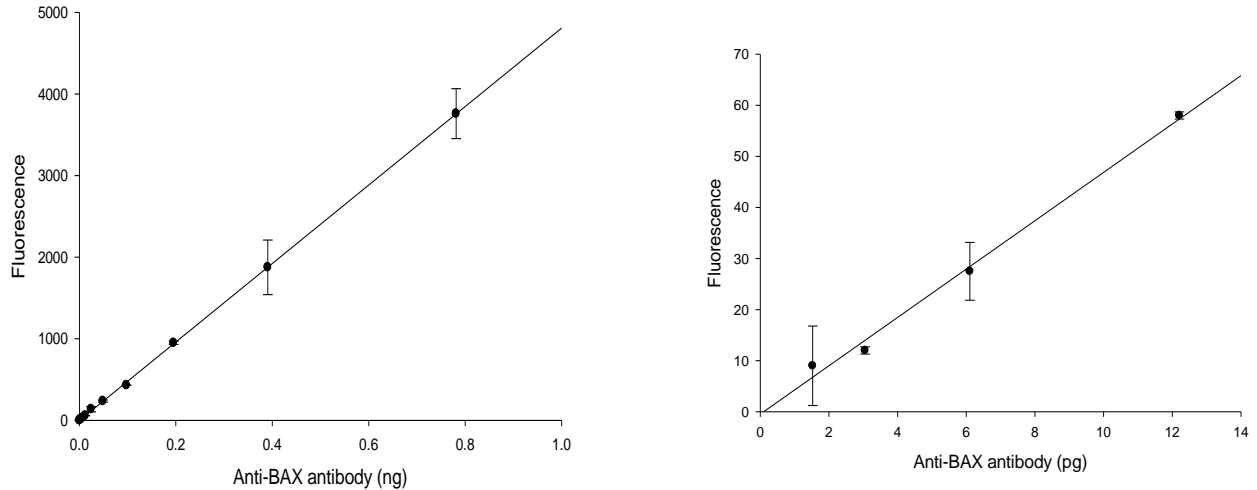


Figure 1. SensoLyte<sup>®</sup> FDP alkaline phosphatase ELISA assay kit was used to detect rabbit anti-BAX antibody.

The assay can detect as low as 1 pg (see right panel). The wells were coated with BAX-BSA. After blocking and washing, serially diluted rabbit anti-BAX antibody was then added into the wells. The wells were washed, and AP-conjugated goat anti-rabbit IgG secondary antibody (1:2000 dilution) was added. After incubation with secondary antibody the wells were washed and FDP reaction mixture was added. In 6 minutes the fluorescence was read by a fluorescence microplate reader at Ex/Em= $485\pm 20$  nm/ $528\pm 20$  nm.

---

## Appendix: General ELISA protocol

### 1. Required buffers:

1. Coating buffer: 1.59 g of  $\text{Na}_2\text{CO}_3$  and 2.93 g of  $\text{NaHCO}_3$  in 1 L of deionized  $\text{H}_2\text{O}$ , pH is 9.6 without adjustment.
2. Tris-buffered saline (TBS): 8.76 g of  $\text{NaCl}$ , 12.1 g of Tris in 800 ml of deionized  $\text{H}_2\text{O}$ . Adjust the pH to 7.4 with  $\text{HCl}$ . Add  $\text{H}_2\text{O}$  to 1L.
3. Blocking buffer: add 10 g of bovine serum albumin (BSA) and 0.2 mL of Tween<sup>®</sup>-20 into 1 L of TBS.
4. EIA buffer: add 1 g of bovine serum albumin (BSA) and 0.2 mL Tween<sup>®</sup>-20 into 1 L of TBS.
5. Wash buffer: add 0.2 mL of Tween<sup>®</sup>-20 into 1 L of TBS.

### 2. ELISA procedures:

1. Coating: Add 100  $\mu\text{L}$ /well of protein to the 96-well plate at a concentration of 10  $\mu\text{g}/\text{mL}$  in coating buffer. Incubate the plate at 4°C overnight.
2. Blocking: Discard the solution. Add 200  $\mu\text{L}$  of blocking buffer and incubate 30 min to 1 h at room temperature. Discard the blocking reagent and dry the plate under vacuum. You can store the plate at 4°C for future use.
3. Washing: Wash the plate with 200  $\mu\text{L}$  of wash buffer per well three to five times. Soak the plate during the last wash step for 5 minutes. Pad dry on paper towel.
4. Add sample: Add 50-100  $\mu\text{L}$ /well of sample to be tested and incubate at room temperature for 1 or more hours on a plate shaker. The sample can be diluted in EIA buffer or other appropriate buffer before adding to the plate.
5. Washing: Repeat Step 3.
6. Add enzyme-conjugated secondary antibody: Dilute alkaline phosphatase conjugated secondary antibody in EIA buffer to an appropriate concentration (1:1,000 – 1:10,000). Alkaline phosphatase-conjugated goat anti rabbit IgG (Component F) is provided in the kit. Add 100  $\mu\text{L}$ /well of diluted secondary antibody and incubate at room temperature for 30 min to 1 h on a plate shaker.
7. Washing: Repeat Step 3.
8. Detect by adding substrate: The plate is now ready for the FDP detection (refer to Protocol).

---

## References

1. Huang, Z., Olson, N.A., You, W., Haugland, R.P. *J.Immunol.Methods* 149, 261-266 (1992).