

## Myeloperoxidase (MPO) Activity Assay Kit

Catalog Number EEA016 (96 tests)

Rev 3.0

For safety and biohazard guidelines, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

### Product description

This kit can be used to detect myeloperoxidase (MPO) activity in serum, plasma, milk, and animal tissue. Myeloperoxidase is a heme-containing cationic glycoprotein that belongs to the heme peroxidase family in mammals. MPO is a dimer formed by polymerization of two subunits. Each subunit contains a heavy chain and a light chain. MPO is abundant in the azurophilic granules of polymorphonuclear leukocytes (PMNLs) and a small number in monocytes and macrophages. Studies have shown that MPO plays an important role in the generation of oxidants and host defense in neutrophils, and is closely related to the pathogenesis of many diseases, including cardiovascular disease, lung injury, and cancer.

Myeloperoxidase reduces hydrogen peroxide to a complex. The complex reacts with o-dianisidine (as hydrogen donor) to produce a yellow product which has a maximum absorption peak at 460 nm. The activity of MPO can be calculated indirectly by measuring the OD value at 460 nm.

## Contents and storage

Kit and components are shipped at 2-8°C. An unopened kit can be stored at 2-8°C for 12 months.

| Components      | Quantity (96 tests) |
|-----------------|---------------------|
| Buffer Solution | 20 mL               |
| Powder A        | Powder × 2 vials    |
| Powder B        | Powder × 2 vials    |
| Saline Solution | 6 mL                |
| Clarificant     | 1.2 mL × 2 vials    |
| Powder C        | Powder × 2 vials    |
| Substrate       | 0.1 mL              |
| Acid Reagent    | 1 mL                |
| Microplate      | 1 plate             |
| Plate Sealer    | 2 pieces            |

## Required materials

- Distilled or deionized water
- Normal saline (0.9% NaCl), PBS (0.01 M, pH 7.4)
- Microtiter plate reader with software capable of measurement at or near 460 nm
- Calibrated adjustable precision pipettes and glass or plastic tubes for diluting solution
- Incubator capable of maintaining 37 °C/ 60 °C .

## Procedural guidelines

---

**IMPORTANT!** Reagents are lot-specific. Do not mix or interchange different reagent lots from various kit lots.

---

Allow reagents to reach room temperature before use. Mix to redissolve any precipitated salts.

### Sample preparation guidelines

**Serum and plasma samples:** Detect the sample directly. If the sample is turbid, centrifuge at 10000 g for 10 min at 4 °C, then take the supernatant for detection.

**Milk samples:** Collect fresh milk sample, centrifuge at 10000 g for 10 min at 4 °C. Discard the upper white liquid and preserve the middle layer liquid on ice for detection.

**Tissue samples:**

- Take 0.02-1 g fresh tissue to wash with PBS(0.01 M, pH 7.4) at 2-8 °C to remove blood cells.
- Absorb the water with filter paper and weigh.
- Homogenize at the ratio of the volume of powder A application solution (mL): the weight of the tissue (g) =19:1. Mechanical homogenate the sample in ice water bath.
- Don't centrifuge, preserve the sample on ice for detection.

## Prepare samples

It is recommended to take 2~3 samples with expected large difference to do a pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (19.42-893.31 U/L).

The recommended dilution factor for different samples is as follows (for reference only):

| Sample type                     | Dilution factor |
|---------------------------------|-----------------|
| Human serum                     | 1               |
| Human plasma                    | 1               |
| Human milk                      | 1               |
| Cell culture supernatant        | 1               |
| Rat serum                       | 1               |
| Rat plasma                      | 1               |
| 5% Rat kidney tissue homogenate | 1               |
| 5% Rat spleen tissue homogenate | 1               |

Note: The diluent is powder A application solution.

## Preparation of buffer solution application solution

Mix 1 part buffer solution with 9 parts distilled water. The prepared solution can be stored at 2-8 °C for 1 month.

## Preparation of powder A application solution

Add vial of powder A with 60 mL buffer solution application solution, and heat at 37 °C to dissolve, store at 2-8 °C for 2 weeks.

## Preparation of powder B application solution

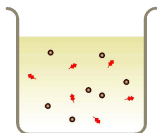
Dissolve vial of powder B with 3 mL Saline Solution. The prepared solution can be store at 2-8 °C for 2 weeks.

## Preparation of chromogenic agent

Dissolve vial of powder C with 12.5 mL buffer application solution, then add 12.5  $\mu$ L of Substrate solution. Mix fully and store at 2-8 °C with shading light.

Note: If clarificant is frozen, shake in 37°C water bath to redissolve fully before use.

## Assay Protocol

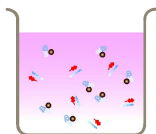


### 1. Sample pretreatment

- Tissue sample: Take 90  $\mu\text{L}$  of tissue homogenate and add 10  $\mu\text{L}$  of powder B application solution, mix fully and incubate at 37  $^{\circ}\text{C}$  for 15 min.
- Serum/plasma: Take 45  $\mu\text{L}$  of sample and add 45  $\mu\text{L}$  of powder A application solution, mix fully, then add 10  $\mu\text{L}$  of powder B application solution and incubate at 37  $^{\circ}\text{C}$  for 15 min.

### 2. Add sample and control

- Sample tube: add 20  $\mu\text{L}$  of sample, 20  $\mu\text{L}$  of clarificant, and 350  $\mu\text{L}$  of chromogenic agent into 1.5 mL microcentrifuge tubes.
- Control well: add 350  $\mu\text{L}$  of distilled water, 20  $\mu\text{L}$  of sample, and 20  $\mu\text{L}$  of clarificant into 1.5 mL microcentrifuge tubes.



### 3. Add substrate

- Oscillate fully with a vortex mixer and incubate for 30 min at 37  $^{\circ}\text{C}$ .
- Add 5  $\mu\text{L}$  of acid reagent, oscillate fully with a vortex mixer and incubate for 10 min at 60  $^{\circ}\text{C}$ .
- Centrifuge the tubes at 3000 g for 10 min and take 300  $\mu\text{L}$  of supernatant for measuring the OD value.
- Measure the OD value with microplate reader at 460 nm.



Target



Horseradish  
peroxidase



Substrate



Enzyme

## Calculation

### Serum (plasma) and other liquid sample:

Unit definition: the amount of MPO in 1 L of sample that catalyze decomposition of 1  $\text{H}_2\text{O}_2$  at 37 °C for 30 min is defined as 1 unit.

$$\text{MPO activity (U/L)} = \frac{\Delta A}{11.3^* \times b} \times V_{\text{Total}} \div \left( \frac{V_{\text{Sample}}}{V_1} \times V_2 \right) \times 1000 \times f = \frac{0.175 \times 1000 \times \Delta A}{V_{\text{Sample}}} \times f$$

### Tissue sample:

Unit definition: the amount of MPO in 1 g wet weight of tissue that catalyze decomposition of 1  $\mu\text{mol H}_2\text{O}_2$  at 37 °C for 30 min is defined as 1 unit.

$$\text{MPO activity (U/g wet weight)} = \frac{\Delta A}{11.3^* \times b} \times V_{\text{Total}} \div \left( \frac{m}{V_3} \times V_2 \times 0.9 \right) = \frac{1.942 \times V_3 \times \Delta A}{m} \times f$$

[Note]

$\Delta A$ :  $\text{OD}_{\text{sample}} - \text{OD}_{\text{control}}$ ;

11.3\*: Constant;

b: Optical path, 1 cm;

$V_{\text{Total}}$ : The total volume of reaction system, 0.395 mL;

$V_{\text{Sample}}$ : The volume of sample added in sample pretreatment step for serum (plasma) and milk sample, 0.045 mL;

$V_1$ : The total volume in sample pretreatment step,  $0.045 + 0.045 + 0.01 = 0.1 \text{ mL}$  or  $0.09 + 0.01 = 0.1 \text{ mL}$ ;

$V_2$ : The volume of sample added to reaction system, 0.02 mL;

$V_3$ : The volume of powder A application solution added into tissue sample in sample preparation step;

1000: 1 L=1000 mL

m: Wet weight of sample, g;

0.9: The ratio of sample volume and total volume in sample pretreatment step,  $0.09 \text{ mL} / 0.1 \text{ mL} = 0.9$ ;

f: Dilution factor of sample before tested

**To easy calculate the test results, refer to the calculation file available on the webpage.**

### Example analysis

For human serum, take 45 µL of human serum, and carry the assay according to the operation steps. The results are as follows: the average OD value of the sample is 0.116, the average OD value of the control is 0.061, and the calculation result is:

$$\text{MPO activity (U/L)} = \frac{0.116-0.061}{0.045} \times 0.175 \times 1000 = 213.89 \text{ U/L}$$

### Performance characteristics

#### ■ Intra-assay Precision

Three human serum samples were assayed in replicates of 20 to determine precision within an assay.

| Parameters | Sample 1 | Sample 2 | Sample 3 |
|------------|----------|----------|----------|
| Mean (U/L) | 36.50    | 364.50   | 579.60   |
| %CV        | 5.8      | 5.4      | 5.0      |

CV = Coefficient of Variation

#### ■ Inter-assay Precision

Three human serum samples were assayed 20 times in duplicate by three operators to determine precision between assays.

| Parameters | Sample 1 | Sample 2 | Sample 3 |
|------------|----------|----------|----------|
| Mean (U/L) | 36.50    | 364.50   | 579.60   |
| %CV        | 7.2      | 7.5      | 7.2      |

CV = Coefficient of Variation

### ▪ Expected values

This assay was tested with human serum, and plasma samples at dilutions from 1:10 to 1:60 in Assay Buffer

| Sample Type  | Range (U/L) | Average (U/L) |
|--------------|-------------|---------------|
| Rat serum    | 400-500     | 442.5         |
| Human plasma | 141-170     | 153.6         |

### ▪ Recovery

Take three samples of high concentration, middle concentration and low concentration to test the samples of each concentration for 6 times parallelly to get the average recovery rate of 104%.

|                      | Sample 1<br>(low conc.) | Sample 2<br>(middle conc.) | Sample 3<br>(high conc.) |
|----------------------|-------------------------|----------------------------|--------------------------|
| Expected Conc. (U/L) | 95.6                    | 342.5                      | 678.2                    |
| Observed Conc. (U/L) | 95.6                    | 363.1                      | 718.9                    |
| Recovery rate (%)    | 100                     | 106                        | 106                      |



▪ Recommended Plate Set Up

|  | 1  | 2   | 3   | 4    | 5   | 6    | 7   | 8    | 9   | 10   | 11  | 12   |
|--|----|-----|-----|------|-----|------|-----|------|-----|------|-----|------|
| A  | S1 | S1' | S9  | S9'  | S17 | S17' | S25 | S25' | S33 | S33' | S41 | S41' |
| B  | S2 | S2' | S10 | S10' | S18 | S18' | S26 | S26' | S34 | S34' | S42 | S42' |
| C  | S3 | S3' | S11 | S11' | S19 | S19' | S27 | S27' | S35 | S35' | S43 | S43' |
| D  | S4 | S4' | S12 | S12' | S20 | S20' | S28 | S28' | S36 | S36' | S44 | S44' |
| E  | S5 | S5' | S13 | S13' | S21 | S21' | S29 | S29' | S37 | S37' | S45 | S45' |
| F  | S6 | S6' | S14 | S14' | S22 | S22' | S30 | S30' | S38 | S38' | S46 | S46' |
| G  | S7 | S7' | S15 | S15' | S23 | S23' | S31 | S31' | S39 | S39' | S47 | S47' |
| H  | S8 | S8' | S16 | S16' | S24 | S24' | S32 | S32' | S40 | S40' | S48 | S48' |
| [Note]: S1-S48, sample wells; S1'-S48', control wells. |    |     |     |      |     |      |     |      |     |      |     |      |

## ■ Sensitivity

The analytical sensitivity of the assay is 19.42 U/L. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 20 times, and calculating the corresponding concentration.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at

**[www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html)**. If you have any questions, please contact Life Technologies at

**[www.thermofisher.com/support](http://www.thermofisher.com/support)**.

Corporate entity: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

© 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.