

EnzyFluo™ Myeloperoxidase Assay Kit (EMPO-100)

Quantitative Fluorimetric Determination of Myeloperoxidase Peroxidation

DESCRIPTION

MYELOPEROXIDASE (MPO; EC 1.11.2.2) is a peroxidase enzyme and can be found in neutrophil, monocytes, and some soft tissue macrophages. MPO has an ability to use chloride as a cosubstrate with hydrogen peroxide to generate hypochlorous acid, a powerful antimicrobial agent produced by neutrophils. However, an excessive production of hypochlorous acid can lead to oxidative stress and tissues damage. Inflammation may also result when MPO oxidizes various substances such as phenols and anilines. Studies show that increased MPO levels may increase the risk of myocardial infarction and cardiovascular disease.

BioAssay Systems' myeloperoxidase (MPO) assay kit is based on the MPO enzyme reaction with hydrogen peroxide (H₂O₂) which oxidizes the dye reagent to a highly fluorescent product. The fluorescence intensity of this product, measured at $\lambda_{\text{ex/em}} = 530/585$ nm, is proportional to the total peroxidation activity in the sample. The provided MPO inhibitor is used to suppress peroxidase activity due to MPO in order to differentiate other peroxidase activities that may be present in the samples.

KEY FEATURES

Fast and sensitive. Linear detection range (20 μ L sample): 0.0025 to 2 U/L for 10 min reaction at 25°C.

Convenient and high-throughput. Homogeneous "mix-incubate-measure" type assay. Can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

APPLICATIONS

MPO peroxidation activity determination in biological samples (e.g. cell lysates, tissues, etc.)

KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

Assay Buffer:	10 mL	Resorufin:	1.5 mL
20x MPO Inhibitor:	120 μ L	Dye Reagent:	120 μ L
3% Stabilized H₂O₂:	100 μ L		

Storage conditions. The kit is shipped on ice. Store all components at -20°C upon receiving. Shelf life: 6 months after receipt.

Precautions: reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

PROCEDURES

This assay is based on a kinetic reaction. To ensure identical incubation time, addition of Working Reagent to samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended. Assays can be executed at any desired temperature (e.g. 25°C or 37°C).

Sample Preparation:

Tissue: prior to dissection, rinse tissue in phosphate buffered saline (pH 7.4) to remove blood. Homogenize tissue (50 mg) with a Dounce homogenizer in ~200 μ L cold 20 mM PBS, pH 7.4. Freeze the homogenized tissue at -80°C to lyse the cells. After freezing, thaw and centrifuge samples at 14,000xg for 20 min at 4°C. Remove supernatant for assay.

Cell Lysate: collect cells by centrifugation at 2,000xg for 5 min at 4°C. For adherent cells, do not harvest cells using proteolytic enzymes; rather use a rubber policeman. Homogenize or sonicate cells in an appropriate volume of cold buffer containing 50 mM potassium phosphate (pH 7.5). Centrifuge at 14,000xg for 10 min at 4°C. Remove supernatant for assay.

All samples can be stored at -20 to -80°C for at least one month.

Reagent Preparation:

Bring all reagents to room temperature prior to assay. Briefly centrifuge tubes before use.

Each sample requires 20 μ L of 1x MPO Inhibitor. Prior to assay, Prepare enough 1x MPO Inhibitor by diluting the provided 20x MPO Inhibitor 20-fold in dH₂O.

Assay Procedure:

- Prepare 250 μ L 30 μ M Resorufin Premix by mixing 15 μ L provided Resorufin and 235 μ L water.
Transfer 100 μ L water and 100 μ L 30 μ M Resorufin into two separate wells of a black flat-bottom 96-well plate.
- For each sample prepare 2 parallel wells. Add 20 μ L of samples to each wells. Add 20 μ L of 1x MPO inhibitor to one of each sample's wells and add 20 μ L assay buffer to the other well. Incubate samples at room temperature for 10 min.
- Prepare 0.07% H₂O₂ by mixing 4.7 μ L 3% H₂O₂ with 195.3 μ L dH₂O. Then to 0.007% H₂O₂ by mixing 60 μ L 0.07% H₂O₂ with 540 μ L dH₂O. Use the 0.007% H₂O₂ within one hour.

Prepare enough Working Reagent (WR) for all reaction wells by mixing, for each 96-well assay, 60 μ L Assay Buffer, 1 μ L 0.007% H₂O₂ and 1 μ L Dye Reagent. Add 60 μ L WR to all sample and inhibitor wells. Tap plate briefly to mix.

- Read fluorescence $\lambda_{\text{ex/em}} = 530/585$ nm at 0 min and 10 min at room temperature.

CALCULATION

The MPO activity in a sample is computed as follows:

$$\text{MPO Activity} = \frac{\Delta R_{\text{SAMPLE}} - \Delta R_{\text{INB}}}{R_{\text{RESORUFIN}} - R_{\text{H}_2\text{O}}} \times \frac{[\text{Resorufin}] (\mu\text{M})}{t (\text{min})} \times \frac{\text{Reaction Vol} (\mu\text{L})}{\text{Sample Vol} (\mu\text{L})} \times n$$

$$= \frac{\Delta R_{\text{SAMPLE}} - \Delta R_{\text{INB}}}{R_{\text{RESORUFIN}} - R_{\text{H}_2\text{O}}} \times 15 \times n \quad (\text{U/L})$$

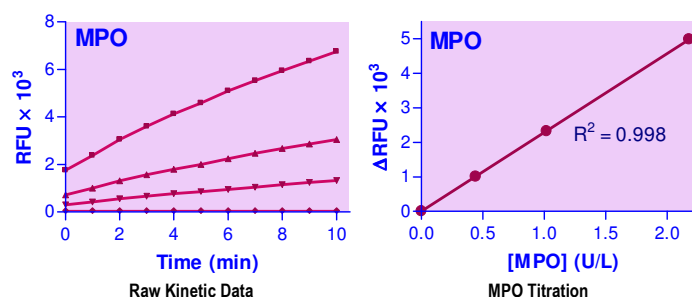
where R_{SAMPLE} , R_{INB} , $R_{\text{RESORUFIN}}$ and $R_{\text{H}_2\text{O}}$ are fluorescence readings of the Sample, Sample Inhibitor, Resorufin and Water wells, respectively. $\Delta R_{\text{SAMPLE}} = R_{\text{SAMPLE},10\text{min}} - R_{\text{SAMPLE},0\text{min}}$ and $\Delta R_{\text{INB}} = R_{\text{INB},10\text{min}} - R_{\text{INB},0\text{min}}$. n is the sample dilution factor. [Resorufin] = 30 μ M, Reaction Vol = 100 μ L, Sample Vol = 20 μ L, Reaction time (t) = 10 min.

Notes: if ΔR_{SAMPLE} values are higher than that of the $R_{\text{RESORUFIN}}$, dilute sample in Assay Buffer and repeat the assay. Multiply the results by the dilution factor, n .

Unit definition: one unit of enzyme will catalyze the formation of 1 μ mole resorufin per min under the assay conditions.

MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, black 96-well plates (e.g. Greiner Bio-One, cat# 655900) and plate reader capable of measuring fluorescence at $\lambda_{\text{ex/em}} = 530/585$ nm.



PUBLICATIONS

- Son, A., et al (2019). Ca²⁺ influx channel inhibitor saraf protects mice from acute pancreatitis. *Gastroenterology*, 157(6), 1660-1672.e2.
- Hashim, A. A., Helmy, M. M., & Mouneir, S. M. (2018). Cysteinyl leukotrienes predominantly mediate cisplatin-induced acute renal damage in male rats. *Journal of Physiology and Pharmacology*, 69(5), 779-787.
- Helmy, M. M., Hashim, A. A., & Mouneir, S. M. (2018). Zileuton alleviates acute cisplatin nephrotoxicity: Inhibition of lipoxygenase pathway favorably modulates the renal oxidative/inflammatory/caspase-3 axis. *Prostaglandins & other lipid mediators*, 135, 1-10.

