

ab283410 – Myeloperoxidase Inhibitor Screening Kit (Fluorometric)

For the screening of potential Myeloperoxidase inhibitors.

For research use only - not intended for diagnostic use.

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab283410>

Storage and Stability

On receipt entire assay kit should be stored at -20°C, protected from light. Upon opening, use kit within 6 months.

Materials Supplied

Item	Quantity	Storage Condition
Assay Buffer LVI/Myeloperoxidase Assay Buffer	50 ml	-20°C
MPO Chlorination Probe/Myeloperoxidase Chlorination Substrate	200 µl	-20°C
OxiRed Probe/Myeloperoxidase Peroxidation Substrate	200 µl	-20°C
MPO Enzyme/Myeloperoxidase Enzyme	1 vial	-20°C
MPO Inhibitor Control/Inhibitor Control	100 µl	-20°C
Hydrogen Peroxide Solution II/Hydrogen Peroxide	50 µl	-20°C

Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Multi-well spectrophotometer
- 96-well white plate with flat bottom

Reagent Preparation

- Before using the kit, spin the tubes prior to opening.

Assay Buffer LVI/Myeloperoxidase Assay Buffer: Bring the Assay Buffer LVI/Myeloperoxidase Assay Buffer to room temperature (RT) before use. Can store at -20°C or +4°C.

MPO Chlorination Probe/Myeloperoxidase Chlorination and OxiRed Probe/Peroxidation Substrate: Aliquot and store at -20°C. Bring to room temperature before use.

MPO Enzyme/Myeloperoxidase Enzyme: Reconstitute with 1.1 ml Assay Buffer LVI/MPO Assay Buffer. Aliquot and store at -70°C. Avoid freeze/thaw. Keep on ice while in use. Use within two months.

MPO Inhibitor Control/Inhibitor Control: Bring to RT before use.

Myeloperoxidase Chlorination Screening Protocol

Screening Compounds, Inhibitor Control, and Enzyme Control Preparation:

1. Dissolve candidate inhibitors into an appropriate solvent at 100X the highest final concentration to be tested.
 2. Dilute to 2X desired test concentration with Assay Buffer LVI/MPO Assay Buffer.
 3. Add 45 µl diluted candidate inhibitor or Assay Buffer LVI/MPO Assay Buffer into desired wells, as Sample Screen [S], or Enzyme Control [EC] (no inhibitor).
 4. For Inhibitor Control (IC), dilute MPO Inhibitor Control/Inhibitor Control 50 times by adding 5 µl MPO Inhibitor Control/Inhibitor Control to 245 µl Assay Buffer LVI/MPO Assay Buffer.
 5. Add 45 µl of diluted MPO Inhibitor Control/Inhibitor Control into desired well(s).
- Δ Note:** *Diluted MPO Inhibitor Control/Inhibitor Control is stable for 4 hours at 4°C.*

Enzyme Solution Preparation:

1. Add 5 µl MPO Enzyme/Myeloperoxidase Enzyme into Sample, Enzyme Control and Inhibitor Control wells.

Substrate Solution Preparation:

1. Prepare 5 mM Hydrogen Peroxide solution by adding 1 µl Hydrogen Peroxide Solution II/Hydrogen Peroxide to 175 µl Assay Buffer LVI/MPO Assay Buffer. Mix. Make enough reagents for the number of assays to be performed. For each well, prepare 50 µl of Substrate solution containing:

	Volume
Assay Buffer LVI/Myeloperoxidase Assay Buffer	46 µl
MPO Chlorination Probe/Myeloperoxidase Chlorination Substrate	2 µl
Diluted Hydrogen Peroxide Solution II/Hydrogen Peroxide	2 µl

2. Mix and add 50 µl of Substrate solution into each well. Mix well. Incubate for 10 minutes at room temperature with gentle shaking, protected from light.

Δ Note: *Do not store diluted Hydrogen Peroxide solution.*

Measurement:

1. Measure fluorescence (Ex/Em = 480/520 nm).

Myeloperoxidase Peroxidation Screening Protocol

Screening Compounds, Inhibitor Control, and Enzyme Control Preparation:

1. Dissolve candidate inhibitors into an appropriate solvent at 100X the highest final concentration to be tested.
2. Dilute to 2X desired test concentration with Assay Buffer LVI/MPO Assay Buffer.
3. Add 45 µl diluted candidate inhibitor or Assay Buffer LVI/MPO Assay Buffer into desired wells, as Sample Screen [S], or Enzyme Control [EC] (no inhibitor).
4. For Inhibitor Control (IC), dilute MPO Inhibitor Control/Inhibitor Control 50 times by adding 5 µl MPO Inhibitor Control/Inhibitor Control to 245 µl Assay Buffer LVI/MPO Assay Buffer.
5. Add 45 µl of diluted MPO Inhibitor Control/Inhibitor Control into desired well(s).

Δ Note: *Diluted MPO Inhibitor Control/Inhibitor Control is stable for 4 hours at 4°C.*

Enzyme Solution Preparation:

1. Mix enough reagents for the number of assays to be performed. For each well, prepare 5 µl of MPO enzyme solution 5 µl MPO Enzyme/Myeloperoxidase Enzyme into Sample, Enzyme Control, and Inhibitor Control wells.

	Volume
Assay Buffer LVI/Myeloperoxidase Assay Buffer	2.5 µl
MPO Enzyme/Myeloperoxidase Enzyme	2.5 µl

2. Mix and add 5 µl of MPO Enzyme Solution into Sample, Enzyme Control, and Inhibitor Control wells.

Substrate Solution Preparation:

1. Prepare 5 mM Hydrogen Peroxide solution by adding 1 µl Hydrogen Peroxide Solution II/Hydrogen Peroxide to 175 µl Assay Buffer LVI/MPO Assay Buffer. Mix. Make enough reagents for the number of assays to be performed. For each well, prepare 50 µl of Substrate solution containing:

	Volume
Assay Buffer LVI/Myeloperoxidase Assay Buffer	46 µl
OxiRed Probe/Myeloperoxidase Peroxidation Substrate	2 µl
Diluted Hydrogen Peroxide Solution II/Hydrogen Peroxide	2 µl

2. Mix and add 50 µl of Substrate solution into each well. Mix well. Incubate for 5 minutes at room temperature with gentle shaking, protected from light.

Measurement:

1. Measure fluorescence (Ex/Em = 535/587 nm).

Calculations:

1. Set the RFU of Enzyme Control (EC) as 100%, and calculate the relative inhibition (%) of the candidate inhibitors as:

$$\% \text{ Inhibitor} = \frac{RFU (EC) - RFU (S)}{RFU (EC)} \times 100$$

Where RFU (EC) is the fluorescence reading of Enzyme Control and RFU (S) is the fluorescence reading of Sample Screen.

Technical Support

Copyright © 2023 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print. For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)