

Product datasheet

Human Myeloperoxidase ELISA Kit (MPO) ab119605

★★★★☆ 1 Abreviews 5 References 1 Image

Overview

| | |
|----------------------------|---|
| Product name | Human Myeloperoxidase ELISA Kit (MPO) |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant, Saliva, Urine, Serum, Cell Lysate, Heparin Plasma, EDTA Plasma, Tissue Homogenate |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 10 pg/ml |
| Range | 312 pg/ml - 20000 pg/ml |
| Assay duration | Multiple steps standard assay |
| Species reactivity | Reacts with: Human |
| Product overview | <p>Abcam's Human Myeloperoxidase (MPO) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Myeloperoxidase in cell culture supernatants, cell lysate, tissue homogenates, serum, plasma (heparin, EDTA), saliva and urine.</p> <p>A Myeloperoxidase specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat specific for Myeloperoxidase is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human Myeloperoxidase amount of sample captured in plate.</p> <p>Get results in 90 minutes with Human Myeloperoxidase ELISA Kit (ab195212) from our SimpleStep ELISA® range.</p> |
| Tested applications | Suitable for: Sandwich ELISA |
| Platform | Microplate |

Properties

Storage instructions Store at -20°C. Please refer to protocols.

| Components | Identifier | 1 x 96 tests |
|---|--------------|--------------|
| ABC Diluent Buffer | Blue Cap | 1 x 12ml |
| Antibody Diluent Buffer | Green Cap | 1 x 12ml |
| Anti-Human Myeloperoxidase antibody Microplate (12 x 8 wells) | | 1 unit |
| Avidin-Biotin-Peroxidase Complex (ABC) | | 1 x 130µl |
| Biotinylated anti-Human Myeloperoxidase antibody | | 1 x 130µl |
| Lyophilized recombinant Human Myeloperoxidase standard | | 2 x 20ng |
| Plate Seal | | 4 units |
| Sample Diluent Buffer | Green Cap | 1 x 30ml |
| TMB Color Developing Agent | Amber Bottle | 1 x 10ml |
| TMB Stop Solution | Yellow Cap | 1 x 10ml |

| | |
|-------------------------------|---|
| Function | Part of the host defense system of polymorphonuclear leukocytes. It is responsible for microbicidal activity against a wide range of organisms. In the stimulated PMN, MPO catalyzes the production of hypohalous acids, primarily hypochlorous acid in physiologic situations, and other toxic intermediates that greatly enhance PMN microbicidal activity. |
| Involvement in disease | Defects in MPO are the cause of myeloperoxidase deficiency (MPD) [MIM:254600]. MPD is an autosomal recessive defect that results in disseminated candidiasis. |
| Sequence similarities | Belongs to the peroxidase family. XPO subfamily. |
| Cellular localization | Lysosome. |

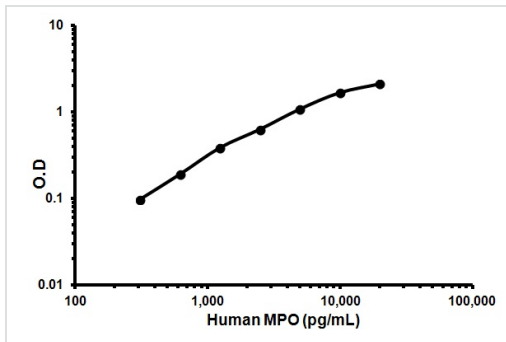
Applications

Our [Abpromise guarantee](#) covers the use of **ab119605** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| Application | Abreviews | Notes |
|----------------|-----------|--|
| Sandwich ELISA | | Use at an assay dependent concentration. |

Images



Representative Standard Curve using ab119605.

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors