

Product datasheet

Human Myeloperoxidase ELISA Kit (MPO) ab119605

★★★★☆ 1 Abreviews 4 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	

Abcam's Human Myeloperoxidase (MPO) *in vitro* 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.

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.

Get results in 90 minutes with Human Myeloperoxidase ELISA Kit ([ab195212](#)) from our SimpleStep ELISA® range.

Tested applications	Suitable for: Sandwich ELISA
Platform	Microplate

Properties

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

Components	1 x 96 tests
ABC Diluent Buffer	1 x 12ml
Antibody Diluent Buffer	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	1 x 30ml
TMB Color Developing Agent	1 x 10ml
TMB Stop Solution	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.

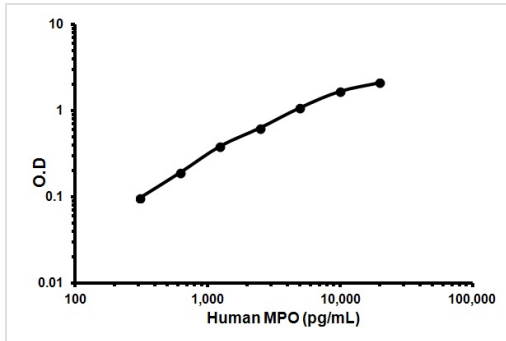
Applications

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