

Human Myeloperoxidase ELISA Kit (MPO) ab119605

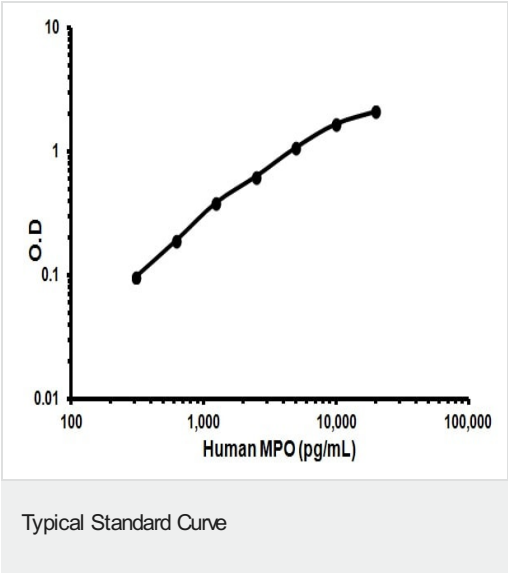
★★★★★ [1 Abreviews](#) [29 References](#) [1 图像](#)

概述	
产品名称	人Myeloperoxidase ELISA试剂盒(MPO)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Saliva, Urine, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate
检测类型	Sandwich (quantitative)
灵敏度	< 10 pg/ml
范围	312 pg/ml - 20000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<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>
平台	Microplate
性能	
存放说明	Store at -20°C. Please refer to protocols.

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Myeloperoxidase antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human Myeloperoxidase antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Myeloperoxidase standard		2 x 20ng	2 x 20ng
Plate Seal		4 units	4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

功能	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.
疾病相关	Defects in MPO are the cause of myeloperoxidase deficiency (MPD) [MIM:254600]. MPD is an autosomal recessive defect that results in disseminated candidiasis.
序列相似性	Belongs to the peroxidase family. XPO subfamily.
细胞定位	Lysosome.

图片



Representative Standard Curve using ab119605.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

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.cn/abpromise> or contact our technical team.

Terms and conditions

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