

Mouse MPO ELISA Kit (Myeloperoxidase) ab155458

★★★★★ [2 Abreviews](#) [23 References](#) [2 Images](#)

Overview

Product name	Mouse MPO ELISA Kit (Myeloperoxidase)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Plasma, Cell Lysate				
Assay type	Sandwich (quantitative)				
Sensitivity	< 0.6 ng/ml				
Range	0.61 ng/ml - 150 ng/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Cell culture supernatant		105.5		96% - 112%
	Plasma		86.89		68% - 105%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's MPO (Myeloperoxidase) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse MPO (Myeloperoxidase) in plasma and cell culture supernatants.				

This assay employs an antibody specific for mouse MPO coated on a 96-well plate. Standards and samples are pipetted into the wells and MPO present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse MPO antibody is added.

After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of MPO bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450nm.

We have not been able to detect the endogenous Mouse MPO (Myeloperoxidase) in normal serum with ab155458, only in serum spiked with Mouse MPO (Myeloperoxidase).

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-mouse MPO	2 vials
MPO Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MPO Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

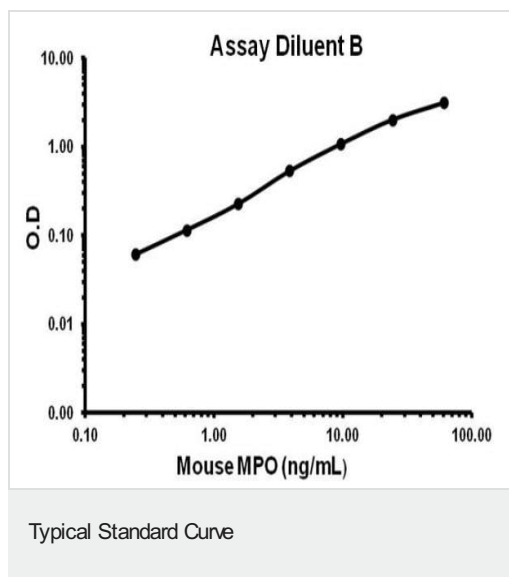
Function Part of the host defense system of polymorphonuclear leukocytes. It is responsible for microbicidal activity against a wide range of organisms. In the stimulated PMN, MPO catalyzes the production of hypohalous acids, primarily hypochlorous acid in physiologic situations, and other toxic intermediates 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.

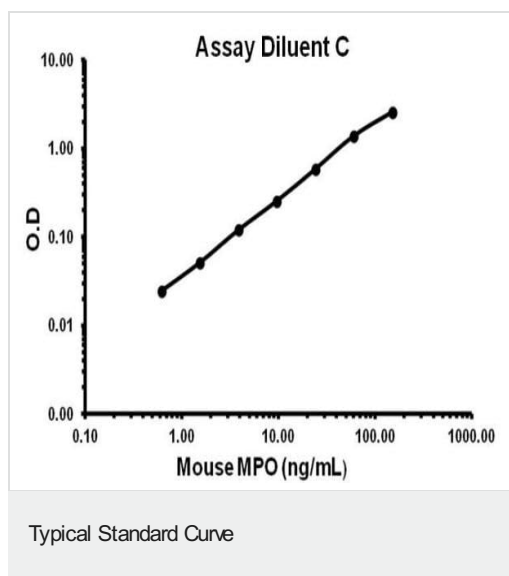
Sequence similarities Belongs to the peroxidase family. XPO subfamily.

Cellular localization Lysosome.

Images



Representative standard curve using ab155458 - Assay Diluent B.



Representative standard curve using ab155458 - Assay Diluent C.

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.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