

Myeloperoxidase (human), ELISA kit

The Myeloperoxidase (human), ELISA kit is a colorimetric immunometric enzyme immunoassay kit

The Myeloperoxidase (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. This sensitive kit lets you quantitate as little as 19 pg/mL MPO in a wide variety of sample matrices that have been validated in-house. Convenient ready-to-use liquid color-coded reagents reduce error and save time in the lab.

Myeloperoxidase (MPO) is synthesized as a single chain precursor, which is then processed into a tetramer consisting of 2 light chains and 2 heavy chains. In response to microbial invasion, MPO is released from the cytoplasmic granules of neutrophils into the extracellular space and phagosome, where it generates hypohalous acids that have a strong microbicidal effect. Myeloperoxidase can be used as a marker of airway inflammation caused either by disease, such as asthma, or by environmental irritants.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

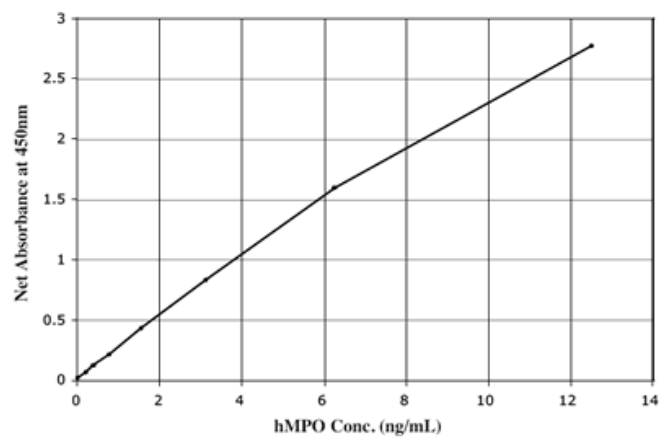
ADI-900-115

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - detect as little as 19 pg/ml MPO
- Reliable - excellent reproducibility and parallelism
- Convenient - ready-to-use liquid color-coded reagents reduce error and save time in the lab
- Higher throughput - results for up to 39 samples in duplicate in just 3 hours



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MPO

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human Myeloperoxidase of culture supernatants, nasal lavage, Lithium-heparin plasma, sputum supernatant, and urine.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Conjugate, Antibody, Assay buffer 13, Assay buffer 31 concentrate, Wash buffer concentrate, Standard, TMB Substrate, Stop solution 2

Crossreactivity Human Myeloperoxidase (100%), human Erythropoietin (1.1%), bovine Lactoperoxidase (0.40%), human Elastase (0.20%), human Lactoferrin (0.20%), sheep COX-2 (0.10%) and <0.10%: human Thyroid Peroxidase, human Eosinophil-derived Neurotoxin

Sensitivity Assay Buffer 13: 0.028ng/mL Assay Buffer 31: 0.019ng/mL (range 0.195 – 12.5 ng/ml)

Species Reactivity Human

UniProt ID P05164

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024