

Myeloperoxidase fluorometric detection kit

The Myeloperoxidase fluorometric detection kit is a one-step no wash homogeneous activity assay that is adaptable to HTS applications. This kit offers highly sensitive fluorescent detection and the ability to monitor multiple time points to follow the kinetics of MPO activity.

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: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

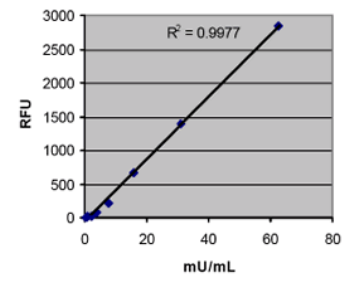
ADI-907-029	500 tests
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

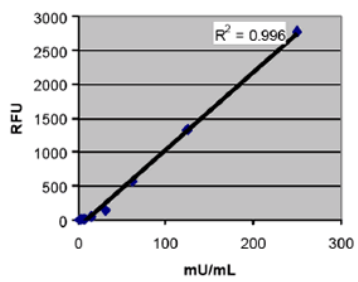
- Fast homogenous assay
- Suitable for High Throughput Screening
- Monitor multiple time points to follow kinetics
- Results in 30 to 60 minutes

MPO without Catalase inhibitor

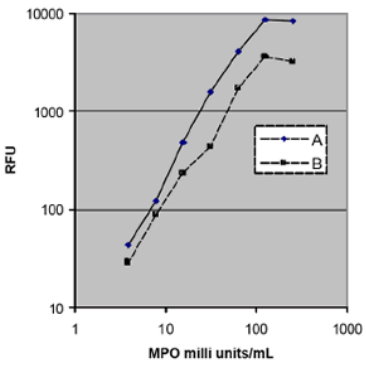


mU/mL	MPO RFU
62.5	2838
31.25	1401
15.625	668
7.8125	207
3.90625	71
1.953125	27
0.976563	14
0.488281	6
0.244141	4

MPO + Catalase inhibitor (20mM)



mU/mL	MPO + Inhib RFU
250	2783
125	1401
62.5	567
31.25	143
15.625	46
7.8125	12
3.90625	6
1.953125	2



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MPO

Application Fluorescence microscopy

Application Notes For the measurement of Myeloperoxidase activity in cells and tissue from any species.

Contents Assay buffer concentrate, Detection Reagent, Hydrogen peroxide, Myeloperoxidase enzyme

Sensitivity 0.00195 units/ml

Species Reactivity Species independent

UniProt ID P05164



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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