

Myeloperoxidase Activity Assay

A plate based colorimetric assay to measure the enzymatic activity of MPO in a wide range of sample types.

Myeloperoxidase (MPO) is an oxidative enzyme abundantly expressed in neutrophils and is involved in the immune response and inflammation. Additionally, MPO can also oxidize lipids, proteins and DNA, leading to oxidative damage in chronic inflammatory conditions such as atherosclerosis, COPD and neurodegenerative disease. AkrivisBio's MPO Assay is a simple sensitive means of measuring MPO enzyme activity below 0.05 mU per well, which can be indicative of neutrophil activation and oxidative stress. In the assay, Myeloperoxidase produces HOCl (hypochlorous acid) from hydrogen peroxide and chloride. HOCl reacts with taurine forming taurine chloramine, which then reacts with TNB, the reduced form of DTNB, Ellman's Reagent, bleaching out the color causing a decrease of absorbance measured at OD 412 nm. The absorbance is inversely proportional to the amount of MPO enzyme.

Ordering Information

[Order Online »](#)

AKR-MA-0135	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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