

Myeloperoxidase (MPO) Activity Assay Kit (Colorimetric) ab105136

★★★★★ [3 Abreviews](#) [67 References](#) [3 图像](#)

概述

产品名称	Myeloperoxidase (MPO) Activity Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Urine, Serum, Other biological fluids, Tissue Homogenate
检测类型	Enzyme activity
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	<p>Myeloperoxidase (MPO) Activity Assay Kit ab105136 is a rapid, simple, sensitive, and reliable colorimetric assay suitable for use as a high throughput MPO activity assay.</p> <p>In the MPO assay protocol, myeloperoxidase produces HClO from H₂O₂ and Cl⁻. The HClO reacts with taurine to generate the taurine chloramine, which subsequently reacts with the DTNB probe to eliminate color (absorbance at 412 nm). The absorbance is inversely proportional to the amount of MPO enzyme.</p> <p>This MPO assay kit can be used to detect myeloperoxidase as low as 0.05 mU per well.</p> <p>MPO assay protocol summary:</p> <ul style="list-style-type: none"> - add samples to wells (and assay buffer only to wells for standards) - add reaction mix to sample wells and incubate for 30 min to 2 hrs - add stop mix and incubate for 10 min - add TNB reagent/standard to all wells and incubate for 5-10 min - analyze with a microplate reader
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K744 Myeloperoxidase (MPO) Colorimetric Activity Assay Kit. K744-100 is the same size as the 100 test size of ab105136.</p> <p>Myeloperoxidase (MPO) is a peroxidase enzyme (EC 1.11.1.7) most abundantly expressed in neutrophil granulocytes. It is a lysosomal protein stored in the azurophilic granules of the neutrophil. Myeloperoxidase contains a heme pigment which causes its green color in secretions rich in neutrophils, such as pus and some forms of mucus. Myeloperoxidase catalyzes the production of hypochlorous acid (HClO) from hydrogen peroxide (H₂O₂) and chloride anion (Cl⁻, or the equivalent from a non-chlorine halide). Myeloperoxidase also oxidizes tyrosine to a tyrosyl</p>

radical using hydrogen peroxide as an oxidizing agent.

平台

Microplate reader

性能

存放说明

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

组件	100 tests
DTNB Probe	1 x 50µl
MPO Assay Buffer II	1 x 25ml
MPO Positive Control	1 vial
Hydrogen Peroxide Solution II	1 x 50µl
Stop Mix	1 x 20µl
TCEP	1 x 50µl

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.

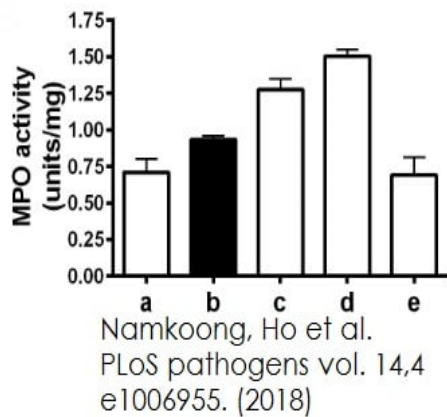
序列相似性

Belongs to the peroxidase family. XPO subfamily.

细胞定位

Lysosome.

图片

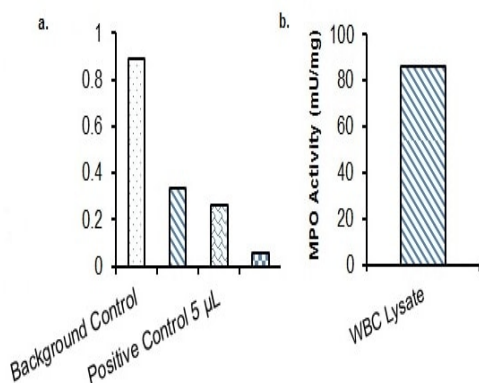


MPO activity in indicated cells was measured using the commercially available MPO Activity Assay Kit (n = 3).

Functional Studies - Myeloperoxidase (MPO)

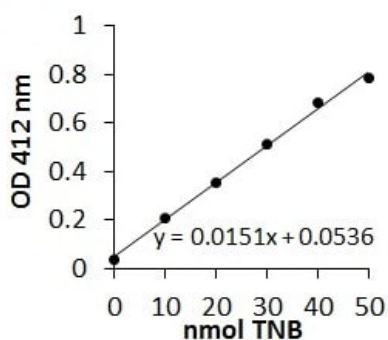
Activity Assay Kit (Colorimetric) (ab105136)

Namkoong, Ho et al., PLoS pathogens vol. 14,4 e1006955., S4 fig, doi:10.1371/journal.ppat.1006955



a. Measurement of MPO activity using WBC lysate (3 µg) and MPO Positive Control (5 and 10 µL). b. MPO specific activity in WBC lysate.

MPO activity



Typical standard curve for ab105136.

Typical Standard Curve

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