

## Product datasheet

# Human Macrophage Stimulating Protein ELISA Kit ab100612

## 2 图像

### 概述

产品名称	人Macrophage Stimulating蛋白ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 8 pg/ml
范围	8.23 pg/ml - 2000 pg/ml
回收率	95 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	97.25	86% - 106%
Serum	94.48	83% - 104%
Plasma	95.29	82% - 105%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's Macrophage Stimulating Protein Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Macrophage Stimulating Protein in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Macrophage Stimulating Protein coated on a 96-well plate. Standards and samples are pipetted into the wells and Macrophage Stimulating Protein present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Macrophage Stimulating Protein 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 Macrophage Stimulating Protein bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**说明** Optimization may be required with urine samples.

**经测试应用** 适用于: Sandwich ELISA

平台 Microplate

## 性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Macrophage Stimulating Protein	2 vials
Macrophage Stimulating Protein Microplate (12 x 8 wells)	1 unit
Recombinant Human Macrophage Stimulating Protein Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**相关性** Macrophage Stimulating Protein contains four kringle domains and a serine protease domain, similar to that found in hepatic growth factor. Despite the presence of the serine protease domain, the encoded protein may not have any proteolytic activity. The receptor for this protein is RON tyrosine kinase, which upon activation stimulates ciliary motility of ciliated epithelial lung cells. This protein is secreted and cleaved to form an alpha chain and a beta chain bridged by disulfide bonds.

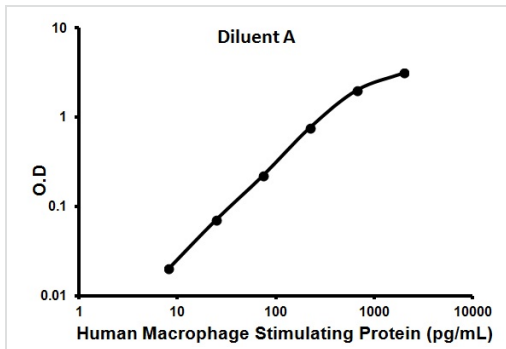
## 应用

Our [Abpromise guarantee](#) covers the use of **ab100612** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

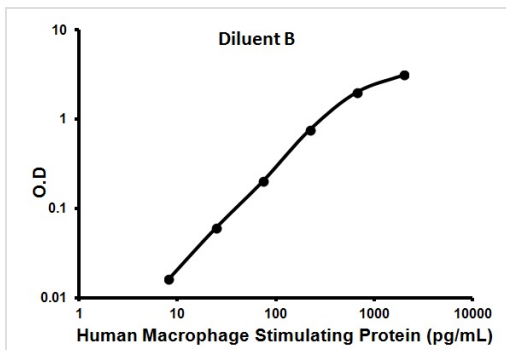
应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

## 图片



Representative Standard Curve using ab100612.

Typical Standard Curve



Representative Standard Curve using ab100612.

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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