

## Product datasheet

# Human Anti-Respiratory syncytial virus IgM ELISA Kit (RSV) ab108766

### 概述

**产品名称** 人Anti-Respiratory syncytial virus IgM ELISA试剂盒(RSV)

**检测方法** Colorimetric

### 精确度

批次内

样品	n	Mean	SD	CV%
Pos.Serum	24			2.6%

批次间

样品	n	Mean	SD	CV%
Pos.Serum	12			5.4%

**样品类型** Serum, Plasma

**检测类型** Indirect

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's anti-Respiratory syncytial virus (RSV) IgM Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Respiratory syncytial virus in Human serum and plasma.

A 96-well plate has been precoated with Respiratory syncytial virus antigens to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidase (HRP) labelled anti-Human IgM conjugate is added to the wells, which binds to the immobilized Respiratory syncytial virus -specific antibodies. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Respiratory syncytial virus IgM sample captured in plate.

**经测试应用** 适用于: Indirect ELISA

**平台** Microplate

### 性能

## 存放说明

Store at +4°C. Please refer to protocols.

组件	标识符	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Anti-Respiratory syncytial virus (IgM) Coated Microplate (12 x 8 wells)		1 unit
IgM Sample Diluent	Colored Green; White cap	1 x 100ml
Respiratory syncytial virus anti-IgM HRP Conjugate	Colored Red; Black cap	1 x 20ml
Respiratory syncytial virus IgM Cut-off Control	Colored yellow; Green cap	1 x 3ml
Respiratory syncytial virus IgM Negative Control	Colored yellow; Blue cap	1 x 2ml
Respiratory syncytial virus IgM Positive Control	Colored yellow; Red cap	1 x 2ml
Stop Solution		1 x 15ml
TMB Substrate Solution	Yellow cap	1 x 15ml

## 应用

Our [Abpromise guarantee](#) covers the use of **ab108766** 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
Indirect ELISA		Use at an assay dependent concentration.

**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 <http://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