

## Product datasheet

# Human Anti-Influenza virus A IgA ELISA Kit ab108743

### 概述

**产品名称** 人Anti-Influenza virus A IgA ELISA试剂盒

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
Pos. Serum	8			2.1%

批次间

样品	n	Mean	SD	CV%
Pos. Serum	19			3.7%

**样品类型** Serum, Heparin Plasma, Citrate 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-Influenza virus A IgA Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA class antibodies against Influenza virus A in Human serum and plasma.

A 96-well plate has been precoated with Influenza virus A 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 IgA conjugate is added to the wells, which binds to the immobilized Influenza virus A-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 Influenza virus A IgA 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
Cover foil		1 unit
IgA Sample Diluent	Yellow, White cap	1 x 100ml
Influenza virus A (IgA) Coated Microplate (12 x 8 wells)		1 unit
Influenza virus A anti-IgA HRP Conjugate	Violet, Black cap	1 x 20ml
Influenza Virus A IgA Cut-off Control	Yellow, Green cap	1 x 3ml
Influenza Virus A IgA Negative Control	Yellow, Blue cap	1 x 2ml
Influenza Virus A IgA Positive Control	Yellow, Red cap	1 x 2ml
Stop Solution		1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

## 应用

Our [Abpromise guarantee](#) covers the use of **ab108743** 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 dilution.

**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