

# Human Anti-Influenza virus A IgG ELISA Kit ab108745

2 References 1 Image

Overview

Product name Human Anti-Influenza virus A IgG ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	20			4.5%
	Pos.Serum	24			3.3%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	12			3.2%
	Pos.Serum	12			2.1%

Sample type Serum, Hep Plasma, C'it plasma

Assay type Indirect

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

**Product overview**

Abcam's anti-Influenza virus A IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG 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 IgG 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 IgG sample captured in plate.

Platform Microplate

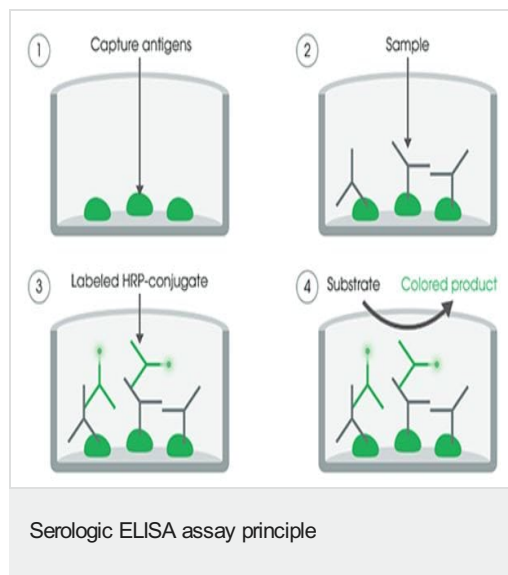
## Properties

### Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
IgG Sample Diluent	Yellow, White cap	1 x 100ml
Influenza virus A (IgG) Coated Microplate (12 x 8 wells)		1 unit
Influenza virus A anti-IgG HRP Conjugate	Blue, Black cap	1 x 20ml
Influenza virus A IgG Cut-off Control	Yellow, Green cap	1 x 3ml
Influenza virus A IgG Negative Control	Yellow, Blue cap	1 x 2ml
Influenza virus A IgG Positive Control	Yellow, Red cap	1 x 2ml
Stop Solution	red cap	1 x 15ml
TMB Substrate Solution	Yellow cap	1 x 15ml

## Images



Specific antigens are coated on the 96-well plate, controls or test samples are added to the well and incubated. The wells are washed to remove any unbound Human anti-antigen antibodies (Ig). A horseradish peroxidase (HRP) labelled anti-Human Ig conjugate is added to the wells. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of yellow coloration is directly proportional to the amount of Human anti-antigen Ig captured on the plate.

**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.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors