

Human Anti-Parainfluenza virus 1,2,3 IgG ELISA Kit

ab108758

1 Image

Overview

Product name Human Anti-Parainfluenza virus 1,2,3 IgG ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	8			6.8%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	3			9.8%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Indirect

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's anti-Parainfluenza virus 1, 2, 3 IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Parainfluenza virus 1, 2, 3 in Human serum and plasma.

A 96-well plate has been precoated with Parainfluenza virus 1, 2, 3 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 Parainfluenza virus 1, 2, 3-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 Parainfluenza virus 1, 2, 3 IgG sample captured in plate.

Platform Microplate

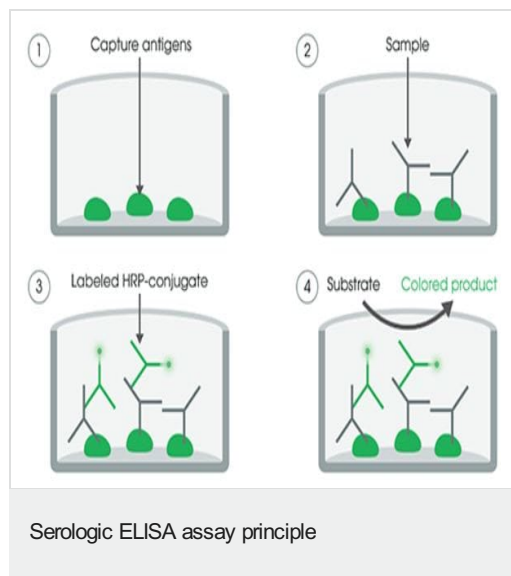
Properties

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	white cap	1 x 100ml
Parainfluenza virus 1, 2, 3 (IgG) Coated Microplate (12 x 8 wells)		1 unit
Parainfluenza virus 1, 2, 3 anti-IgG HRP Conjugate	black cap	1 x 20ml
Parainfluenza virus 1, 2, 3 IgG Cut-off Control	green cap	1 x 3ml
Parainfluenza virus 1, 2, 3 IgG Positive Control	red cap	1 x 2ml
Parainfluenza Virus IgG Negative Control	blue 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