

Human Anti-Parainfluenza virus 1,2,3 IgA ELISA Kit ab108757

1 图像

概述

产品名称 人Anti-Parainfluenza virus 1,2,3 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			2.6%

批次间

样品	n	Mean	SD	CV%
Pos.Serum	24			5.1%

样品类型 Serum, Hep Plasma, Cit 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-Parainfluenza virus 1, 2, 3 IgA Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA 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 IgA 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 IgA sample captured in plate.

平台 Microplate

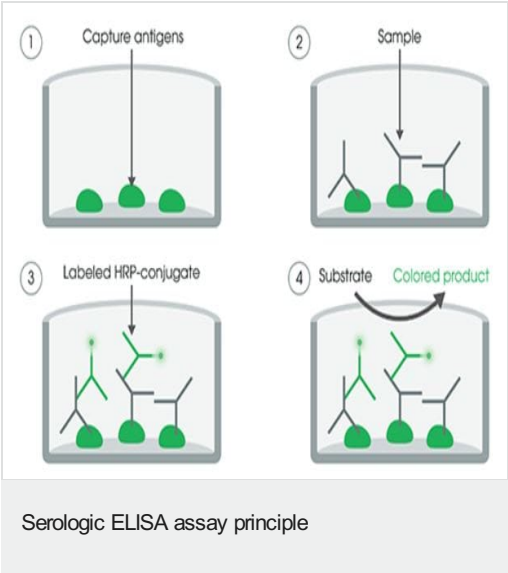
性能

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
Parainfluenza virus 1, 2, 3 (IgA) Coated Microplate (12 x 8 wells)	12 Strips of 8 wells	1 unit
Parainfluenza virus 1, 2, 3 anti-IgA HRP Conjugate	Purple, Black cap	1 x 20ml
Parainfluenza virus 1, 2, 3 IgA Cut-off Control	Yellow, Green cap	1 x 3ml
Parainfluenza virus 1, 2, 3 IgA Negative Control	Yellow, Blue cap	1 x 2ml
Parainfluenza virus 1, 2, 3 IgA Positive Control	Yellow, Red cap	1 x 2ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

图片



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.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