

Human Anti-Chlamydia pneumoniae IgA ELISA Kit ab108719

1 图像

概述

产品名称	人Anti-Chlamydia pneumoniae IgA ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Pos. Serum	20			4.9%
	Pos. Serum	24			8.3%
	批次间				
	样品	n	Mean	SD	CV%
	Pos. Serum	12			9.7%
	Pos. Serum	12			14%
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-Chlamydia pneumoniae IgA Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA class antibodies against Chlamydia pneumoniae in Human serum and plasma.				

A 96-well plate has been precoated with Chlamydia pneumoniae 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 Chlamydia pneumoniae-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 Chlamydia pneumoniae IgA sample captured in plate.

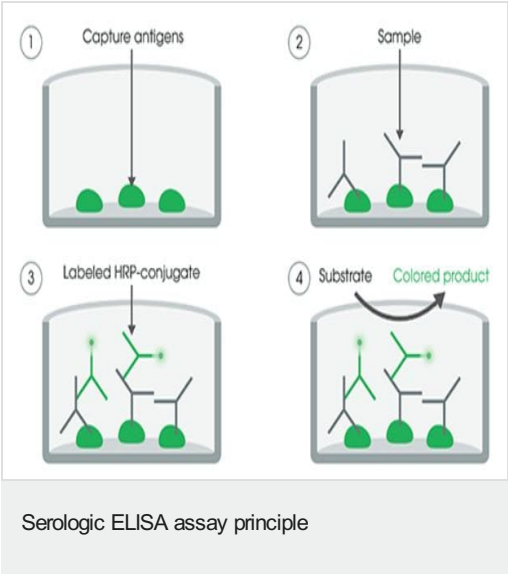
平台	Microplate
----	------------

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
Chlamydia pneumoniae (IgA) Coated Microplate (12 x 8 wells)		1 unit
Chlamydia pneumoniae anti-IgA HRP Conjugate	Colored violet, black cap	1 x 20ml
Chlamydia pneumoniae IgA Cut Off Control	Colored yellow, green cap	1 x 3ml
Chlamydia pneumoniae IgA Negative Control	Colored yellow, blue cap	1 x 2ml
Chlamydia pneumoniae IgA Positive Control	Colored yellow, red cap.	1 x 2ml
IgA Sample Diluent	Colored yellow; white cap	1 x 100ml
Stop Solution	red cap	1 x 15ml
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