

Human Anti-Entamoeba histolytica IgG ELISA Kit ab108734

1 图像

概述

产品名称	人Anti-Entamoeba histolytica IgG ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Pos.Serum	7		2.1%
	批次间			
	样品	n	Mean	SD
	Pos.Serum	6		3.9%
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-Entamoeba histolytica IgG Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Entamoeba histolytica in Human serum and plasma.			

A 96-well plate has been precoated with Entamoeba histolytica antigens to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a Protein A peroxidase labelled anti-Human IgG conjugate is added to the wells, which binds to the immobilized Entamoeba histolytica-specific antibodies. TMB is then catalyzed by the Protein A 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 Entamoeba histolytica IgG sample captured in plate.

平台	Microplate
----	------------

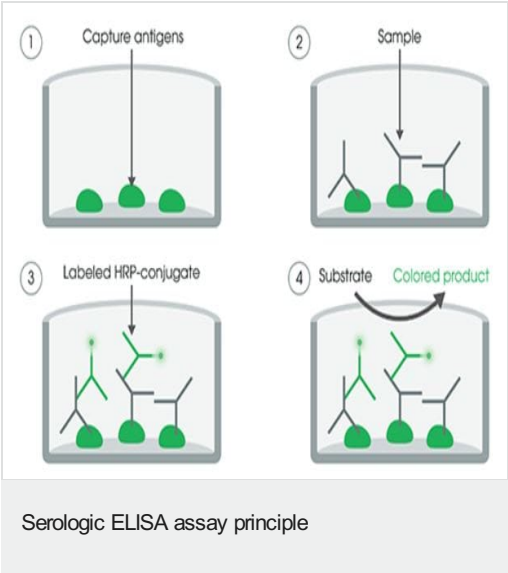
性能

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
Entamoeba histolytica (IgG) Cut-off Control	colored yellow; green cap	1 x 3ml
Entamoeba histolytica (IgG) Negative Control	colored yellow; blue cap	1 x 2ml
Entamoeba histolytica (IgG) Positive Control	colored yellow; red cap	1 x 2ml
Entamoeba histolytica Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Entamoeba histolytica Protein A Conjugate	colored blue; black cap	1 x 20ml
IgG Sample Diluent	colored yellow; white cap	1 x 100ml
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