

Human Anti-Helicobacter pylori IgG ELISA Kit ab108736

5 References 1 图像

概述

产品名称 人Anti-Helicobacter pylori IgG ELISA试剂盒

检测方法 Colorimetric

精确度	批次内				
	样品	n	Mean	SD	CV%
	Pos.Serum	19			6.1%
	Pos.Serum	19			7%

	批次间				
	样品	n	Mean	SD	CV%
	Pos.Serum	13			7.2%
	Pos.Serum	13			3%

样品类型 Serum, Hep Plasma, Cit plasma

检测类型 Indirect

灵敏度 3 U/ml

范围 3 U/ml - 150 U/ml

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's anti-Helicobacter pylori IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of IgG class antibodies against Helicobacter pylori in Human serum and plasma.

A 96-well plate has been precoated with Helicobacter pylori 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 Helicobacter pylori-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 Helicobacter pylori IgG sample

captured in plate.

平台

Microplate

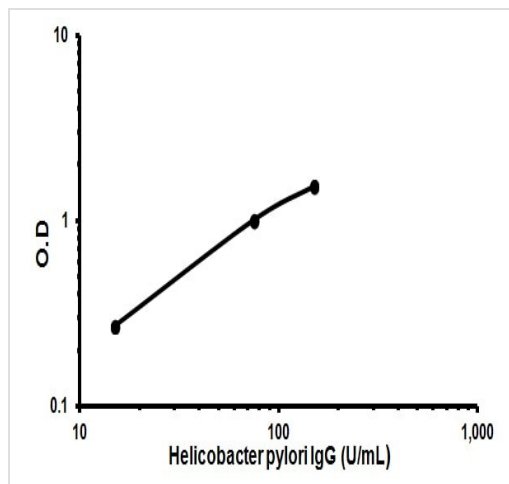
性能

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
Helicobacter pylori (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Helicobacter pylori anti-IgG HRP conjugate	Blue, Black cap	1 x 20ml
Helicobacter pylori IgG Standard A – 0 U/mL	Blue Cap	1 x 2ml
Helicobacter pylori IgG Standard B – 15 U/mL	Green Cap	1 x 2ml
Helicobacter pylori IgG Standard C – 75 U/mL	Yellow cap	1 x 2ml
Helicobacter pylori IgG Standard D – 150 U/mL	Red cap	1 x 2ml
IgG Sample Diluent	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



Representative Standard Curve using ab108736

Typical Standard Curve

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