

Human Anti-Helicobacter pylori IgG ELISA Kit ab108736

5 References 画像数 1

製品の概要

製品名	Human Anti-Helicobacter pylori IgG ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Pos.Serum	19		6.1%
	Pos.Serum	19		7%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Pos.Serum	13		7.2%
	Pos.Serum	13		3%
サンプルの種類	Serum, Hep Plasma, Cit plasma			
アッセイタイプ	Indirect			
検出感度	3 U/ml			
検出範囲	3 U/ml - 150 U/ml			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's anti-Helicobacter pylori IgG Human <i>in vitro</i> 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.

試験プラットフォーム

Microplate

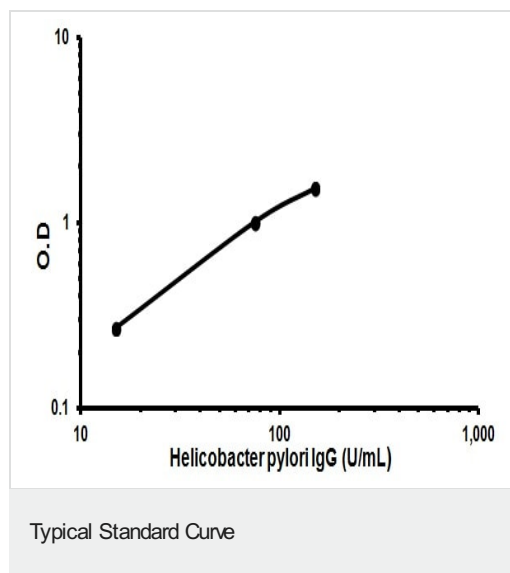
製品の特長

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	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

画像



Representative Standard Curve using ab108736

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors