

Human Anti-Leishmania IgG ELISA Kit ab108749

画像数 1

製品の概要

製品名	Human Anti-Leishmania 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	CV%
Pos.Serum	7			7.4%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Pos.Serum	4			5.9%

サンプルの種類	Serum, Hep Plasma, Cit plasma
アッセイタイプ	Indirect
ステップ	Multiple steps standard assay
種交差性	交差種: Human

製品の概要
Abcam's anti-Leishmania IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Leishmania in Human serum and plasma.

A 96-well plate has been precoated with Leishmania 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 Leishmania-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 Leishmania IgG sample captured in plate.

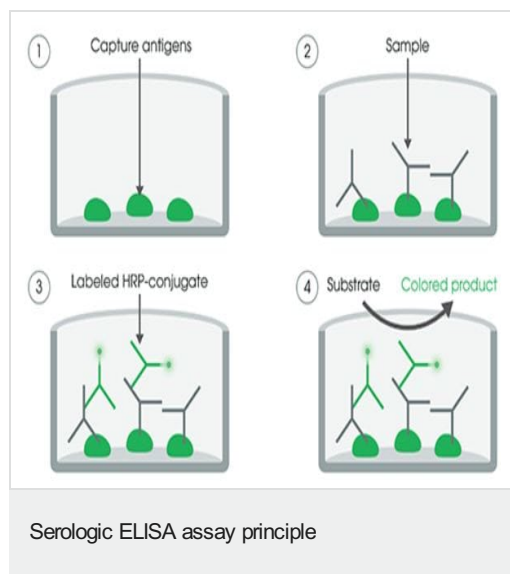
試験プラットフォーム	Microplate
------------	------------

製品の特性

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
IgG Sample Diluent	Yellow, White cap	1 x 100ml
Leishmania (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8	1 unit
Leishmania IgG Cut-off Control	Yellow, Green cap	1 x 3ml
Leishmania IgG Negative Control	Yellow, Blue cap	1 x 2ml
Leishmania IgG Positive Control	Yellow, Red cap	1 x 2ml
Leishmania Protein A Conjugate	Blue, Black cap	1 x 20ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

画像



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.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