

Product datasheet

Human Anti-Candida albicans IgM ELISA Kit ab108716

製品の概要

製品名 Human Anti-Candida albicans IgM 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	6			4.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Pos.Serum	38			5.1%

サンプルの種類 Serum, Heparin Plasma, Citrate 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-Candida albicans IgM Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Candida albicans in Human serum or plasma.

A 96-well plate has been precoated with Candida albicans 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 IgM conjugate is added to the wells, which binds to the immobilized Candida albicans-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 Candida albicans IgM sample captured in plate.

アプリケーション 適用あり: Indirect ELISA

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
Candida albicans (IgM) Coated Microplate (12 x 8 wells)		1 unit
Candida albicans anti-IgM HRP Conjugate	Red Bottle, Black Cap	1 x 20ml
Candida albicans IgM Cut-off Control	Yellow Bottle, Green Cap	1 x 3ml
Candida albicans IgM Negative Control	Yellow Bottle, Blue Cap	1 x 2ml
Candida albicans IgM Positive Control	Yellow Bottle, Red Cap	1 x 2ml
Foil Cover		1 unit
IgM Sample Diluent	Green Bottle, White Cap	1 x 100ml
Stop Solution		1 x 15ml
Strip Holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108716** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Indirect ELISA		Use at an assay dependent concentration.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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