

Product datasheet

Human Anti-Candida albicans IgA ELISA Kit ab108714

概述

产品名称 人Anti-Candida albicans IgA ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Pos.Serum	10			6.6%

批次间

样品	n	Mean	SD	CV%
Pos.Serum	33			9%

样品类型 Serum, 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-Candida albicans IgA Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA class antibodies against Candida albicans in Human serum and 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 IgA 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 IgA sample captured in plate.

经测试应用 适用于: Indirect ELISA

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
Candida albicans (IgA) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Candida albicans anti-IgA HRP Conjugate	Colored violet; Black cap	1 x 20ml
Candida albicans IgA Cut-off Control	Colored yellow; Green cap	1 x 3ml
Candida albicans IgA Negative Control	Colored yellow; Blue cap	1 x 2ml
Candida albicans IgA Positive Control	Colored yellow; Red cap	1 x 2ml
Cover foil		1 unit
IgA Sample Diluent	Colored yellow; White cap	1 x 100ml
Stop Solution		1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

应用

Our [Abpromise guarantee](#) covers the use of **ab108714** in the following tested applications.

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

应用	Ab评论	说明
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 <http://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