

Product datasheet

Human Anti-Respiratory syncytial virus IgG ELISA Kit (RSV) ab108765

1 References

製品の概要

製品名 Human Anti-Respiratory syncytial virus IgG ELISA Kit (RSV)

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	24			8.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Pos.Serum	12			6%

サンプルの種類 Serum, 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-Respiratory syncytial virus (RSV) IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Respiratory syncytial virus in Human serum and plasma.

A 96-well plate has been precoated with Respiratory syncytial virus 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 Respiratory syncytial virus -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 Respiratory syncytial virus IgG 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
Anti-Respiratory syncytial virus (IgG) Coated Microplate (12 x 8 wells)		1 unit
Cover foil		1 unit
IgG Sample Diluent	yellow buffer, white cap	1 x 100ml
Respiratory syncytial virus anti-IgG HRP Conjugate	blue solution, black cap	1 x 20ml
Respiratory syncytial virus IgG Cut-off Control	yellow solution, green cap	1 x 3ml
Respiratory syncytial virus IgG Negative Control	yellow solution, blue cap	1 x 2ml
Respiratory syncytial virus IgG Positive Control	yellow solution, red cap	1 x 2ml
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 **ab108765** 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 <http://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