

## Product datasheet

# Human Anti-Chlamydia trachomatis IgG ELISA Kit ab108720

### Overview

**Product name** Human Anti-Chlamydia trachomatis IgG ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay				
Sample	n	Mean	SD	CV%
Pos. Serum	24			3.7%

Inter-assay				
Sample	n	Mean	SD	CV%
Pos. Serum	12			2.8%

**Sample type** Serum, Heparin Plasma, Citrate Plasma

**Assay type** Indirect

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's anti-Chlamydia trachomatis IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Chlamydia trachomatis in Human serum and plasma.

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

**Tested applications** **Suitable for:** Indirect ELISA

**Platform** Microplate

### Properties

## Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Chlamydia trachomatis (IgG) Coated Microplate (12 x 8 wells)		1 unit
Chlamydia trachomatis anti-IgG HRP Conjugate	Black cap	1 x 20ml
Chlamydia trachomatis IgG Cut-off Control	Green cap	1 x 3ml
Chlamydia trachomatis IgG Negative Control	Blue cap	1 x 2ml
Chlamydia trachomatis IgG Positive Control	Red cap	1 x 2ml
Cover foil		1 unit
IgG Sample Diluent	White cap	1 x 100ml
Stop Solution		1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

## Applications

Our [Abpromise guarantee](#) covers the use of **ab108720** in the following tested applications.

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

Application	Abreviews	Notes
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.com/abpromise> or contact our technical team.

## Terms and conditions

---

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