

Coxiella burnetii Phase 1 IgG ELISA Kit (Q-Fever)
ab178638

1 Image

Overview

Product name	Coxiella burnetii Phase 1 IgG ELISA Kit (Q-Fever)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Pos Serum	15		5.7%
	Pos serum	15		4.7%
	Inter-assay			
	Sample	n	Mean	SD
	Equ serum	5		2.4%
	Pos serum	6		3.5%
Sample type	Serum, Hep Plasma, Cit plasma			
Assay type	Indirect			
Sensitivity	> 90 %			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's anti-Coxiella burnetii (Q-Fever) Phase 1 IgG Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Coxiella burnetii in Human serum.			

A 96-well plate has been precoated with Coxiella burnetii 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 Coxiella burnetii antigens. 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 Coxiella burnetii IgG sample captured in plate.

Platform

Microplate

Properties**Storage instructions**

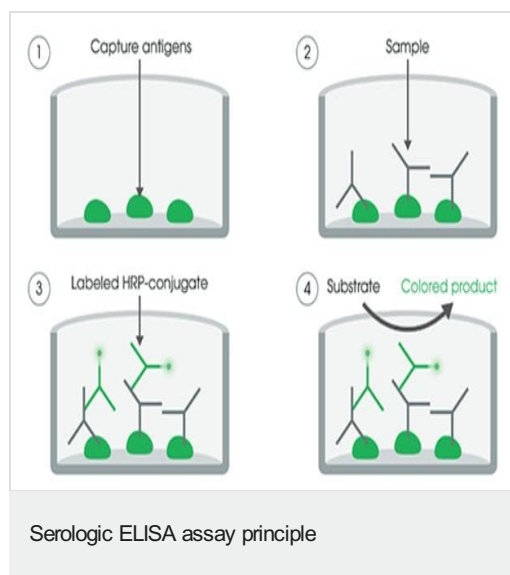
Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover Foil		1 unit
Coxiella burnetii Anti-IgG HRP Conjugate		1 x 20ml
Coxiella burnetii IgG Cut-off Control		1 x 3ml
Coxiella burnetii IgG Negative Control		1 x 2ml
Coxiella burnetii IgG Positive Control		1 x 2ml
Coxiella burnetii Phase 1 (IgG) Coated Microplate (12 x 8 wells)		1 unit
IgG Sample Diluent		1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

Relevance

Coxiella burnetii is an obligate intracellular bacterial pathogen, and is the causative agent of Q fever. The genus Coxiella is morphologically similar to Rickettsia, but with a variety of genetic and physiological differences. C. burnetii is a small Gram-negative bacterium that is highly resistant to environmental stresses such as high temperature, osmotic pressure, and ultraviolet light. These characteristics are attributed to a small cell variant (SCV) form of the organism that is part of a biphasic developmental cycle, including a more metabolically and replicatively active large cell variant (LCV) form. It can survive standard disinfectants, and is resistant to many other environmental changes like those presented in the phagolysosome.

Images



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.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