

Human Anti-Clostridium tetani toxin IgG ELISA Kit
ab108771

1 Image

Overview

Product name Human Anti-Clostridium tetani toxin IgG ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|--------------|----|----------|----|------|
| Weak + Serum | 24 | 0.56U/ml | | 7.1% |
| Pos. Serum | 24 | 1.24U/ml | | 4.9% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|--------------|----|----------|----|-------|
| Weak + Serum | 10 | 0.22U/ml | | 11.3% |
| Pos. Serum | 10 | 0.54U/ml | | 7.7% |

Sample type Serum, Hep Plasma, Cit plasma

Assay type Indirect

Sensitivity < 0.05 U/ml

Range 0.05 U/ml - 1 U/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

A 96-well plate has been precoated with Clostridium tetani toxin 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 Clostridium tetani toxin-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 Clostridium tetani toxin IgG sample captured in plate.

Platform

Microplate

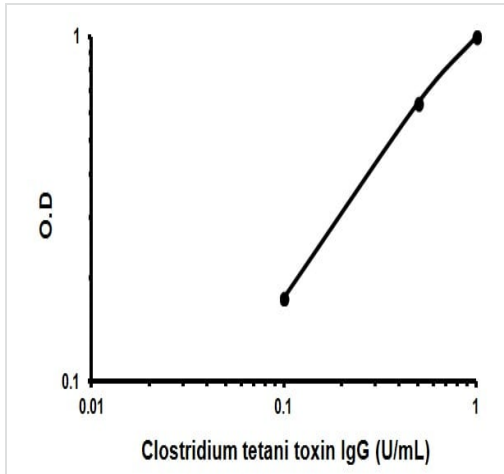
Properties

Storage instructions

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

| Components | Identifier | 1 x 96 tests |
|---|----------------------|--------------|
| 20X Washing Solution | White cap | 1 x 50ml |
| Clostridium tetani toxin (IgG) Coated Microplate (12 x 8 wells) | 12 strips of 8 wells | 1 unit |
| Clostridium tetani toxin anti-IgG HRP Conjugate | Blue, Black cap | 1 x 20ml |
| Clostridium tetani toxin IgG Standard A – 0.0 U/mL | Blue cap | 1 x 2ml |
| Clostridium tetani toxin IgG Standard B – 0.1 U/mL | Green Cap | 1 x 2ml |
| Clostridium tetani toxin IgG Standard C – 0.5 U/mL | Yellow cap | 1 x 2ml |
| Clostridium tetani toxin IgG Standard D – 1.0 U/mL | Red cap | 1 x 2ml |
| Cover foil | | 1 unit |
| IgG Sample Diluent | Yellow, White cap | 1 x 100ml |
| Stop Solution | red cap | 1 x 15ml |
| Strip holder | | 1 unit |
| TMB Substrate Solution | Yellow cap | 1 x 15ml |

Images



Representative Standard Curve using ab108771

Typical Standard Curve

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