

Product datasheet

Human Anti-Corynebacterium diphtheriae toxin IgG ELISA Kit ab108726

[1 Image](#)

Overview

Product name Human Anti-Corynebacterium diphtheriae toxin IgG ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Weak Pos.	24			12.1%
Positive	24			2.1%

Inter-assay

Sample	n	Mean	SD	CV%
Weak Pos	12	0.04U/ml		11.4%
Positive	12	0.16U/ml		2.8%

Sample type Serum, Heparin Plasma, Citrate Plasma

Assay type Indirect

Sensitivity = 0.01 U/ml

Range 0.01 IU/ml - 0.15 IU/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

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

A 96-well plate has been precoated with Corynebacterium diphtheriae 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 Corynebacterium diphtheriae 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 Corynebacterium diphtheriae toxin 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
C. diphtheriae toxin IgG Standard A - 0 U/mL		1 x 2ml
C. diphtheriae toxin IgG Standard B - 0.015 U/mL		1 x 2ml
C. diphtheriae toxin IgG Standard C - 0.075 U/mL		1 x 2ml
C. diphtheriae toxin IgG Standard D - 0.150 U/mL		1 x 2ml
Corynebacterium diphtheriae toxin (IgG) Coated Microplate (12 x 8 wells)		1 unit
Corynebacterium diphtheriae toxin anti-IgG HRP Conjugate	Colored blue, black cap	1 x 20ml
Cover foil		1 unit
IgG 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

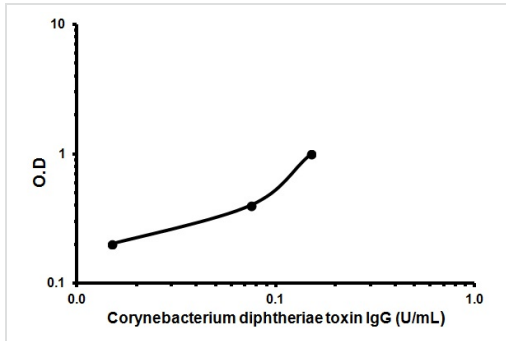
Applications

Our [Abpromise guarantee](#) covers the use of **ab108726** 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.

Images



Representative Standard Curve Using
ab108726

Typical Standard Curve

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