

Human Anti-Corynebacterium diphtheriae toxin IgG ELISA Kit ab108726

1 图像

概述

产品名称	人Anti-Corynebacterium diphtheriae toxin IgG ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Weak Pos.	24			12.1%
	Positive	24			2.1%
	批次间				
	样品	n	Mean	SD	CV%
	Weak Pos	12	0.04U/ml		11.4%
	Positive	12	0.16U/ml		2.8%
样品类型	Serum, Hep Plasma, Cit plasma				
检测类型	Indirect				
灵敏度	= 0.01 U/ml				
范围	0.01 IU/ml - 0.15 IU/ml				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's anti-Corynebacterium diphtheriae toxin IgG Human <i>in vitro</i> 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.

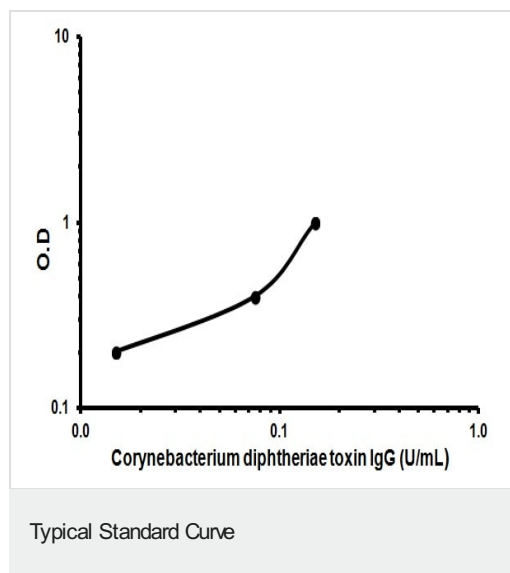
平台 Microplate

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	标识符	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	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

图片



Representative Standard Curve Using ab108726

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.cn/abpromise> or contact our technical team.

Terms and conditions

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