

Human Anti-Corynebacterium diphtheriae toxin IgG ELISA Kit ab108727

画像数 1

製品の概要

製品名	Human Anti-Corynebacterium diphtheriae toxin IgG ELISA Kit
検出方法	Colorimetric
サンプルの種類	Serum, Hep Plasma, Cit plasma
アッセイタイプ	Indirect
検出感度	= 0.01 IU/ml
検出範囲	0.01 IU/ml - 1 IU/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	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 and 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 5S-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.

試験プラットフォーム Microplate

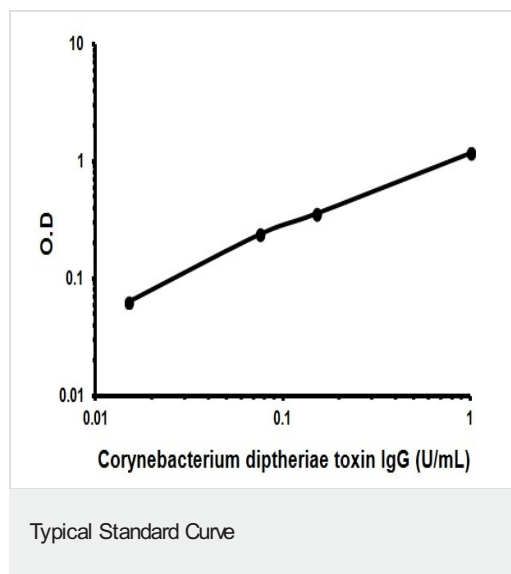
製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml

内容	ラベル	1 x 96 tests
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
Corynebacterium diphtheriae toxin IgG Standard A – 0.000 U/mL		1 x 2ml
Corynebacterium diphtheriae toxin IgG Standard B – 0.015 U/mL		1 x 2ml
Corynebacterium diphtheriae toxin IgG Standard C – 0.075 U/mL		1 x 2ml
Corynebacterium diphtheriae toxin IgG Standard D – 0.150 U/mL		1 x 2ml
Corynebacterium diphtheriae toxin IgG Standard E – 1.000 U/mL		1 x 2ml
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

画像



Representative Standard Curve using ab108727.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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