

CIC-C3d ELISA Kit ab178665

1 Image

Overview

Product name CIC-C3d ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	16			6.1%

Inter-assay

Sample	n	Mean	SD	CV%
Serum				13.9%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (quantitative)

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's CIC-C3d *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of CIC-C3d in Human serum and plasma.

A 96-well plate has been precoated with anti-C3d antibodies. Standards, controls and test samples are added to the wells and incubated. Immune complexes (CIC) in the samples and standards bind to the precoated antibody. Following washing, a horseradish peroxidase (HRP) labelled anti-Human IgG conjugate is added to the wells, which binds to the immobilized CIC-C3d. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of CIC-C3d in the sample and the intensity is measured at 450 nm.

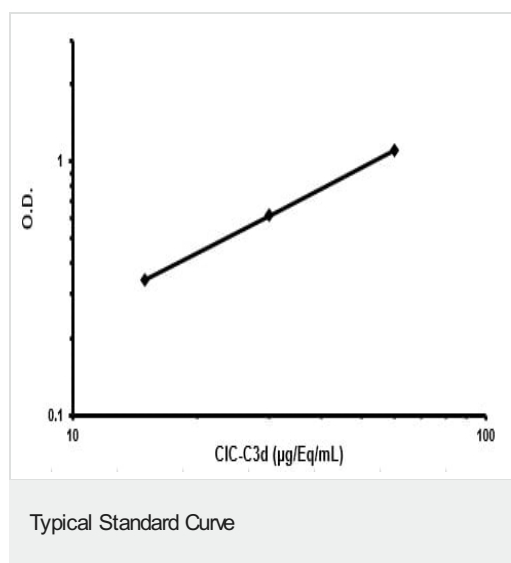
Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
10X Wash Solution	2 x 50ml
Anti-CIC-C3d Coated Microplate (12 x 8 wells)	1 unit
CIC-C3d Anti-IgG HRP Conjugate	1 x 0.5ml
CIC-C3d Negative Control	1 x 1.5ml
CIC-C3d Positive Control	1 x 1.5ml
CIC-C3d Standard 0 - 0 µg Eq/mL	1 x 1.5ml
CIC-C3d Standard 1 - 16 µg Eq/mL	1 x 1.5ml
CIC-C3d Standard 2 - 64 µg Eq/mL	1 x 1.5ml
Conjugate Buffer	1 x 20ml
Cover Foil	1 unit
Incubation Buffer	1 x 50ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Images

Representative standard curve using ab178665.

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