

Human Complement C1r ELISA Kit ab170245

1 Image

Overview

Product name Human Complement C1r ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				5.4%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				9.7%

Sample type Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid

Assay type Sandwich (quantitative)

Sensitivity = 92 pg/ml

Range 0.125 ng/ml - 8 ng/ml

Recovery = 97 %

Assay time 4h 00m

Species reactivity **Reacts with:** Human

Product overview

Abcam's Complement C1r Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Complement C1r concentrations in plasma, serum, saliva, urine, milk and cell culture supernatants.

A Complement C1r specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Complement C1r specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Complement C1r captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Properties

Storage instructions Store at -20°C. Please refer to protocols.

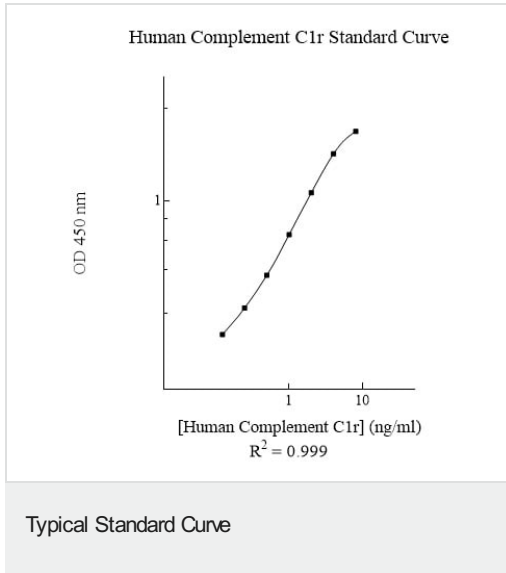
Components	1 x 96 tests
100X Biotinylated Human Complement C1r Antibody	1 x 60µl
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
Chromogen Substrate	1 x 7ml
Complement C1r Microplate (12 x 8 well strips)	1 unit
Complement C1r Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function C1r B chain is a serine protease that combines with C1q and C1s to form C1, the first component of the classical pathway of the complement system.

Sequence similarities Belongs to the peptidase S1 family.
Contains 2 CUB domains.
Contains 1 EGF-like domain.
Contains 1 peptidase S1 domain.
Contains 2 Sushi (CCP/SCR) domains.

Post-translational modifications The iron and 2-oxoglutarate dependent 3-hydroxylation of aspartate and asparagine is (R) stereospecific within EGF domains.

Images



Representative Standard Curve using ab170245

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