

Human Complement C1q ELISA Kit ab170246

15 References 1 Image

Overview

Product name	Human Complement C1q ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Plasma			5.1%
	Inter-assay			
	Sample	n	Mean	SD
	Plasma			9.9%
Sample type	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	> 64 pg/ml			
Range	0.078 ng/ml - 5 ng/ml			
Recovery	= 96 %			
Assay time	4h 00m			
Species reactivity	Reacts with: Human Does not react with: Mouse, Rat, Rabbit, Cow, Pig			
Product overview	Complement C1q Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab170246) is designed for the quantitative measurement of Complement C1q concentrations in plasma, serum, saliva, urine, milk, cerebrospinal fluid and cell culture supernatants.			

A Complement C1q specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Complement C1q 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 C1q captured in plate.

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

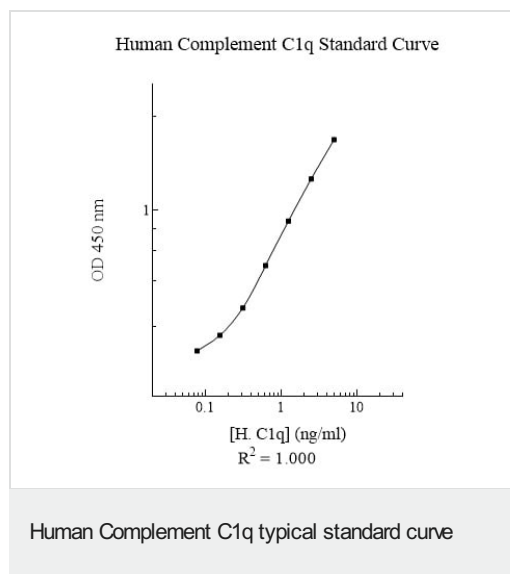
Properties

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Complement C1q Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Complement C1q Microplate (12 x 8 well strips)	1 unit
Complement C1q Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function	C1q associates with the proenzymes C1r and C1s to yield C1, the first component of the serum complement system. The collagen-like regions of C1q interact with the Ca(2+)-dependent C1r(2)C1s(2) proenzyme complex, and efficient activation of C1 takes place on interaction of the globular heads of C1q with the Fc regions of IgG or IgM antibody present in immune complexes.
Involvement in disease	Defects in C1QA are a cause of complement component C1q deficiency (C1QD) [MIM:613652]. A rare defect resulting in C1 deficiency and impaired activation of the complement classical pathway. C1 deficiency generally leads to severe immune complex disease with features of systemic lupus erythematosus and glomerulonephritis.
Sequence similarities	Contains 1 C1q domain. Contains 1 collagen-like domain.
Post-translational modifications	O-linked glycans consist of Glc-Gal disaccharides bound to the oxygen atom of post-translationally added hydroxyl groups.
Cellular localization	Secreted.

Images



Representative Standard Curve using ab170246

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