

Human Complement C2 ELISA Kit ab154132

★★★★★ 1 Abreviews 3 References 1 Image

Overview

Product name	Human Complement C2 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			CV%
				5%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			CV%
				10.8%
Sample type	Milk, Serum, Plasma, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	= 0.0032 µg/ml			
Range	0.006 µg/ml - 0.4 µg/ml			
Recovery	97 %			
Assay time	4h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human Does not react with: Mouse, Rat, Rabbit, Dog, Pig			
Product overview	Abcam's Complement C2 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Complement C2 in plasma, serum, milk, saliva, CSF, cell culture supernatants, cell lysate, and tissue samples.			

A Complement C2 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Complement C2 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 C2 captured in plate.

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

**Platform** Microplate

## 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 N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
70X Biotinylated Human Complement C2 Antibody	1 x 90µl
Chromogen Substrate	1 x 7ml
Complement C2 Microplate (12 x 8 well strips)	1 unit
Complement C2 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

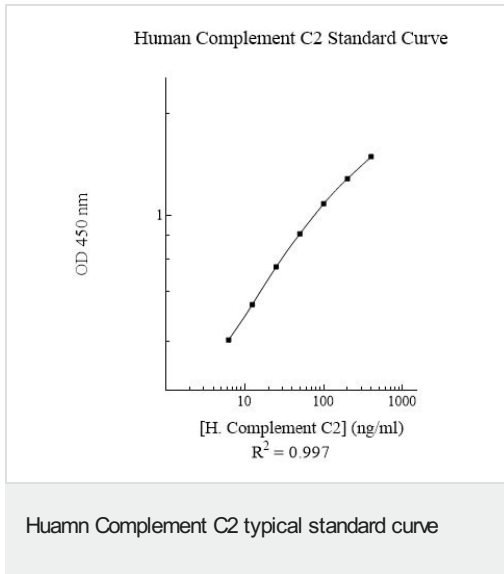
**Function** Component C2 which is part of the classical pathway of the complement system is cleaved by activated factor C1 into two fragments: C2b and C2a. C2a, a serine protease, then combines with complement factor 4b to generate the C3 or C5 convertase.

**Involvement in disease** Defects in C2 are the cause of complement component 2 deficiency (C2D) [MIM:217000]. A deficiency of the complement classical pathway associated with the development of autoimmune disorders, mainly systemic lupus erythematosus. Skin and joint manifestations are common and renal disease is relatively rare. Patients with complement component 2 deficiency are also reported to have recurrent or invasive infections.

**Sequence similarities** Belongs to the peptidase S1 family.  
Contains 1 peptidase S1 domain.  
Contains 3 Sushi (CCP/SCR) domains.  
Contains 1 VWFA domain.

**Cellular localization** Secreted.

## Images



Representative Standard Curve using ab154132

**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