

Human Complement C9 ELISA Kit ab137972

3 References 1 Image

Overview

Product name	Human Complement C9 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			3.2%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			8.4%
Sample type	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	= 0.059 ng/ml			
Range	0.234 ng/ml - 15 ng/ml			
Recovery	= 101 %			
Assay time	4h 0m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Complement C9 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of C9 in Human plasma, serum, saliva, milk, urine, cerebrospinal fluid and cell culture supernatants.			

A Complement C9 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Complement C9 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 C9 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 M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Complement C9 Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Complement C9 Microplate (12 x 8 well strips)	1 unit
Complement C9 Standard	2 vials
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Function** Constituent of the membrane attack complex (MAC) that plays a key role in the innate and adaptive immune response by forming pores in the plasma membrane of target cells. C9 is the pore-forming subunit of the MAC.

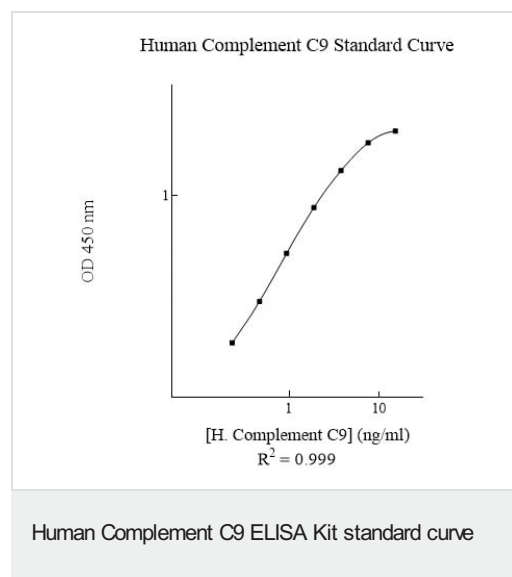
**Tissue specificity** Plasma.

**Involvement in disease** Defects in C9 are a cause of complement component 9 deficiency (C9D) [MIM:613825]. A rare defect of the complement classical pathway associated with susceptibility to severe recurrent infections, predominantly by *Neisseria gonorrhoeae* or *Neisseria meningitidis*.

**Sequence similarities** Belongs to the complement C6/C7/C8/C9 family.  
Contains 1 EGF-like domain.  
Contains 1 LDL-receptor class A domain.  
Contains 1 MACPF domain.  
Contains 1 TSP type-1 domain.

**Post-translational modifications** Thrombin cleaves factor C9 to produce C9a and C9b.  
Phosphorylation sites are present in the extracellular medium.

**Cellular localization** Secreted. Cell membrane. Secreted as soluble monomer. Oligomerizes at target membranes, forming a pre-pore. A conformation change then leads to the formation of a 100 Angstrom diameter pore.



Representative Standard Curve using ab137972

**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