

Human Complement C8 ELISA Kit ab137971

1 Image

Overview

Product name	Human Complement C8 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			4.5%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			8.6%
Sample type	Cell culture supernatant, Saliva, Milk, Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	= 0.3 ng/ml			
Range	0.625 ng/ml - 40 ng/ml			
Recovery	= 103 %			
Assay time	4h 0m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Complement C8 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Complement C8 in Human plasma, serum, saliva, milk, CSF, cell culture supernatants, cell lysates and tissue samples.			

A Complement C8 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Complement C8 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 C8 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

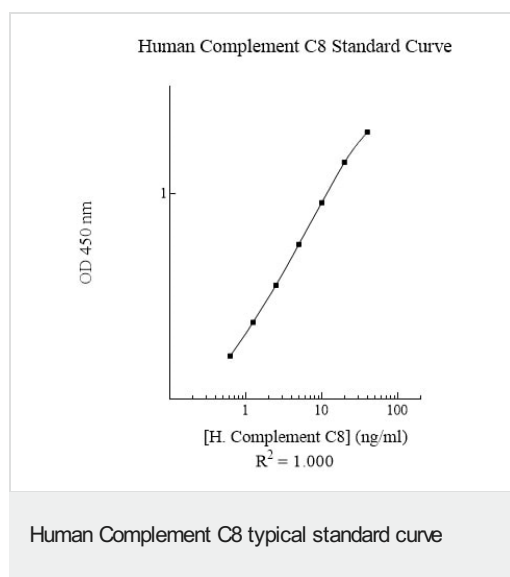
Properties

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

Components	1 x 96 tests
50X Biotinylated Human Complement C8 Antibody	1 vial
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 C8 Microplate (12 x 8 well strips)	1 unit
Complement C8 Standard	1 vial
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. C8A inserts into the target membrane, but does not form pores by itself.
Involvement in disease	Defects in C8A are a cause of complement component 8 deficiency type 1 (C8D1) [MIM:613790]. A rare defect of the complement classical pathway associated with susceptibility to severe recurrent infections, predominantly by <i>Neisseria gonorrhoeae</i> or <i>Neisseria meningitidis</i> .
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 2 TSP type-1 domains.
Cellular localization	Secreted. Cell membrane. Secreted as soluble protein. Inserts into the cell membrane of target cells.

Images



Representative Standard Curve using ab137971

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