

Human Complement C9 ELISA Kit ab137972

3 References [画像数 1](#)

製品の概要

製品名 Human Complement C9 ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall				3.2%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				8.4%

サンプルの種類 Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid

アッセイタイプ Sandwich (quantitative)

検出感度 = 0.059 ng/ml

検出範囲 0.234 ng/ml - 15 ng/ml

添加回収試験 = 101 %

全工程の試験時間 4h 0m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's Complement C9 Human *in vitro* 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.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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

機能

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.

組織特異性

Plasma.

関連疾患

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*.

配列類似性

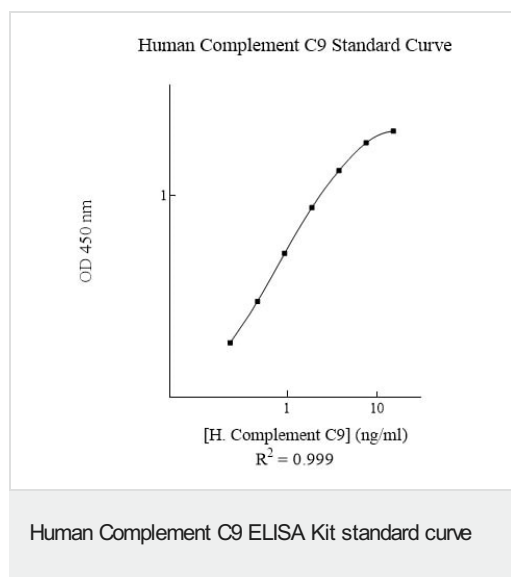
Belongs to the complement C6/C7/C8/C9 family.
Contains 1 EGF-like domain.
Contains 1 LDL-receptor class A domain.
Contains 1 MACPF domain.
Contains 1 TSP type-1 domain.

翻訳後修飾

Thrombin cleaves factor C9 to produce C9a and C9b.
Phosphorylation sites are present in the extracellular medium.

細胞内局在

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors