

Human Complement C1q ELISA Kit ab170246

[15 References](#) [画像数 1](#)

製品の概要

製品名 Human Complement C1q 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%
Plasma				5.1%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Plasma				9.9%

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

アッセイタイプ Sandwich (quantitative)

検出感度 > 64 pg/ml

検出範囲 0.078 ng/ml - 5 ng/ml

添加回収試験 = 96 %

全工程の試験時間 4h 00m

種交差性
交差種: Human
非交差種: Mouse, Rat, Rabbit, Cow, Pig

製品の概要 Complement C1q Human *in vitro* 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.

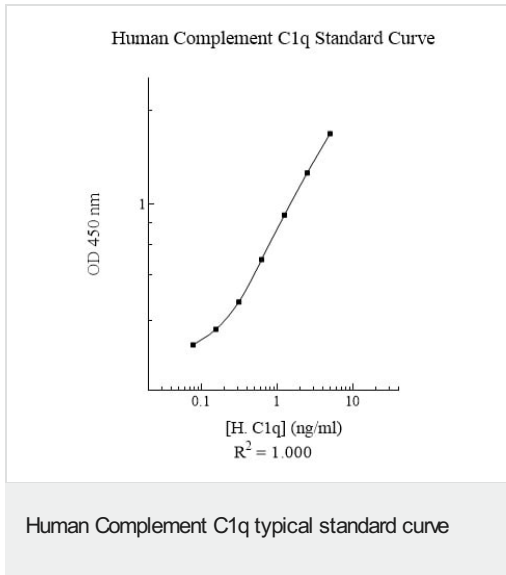
製品の特性

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 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

機能	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.
関連疾患	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.
配列類似性	Contains 1 C1q domain. Contains 1 collagen-like domain.
翻訳後修飾	O-linked glycans consist of Glc-Gal disaccharides bound to the oxygen atom of post-translationally added hydroxyl groups.
細胞内局在	Secreted.

画像



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.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