

Human Complement C4 ELISA Kit ab108824

8 References 画像数 2

製品の概要

| | | | | |
|----------|---|---|-----|-------|
| 製品名 | Human Complement C4 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 |
| | Overall | | | CV% |
| | | | | 5.1% |
| | Inter-Assay (日差再現性) | | | |
| | サンプル | N | 平均値 | SD |
| | Overall | | | CV% |
| | | | | 10.1% |
| サンプルの種類 | Serum, Plasma | | | |
| アッセイタイプ | Competitive | | | |
| 検出感度 | = 0.53 µg/ml | | | |
| 検出範囲 | 160 µg/ml - 480 µg/ml | | | |
| 添加回収試験 | 94 % | | | |
| 全工程の試験時間 | 3h 00m | | | |
| ステップ | Multiple steps standard assay | | | |
| 種交差性 | 交差種: Human | | | |
| 製品の概要 | Complement C4 Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab108834) is designed for the quantitative measurement of Complement C4 levels in plasma and serum. | | | |

A Complement C4 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Complement C4 is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 inversely proportional to the amount of Complement C4 captured in plate.

Reactivity with C4 cleavage products (C4a, C4b, C4d) has not been determined.

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

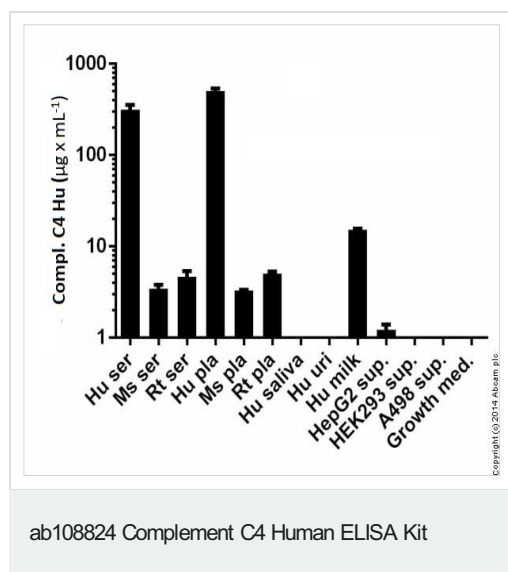
製品の特徴

保存方法

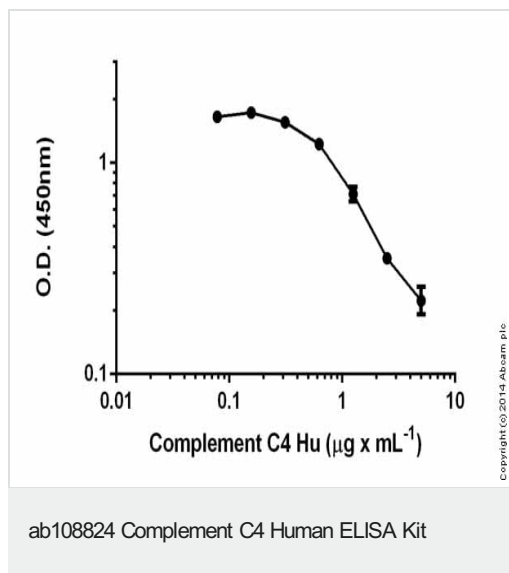
Store at -20°C. Please refer to protocols.

| 内容 | 1 x 96 tests |
|---|--------------|
| 100X Streptavidin-Peroxidase Conjugate | 1 x 80µl |
| 10X Diluent N Concentrate | 1 x 30ml |
| 20X Wash Buffer Concentrate | 1 x 30ml |
| 2X Biotinylated Human Complement C4 (Lyophilized) | 1 vial |
| Chromogen Substrate | 1 x 7ml |
| Complement C4 Microplate (12 x 8 well strips) | 1 unit |
| Complement C4 Standard | 1 vial |
| Sealing Tapes | 3 units |
| Stop Solution | 1 x 11ml |

画像



Complement C4 measured in biological fluids showing quantity (microgram) per mL of tested sample



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