

## Product datasheet

# Human Factor D ELISA Kit ab99969

### 画像数 2

#### 製品の概要

|         |   |
|---------|---|
| 製品名     | Human Factor D ELISA Kit                |
| 検出方法    | Colorimetric                            |
| サンプルの種類 | Cell culture supernatant, Serum, Plasma |
| アッセイタイプ | Sandwich (quantitative)                 |
| 検出感度    | < 4 pg/ml                               |
| 添加回収試験  | < 130 %                                 |

特定サンプルでの回収試験

| サンプルの種類                  | 平均 %  | 測定範囲        |
|--------------------------|-------|-------------|
| Cell culture supernatant | 126.4 | 115% - 133% |
| Serum                    | 130.3 | 118% - 139% |
| Plasma                   | 131.1 | 107% - 140% |

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Human

**製品の概要** Abcam's Human Factor D ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Factor D in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Factor D coated on a 96-well plate. Standards and samples are pipetted into the wells and Factor D present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Factor D antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Factor D bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**特記事項** Optimisation may be required with urine samples.

**アプリケーション** 適用あり: Sandwich ELISA

**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 |
|--|--------------|
| 20X Wash Buffer Concentrate                        | 1 x 25ml     |
| 300X HRP-Streptavidin Concentrate                  | 1 x 200µl    |
| 5X Assay Diluent B                                 | 1 x 15ml     |
| Assay Diluent A                                    | 1 x 30ml     |
| Biotinylated anti-human Adipsin                    | 2 vials      |
| Factor D Microplate (12 strips x 8 wells)          | 1 unit       |
| Recombinant human Factor D Standards (lyophilized) | 2 vials      |
| Stop Solution                                      | 1 x 8ml      |
| TMB One-Step Substrate Reagent                     | 1 x 12ml     |

**機能** Factor D cleaves factor B when the latter is complexed with factor C3b, activating the C3bbb complex, which then becomes the C3 convertase of the alternate pathway. Its function is homologous to that of C1s in the classical pathway.

**関連疾患** Complement factor D deficiency (CFDD) [MIM:613912]: An immunologic disorder characterized by increased susceptibility to bacterial infections, particularly Neisseria infections, due to a defect in the alternative complement pathway. Note=The disease is caused by mutations affecting the gene represented in this entry.

**配列類似性** Belongs to the peptidase S1 family.  
Contains 1 peptidase S1 domain.

**細胞内局在** Secreted.

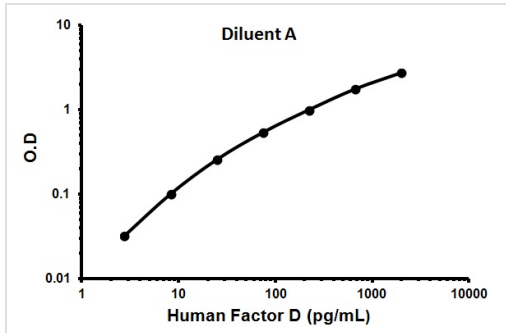
## アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab99969** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

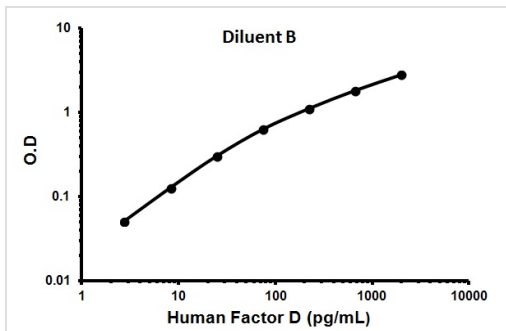
| アプリケーション       | Abreviews | 特記事項                                     |
|----------------|-----------|--|
| Sandwich ELISA |           | Use at an assay dependent concentration. |

## 画像



Representative Standard Curve using ab99969.

Typical Standard Curve



Representative Standard Curve using ab99969.

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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