

# 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	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.

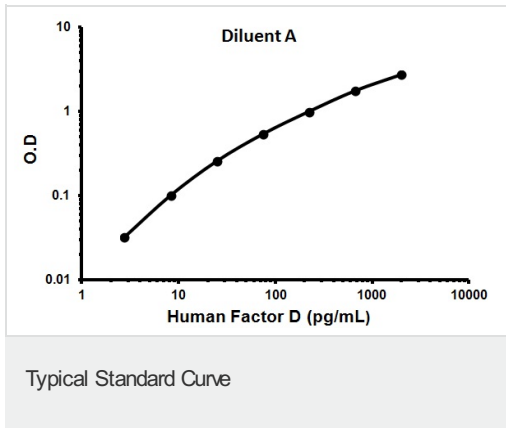
## アプリケーション

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

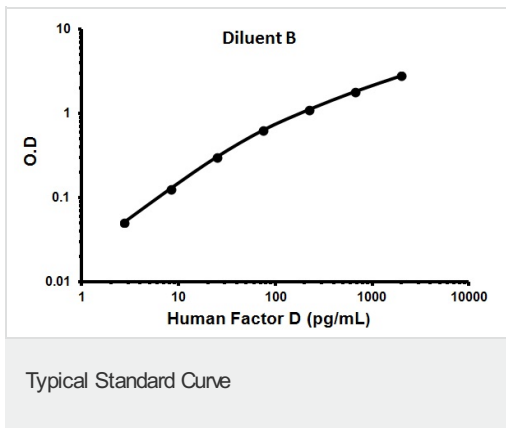
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

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.



Representative Standard Curve using ab99969.

**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 <http://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