

Human Factor D ELISA Kit ab99969

[3 References](#) [2 Images](#)

Overview

Product name	Human Factor D ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 4 pg/ml
Recovery	< 130 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	126.4	115% - 133%
Serum	130.3	118% - 139%
Plasma	131.1	107% - 140%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview 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.

Notes Optimisation may be required with urine samples.

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	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

Function

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.

Involvement in disease

Defects in CFD are the cause of complement factor D deficiency (CFD deficiency) [MIM:134350]. CFD deficiency predisposes to invasive meningococcal disease.

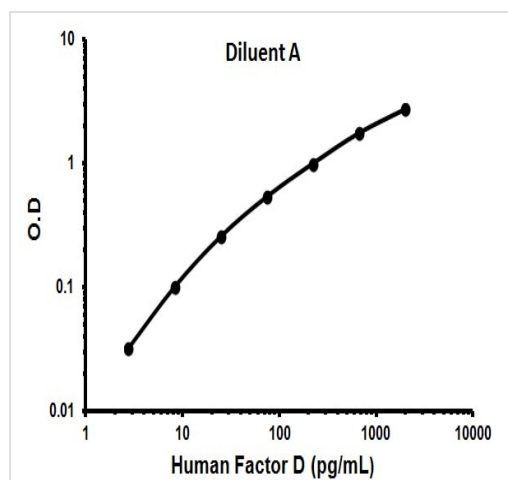
Sequence similarities

Belongs to the peptidase S1 family.
Contains 1 peptidase S1 domain.

Cellular localization

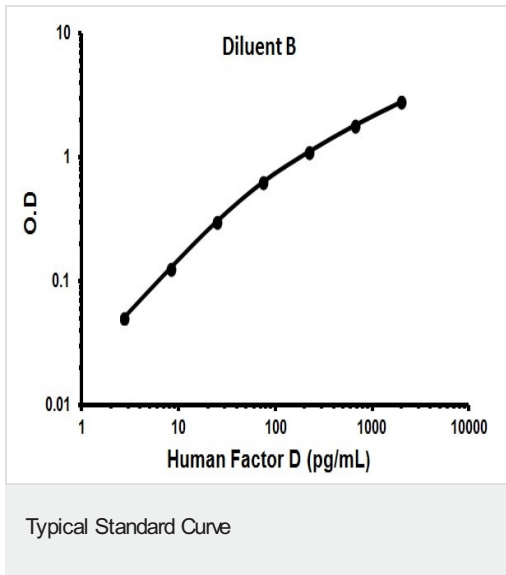
Secreted.

Images



Representative Standard Curve using ab99969.

Typical Standard Curve



Representative Standard Curve using ab99969.

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.com/abpromise> or contact our technical team.

Terms and conditions

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