

Mouse Factor D ELISA Kit ab170244

1 Image

Overview

Product name	Mouse Factor D ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			4.9%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			11%
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Tissue, Cell Lysate			
Assay type	Sandwich (quantitative)			
Sensitivity	19 pg/ml			
Range	50 pg/ml - 500 pg/ml			
Recovery	97 %			
Assay time	4h 00m			
Species reactivity	Reacts with: Mouse			
Product overview	Abcam's Factor D Mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Factor D in cell culture supernatants, urine, plasma and serum.			

A Factor D specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Factor D 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 Factor D captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
10X Diluent M Concentrate	1 x 30ml
50X Biotinylated Mouse Factor D Antibody	1 x 60µl
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
Chromogen Substrate	1 x 7ml
Factor D Microplate (12 x 8 well strips)	1 unit
Factor D Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

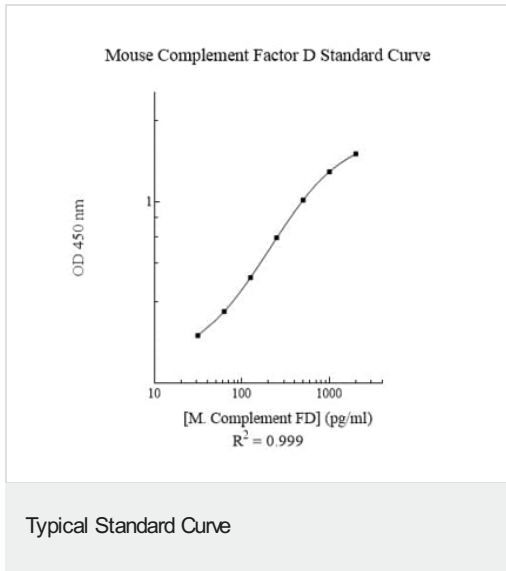
**Function** Factor D cleaves factor B when the latter is complexed with factor C3b, activating the C3bbb complex, which then becomes the C3 convertase of the alternate pathway. Its function is homologous to that of C1s in the classical pathway.

**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 ab170244

**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