

Human Ficolin-2 ELISA Kit ab213778

[1 References](#) [1 Image](#)

Overview

Product name Human Ficolin-2 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	0.977ng/ml	0.044	= 4.6%
2	16	6.182ng/ml	0.497	= 7.3%
3	16	22.201ng/ml	1.354	= 6.1%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	0.979ng/ml	0.052	= 5.4%
2	24	6.528ng/ml	0.509	= 7.8%
3	24	20.934ng/ml	1.611	= 7.7%

Sample type Cell culture supernatant, Serum, Hep Plasma

Assay type Sandwich (quantitative)

Sensitivity < 20 pg/ml

Range 0.78 ng/ml - 50 ng/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

The Human Ficolin-2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213778) is designed for the quantitative measurement of Human Ficolin-2 in cell culture supernatants, serum and plasma (heparin).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for Ficolin-2 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: L26-A313) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Ficolin-2 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human Ficolin-2 amount of sample captured in plate.

Notes Ficolin-2, which was initially identified as L-ficolin, is a protein that in humans is encoded by the FCN2 gene. The product of this gene belongs to the ficolin family of proteins. This family is characterized by the presence of a leader peptide, a short N-terminal segment, followed by a collagen-like region, and a C-terminal fibrinogen-like domain. It is mapped to 9q34.3. This gene is predominantly expressed in the liver, and has been shown to have carbohydrate binding and opsonic activities. Serum concentration of ficolin-2 varies considerably in healthy individuals. FCN2 promoter polymorphisms were associated with marked changes in the ficolin-2 serum concentration.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

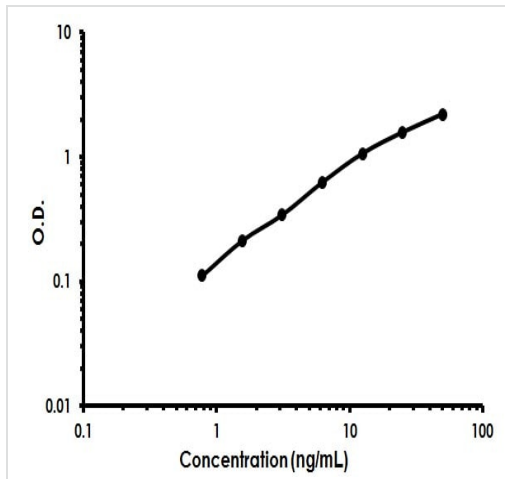
Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Ficolin-2 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human Ficolin-2 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Ficolin-2 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function May function in innate immunity through activation of the lectin complement pathway. Calcium-dependent and GlcNAc-binding lectin. Enhances phagocytosis of *S.typhimurium* by neutrophils, suggesting an opsonic effect via the collagen region.

Tissue specificity Expressed by the liver and secreted in plasma.

Sequence similarities	Belongs to the ficolin lectin family. Contains 1 collagen-like domain. Contains 1 fibrinogen C-terminal domain.
Domain	The fibrinogen-like domain (FBG) contains potential calcium-binding sites that may be involved in carbohydrate binding.
Cellular localization	Secreted.

Images



Human Ficolin-2 ELISA Kit (ab213778) Standard Curve.

Human Ficolin-2 ELISA Kit (ab213778) Standard Curve.

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