

Human alpha Fetoprotein ELISA Kit (AFP) ab108838

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

Overview

Product name Human alpha Fetoprotein ELISA Kit (AFP)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				5.7%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				10.5%

Sample type Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate

Assay type Sandwich (quantitative)

Sensitivity = 0.5 ng/ml

Range 0.625 ng/ml - 40 ng/ml

Recovery 98 %

Assay time 4h 0m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's alpha Fetoprotein (AFP) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of alpha Fetoprotein levels in plasma, serum, cell lysate and culture supernatants, and tissue samples.

An alpha Fetoprotein specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an alpha Fetoprotein specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 alpha Fetoprotein captured in plate.

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

Platform Microplate

Properties

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human alpha Fetoprotein Antibody	1 x 120µl
alpha Fetoprotein Microplate (12 x 8 well strips)	1 unit
alpha Fetoprotein Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function Binds copper, nickel, and fatty acids as well as, and bilirubin less well than, serum albumin. Only a small percentage (less than 2%) of the human AFP shows estrogen-binding properties.

Tissue specificity Plasma. Synthesized by the fetal liver and yolk sac.

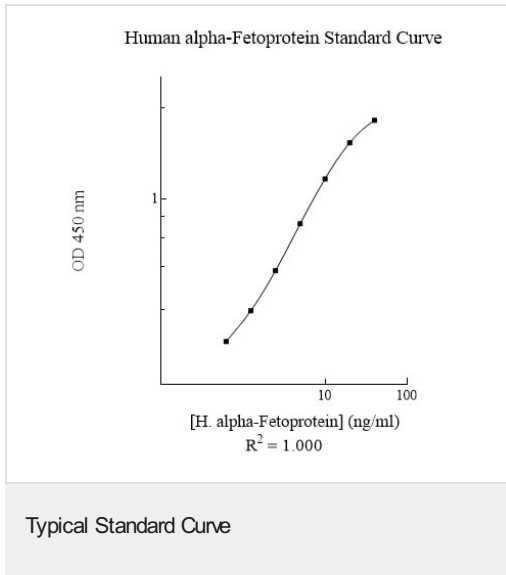
Sequence similarities Belongs to the ALB/AFP/VDB family.
Contains 3 albumin domains.

Developmental stage Occurs in the plasma of fetuses more than 4 weeks old, reaches the highest levels during the 12th-16th week of gestation, and drops to trace amounts after birth. The serum level in adults is usually less than 40 ng/ml. AFP occurs also at high levels in the plasma and ascitic fluid of adults with hepatoma.

Post-translational modifications Independent studies suggest heterogeneity of the N-terminal sequence of the mature protein and of the cleavage site of the signal sequence.
Sulfated.

Cellular localization Secreted.

Images



Representative Standard Curve using ab108838.

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