

Product datasheet

Human alpha Fetoprotein ELISA Kit ab108631

1 Image

Overview

Product name	Human alpha Fetoprotein ELISA Kit
Detection method	Colorimetric
Sample type	Serum
Assay type	Sandwich (quantitative)
Sensitivity	= 2 ng/ml
Range	2 ng/ml - 300 ng/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human

Product overview

Abcam's alpha Fetoprotein (AFP) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of alpha Fetoprotein in Human serum.

A 96-well plate has been precoated with anti-alpha Fetoprotein antibodies. Samples and standards are added to the wells, where alpha Fetoprotein in the sample and standards binds to the precoated antibody. After incubation and washing, added Anti-alpha Fetoprotein-HRP conjugate binds to the antibody-alpha Fetoprotein complex. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of alpha Fetoprotein in the sample and the intensity is measured at 450 nm.

Platform Microplate

Properties

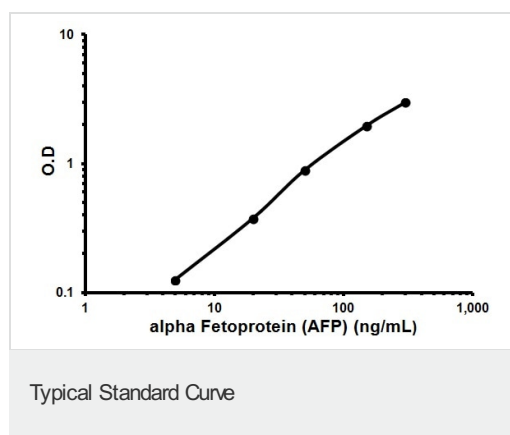
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
alpha Fetoprotein Standard - 0 ng/mL	1 x 0.5ml
alpha Fetoprotein Standard - 150 ng/mL	1 x 0.5ml
alpha Fetoprotein Standard - 20 ng/mL	1 x 0.5ml

Components	1 x 96 tests
alpha Fetoprotein Standard - 300 ng/mL	1 x 0.5ml
alpha Fetoprotein Standard - 5 ng/mL	1 x 0.5ml
alpha Fetoprotein Standard - 50 ng/mL	1 x 0.5ml
Anti-alpha Fetoprotein Coated Microplate (12 x 8 wells)	1 unit
Anti-alpha Fetoprotein HRP Conjugate	1 x 18ml
Stop Solution	1 x 11ml
TMB Reagent	1 x 11ml
Zero Buffer	1 x 13ml

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 ab108631

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