

Human alpha Fetoprotein ELISA Kit (AFP) ab108838

[11 References](#) [画像数 1](#)

製品の概要

製品名 Human alpha Fetoprotein ELISA Kit (AFP)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall				5.7%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				10.5%

サンプルの種類 Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate

アッセイタイプ Sandwich (quantitative)

検出感度 = 0.5 ng/ml

検出範囲 0.625 ng/ml - 40 ng/ml

添加回収試験 98 %

全工程の試験時間 4h 0m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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

機能

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.

組織特異性

Plasma. Synthesized by the fetal liver and yolk sac.

配列類似性

Belongs to the ALB/AFP/VDB family.

Contains 3 albumin domains.

発生段階

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.

翻訳後修飾

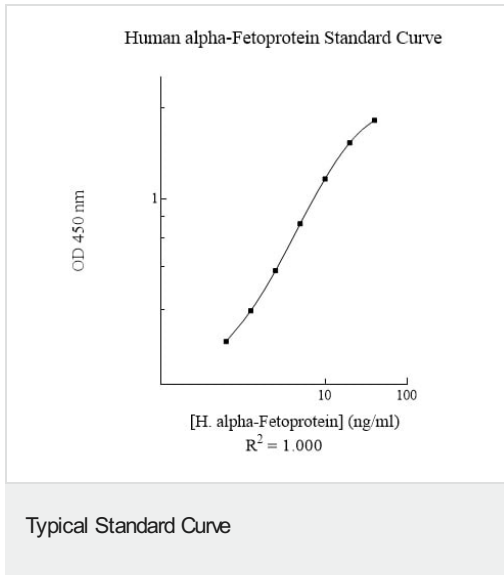
Independent studies suggest heterogeneity of the N-terminal sequence of the mature protein and of the cleavage site of the signal sequence.

Sulfated.

細胞内局在

Secreted.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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