

Human alpha Fetoprotein ELISA Kit (AFP) ab108838

11 References 1 图像

概述

产品名称

检测方法

精确度

样品类型

检测类型

灵敏度

范围

回收率

检测时间

实验步骤

种属反应性

产品概述

人alpha Fetoprotein ELISA试剂盒(AFP)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				5.7%

批次间

样品	n	Mean	SD	CV%
Overall				10.5%

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

Sandwich (quantitative)

= 0.5 ng/ml

0.625 ng/ml - 40 ng/ml

98 %

4h 0m

Multiple steps standard assay

与反应: Human

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.

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

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

功能 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.

组织特异性 Plasma. Synthesized by the fetal liver and yolk sac.

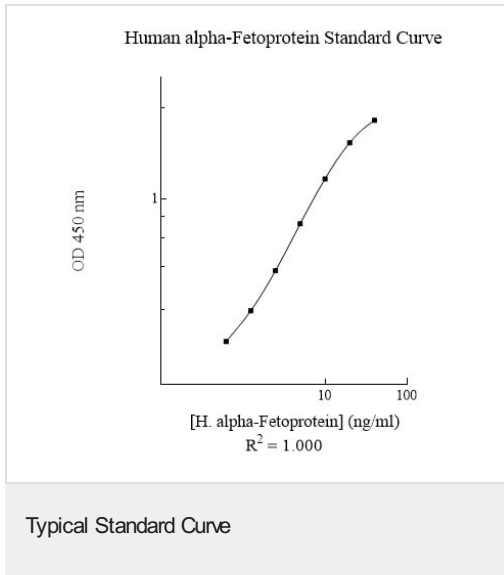
序列相似性 Belongs to the ALB/AFP/VDB family.
Contains 3 albumin domains.

发展阶段 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.

翻译后修饰 Independent studies suggest heterogeneity of the N-terminal sequence of the mature protein and of the cleavage site of the signal sequence.
Sulfated.

细胞定位 Secreted.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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