

Human Transferrin ELISA Kit ab108911

15 References 1 图像

概述

产品名称	人Transferrin ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall			4.2%
	批次间			
	样品	n	Mean	SD
	Overall			10.1%
样品类型	Serum, Plasma			
检测类型	Competitive			
灵敏度	= 89 ng/ml			
范围	0.4 µg/ml - 6.25 µg/ml			
回收率	96 %			
检测时间	2h 00m			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Abcam's Transferrin Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Transferrin in plasma and serum samples.			

A Transferrin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Transferrin 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 inversely proportional to the amount of Transferrin captured in plate.

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	1 x 30ml
6X Biotinylated Human Transferrin (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
Human Transferrin Standard (Lyophilized)	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml
Transferrin Microplate (12 x 8 well strips)	1 unit

功能 Transferrins are iron binding transport proteins which can bind two Fe(3+) ions in association with the binding of an anion, usually bicarbonate. It is responsible for the transport of iron from sites of absorption and heme degradation to those of storage and utilization. Serum transferrin may also have a further role in stimulating cell proliferation.

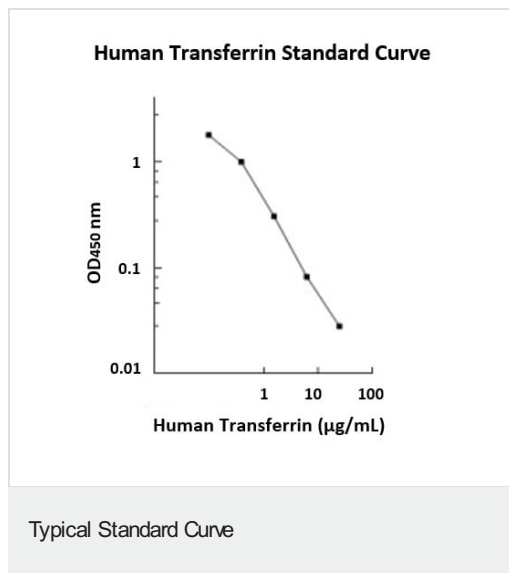
组织特异性 Expressed by the liver and secreted in plasma.

疾病相关 Defects in TF are the cause of atransferrinemia (ATRAF) [MIM:209300]. Atransferrinemia is rare autosomal recessive disorder characterized by iron overload and hypochromic anemia.

序列相似性 Belongs to the transferrin family.
Contains 2 transferrin-like domains.

细胞定位 Secreted.

图片



Representative Standard Curve using ab108911

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