



Product datasheet

Human Transferrin ELISA Kit ab108902

★★★★★ [1 Abreviews](#) [4 References](#) [画像数 1](#)

製品の概要

製品名	Human Transferrin ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				4.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.6%
サンプルの種類	Cell culture supernatant, Saliva, Milk, Tissue, Cell Lysate			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 0.78 ng/ml			
検出範囲	6.25 ng/ml - 50 ng/ml			
添加回収試験	97 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Tranferrin Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Transferrin in milk, urine, saliva, CSF, cell culture, cell lysate, and tissue samples.			

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

Get results in 90 minutes with Human Transferrin ELISA Kit ([ab187391](#)) from our SimpleStep ELISA® range.

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 Transferrin Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
Transferrin Microplate (12 x 8 well strips)	1 unit
Transferrin Standard	1 vial

機能

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.

組織特異性

Expressed by the liver and secreted in plasma.

関連疾患

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

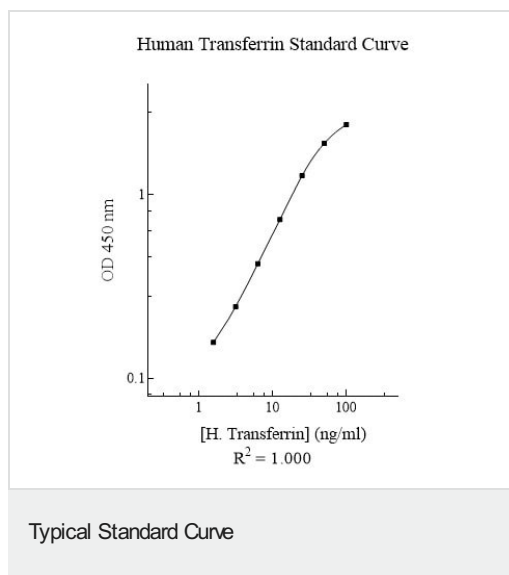
配列類似性

Belongs to the transferrin family.
Contains 2 transferrin-like domains.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab108902.

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