

Human Lactoferrin ELISA Kit ab200015

重组 SimpleStep ELISA

★★★★★ [1 Abreviews](#) [10 References](#) [10 图像](#)

概述

产品名称 人Lactoferrin ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Human Serum	8			5.1%

批次间

样品	n	Mean	SD	CV%
Human Serum	3			5.4%

样品类型 Saliva, Milk, Urine, Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 48.8 pg/ml

范围 156.3 pg/ml - 10000 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Saliva	120	113% - 126%
Milk	109	99% - 118%
Urine	99	99% - 100%
Serum	102	100% - 105%
Cell culture media	86	86% - 86%
Hep Plasma	106	105% - 110%

样品类型	平均%	范围
EDTA Plasma	108	105% - 110%
Cit plasma	110	109% - 111%

检测时间

1h 30m

实验步骤

One step assay

种属反应性

与反应: Human

不与反应: Cow

产品概述

As of October 11 2019, Human Lactoferrin ELISA kit has been re-developed. We have identified new recombinant monoclonal antibodies to provide improved performance and consistency.

Human Lactoferrin ELISA Kit (ab200015) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Lactoferrin protein in edta plasma, hep plasma, milk, saliva, serum, urine, cell culture media, and cit plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Lactoferrin with 48.8 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

ASSAY SPECIFICITY

This kit recognizes native human Lactoferrin protein in serum, plasma, milk, urine, saliva, and cell culture supernatant samples only.

Cell and tissue extract samples have not been tested with this kit.

CROSS REACTIVITY

Recombinant human Transferrin protein was prepared at within the working range of the Lactoferrin standard curve and assayed for cross reactivity. No cross-reactivity was observed.

INTERFERENCE

Recombinant human Transferrin protein was prepared at 50 ng/mL and 2 ng/mL and assayed for interference in the presence 2 ng/mL human Lactoferrin. No Interference was observed.

SPECIES REACTIVITY

This kit recognizes human Lactoferrin protein.

Other species reactivity was determined by measuring 0.5% serum samples of various species, interpolating the protein concentrations from the human standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in human serum assayed at the same dilution.

Reactivity < 3% was determined for the following species: Mouse, Rat, Cow, Monkey

说明

Transferrins are iron binding transport proteins which can bind two Fe^{3+} ions in association with the binding of an anion, usually bicarbonate. Lactoferrin is a member of the transferrin family and is a major iron-binding and multifunctional protein found in exocrine fluids such as breast milk and mucosal secretions. High levels of Lactoferrin are found in saliva and tears, intermediate levels in serum and plasma, and low levels in urine. The canonical sequence of Lactoferrin (known as isoform 1) is 710 amino acids in length and runs at 78.2 kDa. Isoform Delta-Lactoferrin (or DeltaLF) is missing amino acids 1 – 44 of the canonical sequence. Mouse and rat Lactoferrin are 71% and 62% identical to human Lactoferrin, respectively.

Lactoferrin possesses antibacterial, antiviral, anti-inflammatory, and immunoregulatory properties. The presence of Lactoferrin in neutrophils and the release of Lactoferrin during inflammation suggest that Lactoferrin is involved in phagocytic killing and immune responses. Additionally, Lactoferrin mediates bacteriostasis through its ability to sequester free iron and inhibit microbial growth. Lactoferrin also directs bactericidal properties leading to the release of lipopolysaccharides from the bacterial outer membrane. Furthermore, Lactoferrin may play a role in the regulation of bone growth because it has anabolic, differentiating, and anti-apoptotic effects on osteoblasts and can also inhibit osteoclastogenesis. Salivary Lactoferrin plays an important role in protecting the oral cavity against several oral pathogens including *Streptococcus mutans*. Additionally, increased levels of Lactoferrin in fecal samples is used as a marker of active inflammatory bowel disease (IBD).

平台

Microplate (12 x 8 well strips)

性能

存放说明

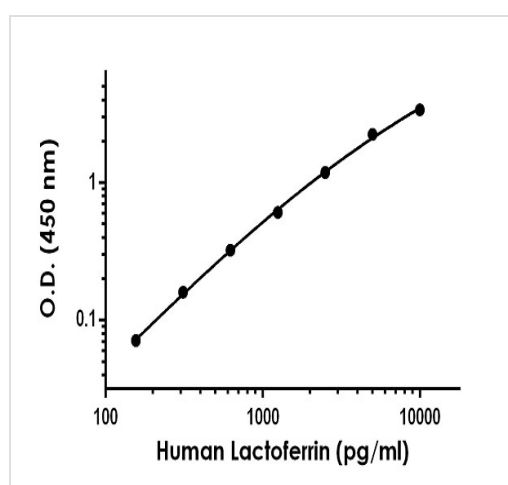
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests	10 x 96 tests
10X Human Lactoferrin Capture Antibody	1 x 600µl	1 x 6000µl
10X Human Lactoferrin Detector Antibody	1 x 600µl	1 x 6000µl

组件	1 x 96 tests	10 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml
Antibody Diluent 5BI	1 x 6ml	10 x 6ml
Human Lactoferrin Lyophilized Purified Protein	2 vials	2 x 10 vials
Plate Seals	1 unit	1 x 10 units
Sample Diluent NS (ab193972)	1 x 50ml	2 x 250ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 x 10 units
Stop Solution	1 x 12ml	1 x 120ml
TMB Development Solution	1 x 12ml	1 x 120ml

功能	<p>Transferrins are iron binding transport proteins which can bind two Fe(3+) ions in association with the binding of an anion, usually bicarbonate.</p> <p>Lactotransferrin has antimicrobial activity which depends on the extracellular cation concentration. Lactoferroxins A, B and C have opioid antagonist activity. Lactoferroxin A shows preference for mu-receptors, while lactoferroxin B and C have somewhat higher degrees of preference for kappa-receptors than for mu-receptors.</p> <p>The lactotransferrin transferrin-like domain 1 functions as a serine protease of the peptidase S60 family that cuts arginine rich regions. This function contributes to the antimicrobial activity.</p>
序列相似性	<p>Belongs to the transferrin family.</p> <p>Contains 2 transferrin-like domains.</p>
细胞定位	<p>Secreted.</p>

图片



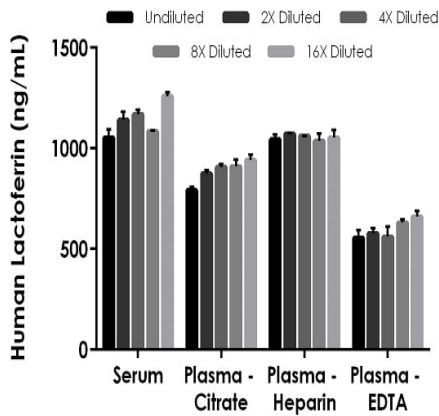
Example of human Lactoferrin standard curve in Sample Diluent NS.

The Lactoferrin standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Standard Curve Measurements			
Concentration (pg/ml)	O.D 450 nm		Mean O.D
	1	2	
0	0.165	0.151	0.158
156.3	0.225	0.233	0.229
312.5	0.317	0.319	0.318
625	0.485	0.479	0.482
1,250	0.762	0.773	0.767
2,500	1.361	1.329	1.345
5,000	2.421	2.384	2.402
10,000	3.538	3.552	3.545

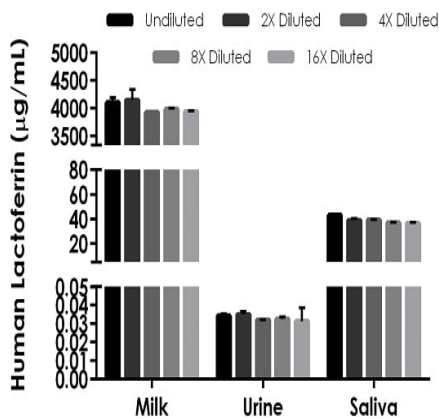
Standard curve

Example of human Lactoferrin standard curve in Sample Diluent NS. The Lactoferrin standard curve was prepared. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



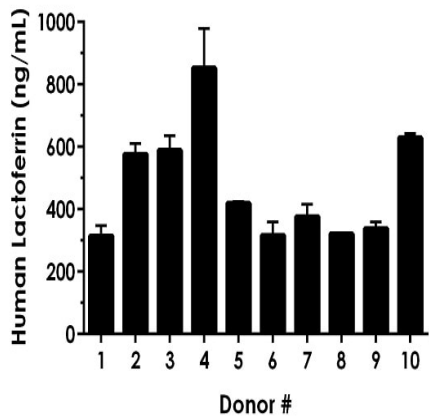
Interpolated concentrations of native Lactoferrin in human serum and plasma samples.

The concentrations of Lactoferrin were measured in duplicates, interpolated from the Lactoferrin standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:200, plasma (citrate) 1:200, plasma (heparin) 1:200, and plasma (EDTA) 1:200. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Lactoferrin concentration was determined to be 1,142 ng/mL in neat serum, 885.1 ng/mL in neat plasma (citrate), 1,054 ng/mL in neat plasma (heparin), and 581.4 ng/mL in neat plasma (EDTA).



Interpolated concentrations of native Lactoferrin in human milk (defatted), urine, and saliva samples.

The concentrations of Lactoferrin were measured in duplicates, interpolated from the Lactoferrin standard curves and corrected for sample dilution. Undiluted samples are as follows: milk 1:500,000, urine 1:20, and saliva 1:10,000. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Lactoferrin concentration was determined to be 4,021.6 µg/mL in neat milk (defatted), 0.033 µg/mL in neat urine, and 39.1 µg/mL in neat saliva.



Serum from ten individual healthy human female donors was measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Lactoferrin concentration was determined to be 473.8 ng/mL with a range of 315.2 – 852.7 ng/mL.

Dilution Factor	Interpolated value	1:200 Human Serum	1:200 Human Plasma (Citrate)	1:200 Human Plasma (Heparin)	1:200 Human Plasma (EDTA)
Undiluted	pg/mL	5,264	3,960	5,223	2,780
	% Expected value	100	100	100	100
2	pg/mL	2,853	2,187	2,680	1,447
	% Expected value	108	110	103	104
4	pg/mL	1,462	1,134	1,327	701.0
	% Expected value	111	114	102	101
8	pg/mL	677.9	568.8	647.7	394.0
	% Expected value	103	115	99	113
16	pg/mL	393.5	294.3	329.1	206.3
	% Expected value	120	119	101	119

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Native Lactoferrin was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

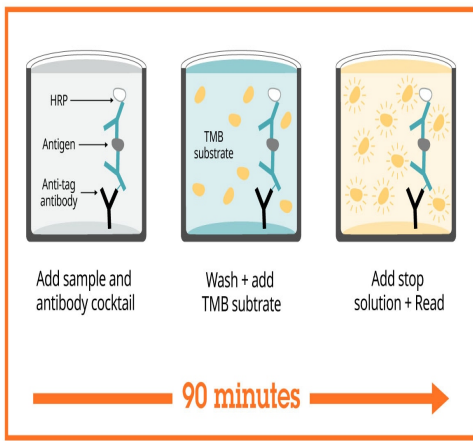
Linearity of dilution.

Dilution Factor	Interpolated value	1:500,000 Human Milk	1:20 Human Urine	1:10,000 Human Saliva	100% Cell Culture Media
Undiluted	pg/mL	8,208	1,714	4,323	1,695
	% Expected value	100	100	100	100
2	pg/mL	4,147	876.8	1,950	893.7
	% Expected value	101	102	90	105
4	pg/mL	1,964	400.9	984.0	476.8
	% Expected value	96	94	91	113
8	pg/mL	996.9	204.0	463.8	252.7
	% Expected value	97	95	86	119
16	pg/mL	492.7	97.99	228.9	102.0
	% Expected value	96	91	85	96

Native Lactoferrin was measured in human milk, urine, and saliva samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

Native purified Lactoferrin protein was spiked into RPMI 1640 cell culture media containing 10% fetal bovine serum and measured in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.





Linearity of dilution.



Sandwich ELISA - Human Lactoferrin ELISA Kit
(ab200015)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Powered by
recombinant antibodies

-  **Research with confidence**
Consistent and reproducible results
-  **Long-term and scalable supply**
Recombinant technology
-  **Success from the first experiment**
Confirmed specificity
-  **Ethical standards compliant**
Animal-free production

Sandwich ELISA - Human Lactoferrin ELISA Kit
(ab200015)

To learn more about the advantages of recombinant antibodies see [here](#).

Get more done with
SimpleStep ELISA

-  **Easy to use**
Single-wash 90-minute protocol
-  **Flexible**
Matched antibody pairs available
-  **Precision antibodies**
High sensitivity, specificity and reproducibility
-  **Scalable**
Now in 10-pack and 384-well formats

Sandwich ELISA - Human Lactoferrin ELISA Kit
(ab200015)

To learn more about the advantages of SimpleStep ELISA[®] kits see [here](#).

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