

Product datasheet

Canine Transferrin ELISA Kit ab157704

1 Image

Overview

Product name	Canine Transferrin ELISA Kit										
Detection method	Colorimetric										
Precision	Intra-assay										
	<table border="1"> <thead> <tr> <th>Sample</th> <th>n</th> <th>Mean</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>< 10%</td> </tr> </tbody> </table>	Sample	n	Mean	SD	CV%	Overall				< 10%
Sample	n	Mean	SD	CV%							
Overall				< 10%							
	Inter-assay										
	<table border="1"> <thead> <tr> <th>Sample</th> <th>n</th> <th>Mean</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>< 10%</td> </tr> </tbody> </table>	Sample	n	Mean	SD	CV%	Overall				< 10%
Sample	n	Mean	SD	CV%							
Overall				< 10%							
Sample type	Serum, Plasma, Other biological fluids										
Assay type	Sandwich (quantitative)										
Sensitivity	4.3054 ng/ml										
Range	6.25 ng/ml - 400 ng/ml										
Recovery	Sample specific recovery										
	<table border="1"> <thead> <tr> <th>Sample type</th> <th>Average %</th> <th>Range</th> </tr> </thead> <tbody> <tr> <td>Serum</td> <td>> 85</td> <td>% - %</td> </tr> </tbody> </table>	Sample type	Average %	Range	Serum	> 85	% - %				
Sample type	Average %	Range									
Serum	> 85	% - %									
Assay duration	Multiple steps standard assay										
Species reactivity	Reacts with: Dog										
Product overview	<p>Abcam's Transferrin Dog ELISA Kit is an <i>in vitro</i> enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of Transferrin levels in serum, tissue extracts and other biological fluids.</p> <p>In this assay the Transferrin present in samples reacts with the anti-Transferrin antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-Transferrin antibodies conjugated with horseradish peroxidase (HRP) are added. These enzyme-labeled antibodies form complexes with the previously bound Transferrin. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of</p>										

bound enzyme varies directly with the concentration of Transferrin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Transferrin in the test sample. The quantity of Transferrin in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X HRP-conjugated anti-dog Transferrin antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chromogen Substrate Solution	1 x 12ml
Dog Transferrin Calibrator (lyophilised)	1 vial
Dog Transferrin ELISA Microplate	1 unit
Stop Solution	1 x 12ml

Function 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.

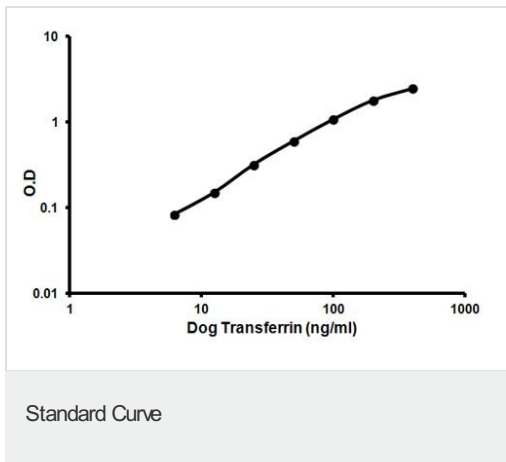
Tissue specificity Expressed by the liver and secreted in plasma.

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

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

Cellular localization Secreted.

Images



Representative standard curve using ab157704 Transferrin Dog ELISA Kit.

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.com/abpromise> or contact our technical team.

Terms and conditions

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