

Product datasheet

Chicken Ovotransferrin ELISA Kit ab157694

1 Image

Overview

Product name Chicken Ovotransferrin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Sample type Serum, Plasma, Other biological fluids

Assay type Sandwich (quantitative)

Sensitivity 0.751 ng/ml

Range 6.25 ng/ml - 400 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	% - %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Chicken

Product overview

Abcam's Ovotransferrin Chicken kit is highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring Ovotransferrin in biological samples of chicken.

In this assay the Ovotransferrin present in samples reacts with the anti-Ovotransferrin antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti-Ovotransferrin antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound Ovotransferrin. 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 bound enzyme varies directly with the concentration of

Ovotransferrin in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of Ovotransferrin in the test sample. The quantity of Ovotransferrin 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-chicken Ovotransferrin antibody	1 x 150µl
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
Chicken Ovotransferrin Calibrator (Lyophilized)	1 vial
Chicken Ovotransferrin ELISA Microplate	1 unit
Chromogen Substrate Solution	1 x 12ml
Stop Solution	1 x 12ml

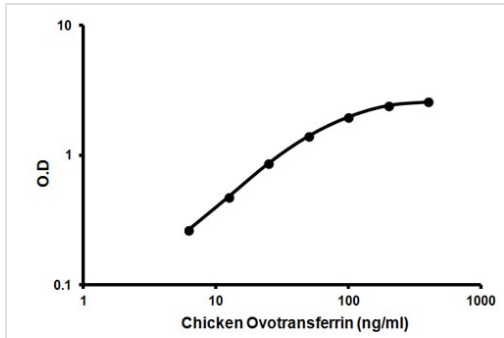
Function Transferrins are iron binding transport proteins which can bind two Fe(3+) ions in association with the binding of an anion, usually bicarbonate. Responsible for the transport of iron from sites of absorption and heme degradation to those of storage and utilization. There are two forms of hen transferrin, ovotransferrin, found in the oviducts and, serum transferrin, secreted by the liver. Serum transferrin may also have a role in stimulating cell proliferation and is regulated by iron levels. Ovotransferrin has a bacteriostatic function and, is not controlled by iron levels.

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

Post-translational modifications Different forms of hen transferrin are distinguished by their carbohydrate composition. Ovotransferrin and embryo serum transferrin but not adult serum transferrin, have bisecting N-acetylglucosamine. Transferrin secreted by embryo hepatocytes in primary culture is marked by the presence of (alpha 1-6) fucosylation of the core N-acetylglucosamine. Serum transferrins also differ in the number of attached neuraminic acid residues. In both embryo forms, sialylation occurs on the Man (alpha 1-3)-linked antennae.

Cellular localization Secreted.

Images



Representative standard curve using ab157694 Ovotransferrin
Chicken ELISA Kit

Standard Curve

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