

Human Ferritin ELISA Kit ab108698

★★★★★ [1 Abreviews](#) [7 References](#) [1 Image](#)

Overview

Product name Human Ferritin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Control	48			< 5.4%

Inter-assay

Sample	n	Mean	SD	CV%
Control	48			< 6.1%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 0.53 ng/ml

Recovery 98.66 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human Ferritin ELISA kit is a competitive ELISA designed for the accurate quantitative measurement of Ferritin (FTL) in serum and plasma.

A 96-well plate has been precoated with anti-Ferritin IgG antibodies. Samples, standards and the Ferritin-HRP conjugate are added to the wells, where Ferritin in the sample and standards binds to the precoated antibody and added Ferritin-HRP conjugate binds to this antibody-AFP complex. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of Ferritin in the sample and the intensity is measured at 450 nm.

Get results in 90 minutes with Human Ferritin ELISA Kit (FLT) ([ab200018](#)) from our SimpleStep ELISA® range.

Platform

Microplate

Properties**Storage instructions**

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

Components	Identifier	1 x 96 tests	1 x 96 tests
10X Washing Solution		1 x 50ml	1 x 50ml
Anti-Ferritin IgG Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit	1 unit
Anti-Ferritin-HRP conjugate		1 x 12ml	1 x 12ml
Cover foil		1 unit	1 unit
Ferritin Control		1 x 1ml	1 x 1ml
Ferritin Standard 0 - 0 ng/mL		1 x 3ml	1 x 3ml
Ferritin Standard 1 - 5 ng/mL		1 x 1ml	1 x 1ml
Ferritin Standard 2 - 20 ng/mL		1 x 1ml	1 x 1ml
Ferritin Standard 3 - 100 ng/mL		1 x 1ml	1 x 1ml
Ferritin Standard 4 - 400 ng/mL		1 x 1ml	1 x 1ml
Ferritin Standard 5 - 800 ng/mL		1 x 1ml	1 x 1ml
Stop Solution		1 x 15ml	1 x 15ml
Strip holder		1 unit	1 unit
TMB Substrate Solution		1 x 15ml	1 x 15ml

Function

Stores iron in a soluble, non-toxic, readily available form. Important for iron homeostasis. Iron is taken up in the ferrous form and deposited as ferric hydroxides after oxidation. Also plays a role in delivery of iron to cells. Mediates iron uptake in capsule cells of the developing kidney.

Involvement in disease

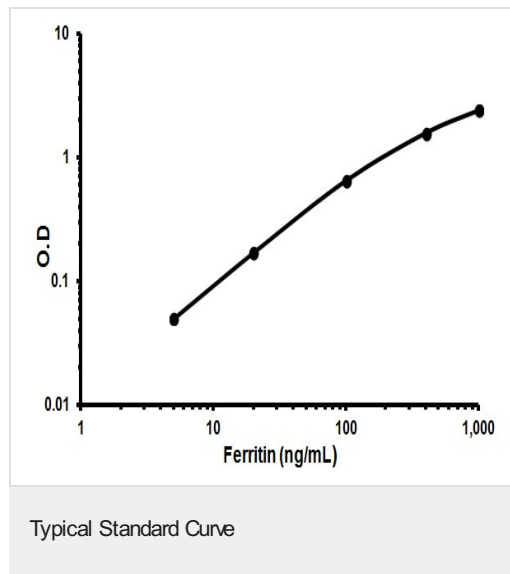
Defects in FTL are the cause of hereditary hyperferritinemia-cataract syndrome (HHCS) [MIM:600886]. It is an autosomal dominant disease characterized by early-onset bilateral cataract. Affected patients have elevated level of circulating ferritin. HHCS is caused by mutations in the iron responsive element (IRE) of the FTL gene.

Defects in FTL are the cause of neurodegeneration with brain iron accumulation type 3 (NBIA3) [MIM:606159]; also known as adult-onset basal ganglia disease. It is a movement disorder with heterogeneous presentations starting in the fourth to sixth decade. It is characterized by a variety of neurological signs including parkinsonism, ataxia, corticospinal signs, mild nonprogressive cognitive deficit and episodic psychosis. It is linked with decreased serum ferritin levels.

Sequence similarities

Belongs to the ferritin family.
Contains 1 ferritin-like diiron domain.

Images



Representative Standard Curve using ab108698

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