

Human Ferritin ELISA Kit ab108698

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

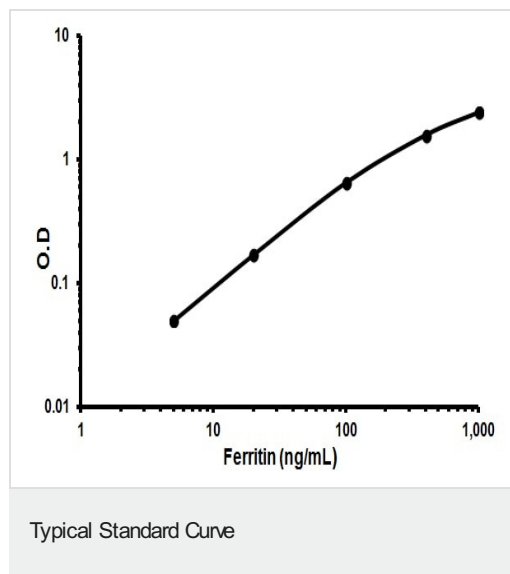
概述

产品名称	人Ferritin ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Control	48		CV% < 5.4%
	批次间			
	样品	n	Mean	SD
	Control	48		CV% < 6.1%
样品类型	Serum, Hep Plasma, Cit plasma			
检测类型	Sandwich (quantitative)			
灵敏度	0.53 ng/ml			
回收率	98.66 %			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	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.

平台	Microplate		
性能			
存放说明	Store at +4°C. Please refer to protocols.		
组 件	标识符	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
功能	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.		
疾病相关	<p>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.</p> <p>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.</p>		
序列相似性	<p>Belongs to the ferritin family.</p> <p>Contains 1 ferritin-like diiron domain.</p>		



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.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