

Product datasheet

Human Frataxin ELISA Kit (10 x 96 well plate) ab201121

SimpleStep ELISA

5 Images

Overview

Product name	Human Frataxin ELISA Kit (10 x 96 well plate)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	HeLa	9			2.3%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	HeLa	3			6.3%
Sample type	Cell culture extracts, Tissue Extracts				
Assay type	Sandwich (quantitative)				
Sensitivity	2.1 pg/ml				
Range	9.88 pg/ml - 800 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		92		86% - 98%
	Cell culture media		97		93% - 101%
	EDTA Plasma		89		88% - 92%
Assay time	1h 30m				
Assay duration	One step assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's Frataxin in vitro SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Frataxin protein in in human cell and tissue extracts.				

The SimpleStep ELISA® employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

Notes

Frataxin is a 17 kDa nuclear-encoded mitochondrial protein. In humans the gene is localized on chromosome 9 and is highly conserved during evolution. The gene is expressed in every cell, although in varying levels in different tissues and during development. The specific function of frataxin is still unknown, but it has been shown to play a role in iron metabolism. Studies have demonstrated that the deletion of the frataxin gene in yeast results in iron accumulation in mitochondria and loss of respiration. Recombinant human frataxin has been shown to bind iron in vitro, and increased mitochondrial iron levels have been observed in patients with Friedreich's Ataxia (FRDA). FRDA is an autosomal recessive, progressive degenerative disease characterized by neurodegeneration and cardiomyopathy it is the most common inherited ataxia.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Microplate

Properties

Storage instructions Please refer to protocols.

Components	10 x 96 tests
10X Human Frataxin Capture Antibody	1 x 6ml
10X Human Frataxin Detector Antibody	1 x 6ml
10X Wash Buffer PT (ab206977)	1 x 200ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 5ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 50ml
Antibody Diluent 4BI	1 x 60ml
Human Frataxin Recombinant Protein	1 x 2ml
Plate Seals	10 units

Components	10 x 96 tests
Sample Diluent NS (ab193972)	1 x 120ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	10 units
Stop Solution	1 x 120ml
TMB Development Solution	1 x 120ml

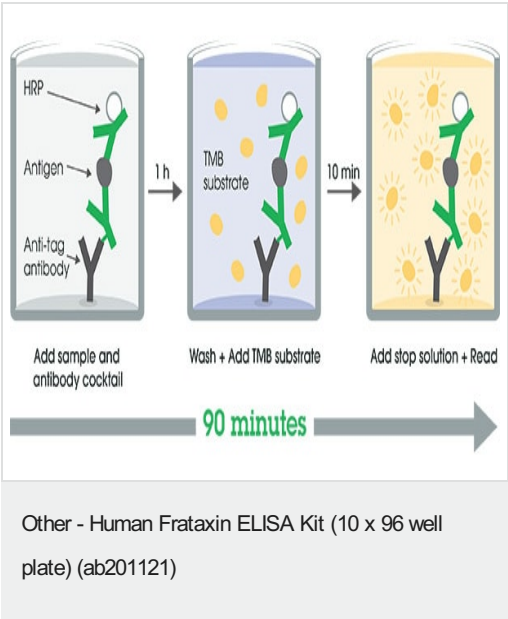
Relevance

Function: Promotes the biosynthesis of heme and assembly and repair of iron-sulfur clusters by delivering Fe2+ to proteins involved in these pathways. May play a role in the protection against iron-catalyzed oxidative stress through its ability to catalyze the oxidation of Fe2+ to Fe3+; the oligomeric form but not the monomeric form has in vitro ferroxidase activity. May be able to store large amounts of iron in the form of a ferrihydrite mineral by oligomerization; however, the physiological relevance is unsure as reports are conflicting and the function has only been shown using heterologous overexpression systems. Modulates the RNA-binding activity of ACO1.

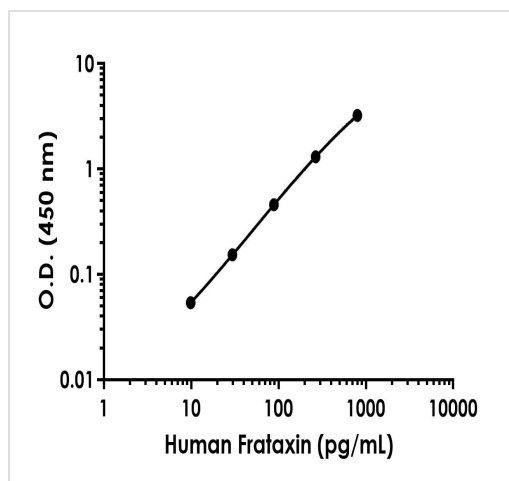
Cellular localization

Cytoplasmic and Mitochondrial

Images

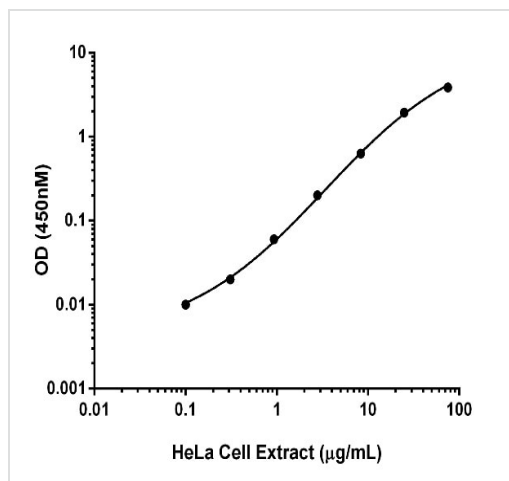


ELISA Protocol Summary



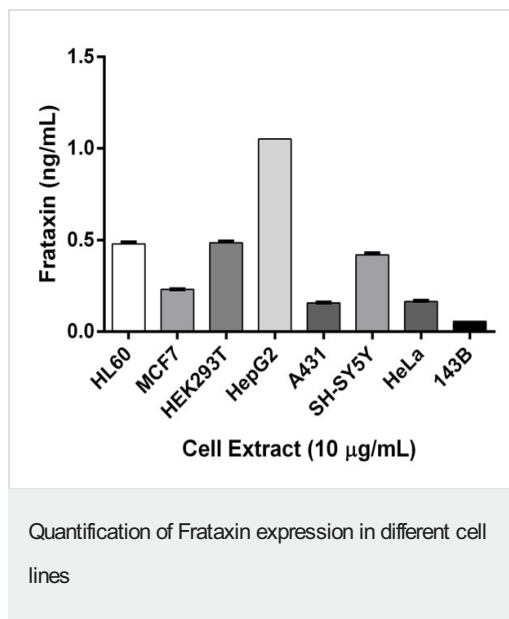
Background-subtracted data values (mean \pm SD) are graphed.

Example of Frataxin standard curve

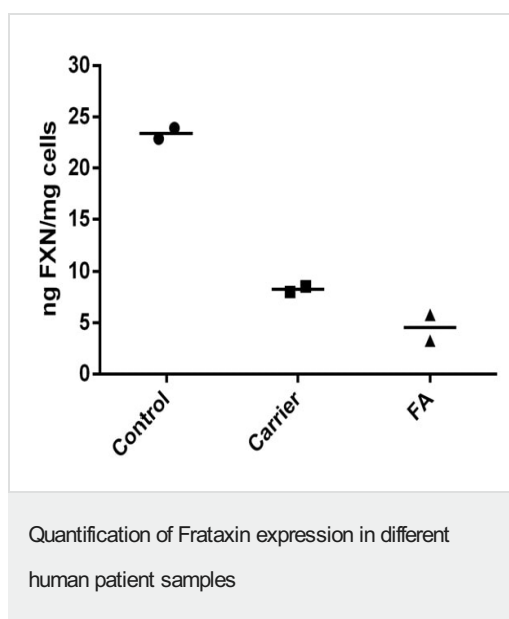


Background subtracted data from duplicate measurements are plotted.

Titration of HeLa cell extract within the working range of the assay



Interpolated values of Frataxin are plotted for the indicated cell lines based on a extract load of 10 µg/mL.



Transformed B lymphocyte cells from Friedreich's Ataxia (FRDA) samples, due to a stable homozygous GAA repeat insertion, are compared to heterozygous carrier B lymphocyte cells and control B lymphocyte cells. B lymphocyte cell extracts were analyzed across a 7-point titration (0.1–100 µg/mL) and frataxin levels were interpolated from the standard curve. Average interpolated values of Frataxin are plotted.

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