

Product datasheet

FRAP Assay Kit (Ferric Reducing Antioxidant Power Assay) ab234626

[3 References](#) [3 Images](#)

Overview

Product name	FRAP Assay Kit (Ferric Reducing Antioxidant Power Assay)
Detection method	Colorimetric
Sample type	Serum, Plasma, Other biological fluids
Assay type	Quantitative
Sensitivity	0.2 mM
Product overview	FRAP Assay Kit ab234626 provides a quick, sensitive and easy way for measuring the antioxidant capacity of various biological samples. FRAP assay stands for Ferric Reducing Antioxidant Power Assay.

The FRAP assay is high-throughput, adaptable and can detect antioxidant capacities as low as 0.2 mM Fe^{2+} equivalents.

The assay measures the antioxidant potential in samples through the reduction of ferric iron (Fe^{3+}) to ferrous iron (Fe^{2+}) by antioxidants present in the samples. Following the reduction of the ferric iron, a blue color develops that is read colorimetrically at 594 nm.

The antioxidant potential of samples is determined using a ferrous iron standard curve and results are expressed as Fe^{2+} equivalents (μM) or FRAP value.

Notes	This product is manufactured by BioVision, an Abcam company and was previously called K515 Ferric Reducing Antioxidant Power (FRAP) Assay Kit (Colorimetric). K515-200 is the same size as the 200 test size of ab234626.
--------------	---

Related products

Review the [oxidative stress marker and assay guide](#) to learn about more assays for oxidative stress.

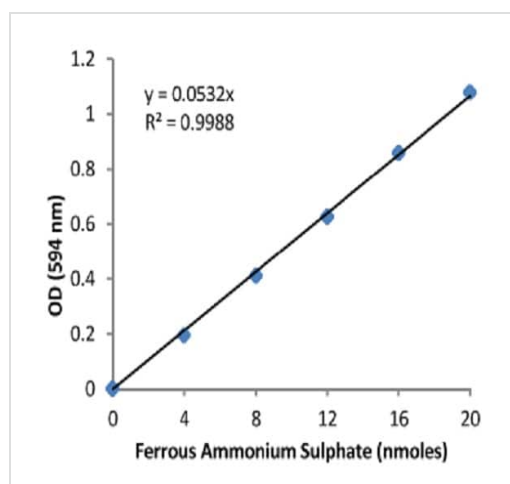
Platform	Microplate reader
-----------------	-------------------

Properties

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

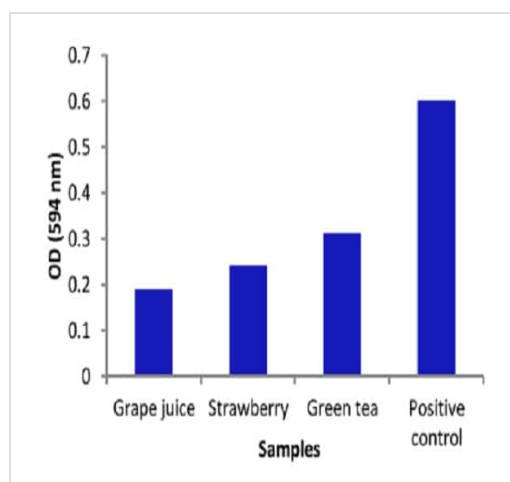
Components	200 tests
Ferric Iron Solution	1 x 4ml
Ferrous Standard	1 x 1ml
FRAP Assay Buffer	1 x 50ml
FRAP Positive Control	1 x 0.1ml
FRAP Probe	1 x 4ml

Images



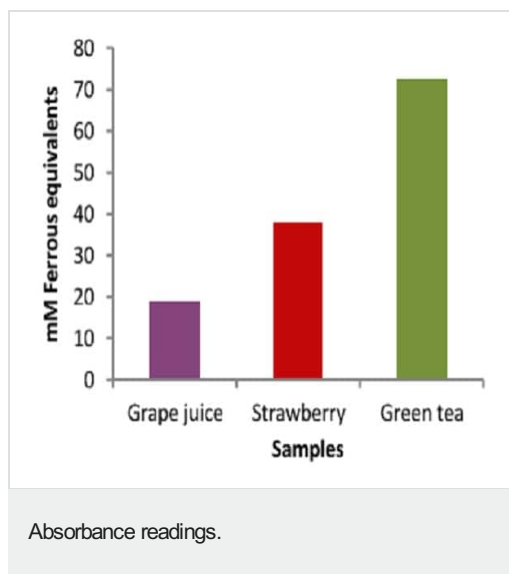
Ferrous standard curve.

Standard Curve



Absorbance readings for positive control and 10 μ l diluted solutions of grape juice (1:50 dilution with dH_2O), strawberry methanolic extract (extract made from 50 mg of freeze-dried strawberries in 5 ml of extraction solvent and final solution diluted 1:80 times with dH_2O) and green tea (brewed for 5 minutes and diluted 1:120 times with dH_2O).

Absorbance readings.



Absorbance readings for mM Ferrous equivalents or FRAP of grape juice, strawberry and green tea.

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