

# FRAP Assay Kit (Ferric Reducing Antioxidant Power Assay) ab234626

**3 References**   **画像数 3**

### 製品の概要

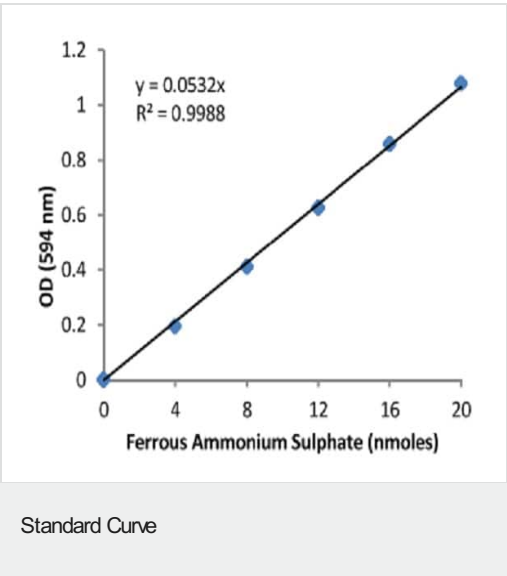
製品名	FRAP Assay Kit (Ferric Reducing Antioxidant Power Assay)
検出方法	Colorimetric
サンプルの種類	Serum, Plasma, Other biological fluids
アッセイタイプ	Quantitative
検出感度	0.2 mM
製品の概要	<p>FRAP Assay Kit ab234626 provides a quick, sensitive and easy way for measuring the antioxidant capacity of various biological samples. FRAP assay stands for Ferric Reducing Antioxidant Power Assay.</p> <p>The FRAP assay is high-throughput, adaptable and can detect antioxidant capacities as low as 0.2 mM <math>\text{Fe}^{2+}</math> equivalents.</p> <p>The assay measures the antioxidant potential in samples through the reduction of ferric iron (<math>\text{Fe}^{3+}</math>) to ferrous iron (<math>\text{Fe}^{2+}</math>) by antioxidants present in the samples. Following the reduction of the ferric iron, a blue color develops that is read colorimetrically at 594 nm.</p> <p>The antioxidant potential of samples is determined using a ferrous iron standard curve and results are expressed as <math>\text{Fe}^{2+}</math> equivalents (<math>\mu\text{M}</math>) or FRAP value.</p>
特記事項	<p>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.</p> <p><b>Related products</b></p> <p>Review the <a href="#">oxidative stress marker and assay guide</a> to learn about more assays for oxidative stress.</p>
試験プラットフォーム	Microplate reader

### 製品の特性

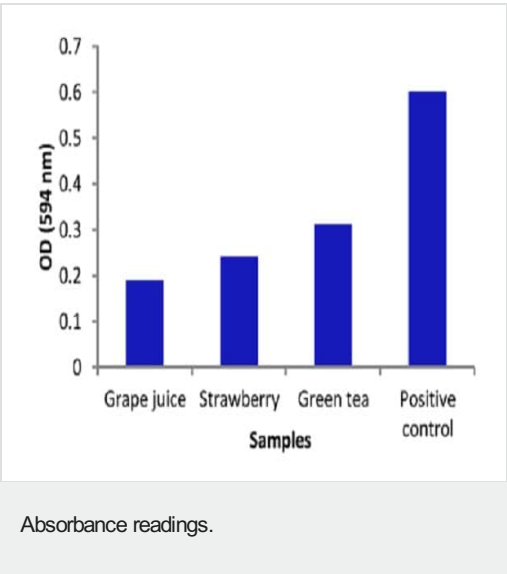
**保存方法**   Store at +4°C. Please refer to protocols.

内容	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

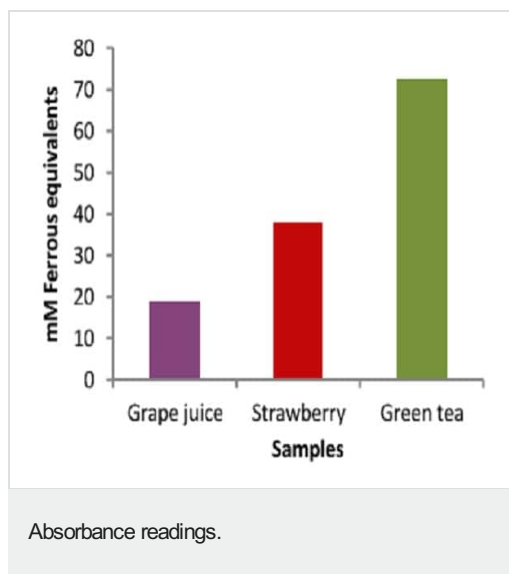
画像



Ferrous standard curve.



Absorbance readings for positive control and 10 µl diluted solutions of grape juice (1:50 dilution with dH<sub>2</sub>O), 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<sub>2</sub>O) and green tea (brewed for 5 minutes and diluted 1:120 times with dH<sub>2</sub>O).



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.co.jp/abpromise> or contact our technical team.

#### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors