

Total Iron-Binding Capacity (TIBC) and Serum Iron Assay Kit ab239715

★★★★★ [3 Abreviews](#) [3 References](#) [2 Images](#)

Overview

Product name	Total Iron-Binding Capacity (TIBC) and Serum Iron Assay Kit
Detection method	Colorimetric
Sample type	Serum, Plasma
Product overview	Total Iron-Binding Capacity (TIBC) and Serum Iron Assay Kit (ab239715) provides a simple, sensitive, and high-throughput adaptable approach to detect physiological Total iron-binding capacity (TIBC) and Serum iron. Those values indicate the requisite iron for transferrin saturation and Serum Iron respectively. Those measurements can be used to detect and monitor transferrin saturation and also iron-deficiency anemia and chronic inflammatory diseases.
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K392 Total Iron-Binding Capacity (TIBC) and Serum Iron Assay Kit (Colorimetric). K392-100 is the same size as the 100 test size of ab239715.</p> <p>The ab239715 assay is not compatible with plasma samples collected with EDTA or citrate anticoagulants. Plasma samples collected in heparin are compatible.</p>
Platform	Microplate reader

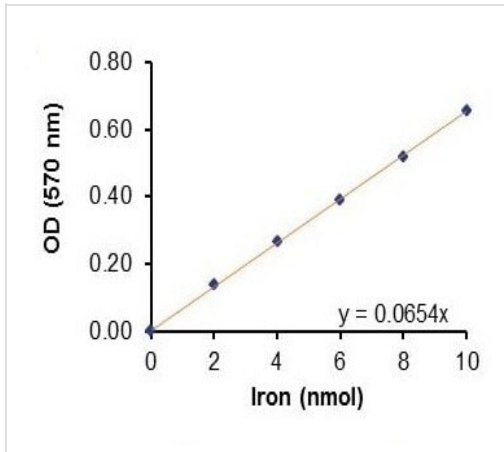
Properties

Storage instructions	Store at -20°C. Please refer to protocols.	
Components	100 tests	2000 tests
Iron Solution	1 x 100µl	20 x 100µl
Iron Standard II	1 x 100µl	20 x 100µl
TIBC Assay Buffer	1 x 25ml	20 x 25ml
TIBC Detector	2 x 1.5ml	40 x 1.5ml
TIBC Developer Solution	1 x 5ml	20 x 5ml

Relevance Iron is essential to nearly all known organisms. It is generally stored in the centre of

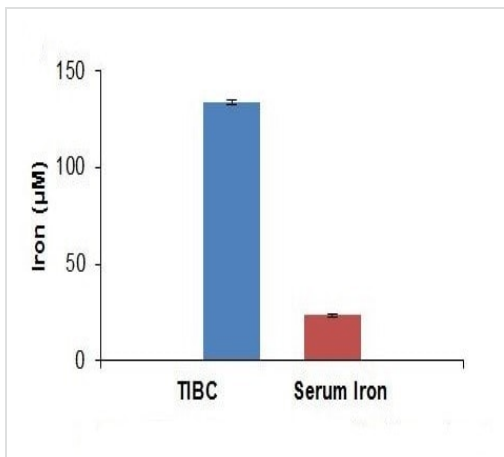
metalloproteins, in the heme complex, and in oxygen carrier proteins. Inorganic iron also contributes to redox reactions in the iron-sulfur clusters of many enzymes, such as nitrogenase and hydrogenase.

Images



Iron standard Curve generated using the kit protocol.

Iron Standard Curve



Serum Iron and TIBC determination of serum. Assays were performed following the kit protocol.

Serum Iron and TIBC determination of serum.

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