

BCA protein assay kit for low concentrations ab207002

14 References 2 图像

概述

产品名称	BCA蛋白assay试剂盒for low concentrations
检测方法	Colorimetric
样品类型	Cell Lysate, Purified protein, Tissue Lysate
产品概述	<p>Abcam's BCA protein assay kit for low concentrations (ab207002) measures total protein concentration of pure proteins, extracts or lysates. It provides an ultra-sensitive method to detect and quantify total protein concentration. The kit has been optimized for determining total protein content in low protein concentration samples (0.5-40 µg/mL), even in the presence of detergents.</p> <p>The assay is based on the reaction of bicinchoninic acid (BCA) with the cuprous cation (Cu<sup>1+</sup>), which is generated by reduction of cupric cation (Cu<sup>2+</sup>) by the protein in alkaline conditions. The Cu<sup>1+</sup>-BCA chelate is a water-soluble complex and exhibits a strong absorbance at 562 nm that is linear over a concentration range of 0.5-40 µg/mL. The kit includes Bovine Serum Albumin (BSA) as a reference protein standard.</p> <p>Support on how to determine protein concentration can be found <a href="#">here</a>.</p>
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K814 Extra Sense™ BCA Protein Assay Kit. K814-2500 is the same size as the 2500 test size of ab207002.</p>
平台	Microplate reader

性能

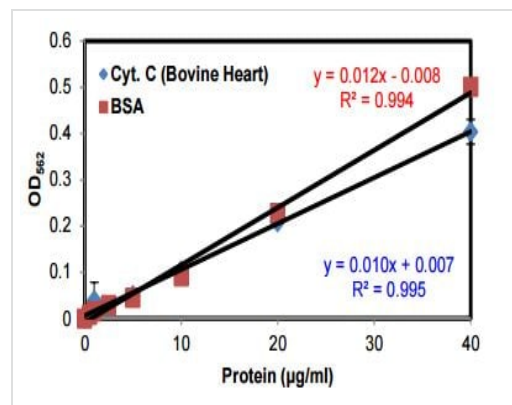
存放说明Store at room temperature. Please refer to protocols.

组件	2500 tests
BCA Reagent A	1 x 190ml
BCA Reagent B	1 x 190ml
BCA Reagent C	1 x 9ml
BSA Standard II	5 x 1ml

## 相关性

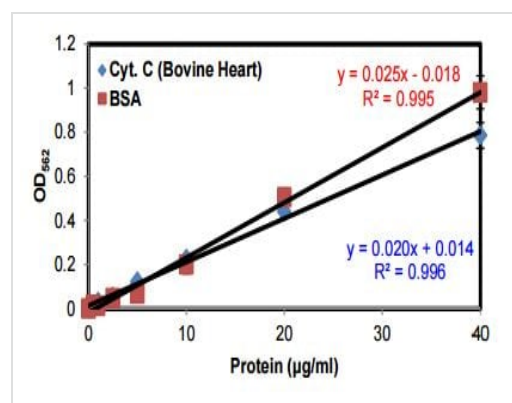
The BCA Protein assay (or Smith assay) is a biochemical assay for determining the total level of protein in a solution. The total protein concentration is determined by measuring the color change of the sample solution from green to purple which is proportional to the protein concentration.

## 图片



Typical standard curve obtained for BSA and Cytochrome C from bovine heart using ab207002 and the microplate procedure (37°C for 2 hours).

Typical absorbance plot obtained for BSA and Cytochrome C from Bovine Heart by using microplate procedure (37°C for 2 hours).



Typical standard curve obtained for BSA and Cytochrome C from bovine heart by using ab207002 and the test tube procedure (60°C for 1 hour).

Typical absorbance plot obtained for BSA and Cytochrome C from Bovine Heart by using test tube procedure (60°C for 1 hour).

**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