

BCA protein assay kit for low concentrations ab207002

[14 References](#) [2 Images](#)

Overview

Product name	BCA protein assay kit for low concentrations
Detection method	Colorimetric
Sample type	Cell Lysate, Purified protein, Tissue Lysate
Product overview	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.

The assay is based on the reaction of bicinchoninic acid (BCA) with the cuprous cation (Cu^{1+}), which is generated by reduction of cupric cation (Cu^{2+}) by the protein in alkaline conditions. The Cu^{1+} -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.

Support on how to determine protein concentration can be found [here](#).

Notes	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.
--------------	----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------

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

Properties

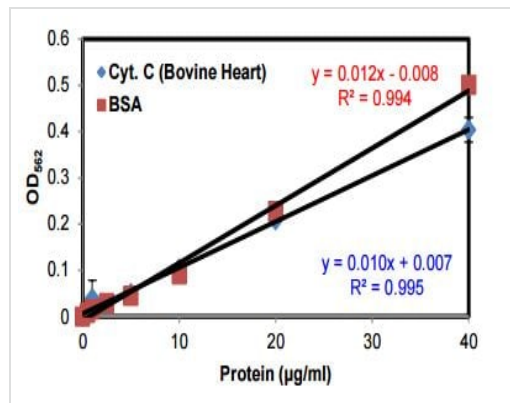
Storage instructions	Store at room temperature. Please refer to protocols.
-----------------------------	-------------------------------------------------------

Components	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

Relevance

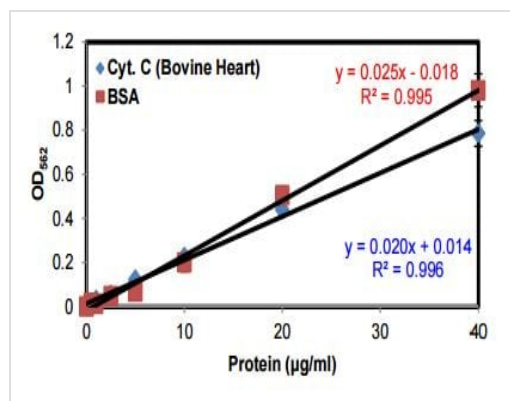
The BCA Protein assay (or Smith assay) is a biochemical assay for determining the total level of protein in a solution. The total protein concentration is determined by measuring the color change of the sample solution from green to purple which is proportional to the protein concentration.

Images



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 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 test tube procedure (60°C for 1 hour).

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).

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