

Product datasheet

BCA protein assay kit reducing agent compatible (test tube) ab207004

3 Images

Overview

Product name	BCA protein assay kit reducing agent compatible (test tube)
Detection method	Colorimetric
Sample type	Cell Lysate, Purified protein, Tissue Lysate
Product overview	Abcam's BCA protein assay kit reducing agent compatible (test tube) (ab207004) measures total protein concentration of pure proteins, extracts or lysates in the presence of reducing agents.

Unlike the majority of kits on the market, ab207004 is compatible with strong reducing agents such as TCEP up to 20 mM, DTT up to 10 mM and β -mercaptoethanol up to 35 mM. This kit is based on the chelation 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 assay is linear over a wide range of protein concentrations (25-2000 $\mu\text{g}/\text{mL}$). The kit also includes Bovine Serum Albumin (BSA) as a reference protein standard. The kit has been formatted for test tube-based assays.

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

Properties

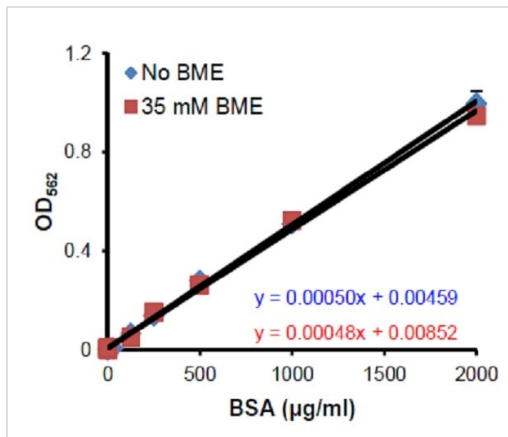
Storage instructions Store at room temperature. Please refer to protocols.

Components	250 tests
BCA Reagent A	1 x 250ml
BCA Reagent B	1 x 20ml
Blocking Reagent	20 vials
Blocking Reagent Buffer	1 x 20ml
BSA Standard	10 x 1ml

Relevance

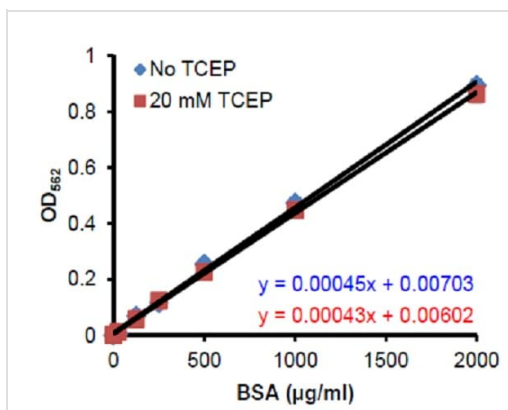
The BCA Protein assay (or Smith assay) is a biochemical assay for determining the total level of protein in a solution. The total protein concentration is determined by measuring the color change of the sample solution from green to purple which is proportional to the protein concentration.

Images



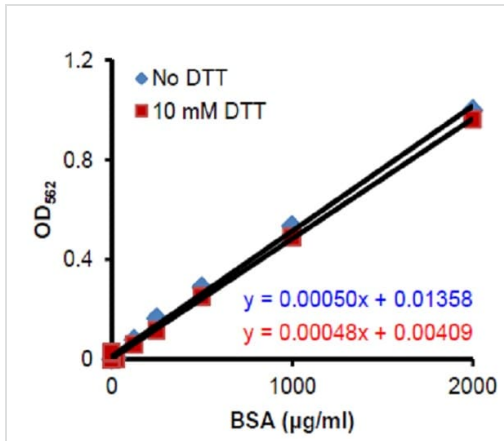
Standard curves for BSA containing blocking reagent in the presence and absence of 35 mM β -mercaptoethanol (BME). Generated using ab207004.

Standard Curves for BSA containing blocking reagent in the presence and absence of 35 mM β -mercaptoethanol (BME)



Standard curves for BSA containing blocking reagent in the presence and absence of 20 mM TCEP. Generated using ab207004.

Standard Curves for BSA containing blocking reagent in the presence and absence of 20 mM TCEP



Standard curves for BSA containing blocking reagent in the presence and absence of 10 mM DTT obtained using ab207004.

Standard Curves for BSA containing blocking reagent in the presence and absence of 10 mM DTT

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