

## BCA Protein Assay Kit

### Table of Contents

|  |   |
|--|---|
| About the Kit .....                      | 2 |
| Description .....                        | 2 |
| Components .....                         | 2 |
| Storage .....                            | 2 |
| BCA Protein Assay .....                  | 3 |
| Preparation of BSA standards .....       | 3 |
| Preparation of BCA working reagent ..... | 3 |
| Assay protocols .....                    | 4 |
| Troubleshooting .....                    | 5 |
| Substance Compatibility .....            | 5 |
| References .....                         | 7 |

© 2011 EMD Chemicals, Inc., an affiliate of Merck KGaA, Darmstadt, Germany. All rights reserved. BugBuster<sup>®</sup>, PopCulture<sup>®</sup>, the Novagen name and logo are registered trademarks of EMD Chemicals, Inc. in the United States and certain other jurisdictions. CytoBuster<sup>™</sup>, Non-Interfering Protein Assay<sup>™</sup>, and Reportasol<sup>™</sup> are trademarks of EMD Chemicals, Inc. Triton<sup>®</sup> is a registered trademark of Rohm & Haas Co. Brij<sup>®</sup> and Tween<sup>®</sup> are registered trademarks of ICI Americas. Zwittergent<sup>®</sup> is a registered trademark of American Hoechst Corporation. BCA Protein Assay Kits are limited for research purposes only.

---

#### USA and Canada

Tel (800) 628-8470  
bioscienceshelp@  
emdchemicals.com

#### Europe

**France**  
Freephone  
0800 126 461

**Germany**  
Freecall  
0800 100 3496

**Ireland**  
Toll Free  
1800 409 445

**United Kingdom**  
Freephone  
0800 622 935

**All other  
European Countries**  
+44 115 943 0840

#### All Other Countries

**Contact Your Local Distributor**  
www.merck4biosciences.com  
bioscienceshelp@  
emdchemicals.com

---

techservice@merckbio.eu

www.merck4biosciences.com

FOR RESEARCH USE ONLY. NOT FOR HUMAN OR DIAGNOSTIC USE.

## About the Kit

BCA Protein Assay Kit

71285-3

### Description

The BCA Protein Assay is based on a biuret reaction, the reduction of  $\text{Cu}^{2+}$  to  $\text{Cu}^{1+}$  by protein in an alkaline solution, and a concentration-dependent detection of the monovalent copper ions ( $\text{Cu}^{1+}$ ) produced. Bicinchoninic acid is a chromogenic reagent that chelates with the reduced copper, producing a purple reaction complex with strong absorbance at 562 nm (1). The kit can be used to determine protein concentrations in the range of 20–2,000  $\mu\text{g}/\text{ml}$  in either the standard assay or microassay size. Kit components are sufficient to complete 500 standard size reactions (50  $\mu\text{l}$  protein sample plus 1 ml reagent) or 2,500 micro-scale reactions (25  $\mu\text{l}$  protein sample plus 200  $\mu\text{l}$  reagent in 96-well plate). For convenient preparation of protein concentration standard curves, a bovine serum albumin (BSA) standard (2 mg/ml) is provided.

The BCA Protein Assay is robust and can be performed in the presence of many chemical compounds and detergents including the Novagen BugBuster<sup>®</sup>, PopCulture<sup>®</sup>, CytoBuster<sup>™</sup>, Reportasol<sup>™</sup> and Insect PopCulture<sup>®</sup> lysis reagents. Some reagents, including chelating agents, strong acids or bases, and reducing agents interfere with the reduction and chelating reactions on which this assay depends.

### Components

- 500 ml BCA Solution (bicinchoninic acid, sodium carbonate, sodium tartrate, and sodium bicarbonate in 0.1 M NaOH, pH 11.25)
- 15 ml 4% Cupric Sulfate
- 3 × 1 ml BSA Standard (2 mg/ml)

### Storage

Store all components at room temperature.

## BCA Protein Assay

### Preparation of BSA standards

Table 1 below gives guidelines for preparing diluted BSA standards. Prepare each standard in a clean vial. To ensure the effects of the buffer are taken into consideration, it is recommended that the dilutions are made in the same buffer as the samples. Deionized water may be used instead of the sample buffer, however, any interference from the buffer will not be corrected for in the BSA standard curve.

**Table 1 Preparation of BSA Standards**

| <b>Dilutions for standard assay</b> |                              |                          |                                |
|-------------------------------------|------------------------------|--------------------------|--------------------------------|
| <b>Tube</b>                         | <b>Volume of BSA</b>         | <b>Volume of diluent</b> | <b>Final BSA concentration</b> |
| 1                                   | 250 µl from 2 mg/ml solution | 250 µl                   | 1,000 µg/ml                    |
| 2                                   | 250 µl from tube 1           | 250 µl                   | 500 µg/ml                      |
| 3                                   | 250 µl from tube 2           | 250 µl                   | 250 µg/ml                      |
| 4                                   | 300 µl from tube 3           | 300 µl                   | 125 µg/ml                      |
| 5                                   | 100 µl from tube 4           | 400 µl                   | 25 µg/ml                       |
| 6                                   | 0                            | 400 µl                   | 0 µg/ml                        |

  

| <b>Dilutions for enhanced assay</b> |                              |                          |                                |
|-------------------------------------|------------------------------|--------------------------|--------------------------------|
| <b>Tube</b>                         | <b>Volume of BSA</b>         | <b>Volume of diluent</b> | <b>Final BSA concentration</b> |
| 1                                   | 100 µl from 2 mg/ml solution | 700 µl                   | 250 µg/ml                      |
| 2                                   | 400 µl from tube 1           | 400 µl                   | 125 µg/ml                      |
| 3                                   | 300 µl from tube 2           | 450 µl                   | 50 µg/ml                       |
| 4                                   | 200 µl from tube 3           | 200 µl                   | 25 µg/ml                       |
| 5                                   | 100 µl from tube 4           | 400 µl                   | 5 µg/ml                        |
| 6                                   | 0                            | 400 µl                   | 0 µg/ml                        |

### Preparation of BCA working reagent

Table 2 below gives guidelines for preparing the BCA working reagent for both the test tube and microtiter assays. Working reagent is prepared by mixing 50 parts BCA solution with 1 part of 4% Cupric Sulfate.

*Note: The volume of BCA working reagent for the test tube assay is based on 1 ml of reagent per reaction. This volume is sufficient if reading the absorbance in a microcuvette (0.7 ml). If the cuvette requires a larger minimum volume, up to 3 ml of reagent may be need per reaction.*

| <b>Table 2 Preparation of BCA working reagent</b> |                        |             |
|---|------------------------|-------------|
|   | <b>For each sample</b> | <b>× 20</b> |
| <b>Test tube assay</b>                            |                        |             |
| BCA Solution                                      | 1 ml                   | 20 ml       |
| 4% Cupric Sulfate                                 | 20 µl                  | 400 µl      |
| <b>Micro-scale assay</b>                          |                        |             |
| BCA Solution                                      | 200 µl                 | 4 ml        |
| 4% Cupric Sulfate                                 | 4 µl                   | 80 µl       |

## Assay protocols

The following procedures describe two assay protocols. The test tube assay requires a larger volume of protein sample (50 µl). However, the effect of interfering substances is minimized because the ratio of BCA working reagent to protein sample is 20:1. The micro-scale assay requires less of the protein sample (25 µl) and offers the convenience of a 96-well plate. However, because the ratio of BCA working reagent to protein sample is 8:1 the effect of interfering substances will be greater.

### Test tube assay

1. Pipette 50 µl each standard or protein sample replicates into labeled test tubes.
2. Add 1.0 ml BCA working reagent. Mix by gentle vortexing.  
*Note: The volume of the reaction may need to be increased to 3.15 ml (3 ml BCA working reagent plus 150 µl standard or protein sample) if the cuvette requires a larger minimum volume for measuring the absorbance.*
3. **Standard assay:** Incubate reactions at 37°C for 30 min or at room temperature for 2–16 h.  
**Enhanced assay:** Incubate reactions at 60°C for 15 min.
4. Allow tubes to cool to room temperature.  
*Note: Color development continues after cooling to room temperature. However, no significant error in protein concentration is observed when all the reactions are read within 10 min.*
5. Add 1 ml water to a clean cuvette and adjust the absorbance reading at 562 nm to zero.
6. Transfer the reactions to clean cuvettes.
7. Measure and record the absorbance (A<sub>562</sub>) of all reactions within 10 min.
8. To obtain the corrected absorbances, subtract the absorbance of the blank standard from the absorbance measurement of all other standard and protein samples.
9. Plot the corrected absorbance versus the known mass of the BSA standards to generate the standard curve.
10. Using the standard curve, interpolate the recorded corrected absorbance reading for the samples assayed which fall within the linear range of the standard curve.
11. Calculate the amount of protein present in the original sample by correcting for the dilution and sample volume.

### Micro-scale assay

1. Pipette 25 µl each standard or protein sample replicate into individual wells of a 96-well plate.
2. Add 200 µl BCA working reagent to each well. Mix on plate shaker for 30 seconds. Cover plate.
3. **Standard assay:** Incubate reactions at 37°C for 30 min or at room temperature for 2–16 h.  
**Enhanced assay:** Incubate reactions at 60°C for 15 min.
4. Cool plate to room temperature.
5. Measure the absorbance at 562 nm on a plate reader.  
*Note: If the plate reader does not have a 562 nm filter, wavelengths from 540–590 nm have been used successfully.*
6. To obtain the corrected absorbances, subtract the absorbance of the blank standard from the absorbance measurement of all other standard and protein samples.
7. Plot the corrected absorbance versus the known mass of the BSA standards to generate the standard curve.  
*Note: If a curve-fitting algorithm associated with the microplate reader is used for preparing the standard curve, a four-parameter (quadratic) or best-fit curve will provide the most accurate results.*
8. Using the standard curve, interpolate the recorded corrected absorbance reading for the samples assayed which fall within the linear range of the standard curve.
9. Calculate the amount of protein present in the original sample by correcting for the dilution and sample volume.

## Troubleshooting

| Symptom   | Possible cause                                   | Solution  |
|---|--|---|
| No color development in any of the reactions  | Sample contains a copper chelating agent         | Dialyze, desalt, or dilute protein sample.<br>Increase copper concentration in BCA working reagent (i.e., use a 50:2 ratio of BCA solution to 4% Cupric Sulfate). |
| All reactions, including blank are dark purple  | Buffer contains a thiol                          | Dialyze or dilute sample.   |
|   | Buffer contains a reducing agent                 | Dialyze or dilute sample.   |
|   | Buffer contains biogenic amines (catecholamines) | Dialyze or dilute sample.   |
| Blank absorbance is OK, but standards and samples absorbance measurements are lower than expected | Absorbance measured at incorrect wavelength      | Measure absorbance at 562 nm. Absorbance may be measured between 540–590 nm, although the sensitivity will be reduced.  |
|   | Sample is in a strong acid or alkaline buffer    | Dialyze, desalt, or dilute sample.  |
| Absorbance measurements of protein samples higher than expected                                   | Sample contains lipids or lipoproteins           | Add SDS to the protein sample (final concentration 2%) to reduce interference from lipids (2).  |
|   | Protein concentration is too high                | Dilute sample.  |

## Substance Compatibility

### Incompatible substances

The following substances have been found to interfere with the BCA Protein Assay Kit at small concentrations and should be avoided as components of the sample buffer. If interfering substances cannot be removed, consider using the Non-Interfering Protein Assay™ Kit (Cat. No. 488250) for protein concentration determination. This table is not comprehensive, and it is not certain whether other substances may interfere with the assay.

|                |                   |            |                  |
|----------------|-------------------|------------|------------------|
| Asorbic acid   | EGTA              | Iron       | Sucrose (impure) |
| Catecholamines | Glycerol (impure) | Lipids     | Tryptophan       |
| Creatinine     | Hydrogen peroxide | Melibiose  | Tyrosine         |
| Cysteine       | Hydrazides        | Phenol Red | Uric acid        |

## Compatible substances

The following table lists substances that are compatible with the BCA Protein Assay Kit and the maximal non-interfering concentration. This table is not comprehensive, and it is not certain whether other substances may interfere with the assay.

| Substance   | Concentration |
|---|---------------|
| ACES, pH 7.8  | 25 mM         |
| Acetone   | 10%           |
| Acetonitrile  | 10%           |
| Ammonium sulfate  | 1.5 M         |
| Aprotinin   | 10 mg/l       |
| Bicine, pH 8.4  | 20 mM         |
| Bis-Tris, pH 6.5  | 33 mM         |
| Borate, pH 8.5  | 50 mM         |
| Brij <sup>®</sup> -35   | 5%            |
| Brij-56, Brij-58  | 1%            |
| BugBuster <sup>®</sup> Protein Extraction Reagent (Cat. No. 70584)  | undiluted     |
| Calcium chloride (in TBS, pH 8.0)                                   | 10 mM         |
| Cesium bicarbonate  | 100 mM        |
| CHAPS   | 5%            |
| CHAPSO  | 5%            |
| CHES, pH 9.0  | 100 mM        |
| Cobalt chloride (in TBS, pH 8.0)                                    | 0.8 mM        |
| CytoBuster <sup>™</sup> Protein Extraction Reagent (Cat. No. 71009) | undiluted     |
| Deoxycholic acid  | 5%            |
| DMF   | 10%           |
| DMSO  | 10%           |
| Dithioerythritol (DTE)  | 1 mM          |
| Dithiothreitol (DTT)  | 0.5 mM        |
| EDTA  | 10 mM         |
| EPPS, pH 8.0  | 100 mM        |
| Ethanol   | 10%           |
| Ferric chloride (in TBS, pH 8.0)                                    | 10 mM         |
| Glucose   | 10 mM         |
| Glycine-HCl, pH 2.8   | 100 mM        |

| Substance  | Concentration |
|--|---------------|
| Glycercol (pure)   | 10%           |
| Guanidine-HCl  | 4 M           |
| HEPES  | 100 mM        |
| Hydrochloric acid  | 100 mM        |
| Imidazole, pH 7.0  | 50 mM         |
| Insect PopCulture <sup>®</sup> Reagent (Cat. No. 71187)    | undiluted     |
| Leupeptin  | 10 mg/l       |
| 2-Mercaptoethanol  | 0.01%         |
| MES, pH 6.1  | 100 mM        |
| Methanol   | 10%           |
| MOPS, pH 7.2   | 100 mM        |
| Nickel chloride (in TBS, pH 8.0)                           | 10 mM         |
| Nonidet P-40 (NP-40)                                       | 5%            |
| Octyl $\beta$ -glucoside                                   | 5%            |
| Octyl $\beta$ -thioglucopyranoside                         | 5%            |
| PIPES, pH 6.8  | 100 mM        |
| PMSF   | 1 mM          |
| PopCulture Reagent (Cat. No. 71092)                        | undiluted     |
| Potassium thiocyanate                                      | 3.0 M         |
| Reportasol <sup>™</sup> Extraction Buffer (Cat. No. 70909) | undiluted     |
| SDS  | 5%            |
| Sodium acetate, pH 4.8                                     | 200 mM        |
| Sodium azide   | 0.2%          |
| Sodium bicarbonate   | 100 mM        |
| Sodium chloride  | 1 M           |
| Sodium citrate, pH 4.8 or pH 6.4                           | 200 mM        |
| Sodium hydroxide   | 100 mM        |
| Sodium phosphate   | 100 mM        |
| Sucrose  | 40%           |

| <b>Substance (continued)</b>               | <b>Concentration</b> |
|--|----------------------|
| TBS (150 mM NaCl, 100 mM Tris-HCl, pH 8.0) | undiluted            |
| Tricine, pH 8.0                            | 25 mM                |
| Triethanolamine, pH 7.8                    | 25 mM                |
| Tris                                       | 250 mM               |
| Tris(hydroxypropyl)phosphine (THP)         | 1 mM                 |
| Triton <sup>®</sup> X-100                  | 5%                   |
| Triton X-114, X-305, X-405                 | 1%                   |
| Tween <sup>®</sup> 20, Tween 60, Tween 80  | 5%                   |
| Zinc chloride (in TBS, pH 8.0)             | 10 mM                |
| Zwittergent <sup>®</sup> 3-14              | 1%                   |

## References

- Smith, P.K., et al. 1985. *Anal. Biochem.* **150**, 76.
- Kessler, R. J. and Fanestil, D. D. 1986. *Anal Biochem.* **159**, 138.
- Wiechelman, K., et al. 1998. *Anal Biochem.* **175**, 231.
- Brown, R., et al. 1989. *Anal Biochem.* **180**, 136.