## **INSTRUCTIONS**

# Micro BCA Protein Assay Kit



23235	0412.6
Number	Description
23235	Micro BCA Protein Assay Kit, sufficient reagents for 480 tube assays or 3200 microplate assays
	Kit Contents:
	Micro BCA Reagent A (MA), 240mL
	Micro BCA Reagent B (MB), 240mL
	Micro BCA Reagent C (MC), 12mL
	<b>Bovine Serum Albumin Standard Ampules, 2mg/mL</b> , 10 × 1mL ampules containing bovine serum albumin (BSA) at 2.0mg/mL in a solution of 0.9% saline and 0.05% sodium azide
	Storage: Upon receipt store product at room temperature. Product shipped at ambient temperature.
	<b>Note</b> : If either Reagent MA or Reagent MB precipitates upon shipping in cold weather or during long- term storage, dissolve precipitates by gently warming and stirring solutions. Discard any reagent that shows discoloration or evidence of microbial contamination.

## **Table of Contents**

Introduction	1
Preparation of Standards and Working Reagent	2
Test Tube Procedure (linear working range of 0.5-20µg/mL)	3
Microplate Procedure (linear working range of 2-40µg/mL)	3
Troubleshooting	3
Additional Information	4
Related Thermo Scientific Products	5
References	5

## Introduction

The Thermo Scientific<sup>TM</sup> Micro BCA<sup>TM</sup> Protein Assay Kit is a detergent-compatible bicinchoninic acid formulation for the colorimetric detection and quantitation of total protein. An adaptation of the Thermo Scientific<sup>TM</sup> BCA Protein Assay Kit (Product No. 23225), the Micro BCA Kit has been optimized for use with dilute protein samples (0.5-20µg/mL). The unique, patented method uses bicinchoninic acid (BCA) as the detection reagent for Cu<sup>+1</sup>, which is formed when Cu<sup>+2</sup> is reduced by protein in an alkaline environment.<sup>1</sup> A purple-colored reaction product is formed by the chelation of two molecules of BCA with one cuprous ion (Cu<sup>+1</sup>). This water-soluble complex exhibits a strong absorbance at 562nm that is linear with increasing protein concentrations.

The macromolecular structure of protein, the number of peptide bonds and the presence of four amino acids (cysteine, cystine, tryptophan and tyrosine) are reported to be responsible for color formation with BCA.<sup>2</sup> Studies with di-, tri- and tetrapeptides suggest that the extent of color formation is caused by more than the mere sum of individual color-producing functional groups.<sup>2</sup>

The Micro BCA Protein Assay Kit uses concentrated reagents and a protocol that utilizes an extended incubation time at an elevated temperature (60°C, Test Tube Procedure only). The result is an extremely sensitive colorimetric protein assay in a test tube or microplate assay format.

**Note:** For peptide sample concentration measurements, use the Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> Quantitative Fluorometric Peptide Assay or the Pierce<sup>TM</sup> Quantitative Colorimetric Peptide Assay (see Related Thermo Scientific Products).



## **Preparation of Standards and Working Reagent**

### A. Preparation of Diluted Albumin (BSA) Standards

Use Table 1 as a guide to prepare a set of protein standards. Dilute the contents of one Albumin (BSA) Standard ampule into several clean vials, preferably using a diluent that is similar to the sample buffer. Each 1mL ampule of 2.0mg/mL Albumin Standard is sufficient to prepare a set of diluted standards such that three replicates of each dilution may be included in the Test Tube Procedure.

Vial	Volume of Diluent	Volume and Source of BSA	Final BSA Concentration
А	4.5mL	0.5mL of Stock	200µg/mL
В	8.0mL	2.0mL of vial A dilution	$40 \mu g/mL$
С	4.0mL	4.0mL of vial B dilution	20µg/mL
D	4.0mL	4.0mL of vial C dilution	$10\mu g/mL$
Е	4.0mL	4.0mL of vial D dilution	5µg/mL
F	4.0mL	4.0mL of vial E dilution	2.5µg/mL
G	4.8mL	3.2mL of vial F dilution	$1 \mu g/mL$
Н	4.0mL	4.0mL of vial G dilution	$0.5 \mu g/mL$
Ι	8.0mL	0	$0\mu g/mL = Blank$

Table 1. Preparation of Diluted Albumin (BSA) Standards

#### B. Preparation of the Micro BCA Working Reagent (WR)

1. Use the following formula to determine the total volume of WR required:

(# standards + # unknowns) × (# replicates) × (volume of WR per sample) = total volume WR required

Example: for the standard Test Tube Procedure with 3 unknowns and 2 replicates of each sample:

 $(9 \text{ standards} + 3 \text{ unknowns}) \times (2 \text{ replicates}) \times (1 \text{ mL}) = 24 \text{ mL WR required (round up to 25 mL)}$ 

**Note:** 1mL of the WR is required for each sample in the Test Tube Procedure, while only  $150\mu$ L of WR is required for each sample in the Microplate Procedure.

Prepare WR by mixing 25 parts of Micro BCA Reagent MA and 24 parts Reagent MB with 1 part of Reagent MC (25:24:1, Reagent MA:MB:MC). For the above example, combine 12.5mL of Reagent MA and 12.0mL Reagent MB with 0.5mL of Reagent MC.

**Note:** When Reagent MC is initially added to Reagents MA and MB, turbidity occurs that quickly disappears upon mixing to yield a clear-green solution. Prepare sufficient volume of WR based on the number of samples to be assayed. The WR is stable for one day when stored in a closed container at room temperature (RT). It is not necessary to protect the solution from light.



## **Procedure Summary (Test Tube Procedure)**



## Test Tube Procedure (linear working range of 0.5-20µg/mL)

- 1. Pipette 1.0mL of each standard and unknown sample replicate into appropriately labeled test tubes.
- 2. Add 1.0mL of the WR to each tube and mix well.
- 3. Cover tubes and incubate at 60°C in a water bath for 1 hour.
- 4. Cool all tubes to room temperature (RT).
- 5. With the spectrophotometer set to 562nm, zero the instrument on a cuvette filled only with water. Subsequently, measure the absorbance of all the samples within 10 minutes.

**Note:** Color development continues even after cooling to RT. However, the rate of development at RT is sufficient low that no significant error is introduced if all absorbance measurements are made within a 10-minute period.

- 6. Subtract the average 562nm absorbance reading of the Blank standard replicates from the 562nm reading of all other individual standard and unknown sample replicates.
- 7. Prepare a standard curve by plotting the average Blank-corrected 562nm reading for each BSA standard vs. its concentration in  $\mu$ g/mL. Use the standard curve to determine the protein concentration of each unknown sample.

## Microplate Procedure (linear working range of 2-40µg/mL)

- 1. Pipette 150µL of each standard or unknown sample replicate into a microplate well (Product No. 15041).
- 2. Add 150µL of the WR to each well and mix plate thoroughly on a plate shaker for 30 seconds.
- 3. Cover plate using Sealing Tape for 96-Well Plates (Product No. 15036) and incubate at 37°C for 2 hours.

**Note:** Limit incubations of microplate to less than or equal to 37°C, otherwise high background and aberrant color development may result. Most polystyrene assay plates deform, leach, and become cloudy at 60°C.

- 4. Cool plate to room temperature (RT).
- 5. Measure the absorbance at or near 562nm on a plate reader.
- 6. Subtract the average 562nm absorbance reading of the Blank standard replicates from the 562nm reading of all other individual standard and unknown sample replicates.
- 7. Prepare a standard curve by plotting the average Blank-corrected 562nm reading for each BSA standard vs. its concentration in  $\mu$ g/mL. Use the standard curve to determine the protein concentration of each unknown sample.

**Note:** If using curve-fitting software, use a best-fit polynomial equation rather than a linear equation for the standard curve. If plotting results by hand, a point-to-point fit is preferable to a linear fit to the standard points.

## Troubleshooting

Problem	Possible Cause	Solution
No color in any tubes	Sample contains a copper chelating agent	Dialyze, desalt, or dilute sample
		Increase copper concentration in working
		reagent (e.g., use more Reagent MC)
Blank absorbance is OK, but	Strong acid or alkaline buffer, alters	Dialyze, desalt, or dilute sample
standards and samples show	working reagent pH	
less color than expected	Color measured at the wrong wavelength	Measure the absorbance at 562nm
Color of samples appears	Protein concentration is too high	Dilute sample
darker than expected	Sample contains lipids or lipoproteins	Add 2% SDS to the sample <sup>3</sup>
All tubes (including blank) are	Buffer contains a reducing agent	Dialyze or dilute sample
dark purple	Buffer contains a thiol	
	Buffer contains biogenic amines	
	(catecholamines)	
Need to measure color at a	Spectrophotometer or plate reader does	Wavelengths between 540nm and 590nm
different wavelength	not have 562nm filter	can be used, but standard curve slope and
		overall assay sensitivity will be decreased.
		See Tech Tip on website

thermofisher.com/techresources



## Additional Information

#### A. Interfering Substances

Certain substances are known to interfere with the Micro BCA Assay including those with reducing potential, chelating agents, and strong acids or bases. Avoid the following substances as components of the sample buffer:

Ascorbic Acid	Hydrogen Peroxide	Iron	Reducing Sugars
Catecholamines	Hydrazides	Lipids	
Cysteine	Impure Glycerol	Phenol Red	Tyrosine
EGTA	Impure Sucrose	Reducing Agents	Uric Acid

Maximum compatible concentrations for many substances in the Test Tube Procedure are listed in Table 2 (see last page). Substances were considered compatible at the indicated concentration if the error in protein concentration estimation caused by the presence of the substance in the sample was less than or equal to 10%. The substances were tested using WR prepared immediately before each experiment. The Blank-corrected 562nm absorbance measurements (for the 10µg/mL BSA standard + substance) were compared to the net 562nm readings of the same standard prepared in 0.9% saline.

#### B. Strategies for Eliminating or Minimizing the Effects of Interfering Substances

The effects of interfering substances in the Micro BCA Protein Assay may be eliminated or overcome by several methods.

- Remove the interfering substance by dialysis or gel filtration (see Related Thermo Scientific Products).
- Dilute the sample until the substance no longer interferes. (This is only effective for relatively concentrated samples.)
- Precipitate proteins with acetone or trichloroacetic acid (TCA).<sup>4</sup>
- Increase the amount of copper in the WR (prepare WR using a greater proportion of Reagent MC; e.g., MA:MB:MC equal to 25:24:2 or 25:24:3), which may eliminate interference by copper chelating agents.
  - **Note:** For greatest accuracy, the protein standards must be treated identically to the sample(s).

#### C. Response Characteristics for Different Proteins

Each of the commonly used total protein assay methods exhibits some degree of varying response toward different proteins (Table 3). These differences relate to amino acid sequence, pI, structure and the presence of certain side chains or prosthetic groups that can dramatically alter the protein's color response. Most protein assay methods use BSA or immunoglobulin (IgG) as the standard against which the concentration of protein in the sample is determined. Pierce Albumin Standard (BSA) (Product No. 23209) provides a consistent standard for protein estimations. Nevertheless, individual proteins, including BSA and IgG, differ slightly in their color responses in the Micro BCA Assay (Figure 1). Therefore, for maximum accuracy use a purified (known concentration) sample of the target protein as the assay standard.



Micro BCA™ Reagent Protein Assay

Table 3. Protein-to-Protein Variation. Absorbance ratios (562nm) for proteins relative to BSA using the Test Tube Procedure.

Ratio = (avg. "test" net Abs.) / (avg. BSA net Abs.)		
Protein Tested	<u>Ratio</u>	
Albumin, bovine serum	1.00	
Aldolase, rabbit muscle	0.80	
$\alpha$ -Chymotrypsinogen, bovine	0.99	
Cytochrome C, horse heart	1.11	
Gamma globulin, bovine	0.95	
IgG, bovine	1.12	
IgG, human	1.03	
IgG, mouse	1.23	
IgG, rabbit	1.12	
IgG, sheep	1.14	
Insulin, bovine pancreas	1.22	
Myoglobin, horse heart	0.92	
Ovalbumin	1.08	
Transferrin, human	0.98	
Average Ratio	1.05	
Standard Deviation	0.12	
Coefficient of Variation	11.4%	

Figure 1. Typical color response curves for BSA and BGG using the Test Tube Procedure.

(815) 968-0747 (815) 968-7316 fax



#### D. Alternative Total Protein Assay Reagents

If interference by a reducing substance or metal-chelating substance contained in the sample cannot be overcome, try the Thermo Scientific<sup>TM</sup> Coomassie Plus<sup>TM</sup> (Bradford) Protein Assay Kit (Product No. 23236), which is less sensitive to such substances.

#### E. Cleaning and Re-using Glassware

Care must be exercised when re-using glassware. The Micro BCA WR is sensitive to metal ions, especially copper ions. All glassware must be cleaned and then given a thorough final rinse with ultrapure water.

## **Related Thermo Scientific Products**

15041	Pierce <sup>TM</sup> 96-Well Plates, 100/pkg
15075	Reagent Reservoirs, 50mL, 200/pkg
15036	Sealing Tape for 96-Well Plates, 100/pkg
23209	Bovine Serum Albumin (BSA) Standard Ampules, 2mg/mL, $10 \times 1mL$
23212	Bovine Gamma Globulin (BGG) Standard Ampules, 2mg/mL, $10 \times 1$ mL
23290	Pierce Quantitative Fluorometric Peptide Assay
23275	Pierce Quantitative Colorimetric Peptide Assay
23236	Coomassie Plus <sup>TM</sup> (Bradford) Protein Assay Kit
89882	Zeba <sup>™</sup> Spin Desalting Columns, 0.5mL
89889	Zeba Spin Desalting Columns, 2mL
69576	Slide-A-Lyzer <sup>TM</sup> MINI Dialysis Units
69576	Slide-A-Lyzer Dialysis Cassettes

#### **Cited References**

- 1. Smith, P.K., et al. (1985). Measurement of protein using bicinchoninic acid. Anal Biochem 150:76-85.
- 2. Wiechelman, *et al.* (1988). Investigation of the bicinchoninic acid protein assay: Identification of the groups responsible for color formation. *Anal Biochem* **175**:231-7.
- 3. Kessler, R. and Fanestil, D. (1986). Interference by lipids in the determination of protein using bicinchoninic acid. Anal Biochem 159:138-42.
- 4. Brown, R., et al. (1989). Protein measurement using bicinchoninic acid: elimination of interfering substances. Anal Biochem 180:136-9.

#### **General References**

Gembitsky, D.S., *et al.* (2004). A prototype antibody microarray platform to monitor changes in protein tyrosine phosphorylation. *Mol Cell Proteomics* **3:**1102-1118.

O'Nuallain, B. and Wetzel, R. (2002). Conformational Abs recognizing a generic amyloid fibril epitope. Proc Natl Acad Sci USA 99:1485-90.

Paratcha, G., et al. (2003). The neural cell adhesion molecule NCAM is an alternative signaling receptor for GDNF family ligands. Cell 113:867-79.

Yang, G., et al. (2005). Activation-induced deaminase cloning, localization, and protein extraction from young Vh-mutant rabbit appendix. Proc Natl Acad Sci USA 102:17083-8.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com/techresources. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2015 Thermo Fisher Scientific Inc. All rights reserved. Triton is a trademark of The Dow Chemical Company. Brij, Tween and Span are trademarks of Croda International Plc. Zwittergent is a trademark of Merck KGaA. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.



### Table 2. Compatible Substance Concentrations in the Micro BCA Protein Assay (see text for details).

Substance	Compatible Concentration	Substance
Salts/Buffers		Deter
ACES, pH 7.8	10mM	Brij™-35
Ammonium sulfate		Brij-56, Brij-58
Bicine, pH 8.4	2mM	CHAPS (CHAPSO)
Bis-Tris, pH 6.5	0.2mM	Deoxycholic acid
Borate (50mM), pH 8.5 (#28384)	1:4 dilution*	Nonidet P-40 (NP-4
B-PER <sup>®</sup> Reagent (#78248)	1:10 dilution*	Octyl β-glucoside
Calcium chloride in TBS, pH 7.2	10mM	Octyl B-thioglucopy
Na-Carbonate/Na-Bicarbonate (0.2 M), pH 9.4 (#28382)	undiluted	SDS Span™ 20
Cesium bicarbonate	100mM	Triton™ X-100
	100mM	Triton X-114
No Citrate $(0.6 \text{ M})$ No Corbonate $(0.1 \text{ M})$	1:600 dilution*	Triton X 205 X 405
$M_{\rm p} = 0.0  (\#28388)$	1.000 ullution	Thion X-303; X-403
	1.600 dilution*	Tween 60
(#28386)	1.600 dilution	
Cobalt oblarida in TPC all 7.2		
Cobait chioride in TBS, pH 7.2		Chelati
EPPS, pH 8.0		EDIA
Ferric chloride in TBS, pH 7.2	0.5mM	EGIA
Glycine	n/a	Sodium citrate, pH
Guanidine•HCl	4M	Reducing & Thio
HEPES, pH 7.5	100mM	N-acetylglucosamin
Imidazole, pH 7.0	12.5mM	Ascorbic acid
MES, pH 6.1	100mM	Cysteine
MES (0.1M), NaCl (0.9%), pH 4.7 (#28390)	1:4 dilution*	Dithioerythritol (DTE
MOPS, pH 7.2	100mM	Dithiothreitol (DTT)
Modified Dulbecco's PBS, pH 7.4 (#28374)	undiluted	Glucose
Nickel chloride in TBS, pH 7.2	0.2mM	2-Mercaptoethanol
PBS; Phosphate (0.1M), NaCl (0.15M), pH 7.2 (#28372)	undiluted	Thimerosal
	100mM	Misc. Reage
PIPES, PH 0.0	1.10 dilution*	Acetone
0.5% DOC 1% NP-40 0.1% SDS pH 8.0	1:10 dilution"	Acetonitrile
		Aprotinin
Sodium acetate, pH 4.8	200mivi	DMF, DMSO
	0.2%	Ethanol
Sodium bicarbonate	100mM	Glycerol (Fresh)
Sodium chloride	1M	Hydrazide
Sodium citrate, pH 4.8 (or pH 6.4)	5mM (16.7mM)	Hydrides (Na <sub>2</sub> BH <sub>4</sub> o
Sodium phosphate	100mM	Hydrochloric Acid
Tricine, pH 8.0	2.5mM	Leupeptin
Triethanolamine, pH 7.8	0.5mM	Methanol
Tris	50mM	Phenol Red
TBS; Tris (25mM), NaCl (0.15M), pH 7.6	1:10 dilution*	PMSF
(#28376)		Sodium Hydroxide
Tris (25mM), Glycine (192mM), pH 8.0	1:10 dilution*	Sucrose
(#28380)		TLCK
Tris (25mM), Glycine (192mM), SDS	undiluted	TPCK
(0.1%), pH 8.3 (#28378)		Urea
Zinc chloride in TBS, pH 7.2	0.5mM	o-Vanadate (sodium

Substance	Compatible Concentration
Detergents**	
Brij™-35	5.0%
Brij-56, Brij-58	1.0%
CHAPS (CHAPSO)	1.0% (5.0%)
Deoxycholic acid	5.0%
Nonidet P-40 (NP-40)	5.0%
Octyl β-glucoside	0.1%
Octyl β-thioglucopyranoside	5.0%
SDS	5.0%
Span™ 20	1.0%
Triton™ X-100	5.0%
Triton X-114	0.05%
Triton X-305, X-405	1.0%
Tween™-20, Tween-80	5.0%
Tween-60	0.5%
Zwittergent™ 3-14	
Chelating agents	
EDTA	0.5mM
EGTA	
Sodium citrate, pH 4.8 (or pH 6.4)	5mM (16.7mM)
Reducing & Thiol-Containing Agents	
N-acetylglucosamine in PBS, pH 7.2	
Ascorbic acid	
Cysteine	
Dithioerythritol (DTE)	
Dithiothreitol (DTT)	
Glucose	1mM
2-Mercaptoethanol	1mM
Thimerosal	
Misc. Reagents & Solvents	
Acetone	1.0%
Acetonitrile	1.0%
Aprotinin	1mg/L
DMF, DMSO	1.0%
Ethanol	1.0%
Glycerol (Fresh)	1.0%
Hydrazide	
Hydrides (Na <sub>2</sub> BH <sub>4</sub> or NaCNBH <sub>3</sub> )	
Hydrochloric Acid	10mM
Leupeptin	10mg/L
Methanol	1.0%
Phenol Red	
PMSF	1mM
Sodium Hydroxide	50mM
Sucrose	4%
TLCK	0.1mg/L
ТРСК	0.1mg/L
Urea	3M
o-Vanadate (sodium salt), in PBS, pH 7.2	1mM
· · · · ·	

\* Diluted with ultrapure water \*\* Detergents were tested using high-purity Thermo Scientific<sup>™</sup> Surfact-Amps<sup>™</sup> Products, which have low peroxide content

--- Dashed-line entry indicates that the material is incompatible with the assay

(815) 968-0747 (815) 968-7316 fax