## **INSTRUCTIONS**



# Pierce Quantitative Colorimetric Peptide Assay, 500 Assays

23275

Number

## **Description**

23275

Pierce Quantitative Colorimetric Peptide Assay, sufficient reagents for 500 microplate assays

**Kit Contents:** 

Colorimetric Peptide Assay Reagent A, 50mLColorimetric Peptide Assay Reagent B,  $2 \times 25mL$ Colorimetric Peptide Assay Reagent C, 2mLPeptide Digest Assay Standard (1mg/mL), 1.5mL

**Storage:** Upon receipt store at 4°C. Product shipped on ice packs.

**Note:** If Colorimetric Peptide Assay Reagent A, B or C precipitates upon shipping in cold weather or during long-term storage, dissolve precipitates by gently warming and stirring solution. Discard any kit reagent that shows discoloration or evidence of microbial contamination.

#### Introduction

The Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> Quantitative Colorimetric Peptide Assay provides modified BCA reagents along with a proprietary chelator and uses the biuret reduction of Cu<sup>+2</sup> to Cu<sup>+1</sup> to analyze peptide mixtures. In the reaction, copper is first reduced by the amide backbone of peptides under alkaline conditions (the biuret reaction), followed by the proprietary chelator coupling with the reduced copper to form a bright red complex with absorbance of 480nm. The signal produced from this reaction is three- to four-fold more sensitive than the Thermo Scientific<sup>TM</sup> Micro BCA<sup>TM</sup> Protein Assay Kit for peptide analysis. The kit also contains a high quality peptide digest reference standard to generate linear standard curves and calibration controls. This colorimetric peptide assay requires a small amount of sample (20µL) and has a working peptide concentration range of 25-1,000µg/mL. The increased sensitivity, low sample assay volume, and included reference standard are essential for accurate, robust measurement of peptide digest samples, specifically for mass spectrometry applications.

#### **Important Product Information**

- The Standard provided with this assay is optimized to produce concentration measurements that are consistent with the average results from tryptic digests of proteins or cell lysates. **This assay is not recommended for individual peptide quantitation.** The peptide hydrophobicity; length; and specific amino acid content, such as the presence of cysteine, tryptophan and tyrosine residues are all known to influence the rate of copper reduction, which in turn affect color formation at 480nm.
- Since the assay is based on copper reduction by amide bonds, full-length protein or partially digested protein will produce a colorimetric assay response that can interfere with results. Ensure complete peptide digestion before analysis. If the sample contains significant amounts of undigested protein, use the Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> Quantitative Fluorometric Peptide Assay (Product No. 23290).
- Reducing and chelating reagents such as DTT and EDTA will interfere with the assay.
- Amine-reactive labeling reagents (e.g., Thermo Scientific<sup>TM</sup> TMT<sup>TM</sup> Isobaric Labeling Reagents or biotinylation reagents) are compatible with this assay.
- This method is not a true end-point method; that is, the copper continues to be reduced and color develops over time. However, the rate of color development is sufficiently slow at room temperature, allowing large numbers of samples to be assayed together with minimal error based on the normal timing of reagent additions.

Thermo Fisher Scientific PO Box 117

(815) 968-0747 or (800) 874-3723

(815) 968-7316 fax

thermoscientific.com/pierce



• This assay is optimized to be performed in a 96-well microplate with a sample volume of  $\leq 20\mu L$  and total assay volume of  $200\mu L$ . If a larger assay volume is desired, e.g., for cuvette-based measurements, use a Sample:Assay Working Reagent ratio of 1:33 (e.g., use  $30\mu L$  of sample and  $970\mu L$  of working reagent to have a 1mL total solution volume). If larger volumes of sample are used, the number of assays per kit will be less than 500.

#### **Additional Material Required**

- Ultrapure water
- Sealing tape (e.g., Sealing Tape for 96-Well Plates, Product No. 15036)
- 96-well microplate (e.g., Thermo Scientific<sup>TM</sup> Pierce<sup>TM</sup> 96-Well Plates, Product No. 15041)
- 96-well microplate reader compatible with absorbance measurement at 480nm (e.g., Thermo Scientific<sup>TM</sup> VarioSkan<sup>TM</sup> Flash Multimode Reader)

#### **Preparation of Standards**

Use the procedure in Table 1 to prepare a dilution series of the Peptide Digest Assay Standard to generate a standard curve. Dilute the Peptide Digest Assay Standard into clean vials, preferably using the same diluent as the sample(s). This will provide sufficient volume to run an 8-point standard curve (from 0-1000µg/mL) in triplicate.

Table 1. Preparation of diluted Peptide Digest Assay Standards.

Centrifuge tubes	<u>Volume of Water</u> (μL)	<u>Volume of Digest</u> (μL)	Final Concentration of Digest (µg/mL)
A	0	150 of Stock	1000
В	75	75 of Vial A	500
C	75	75 of Vial B dilution	250
D	75	75 of Vial C dilution	125
E	75	75 of Vial D dilution	62.5
F	75	75 of Vial E dilution	31.3
G	75	75 of Vial F dilution	15.6
Blank		Use water for blank	0

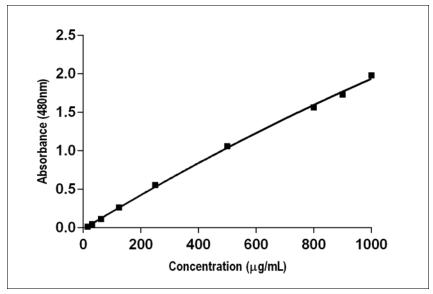


Figure 1. Example standard curve.



## **Preparation of Working Reagent**

1. Use the following formula to determine the total volume of Working Reagent (WR) required:

[(# of standards) + (# of unknowns)]  $\times$  (# of replicates)  $\times$  (0.18mL of WR per sample) = total volume WR required Example: For an 8-point standard curve and 5 unknowns in triplicate:

 $(8 \text{ standards} + 5 \text{ unknowns}) \times (3 \text{ replicates}) \times (0.18\text{mL}) = 7.02\text{mL} \text{ of WR required}$ 

**Note:** For the above example, make at least 7.5mL of WR to ensure sufficient reagent for use.

- 2. Prepare WR by mixing:
  - 50 parts of Colorimetric Peptide Assay Reagent A
  - 48 parts of Colorimetric Peptide Assay Reagent B
  - 2 parts of Colorimetric Peptide Assay Reagent C

For the above example, combine 3.75mL of Peptide Assay Reagent A, 3.6mL of Peptide Assay Reagent B and 0.15mL of Peptide Assay Reagent C.

**Note:** After preparation, the WR can be kept at room temperature for  $\leq 30$  minutes.

#### **Procedure**

- 1. Prepare samples and standards.
- 2. Pipette 20µL of each standard or unknown sample replicate into a well of a 96-well microplate.
- 3. Add 180µL of the WR to each well and mix plate thoroughly on a plate shaker for 30 seconds to 1 minute.
- 4. Cover plate and incubate at 22°C (room temperature) for 30 minutes or 37°C for 15 minutes.

**Note:** Incubation time can be increased to greater than 15 minutes if there is no sufficient color development within 15 minutes at 37°C. Increased incubation time or temperature will increase the assay response; however, prolonged incubation at elevated temperatures may reduce the linearity of the assay response. Follow instrument instructions for determining sample concentrations based on the standard curve. Alternatively, prepare a standard curve by plotting the average blank-corrected 480nm measurement for each standard versus its concentration in µg/mL. Use the standard curve to determine the peptide concentration of each unknown sample based on the average blank-corrected absorbance value of the samples.

**Note:** Assay produces a near-linear response across the recommended concentration ranges. However, if using curve-fitting algorithms associated with a microplate reader, a four-parameter (quadratic) can also be used as an alternative to a linear fit.

## **Troubleshooting**

Problem	Possible Cause	Solution
No color in wells	Sample contained a copper chelating agent	Clean-up sample on a C18 column
		Dilute sample
Blank absorbance measures	Strong acid or alkaline buffer altered WR	Dilute sample
correctly, but standards and	pH	Neutralize pH
samples show less color than expected	Color measured at the wrong wavelength	Measure the absorbance at 480nm
Sample colors appear darker	Peptide concentration was too high	Dilute sample
than expected	Tryptic-digested sample contained undigested protein  Sample contained a reducing agent such as	Confirm digestion efficiency by SDS-PAGE. Optimize digestion protocol Clean-up sample on a C18 column
	DTT or EDTA	The state of the s
Sample colors appear very light,	Peptide sample precipitated	Dissolve the sample in 50% acetonitrile
but standards measure correctly	Peptide bound to tubing	and perform assay at 37°C for
	Peptide had significant secondary structure	30 minutes



## **Compatible Substances**

- Reducing and chelating reagents such as DTT and EDTA will interfere with the assay. Clean samples to remove residual interfering reagents before assaying.
- Other than reducing agents and chelators, the Pierce Quantitative Colorimetric Peptide Assay is compatible with a wide range of standard laboratory chemicals. The maximum concentrations tested and shown to be compatible for many substances are listed in Table 2. Substances were compatible at the indicated concentration if the error in the concentration estimation caused by the presence of the substance was ≤ 15%.

Table 2. Compatible substance concentrations in the Thermo Scientific Pierce Quantitative Colorimetric Peptide Assay.

Substance	<b>Compatible Concentration</b>
Acetone	50%
Acetonitrile	50%
Ammonium bicarbonate	50mM
DMSO	50%
DTT(dithiothreitol)	Not compatible
EDTA	5mM
Formic acid	0.5%
Guanidine	0.25M
Iodoacetamide	1M
Methanol	25%
SDS	1%
Sodium azide	1%
TCEP	Not compatible
TEA acetate	5mM
TEA bicarbonate	5mM
Trifluoroacetic acid	0.5%
Tris	100mM
Urea	1M

**Note:** It is possible to have a substance additive effect even though a single component may be present at a concentration below its listed compatibility. A sample buffer containing a combination of substances could potentially interfere with the assay. Combinations of reagents have not been examined in detail.

#### **Related Thermo Scientific Products**

15041	Pierce 96-Well Plates, 100/pkg
15075	Reagent Reservoirs, 200/pkg
15036	Sealing Tape for 96-Well Plates, 100/pkg
23290	$\textbf{Pierce Quantitative Fluorometric Peptide Assay,} \ 500 \ assays$
28904	Trifluoroacetic acid, $10 \times 1 \text{mL}$
51101	Acetonitrile, 1L
87782	Pierce <sup>TM</sup> C18 Tips, 100μL bed column, 96/pkg
84840	Mass Spec Sample Prep Kit for Cultured Cells, 20-rxn kit



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 <a href="www.thermoscientific.com/pierce">www.thermoscientific.com/pierce</a>. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2014 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.