

# Pierce<sup>®</sup> Iodination Beads

28665 28666 0331.3

Number Description

**Pierce Iodination Beads**, 50 beads, contains sufficient material to perform 10 to 50 reactions using

0.1 to 0.5mg of protein each

**28666** Pierce Iodination Beads, 250 beads, contains sufficient material to perform 25 to 250 reactions using

0.1 to 0.5mg of protein each

Bead diameter: 3.175mm

Oxidative capacity:  $0.55 \pm 0.05 \mu \text{mol/bead}$ 

Storage: Upon receipt store product desiccated at 4°C. Avoid exposing beads to moisture

or reducing agents. Product is shipped at ambient temperature.

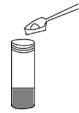
#### Introduction

The Thermo Scientific Pierce Iodination Beads (previously called "IODO-BEADS") consist of the iodination reagent *N*-chloro-benzenesulfonamide (sodium salt) immobilized on nonporous, polystyrene beads. This oxidizing agent was first reported by Markwell<sup>1</sup> as a means for iodinating proteins. Iodination Beads provide for effective iodination of peptides and proteins in 2-15 minutes across a broad pH range.

Radioactive iodine (<sup>131</sup>I or <sup>125</sup>I) can be incorporated into protein either by enzymatic or chemical oxidation. In the chemical oxidation method, Na<sup>125</sup>I or Na<sup>131</sup>I is converted to its corresponding reactive radioactive iodine form. Historically, the oxidizing agent of choice was chloramine-T. However, because of its strong oxidizing properties, chloramine-T may damage proteins and thereby adversely affect their biological function for the intended experiment. Iodination with these beads is milder than chloramine-T, generates sufficient radioactive iodine, and does not require a reduction step, which makes it advantageous for maintaining biological activity of proteins. Furthermore, Iodination Beads allow for quick and complete separation of the reagent from the reaction mixture, making it easy to control the reaction time and to limit direct contact of the oxidant with the protein.

Iodination with Pierce Iodination Beads is compatible with many common buffer components. Detergents such as SDS, NP-40, and Triton® X-100, and denaturants such as urea, and high-salt concentrations (1.0 M NaCl) will not harm the reaction and may improve incorporation of the radioactive iodine by making tyrosines less hindered. Iodinations may be performed in phosphate, Tris, HEPES and other common buffers. The beads are not compatible with reducing agents.

## **Procedure Summary**



**1.** Add beads to an <sup>125</sup>I-buffered solution in the reaction vial. Incubate for 5 minutes at room temperature.



**2.** Add protein in buffer to the reaction vial. Allow reaction to proceed for 2-15 minutes.



**3.** Separate reaction volume from the beads to terminate the iodination reaction.



## **Important Product Information**

**Beads and Protein:** Use at least one bead for each 5-500µg of tyrosine-containing peptide or protein dissolved in 0.1-1mL of Reaction Buffer. Because beads can vary in oxidative capacity, use at least two beads per reaction for best results. The level of protein iodination can be controlled by varying the bead:protein ratio, the time of reaction, and amount of NaI added. If the protein or peptide does not contain tyrosine residues, which are the target of iodination, use SHPP or Sulfo-SHPP (see Related Thermo Scientific Products Section) to add these tyrosyl groups to the molecule.

**Reaction Buffer:** Phosphate-buffered saline (e.g., 0.1M phosphate, 0.15M NaCl; pH 7.2; Product No. 28372) or Tris, pH 5.5-7.5 (pH 6.5 is optimal). Phosphate buffered Saline (PBS) results in higher <sup>125</sup>I incorporation than Tris. Other buffers may be used provided they are free of reducing agents and anti-oxidants. Avoid organic solvents, such as DMSO and DMF, which readily dissolve polystyrene.

**Temperature and time for iodination:** Room temperature is optimal for iodine labeling. However, the reaction may also be performed at 4°C, but with reduced efficiency. Typical reaction time is 2-15 minutes. The optimal reaction time for each protein must be determined empirically.

## **Example Procedure for Protein Iodination**

## A. Additional Materials Required

- Reaction Buffer: see description in Important Product Information above
- Sodium iodide (Na<sup>125</sup>I), carrier-free
- Microcentrifuge tube or glass reaction vial
- Desalting column or dialysis unit to separate labeled protein/peptide from excess unincorporated iodine (see Step 5)

#### B. Procedure

- 1. Immediately before use, wash beads with 500µL of Reaction Buffer per bead. Dry the beads on filter paper. (This wash step removes any loose particles and reagent.)
- 2. Add beads to a solution of carrier-free Na<sup>125</sup>I (approximately 1 mCi per 100μg of protein) diluted with Reaction Buffer and incubate for 5 minutes. A small reaction vial is optimal for this step.
- 3. Dissolve or dilute protein in Reaction Buffer and add to the reaction vessel. Incubate reaction mixture for 2-15 minutes.
  - **Note:** For best results, perform a time study (see Time Study Procedure below) to determine optimal reaction time.
- 4. Stop the reaction by removing the solution from the reaction vessel. Once beads are separated from the solution, the reaction will stop.
  - **Note:** Beads may be washed with a small quantity of Reaction Buffer to recover any protein from the bead surface.
- 5. Thermo Scientific Zeba Spin Desalting Columns (e.g., Product No. 89891) may be used to remove excess Na<sup>125</sup>I or unincorporated <sup>125</sup>I from the iodinated protein (if the protein is larger than 7kDa). Alternatively, use a Thermo Scientific Slide-A-Lyzer Dialysis Cassette (e.g., Product No. 66382) and dialyze the iodinated protein against a buffer suitable for the specific downstream application. For peptides larger than 2kDa, use Slide-A-Lyzer<sup>®</sup> Dialysis Cassette with a 2K MWCO (see web site for complete listing).

## **Example Time Study Procedure**

This is a summary of the method described by Cheng and Rudick.<sup>2</sup>

- 1. Remove  $5\mu L$  of the reaction mixture every 30 seconds, removing the first aliquot when the bead is added (time zero).
- 2. Apply each aliquot to separate squares of nitrocellulose. Support the squares using a non-adsorbent surface.
- 3. Place squares in a vials and count for gamma radiation. The total counts will be approximately the same for each square.
- 4. Prepare Tris-glycine/NaI buffer by titrating 25mM Tris base with 192mM glycine to pH 8.3, add methanol to 20% and sodium iodide to a final concentration of 10mM.
- 5. To remove free  $^{125}$ I, wash squares  $4 \times 5$  minutes with gentle agitation with 50mL of Tris-glycine/NaI buffer.
- 6. Count gamma radiation on the washed squares. Subtract the zero time counts of washed squares from the other counts to determine net radioactivity for each incubation-time aliquot. This provides the radioactivity incorporated into the protein.
- 7. The efficiency of radioiodine incorporation may be measured by dividing the net radioactivity of the washed membranes by the radioactivity determined from the unwashed membranes after sample application.



## **Example Procedure for Cell-Surface Iodination**

#### A. Additional Materials Required

- Phosphate-buffered saline (PBS): 0.1M phosphate, 0.15M NaCl; pH 7.2 (Product No. 28372)
- Na<sup>125</sup>I
- Carrier iodide: 25mM NaI in PBS (optional)

#### B. Procedure

- 1. Wash  $10^5$ - $10^6$  cells 3 times with PBS.
- 2. Prepare Iodination Beads by washing them twice with 0.5mL PBS per bead. Dry the beads on filter paper. (This wash step removes any loose particles and reagent.)

**Note:** For best results, use six beads per  $10^5$ - $10^6$  cells. Specific activity may be controlled by changing the number of beads. Using more than six beads per  $10^5$ - $10^6$  cells will not increase specific activity and may reduce protein recovery.

3. Resuspend cells in 4mL PBS and add  $^{125}$ I (approximately 1mCi per 100 $\mu$ g wet cells), 40 $\mu$ L carrier iodide (optional) and the iodinating beads.

Note: For best results, determine the optimal amount of carrier iodide to add for the specific application.

4. Incubate the samples at room temperature for 2-15 minutes.

**Note:** For best results, perform a time study to determine optimal reaction time. Iodination may be performed at 4°C; however, the efficiency will be reduced.

- 5. Stop the reaction by removing the iodinating beads with forceps.
- 6. Wash cells 3 times with PBS, to remove non-reacted <sup>125</sup>I. Cell-surface proteins may be solubilized and analyzed by SDS-PAGE.

## **Troubleshooting**

Problem	Cause	Solution
Iodination of protein is low	Absence of tyrosines available for labeling	Use Bolton-Hunter Reagent (Product No. 27710) or its water-soluble analog Sulfo-SHPP (Product No. 27712) to modify primary amines on the protein with tyrosine-"like" functional groups
		<b>Note:</b> These reagents may be iodinated first and the protein subsequently reacted with the iodination reagent
	Insufficient surface tyrosines available for labeling	Use high-salt (1.0M NaCl), detergents or denaturants to allow better access to "buried" tyrosines
Protein could not be iodinated	Reducing agents or anti- oxidants are interfering with labeling	Avoid 2-mercaptoethanol, dithiothreitol, sodium borohydride, sodium cyanoborohydride, glycerol or any other compound with reducing properties
Protein is being degraded	The protein is susceptible to oxidation	Allow the bead to oxidize the Na <sup>125</sup> I solution as a separate step from the protein iodination and then transfer the oxidized iodine to the protein solution
		<b>Note:</b> Iodine incorporation is much less efficient than the standard method but may be necessary for easily oxidized proteins
		Perform a time study to determine the optimal reaction time



## **Appendix**

Pierce Iodination Beads are susceptible to inactivation from exposure to reducing agents or moisture. To verify if the beads have sufficient activity, use this simple qualitative test, which was first described by Lee and Griffiths.<sup>3</sup>

- 1. Dissolve 125μg of potassium iodide in 100μL of 0.1M sodium phosphate buffer, pH 7.0.
- 2. Add one unused bead to this solution.

An active bead will turn yellow to brown in less than 15 seconds and the solution will turn yellow a few seconds later. A bead with low activity will require several minutes or longer to produce the color change.

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

27710 Bolton-Hunter Reagent (SHPP), 1g

27712 Water-Soluble Bolton-Hunter Reagent (Sulfo-SHPP), 100mg

**28601** Pierce Iodination Tubes, 10/pkg, glass test tubes coated with Pierce Iodination Reagent

**28600** Pierce Iodination Reagent, 1g

#### **Cited References**

- 1. Markwell, M.A. (1982). A new solid-state reagent to iodinate proteins: conditions for the efficient labeling of antiserum. Anal Biochem 125: 427-32.
- Cheng, Hui and Rudick, M.J. (1991). A membrane blotting method for following the time course of protein radioiodination background references using IODO-BEADS. Anal Biochem 198: 191-193.
- Lee, D.S.C. and Griffiths, B.W. (1984). Comparative studies of IODO-BEADS and chloramine-T methods for the radioiodination of human alphafetoprotein. J Immunol Meth 74(181).

#### **Product References**

Bohnsack, J.F., et al. (1985). Dependence of I-laminin binding to immune complexes in the presence of Clq. Proc Natl Acad Sci USA 82: 3824-8.

Garrison, S., et al. (2003). Surface translocation of pacolus is induced by cell activation and death but is not required for neutrophil migration and function. J Immunol 171: 6795-806.

Holmes, K.L., et al. (1985). Radiolabeling mouse IgG Fc receptor proteins on murine tumor cell lines and spleen cells. *Proc Natl Acad Sci USA* 82:7706-10. Joiner, K.A., et al. (1985). Binding of I-labeled C7 and C9 to E. coli J5 cells. *Proc Natl Acad Sci USA* 82: 4808-12.

Maler, T., Rosenblum, B.B. and Jourdian, G.W. (1985). Locating Phosphomannosyl receptor proteins in 2D-PAGE gels by iodinated protein A. *Proc Natl Acad Sci USA* 82: 8379-83.

Parise, L.V. and Phillips, D.R. (1985). I-fibronogen binding properties of the glycoprotein IIb-IIIa complex. J Biol Chem 260: 10698-707.

Schagat, T.L., et al. (2003). Surfactant protein A differentially regulates peripheral and inflammatory neutrophil chemotaxis. *Am J Physiol Lung Cell Physiol* **284:** L140-7.

Udenfriend, S., et al. (1985). Scintillation proximity radioimmunoassay utilizing I-labeled ligands. Proc Natl Acad Sci USA 82: 8672-6.

Vogel, T., et al. (1985). Lipoprotein receptor binding assays with recombinant I-human apolipoprotein E. Proc Natl Acad Sci USA 82: 8696-700.

Zhang, L., et al. (2004). RNA interference-mediated silencing of the S100A10 gene attenuates plasmin generation and invasiveness of Colo222 colorectal cancer cells. J Biol Chem 279(3): 2053-62.

Triton is a trademark of The Dow Chemical Company.

This product ("Product") is 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") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. 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 the original purchaser of the Product ("Buyer").

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 replacement of or refund for the non-conforming Product(s).

There is no obligation to replace 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.

Current product instructions are available at <a href="https://www.thermoscientific.com/pierce">www.thermoscientific.com/pierce</a>. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2011 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.