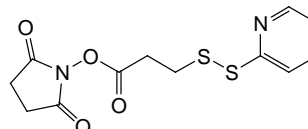


Pierce Premium Grade SPDP

PG82087 PG82088

2543.0

Number	Description
PG82087	Premium Grade SPDP, <i>N</i> -succinimidyl 3-(2-pyridyldithio) propionate, 100mg
PG82088	Premium Grade SPDP, 1g Molecular Weight: 312.37 Spacer Arm Length: 6.8Å



Storage: Upon receipt store product at -20°C protected from moisture. Product is shipped at ambient temperature.

Introduction

Thermo Scientific™ Pierce™ Premium Grade Reagents are high-quality formulations of selected chemical modification reagents, specially characterized for applications where product integrity and risk minimization are critical. Compared to standard grade equivalents, Pierce Premium Grade Reagents provide more clearly defined quality and product support by including: (a) increased analytical testing and product characterization, (b) greater batch-specific information and quality assurance review, (c) extensive lot sample retention, and (d) change control notification.

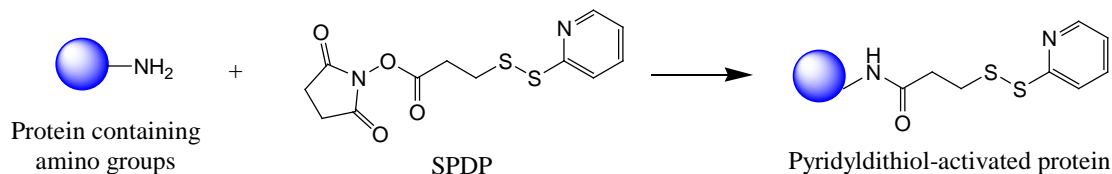
Thermo Scientific™ Pierce™ Premium Grade SPDP Reagent is a unique heterobifunctional crosslinker containing amine- and sulfhydryl-reactive groups. Whether it is used to form amine-to-amine or amine-to-sulfhydryl crosslinks among molecules, Pierce Premium Grade SPDP Reagent produces disulfide-containing linkages that can be cleaved later with reducing agents such as dithiothreitol (DTT).

The amine-reactive portion of SPDP is the *N*-hydroxysuccinimide (NHS) ester. Reactions are most commonly performed in phosphate, carbonate/bicarbonate or borate buffers at pH 7-8. Other buffers can be used provided they do not contain primary amines, thiols or disulfide reducing reagents (see discussion of the sulfhydryl reactivity, below). The rate of reaction and degradation by hydrolysis increases with increasing pH; for example, the half-life of the NHS ester is several hours at pH 7 and less than 10 minutes at pH 9. NHS-ester reagents like SPDP have limited aqueous solubility and must be dissolved in organic solvent before adding them to a reaction mixture. The sulfhydryl-reactive portion of SPDP is the 2-pyridyldithio group, which reacts optimally with sulfhydryls between pH 7 and 8.¹ The reaction results in displacement of a pyridine-2-thione group, the concentration of which can be determined by measuring the absorbance at 343nm (see Additional Information section). Reaction buffers must be free of thiols and disulfide reducing agents until quenching or reduction of the 2-pyridyl disulfide is desired.

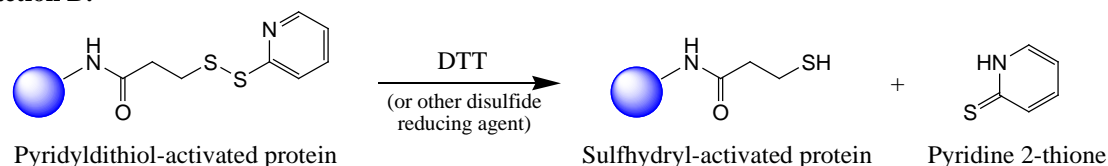
Two basic strategies can be used to form cleavable crosslinks between proteins with SPDP, depending whether one or neither protein already possesses sulfhydryl groups (-SH) in addition to primary amines (Figure 1). Both conjugation methods result in crosslinks that contain a disulfide bond in the spacer arm, which can be cleaved by reduction with dithiothreitol (DTT) or another reducing agent. In most cases, crosslinks created using SPDP reagent can be cleaved with 25mM DTT at pH 4.5 without reducing native protein disulfide bonds. However, when preservation of native disulfide bonds is not a concern, cleavage with DTT can be performed most efficiently at pH 7-9.¹

Crosslinking experiments with SPDP are not limited to those involving proteins. Any of a variety of molecules with primary amines (-NH₂) and sulfhydryl groups can be modified or crosslinked using Pierce Premium Grade SPDP Reagent (see Additional Information section).

Reaction A:



Reaction B:



Reaction C:

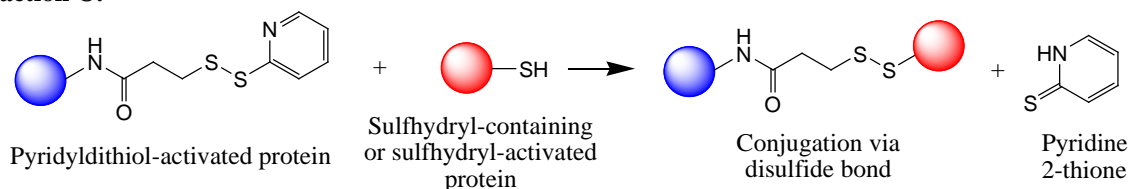


Figure 1. Two protein-to-protein conjugation strategies involving SPDP. If one protein already contains sulfhydryl groups, then only one protein must be modified by SPDP Reagent and only Reactions A and C are used. If neither protein contains available sulfhydryl groups, then both proteins are modified separately with SPDP Reagent (Reaction A), one protein is treated with reducing agent to expose sulfhydryl groups (Reaction B), and finally the two proteins are conjugated (Reaction C).

Procedure for Conjugating Two Proteins, One of Which Contains Sulfhydryls

To conjugate between primary amines on one protein and sulfhydryls (-SH) already present on a second protein, only the first protein must be modified with SPDP.² Then, after desalting the SPDP-modified protein to remove reaction by-products and excess non-reacted reagent, the sulfhydryl-containing protein is added to make the final conjugate (Figure 1, Reactions A and C).

A. Additional Materials Required

- Water-miscible organic solvent such as dimethylsulfoxide (DMSO) or dimethylformamide (DMF)
- Phosphate-buffered saline with EDTA (PBS-EDTA): 100mM sodium phosphate, 150mM NaCl, 1mM EDTA, 0.02% sodium azide, pH 7.5
- Desalting column (e.g., Thermo Scientific™ Zeba™ Spin Desalting Columns, Product No. 89891) sufficient to process 1-2mL samples
- Proteins to be conjugated: In the procedure, IgG and β -galactosidase represent the amine-containing and sulfhydryl-containing proteins, respectively.

B. Procedure (For larger-scale coupling, scale accordingly)

1. Equilibrate the vial of Pierce Premium Grade SPDP Reagent to room temperature before opening.
2. Immediately before use, prepare a 20mM solution of Pierce Premium Grade SPDP Reagent:
 - For example, dissolve 2mg reagent in 320 μ L of DMSO or DMF.
3. Add 25 μ L of the 20mM SPDP solution to 2-5mg IgG dissolved in 1.0mL of PBS-EDTA.
4. Incubate for 30 minutes at room temperature.

5. Equilibrate a desalting column with PBS-EDTA, and buffer exchange the SPDP-modified IgG to remove reaction by-products and excess non-reacted SPDP.
6. Add 1-3mg of β -galactosidase (~1-3 moles β -gal per mole of IgG) to the IgG solution and incubate the reaction mixture overnight at room temperature.
7. The proteins (IgG and β -galactosidase in this example) are now conjugated.

Procedure for Conjugating Proteins, Neither of Which Contains Sulfhydryls

To conjugate between primary amines of both proteins, each protein must be separately modified with SPDP.³ The pyridine-2-thione group is then removed from one of the modified proteins using a reducing agent. Finally, after removing the reducing agent using a desalting column, the resulting sulfhydryl-modified protein and the SPDP-modified protein are incubated together to make the final conjugate (Figure 1, Reactions A, B and C).

A. Additional Materials Required

- Water-miscible organic solvent such as dimethylsulfoxide (DMSO) or dimethylformamide (DMF)
- Phosphate-buffered saline with EDTA (PBS-EDTA): 20mM sodium phosphate, 150mM NaCl, 1mM EDTA, 0.02% sodium azide, pH 7.5. **Note:** If acetate buffer will be used in Section C, do not exceed 20mM sodium phosphate in this buffer. Otherwise 20-100mM sodium phosphate can be used.
- Acetate buffer: 100mM sodium acetate buffer, 100mM NaCl, pH 4.5
- Desalting columns (e.g., Thermo Scientific™ Zeba™ Spin Desalting Columns, Product No. 89891) sufficient to process 1-2mL samples
- Dithiothreitol (DTT), (e.g., Thermo Scientific, Product No. 20290)
- Two proteins to be conjugated

B. SPDP-modify Both Proteins in Separate Reactions

Note: Perform this section of the procedure separately for both proteins.

1. Equilibrate the vial of Pierce Premium Grade SPDP Reagent to room temperature before opening.
2. Immediately before use, prepare a 20mM solution of Pierce Premium Grade SPDP Reagent:
 - For example, dissolve 2mg reagent in 320 μ L of DMSO or DMF.
3. Add 25 μ L of the 20mM SPDP solution to 2-5mg protein dissolved in 1.0mL of PBS-EDTA.
4. Incubate for 30-60 minutes at room temperature.
5. Equilibrate two desalting columns with PBS-EDTA and buffer exchange the SPDP-modified proteins to remove reaction by-products and excess non-reacted SPDP Reagent. Alternatively, the samples can be dialyzed using a Thermo Scientific™ Slide-A-Lyzer™ Dialysis Cassette and concentrated using Slide-A-Lyzer Concentrating Solution.
6. At this point, both proteins are SPDP-modified and one of them must be reduced (Section C) before conjugation.

C. Reduce SPDP Disulfides in One Protein

1. Choose only one of the two SPDP-modified proteins for this section of the procedure. Use the protein whose function or activity is least dependent on native disulfide bonds.
2. Dissolve 23mg DTT in 1mL acetate buffer or PBS-EDTA (makes 150mM solution). Use acetate buffer to avoid reducing native disulfide bonds in the protein.
3. Add 0.5mL DTT solution per 1mL of SPDP-modified protein (results in 50mM DTT).
4. Incubate for 30 minutes.
5. Equilibrate a desalting column with PBS-EDTA and desalt the protein (now sulfhydryl-modified) to remove the DTT.

D. Conjugate the Activated Proteins

1. Mix the SPDP-modified and sulfhydryl-modified proteins and incubate for 18 hours at room temperature or 4°C.
2. To isolate conjugate from the two unconjugated proteins, use an appropriate size-exclusion chromatography method.

Additional Information

A. Please visit our website for additional information on this product including the following items:

- Tech Tip #1: Attach a protein onto glass, silica or quartz surface using a cleavable crosslinker
- Tech Tip #26: Prepare a reversible and measureable sulfhydryl-reactive affinity column

B. Protocol For Pyridine-2-Thione Assay to Determine Level of SPDP-modification

1. Dilute 100µL of SPDP-modified and desalted protein to 1mL with PBS.
2. Measure and record the absorbance at 343nm of the protein sample compared to PBS-EDTA blank (test in triplicate).
3. Add 10µL of 15mg/mL DTT to the 1mL SPDP-modified protein sample and mix.
4. After exactly 15 minutes, measure and record the absorbance at 343nm of the reduced sample.
5. Calculate the change in absorbance: $\Delta A_{343} = (\text{Ave. } A_{343} \text{ after DTT}) - (\text{Ave. } A_{343} \text{ before DTT})$
6. Calculate molar ratio of SPDP to protein using the following equation:

$$\frac{\Delta A}{8080} \times \frac{\text{MW of Protein}}{\text{mg/mL of Protein}} = \text{moles of SPDP per mole of Protein}$$

Where the value 8080 reflects the extinction coefficient for pyridine-2-thione at 343 nm: $8.08 \times 10^3 \text{ M}^{-1}\text{cm}^{-1}$.

Related Thermo Scientific Products

21650	Pierce Sulfo-LC-SPDP , sulfosuccinimidyl 6-[3'-(2-pyridyldithio)-propionamido] hexanoate, 50mg
21651	Pierce LC-SPDP , succinimidyl 6-[3'-(2-pyridyldithio)-propionamido] hexanoate, 50mg
21857	Pierce SPDP , <i>N</i> -succinimidyl 3-(2-pyridyldithio) propionate, 50mg
89889	Zeba Spin Desalting Columns , 2mL
89891	Zeba Spin Desalting Columns , 5mL
89893	Zeba Spin Desalting Columns , 10mL
20290	DTT (Dithiothreitol) , 5g
85190	Dimethylsulfoxide (DMSO) , 50mL
20673	Dimethylformamide (DMF) , 50mL

Cited References

1. Carlsson, J., *et al.* (1978). Protein thiolation and reversible protein-protein conjugation. *Biochem J* **173**:723-7.
2. Cumber, A. J., *et al.* (1985). Preparation of antibody-toxin conjugates. *Meth Enzymol* **112**:207-25.
3. Neurath, A. R. and Strick, N. (1981). Enzyme-linked fluorescence immunoassays using β -galactosidase and antibodies covalently bound to polystyrene plates. *J Virol Meth* **3**:155-65.

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

© 2013 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.