

DSS and BS³ Crosslinkers

MAN0011240 Rev. B.0 Pub. Part No. 2160418

21555 21580 A39266 21586 21655 A39267

Formula: $C_{16}H_{18}N_2Na_2O_{14}S_2$

Number	Description	
21555	DSS (disuccinimidyl suberate), 1g	
21655	DSS , 50mg	~ 0 0
A39267	DSS, No-Weigh Format, $10 \times 2mg$	NO N
	Molecular Weight: 368.35 Spacer Arm: 11.4Å Formula: C ₁₆ H ₂₀ N ₂ O ₈	
21586	BS ³ (bis[sulfosuccinimidyl] suberate), 1g	
21580	BS ³ , 50mg	Na ⁺ O O
A39266	BS ³ , No-Weigh Format, $10 \times 2mg$	0-\$ N 0 N 0 0
	Molecular Weight: 572.43 Spacer Arm: 11.4Å	O O Na ⁺

Storage: Upon receipt store product desiccated at 4°C. Products are shipped at ambient temperature.

Note: Product labels have been provided for your convenience. Please label the vials using one of the labels provided in the Al foil pouch to avoid any confusion as you work with this No-Weigh reagent.

Introduction

Thermo ScientificTM DSS is a water-insoluble, homobifunctional *N*-hydroxysuccinimide ester (NHS ester), and BS³ is its water-soluble analog. NHS esters react efficiently with primary amino groups (-NH₂) in pH 7 to 9 buffers to form stable amide bonds. The reaction results in the release of *N*-hydroxysuccinimide. Proteins, including antibodies, generally have several primary amines in the side chain of lysine (K) residues and the N-terminus of each polypeptide that are available as targets for NHS-ester reagents.

The water-soluble and insoluble forms of NHS-esters have essentially identical reactivity toward primary amines. BS³ is supplied as a sodium salt and is water-soluble up to 100mM. DSS is hydrophobic and dissolved in an organic solvent such as DMSO or DMF then added to the aqueous reaction mixture. DSS does not possess a charged group and is lipophilic and membrane-permeable, which makes it useful for intracellular and intramembrane conjugations. Water-soluble BS³ possess a charged group and is useful for cell-surface protein crosslinking.

Important Product Information

- DSS and BS³ are moisture-sensitive. To avoid moisture condensation onto the product, vial must be equilibrated to room temperature before opening.
- Prepare these crosslinkers immediately before use. The NHS-ester moiety readily hydrolyzes and becomes non-reactive; therefore, do not prepare stock solutions for storage. Discard any unused reconstituted crosslinker.
- Thermo ScientificTM No-WeighTM tube handling: Each of the 10 microtubes contains 2mg of crosslinker. Uncap the vial immediately before use and add DMSO or DMF to the DSS or aqueous buffer to the BS³. Store the unused vials in the foil pouch provided. The maximum useable volume of the vial is 800μL.



- Crosslinking proteins with biological activity (i.e., enzymes, antibodies etc.) can result in activity loss upon conjugation possibly caused by conformational changes of the molecule. Activity loss also may occur when the crosslinker modifies lysine groups involved in binding substrate or antigen. Adjusting the molar ratios of reagent to the target may overcome activity loss. Alternatively, use a crosslinker that targets a different functional group.
- Hydrolysis of the NHS ester is a competing reaction and increases with increasing pH. Hydrolysis occurs more readily in dilute protein or peptide solutions. In concentrated protein solutions, the acylation reaction is favored.

Procedure for Crosslinking Proteins

The following protocol is an example application for this product. Specific applications will require optimization.

A. Materials Required

- Dry dimethyl sulfoxide (DMSO) or dimethyl formamide (DMF) for use with DSS
- Conjugation Buffer: Use a non-amine-containing buffer at pH 7 to 9, such as 100mM sodium phosphate, 0.15M NaCl (Product No. 28372); 20mM HEPES; 100mM carbonate/bicarbonate; or 50mM borate
- Quenching Buffer: 1M Tris•HCl, pH 7.5 (1M glycine or lysine also may be used) non-reacted reagent can be removed by dialysis or gel filtration.

B. Procedure

- 1. Prepare protein in Conjugation Buffer.
- 2. Prepare crosslinker immediately before use. Dissolve BS³ first in water or 20mM sodium phosphate buffer, as more concentrated buffer salt may interfere with initial solubility of the reagent; once the BS³ is dissolved, the solution can be diluted or added to more concentrated buffer solutions without adversely affecting its solubility. Prepare DSS by dissolving in DMSO or DMF. Examples for preparations are as follows:
 - **No-Weigh:** Each No-Weigh tube contains 2mg of crosslinker.
 - **Powder:** Weigh 2mg of crosslinker into a microcentrifuge tube.

Solvent volume to add to 2mg DSS	Buffer volume to add to 2mg BS ³	Crosslinker Concentration
432μL*	277μL*	12.5mM
216μL	140μL	25mM
108μL	70μL	50mM
54μL	35μL	100mM

^{*}The maximum volume that can be added to the No-Weigh vial is $800\mu L. \label{eq:local_local_problem}$

- 3. Add crosslinker to the protein sample. If the protein concentration is greater than 5mg/mL, use a 10-fold molar excess of the crosslinker. For samples < 5mg/mL, use a 20- to 50-fold molar excess of the crosslinker. Use a final concentration of crosslinker at 0.25-5mM.
- 4. Incubate the reaction mixture at room temperature for 30 minutes or on ice for 2 hours.
- 5. Quench the reaction using by adding Quenching Buffer to a final concentration of 20-50mM Tris. Alternatively, remove the non-reacted reagent by dialysis or desalting.
- 6. Incubate the quenching reaction at room temperature for 15 minutes.

Procedure for Intra- and Extracellular Crosslinking

Crosslinking may be performed on cells in suspension or on adherent cells in culture plates. In the latter situation, diffusion of the crosslinking reagent to all surfaces of the cells will be limited and will occur predominately on the exposed surface. Culture media must be washed from the cells otherwise amine-containing components will quench the reaction. Using a more concentrated cell suspension is most effective as less reagent will be required in the reaction. Generally, a final concentration of 1-5mM reagent is effective. NHS reactions occur more rapidly with increasing pH; therefore, pH 8.0 is used in the following example so the reaction can be completed quickly.

Note: Use membrane-insoluble BS³ for crosslinking molecules on the cell surface. Use DSS when crosslinking within the cell is required.



A. Materials Required

- Crosslinker Solution: Immediately before use, dissolve the DSS in dry DMSO at 10-25mM. Dissolve BS³ in water or buffer. BS³ may be added directly to the cells to decrease the extent of hydrolysis.
- Phosphate-buffered saline (PBS): 20mM sodium phosphate, 0.15M NaCl; pH 8. HEPES, bicarbonate/carbonate or borate buffers between pH 7 and 9 may be used as alternative buffers.
- Quench Solution: 1M Tris, pH 7.5 (Tris or glycine will quench the reaction.)

B. Procedure

- 1. Suspend cells at $\sim 25 \times 10^6$ cells/mL in PBS (pH 8.0).
- 2. Wash cells three times with ice-cold PBS (pH 8.0) to remove amine-containing culture media and proteins from the cells.

Note: For cell-surface interaction studies, add ligands to the cells and incubate for 1 hour at 4°C.

- 3. Add the DSS or BS³ solution to a final concentration of 1-5mM.
- 4. Incubate the reaction mixture for 30 minutes at room temperature.

Note: Performing this incubation at 4°C may reduce active internalization of BS³.

- 5. Add the Quench Solution to a final concentration of 10-20mM Tris.
- 6. Incubate the quenching reaction for 15 minutes at room temperature.

Related Thermo Scientific Products

26166	Bioconjugate Techniques, softcover	
66382	Slide-A-Lyzer TM Dialysis Cassette Kit, 10K MWCO, 3mL, 10 cassette kit	
66807	Slide-A-Lyzer Dialysis Cassette Kit, 10K MWCO, 12mL, 8 cassette kit	
22585	DSP (dithiobis[succinimidylpropionate]), 1g	
21578	DTSSP (3,3'-dithiobis[sulfosuccinimidylpropionate]), 50mg	
28372	BupH™ Phosphate Buffered Saline Packs, 40 pack	

General References

Cox, G.W., et al. (1990). Characterization of IL-2 receptor expression and function on murine macrophages. J Immunol 145:1719-26.

Knoller, S., et al. (1991). The membrane-associated component of the amphiphile-activated, cytosol-dependent superoxide-forming NADPH oxidase of macrophages is identical to cytochrome b559. J Biol Chem 266:2795-2804.

Partis, M.D., et al. (1983). Cross-linking of protein by ω-maleimido alkanoyl N-hydroxysuccinimido esters. J Prot Chem 2(3):263-77.

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

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. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2018 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.