

MBS and Sulfo-MBS

22311 22312

0438.2

Number Description

MBS (*m*-maleimidobenzoyl-*N*-hydoxysuccinimide ester), 50mg

Molecular Weight: 314.25 Spacer Arm Length: 7.3Å

CAS #: 58626-38-3

Storage: Upon receipt store desiccated at 4°C. Product is shipped at ambient temperature.

Sulfo-MBS (*m*-maleimidobenzoyl-*N*-hydoxysuccinimide ester), 50mg

Molecular Weight: 416.30 Spacer Arm Length: 7.3Å

Storage: Upon receipt store desiccated at -20° C. Product is shipped at ambient temperature.

Introduction

Thermo Scientific MBS and its water-soluble analog Sulfo-MBS are heterobifunctional crosslinkers that contain *N*-hydroxysuccinimide (NHS) ester and maleimide groups that allow covalent conjugation of amine- and sulfhydryl-containing molecules. NHS esters react with primary amines at pH 7-9 to form amide bonds, while maleimides react with sulfhydryl groups at pH 6.5-7.5 to form stable thioether bonds. In aqueous solutions, hydrolytic degradation of the NHS ester is a competing reaction whose rate increases with pH. The maleimide group is more stable than the NHS-ester group but will slowly hydrolyze and also lose its reaction specificity for sulfhydryls at pH values > 7.5. For these reasons, conjugation experiments involving these crosslinkers are usually performed at pH 7.2-7.5, with the NHS-ester (amine-targeted) reaction being accomplished before or simultaneously with the maleimide (sulfhydryl-targeted) reaction.

MBS and Sulfo-MBS are often used to prepare antibody-enzyme and hapten-carrier protein conjugates in a two-step reaction scheme. First, the amine-containing protein is reacted with a several-fold molar excess of the cross-linker, followed by removal of excess (nonreacted) reagent by desalting or dialysis; finally, the sulfhydryl-containing molecule is added to react with the maleimide groups already attached to the first protein.

Sulfo-MBS is soluble in water and many other aqueous buffers to approximately 10mM, although solubility decreases with increasing salt concentration. MBS is not directly water-soluble and must be dissolved first in an organic solvent such as dimethylsulfoxide (DMSO) or dimethylformamide (DMF); subsequent dilution into aqueous reaction buffer is generally possible, and most protein reactants will remain soluble if the final concentration of organic solvent is less than 10%.

Important Product Information

- MBS and Sulfo-MBS crosslinkers are moisture-sensitive. Store vial of reagent in desiccant at the specified temperature.
 Equilibrate vial to room temperature before opening to avoid moisture condensation inside the container. Dissolve needed amount of reagent and use it immediately before hydrolysis occurs. Discard any unused reconstituted reagent. Do not attempt to make and store stock solutions.
- Avoid buffers containing primary amines (e.g., Tris or glycine) and sulfhydryls during conjugation because they will
 compete with the intended reaction. If necessary, dialyze or desalt samples into an appropriate buffer such as phosphate
 buffered saline (PBS).



• Molecules to be reacted with the maleimide moiety must have free (reduced) sulfhydryls. Reduce peptide disulfide bonds with Thermo Scientific Immobilized TCEP Disulfide Reducing Gel (Product No. 77712). Reduce disulfide bonds in high molecular weight proteins using 5mM TCEP (1:100 dilution of Thermo Scientific Bond-Breaker TCEP Solution, Product No. 77720) for 30 minutes at room temperature, followed by two passes through an appropriate desalting column (e.g., Thermo Scientific Zeba Spin Desalting Columns). Be aware that proteins (e.g., antibodies) may be inactivated by complete reduction of disulfide bonds they contain. Selective reduction of hinge-region disulfide bonds in IgG may be accomplished with 2-Mercaptoethylamine•HCl (2-MEA, Product No. 20408). Sulfhydryls may be added to molecules using *N*-succinimidyl *S*-acetylthioacetate (SATA, Product No. 26102) or 2-iminothiolane•HCl (Traut's Reagent, Product No. 26101), which modify primary amines.

Procedure for Two-step Protein Crosslinking

Generally, a 10- to 50-fold molar excess of crosslinker over the amount of amine-containing protein results in sufficient maleimide activation to enable several sulfhydryl-containing proteins to be conjugated to each amine-containing protein. More dilute protein solutions require greater fold molar excess of reagent to achieve the same level of activation. Empirical testing is necessary to determine activation levels and final conjugation ratios that are optimal for the intended application.

A. Material Preparation

- Conjugation Buffer: Phosphate buffered saline (PBS, pH 7.2; e.g., Product No. 28372) or other amine- and sulfhydryl-free buffer at pH 6.5-7.5 (see Important Product Information) adding EDTA to 1-5mM helps to chelate divalent metals, thereby preventing disulfide formation in the sulfhydryl-containing protein
- Desalting column to separate modified protein from excess cross-linker and reaction byproducts (e.g., ZebaTM Spin Desalting Columns)
- Amine-containing protein (Protein-NH₂) and sulfhydryl-containing protein (Protein-SH) to be conjugated

B. Protocol

Note: For best results, ensure that Protein-SH is prepared (see Important Product Information) and ready to combine with Protein-NH₂ in step 5.

- 1. Dissolve Protein-NH₂ in Conjugation Buffer at 0.1mM (e.g., 5mg in 1mL for a 50kDa protein).
- 2. Add cross-linker to dissolved Protein-NH₂ at 1 mM final (= 10-fold molar excess):
 - For Sulfo-MBS, add 0.416mg per milliliter of Protein-NH₂ solution or dissolve 4.16mg Sulfo-MBS in 1mL Conjugation Buffer (makes 10mM temporary stock) and immediately add 100μL/mL of Protein-NH₂ solution.
 - For MBS, dissolve 3.14mg MBS in 1mL DMSO (makes 10mM) and then add 100µL/mL of Protein-NH₂ solution.
- 3. Incubate reaction mixture for 30 minutes at room temperature or 2 hours at 4°C.
- 4. Remove excess cross-linker using a desalting column equilibrated with Conjugation Buffer.

Note: Follow the desalting column product instructions to determine which fractions contain Protein-NH₂. Alternatively, locate the protein by measuring for fractions having peak absorbance at 280nm; however, be aware that the NHS-ester leaving group also absorbs strongly at 280nm.

- 5. Combine and mix Protein-SH and desalted Protein-NH₂ in a molar ratio corresponding to that desired for the final conjugate and consistent with the relative number of sulfhydryl and activated amines that exist on the two proteins.
- 6. Incubate the reaction mixture at room temperature for 30 minutes or 2 hours at 4°C.

Note: Generally, there is no harm in allowing the reaction to proceed for several hours or overnight; however, the reaction will normally be complete in the specified time. To stop the conjugation reaction before completion, add buffer containing reduced cysteine at a concentration several times greater than the sulfhydryls of Protein-SH.

Note: Conjugation efficiency may be estimated by electrophoresis separation and subsequent protein staining.



Related Thermo Scientific Products

Table 1. Noncleavable NHS/Maleimide crosslinkers.

Cross-linker Name	Spacer Arm Length (Å)	Spacer Arm Composition (between ester and maleimide)	Product No. (NHS)	Product No. (Sulfo-NHS)
AMAS	4.4	Alkane	22295	NA
BMPS	5.9	Alkane	22298	NA
GMBS	7.3	Alkane	22309	22324
MBS	7.3	Aromatic	22311	22312
SMCC	8.3	Cyclohexane	22360	22322
EMCS	9.4	Alkane	22308	22307
SMPB	11.6	Alkane/Aromatic	22416	22317
SMPH	14.2	Alkane/Amide	22363	NA
LC-SMCC	16.2	Alkane/Amide/Cyclohexane	22362	NA
KMUS	16.3	Alkane	NA	21111

Product References

- 1. Kitagawa, T. and Aikawa, T. (1976). Enzyme coupled immunoassay of insulin using a novel coupling reagent. J. Biochem. 79: 233-6.
- 2. Kitagawa, T., et al. (1978). Enzyme immunoassay of viomycin. J Biochem 83: 1493-1501.
- 3. Liu, F-T., et al. (1979). New procedures for preparation and isolation of conjugates of proteins and a synthetic copolymer of D-amino acids and immunochemical characterization of such conjugates. Biochemistry 18: 690-7.
- O'Sullivan, M.J., et al. (1979). Comparison of two methods of preparing enzyme-antibody conjugates: application of these conjugates for enzyme immunoassay. Anal Biochem 100: 100-8.
- 5. Youle, R.J. and Nevelle, Jr., D.M. (1980). Anti-Thy 1.2 monoclonal antibody linked to ricin is a potent cell-type-specific toxin. *Proc Natl Acad Sci USA* 77(9): 5483-6.
- Lerner, R.A., et al. (1981). Chemically synthesized peptides predicted from the nucleotide sequence of the hepatitis B virus genome elicit antibodies reactive with the native envelope protein of Dane particles. Proc Natl Acad Sci USA 78(6): 3403-7.
- Tateishi, K., et al. (1981). A novel immunization procedure for production of anti-cholecystokinin specific antiserum of low cross-reactivity. J Immunol Meth 47: 249-58.
- 8. Hasimura, E., et al. (1982) Production of rabbit antibody specific for amino-terminal residues of cholecystokinin octapeptide (CCK-8) by selective suppression fo cross-reactive antibody response. J Immunol Meth 55: 375-87.
- 9. T. Kitagawa, T. Kawasaki, and H. Munechika (1982). Enzyme immunoassay of blasticidin S with high sensitivity: a new and convenient method for preparation of immunogenic (hapten-protein) conjugates. *J Biochem (Tokyo)* **92:** 585-90.
- 10. Freytag, J.W., Dickinson, J.C. and Tseng, S.Y. (1984). A highly sensitive affinity-column-mediated immunometric assay, as exemplified by digoxin. *Clin Chem* **30(3)**: 417-20.
- 11. Niman, H.L., et al. (1985). Anti-peptide antibodies detect oncogene-related proteins in urine. Proc Natl Acad Sci USA 82: 7924-8.
- 12. Dell'Arciprete, L, et al. (1988). A C terminus cysteine of diptheria toxin B chain involved in immunotoxin cell penetration and cytotoxicity. J Immunol 140: 2466-71.
- 13. Aithal, H.N., *et al.* (1988). An alternate method utilizing small quantities of ligand for affinity purification of monospecific antibodies. *J Immunol Meth* **112:** 63-70.
- 14. Chamberlain, N.R., et al. (1989). Acylation of the 47-kilodalton major membrane immunogen of Treponema pallidum determines its hyrophobicity. *Infec Immunity* **57(9)**: 2878-85.
- Edwards, R.J., et al. (1989). Cross-reaction of antibodies to coupling groups used in the production of anti-peptide antibodies. J Immunol Meth 117: 215-20.
- 16. Peeters, J.M., et al. (1989). Comparison of four bifunctional reagents for coupling peptides to proteins and the effect of the three moieties on the immunogenicity of the conjugates. J Immunol Meth 120: 133-43.
- 17. Miller, M.D., et al. (1989). A novel polypeptide secreted by activated human T lymphocytes. J Immunol 143(9): 2907-16.
- 18. Swanson, S.J., *et al.* (1991). A synthetic peptide corresponding to the phosphorylcholine (PC)-binding region of human C-reactive protein possesses the TEPC-15 myeloma PC-idiotype. *J Immunol* **146(5)**: 1596-1601.
- 19. Myers, D.E., et al. (1989). The effects of aromatic and aliphatic maleimide crosslinkers on anti-CD5 ricin immunotoxins. J Immunol Meth 121:129-42.
- 20. Shoji, H., et al. (1992). Cross-linking of proteins in acetylcholine receptor-rich membranes from *Torpedo californica*: relation of 43-kD protein and *Torpedo* dystrophin to acetylcholine receptor. *Biochemistry International* **28(6)**: 1071-77.
- 21. Hicks, G.R. and Raikhel, N.V. (1995). Nuclear localization signal binding proteins in higher plant nuclei. Proc Natl Acad Sci USA 92:734-8.
- 22. Capala, J., et al. (1996) Boronated epidermal growth factor as a potential targeting agent for boron neutron capture therapy of brain tumors. Bioconjugate Chemistry 7: 7-15.
- 23. Lagriffoul, A., et al. (1996). Secretion of protease nexin-1 by C6 glioma cells is under the control of a heterotrimeric G protein G_{o1}. J Biol Chem **271(49)**: 31508-16.
- 24. Zhang, S., et al. (1996). Augmenting the immunogenicity of synthetic MUC1 peptide vaccines in mice. Cancer Res 56: 3315-9.



- 25. Laird, V. and High, S. (1997). Discrete cross-linking products identified during membrane protein synthesis. J Biol Chem 272(3): 1983-9.
- 26. Sabbatini, A.R.M., *et al.* (1999). Presence in human skeletal muscle of an AMP deaminase-associated protein that reacts with an antibody to human plasma histidine-proline-rich glycoprotein. *J Histochem Cytochem* **47:** 255-60.
- 27. Adessi, C., et al. (2000). Solid phase DNA amplification: characterization of primer attachment and amplification mechanisms. *Nucl Acid Res* **28(20)**:e87.
- 28. Chen, Z., et al. (2003). Spatial and dynamic interactions between phospholamban and the canine cardiac Ca²⁺ pump revealed with use of heterobifunctional cross-linking agents. J Biol Chem 278(48): 48348-56.
- 29. Keding, S.J. and Danishefsky, S.J. (2004). Prospects for total synthesis: a vision for a totally synthetic vaccine targeting epithelial tumors. *Proc Natl Acad Sci USA* **101:** 11937-42.
- 30. Ostrowski, M., et al. (2002). Identification of neutralizing and nonneutralizing epitopes in the porcine reproductive and repiratory syndrome virus GP5 ectodomain. J Virol 76(9): 4241-50.
- 31. Hofmann, S., et al. (2003). Wolfram syndrome: structural and functional analyses of mutant and wild-type wolframin, the WFS1 gene product. Hum Mol Genet 12(16): 2003-12.
- 32. Heit, A., et al. (2003). Cutting edge: toll-like receptor 9 expression is not required for CpG DNA-aided cross-presentation of DNA-conjugated antigen but essential for cross-priming of CD8 T cells. *J Immunol* 170: 2802-5.
- 33. Heit, A., et al. (2004). CpG-DNA Aided cross-priming by cross-presenting B cells. J Immunol 172: 1501-7.
- 34. Wick, M.J., et al. (2003). Mouse 3-phosphoinositide-dependent protein kinase-1 undergoes dimerization and trans-phosphorylation in the activation loop. J. Biol. Chem. 278(44): 42913-9
- 35. Swiderska, A., et al. (2001). Inhibition of the Agrobacterium tumefaciens TraR Quorum-sensing regulator. J Biol Chem 276(52): 49449-58.
- 36. Mullick, J., *et al.* (2001). Physical interaction and functional synergy between glucocorticoid receptor and Ets2 proteins for transcription activation of the rat cytochrome P-450c27 promoter. *J Biol Chem* **276(21)**: 18007-17.
- 37. Dong, M., et al. (2004). Spatial approximation between the amino terminus of a peptide agonist and the top of the sixth transmembrane segment of the secretin receptor. *J Biol Chem* **279(4)**: 2894-903.

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 www.thermoscientific.com/pierce. 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.