

Coupling of Oligosaccharides to Thermo Scientific Nunc CovaLink NH Modules

Key Words

Thermo Scientific™ Nunc™ CovaLink™ NH Modules, solid phase, ELISA, activation with NHS/EDC, carbohydrates, oligosaccharides, binding of activated lactose.

Goal

The goal of this application note is to describe the use of carbohydrates covalently immobilized and detected on the Nunc CovaLink NH Modules. This is exemplified by the Erythrina corallodendron – lactose interaction 5.

The interest in carbohydrates has increased. One of the main reasons for this is due to the results obtained through research in cell communication, e.g. between leucocytes and in bacterial adhesion.

From that point of view carbohydrates are remarkable information molecules compared to peptides and nucleotides.

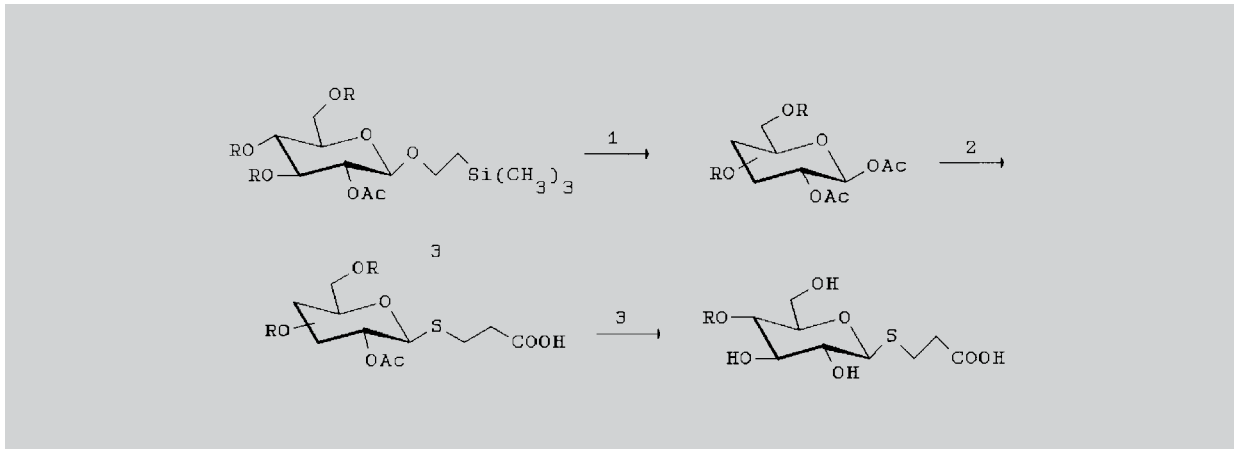
If, for instance, two identical monosaccharides are combined, 11 different disaccharides can be built. Whereas, if two identical aminoacids are combined, only one dipeptide can be formed, and with four different monosaccharides 35,560 unique tetrasaccharides can be created. It is no wonder then that carbohydrates by nature have been assigned a role in communication within living creatures ^{1,2}.

The diversity of polysaccharides has, however, been a problem in the search for biologically relevant carbohydrate structures for use in the struggle against infection and cancer. Not only is it difficult to collect enough material for examination, it has also been difficult to perform carbohydrate chemistry.



However, synthetically made oligosaccharides have become a potential tool in the search for active structures. The technology now has evolved to a level comparable to that employed in peptide and nucleotide synthesis, hence a wide range of synthetic saccharides are now commercially available.

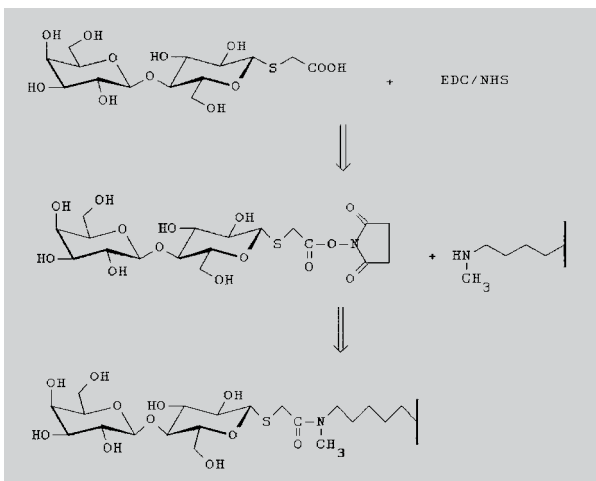
Synthesis of oligosaccharide utilizing the 2-(Tri-methylsilyl)-ethyl (TMSEt) group for anomeric protection.



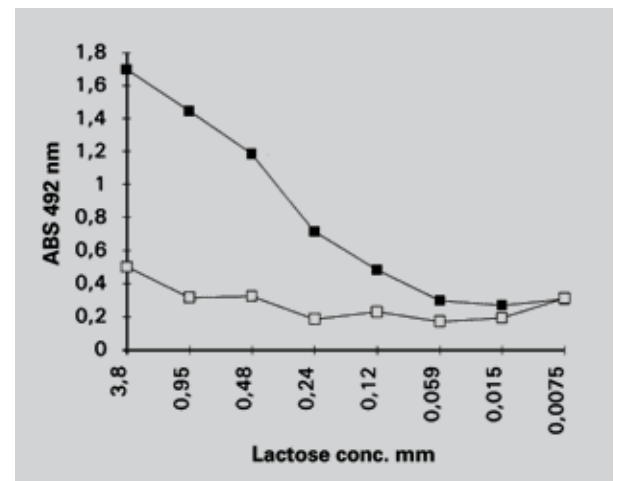
R = acryl group or »protected« sugar residues
A = acetate group

- 1) Typically 3 equivalents of acetic anhydride and 1.1 equivalents of boron trifluoride-etherate in chloroform or toluene at room temperature for 1 hour. Purification by column chromatography³.
- 2) Mercaptopropionic acid (2.5-3 equivalents) and boron trifluoride-etherate (1.2 equivalents) in methylene chloride at 0°C⁴. Purification by column chromatography.
- 3) Deacetylation using sodium methoxide in methanol. Neutralization and chromatographic purification normally produces the deblocked sugar in quantitative yield.

Coupling to Nunc Covalink NH



Binding of activated lactose on Nunc Covalink NH



Coupling of sugar to Nunc Covalink NH and detection of binding exemplified by the Erythrina corallodendron – lactose interaction 5.

Lactose activated (A-Lac) according to the method described above was used:

1. A 1:2 dilution of A-Lac was made in Nunc Covalink NH wells. 100 μ L 7.8 mM A-Lac in H₂O were added to wells in column 1. 50 μ L H₂O were added to the rest of the wells, 50 μ L were transferred from column 1 to column 2, etc. After mixing, 50 μ L were disposed of from column 12. Finally NHS/EDC in aqueous solution (25 μ g each/well) was added to all wells (50 μ L/well).
2. The plate was incubated on shaker at RT for 1.5 hour.
3. Blocked using PBS + 0.5% BSA overnight at 4°C (200 μ L/well).
4. Rinsed with PBS + 0.5% BSA + 0.05% Triton X-100, 3 times 200 μ L/well.
5. Biotinylated Erythrina corallodendron (10 μ g/mL PBS + 0.5% BSA) corresponding to 0.36 μ m regarding binding sites (100 μ L/well).
6. Incubated for 2 hours at RT.
7. Rinsed using CovaBuffer (PBS + 2 M NaCl + 65 mM MgSO₄·2 H₂O + 0.05% Tween 20) 3 times 200 μ L/well.
8. Horseradish peroxidase conjugated Avidine added (1:1000 v/v in PBS + 0.5% BSA) (100 μ L/well).
9. Incubated for 1 hour at RT.
10. Rinsed using CovaBuffer 3 times 200 μ L/well.
11. Substrate (H₂O₂, OPD; 100 μ L/well) 5 min. at RT; the reaction stopped by adding 1M H₂SO₄ (100 μ L/well).
12. The results read at 492 nm.

References

- 1: Nathan Sharon and Halina Lis. Carbohydrates in Cell Recognition. Scientific American, January 1993, 74-81.
- 2: Nathan Sharon and Halina Lis. Lectins as Cell Recognition Molecules. Science, 246, (1989), 227-234.
- 3: Karl Jansson, Stefan Ahlfors, Torbjörn Frejd, Jan Kihlberg, Göran Magnusson, Jan Dahmén, Ghazi Noorl and Kristina Stenvall. 2-(Trimethylsilyl)ethyl Glycosides. 3. Synthesis, Anomeric Deblocking, and Transformation into 1,2-trans 1-O-acyl Sugars. J. Org. Chem., 53, (1988), 5629-5647.
- 4: Mikael Elofsson, Björn Walse and Jan Kihlberg. Building Blocks for Glycopeptide Synthesis: Glycosylation of 3-Mercaptopropionic Acid and Fmoc Amino Acids with Unprotected Carboxyl Groups. Tetrahedron Lett., 32, (1991) 7613-7616.
- 5: N. Gilboa-Garber and L. Mizrahi. A New Mitogenic D-galactosephilic Lectin Isolated from Seeds of the Coral-tree Erythrina Corallodendron. Comparison with Glycine Max (Soybean) and Pseudomonas Aeruginosa Lectins. Can. J. Biochem., 59, (1981) 315-320.

thermoscientific.com/oemdiagnostics

© 2014 Thermo Fisher Scientific Inc. All rights reserved. "Triton" is a registered trademark of Dow Chemical Company, and "Tween" is a registered trademark of Uniqema Americas. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

ANZ: Australia: 1300 735 292, New Zealand: 0800 933 966; **Asia:** China Toll-free: 800-810-5118 or 400-650-5118; India: +91 22 6716 2200, India Toll-free: 1 800 22 8374; Japan: +81-3-5826-1616; Other Asian countries: 65 68729717
Europe: Austria: +43 1 801 40 0; Belgium: +32 2 482 30 30; Denmark: +45 4631 2000; France: +33 2 2803 2180; Germany: +49 6184 90 6000, Germany Toll-free: 0800 1-536 376; Italy: +39 02 95059 554; Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 10200; Russia/CIS: +7 (812) 703 42 15; Spain/Portugal: +34 93 223 09 18; Switzerland: +41 44 454 12 22; UK/Ireland: +44 870 609 9203
North America: USA/Canada +1 585 586 8800; USA Toll-free: 800 625 4327
South America: USA sales support: +1 585 899 7198 **Countries not listed:** +49 6184 90 6000 or +33 2 2803 2000

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand