



In Situ Studies of a Fungal Polysaccharide Using MAC Mode AFM

Application Note

Theresa M. McIntire¹ and Judy Y. Zhu²

¹ Department of Chemistry, University of California, Irvine

² Molecular Imaging (Agilent Technologies)

Introduction

In atomic force microscopy (AFM), a sharp tip is scanned over a sample while maintaining a constant interatomic force between atoms on the end of the tip and atoms on the surface of interest (1). Early implementation of AFM employed the contact mode of operation in which the repulsive force experienced by the tip was measured by recording the cantilever deflection. Under ambient conditions, these repulsive forces range from a few tens to hundreds of nanonewtons. The forces between the AFM tip and the sample cause pressures large enough to induce distortion, remove portions of the sample from the substrate, and even damage the tip. By scanning in a liquid, capillary forces are reduced, which decreases the forces between the tip and sample by orders of magnitude. Other forces such as contact pressures, however, can still lead to dulling of the AFM tip and deformation of soft biological samples.

An effective method to reduce destructive lateral forces and to enhance sensitivity and reproducibility involves dynamic oscillation of the AFM tip while imaging. Agilent's MAC Mode AFM was developed (2) to reduce damage-causing vertical interactions

between the sample and the AFM tip. In MAC Mode, a cantilever coated with a paramagnetic film is oscillated directly by a solenoid located underneath or above the sample. The oscillation amplitude that is utilized to drive the cantilever in MAC Mode is better controlled and can be much smaller than that used in other dynamic-force AFM modes (3). Compared to other AFM imaging modes, DNA molecules resolution has been greatly improved using MAC Mode AFM (4).

In this application note, we demonstrate the ability to image a (1→6)-branched (1→3)-linked-β-D-glucan polysaccharide of fungal origin, in liquid, using MAC Mode AFM. These polysaccharides are widely distributed in plants, fungi, bacteria, and yeast, and serve as structural components of cell walls or as reserve polysaccharides (5). Scleroglucan (6), also known as schizophyllan (7) or lentinan, is a fungal polysaccharide with immunostimulatory properties (8, 9). The rod-like Scleroglucan triple helix structure can be dissociated or denatured in dilute solution into random coil chains (10) and subsequently renatured as linear and circular forms (11), depending on annealing temperature and polymer molecular weight (11-15).



Agilent Technologies

Figure 1a Computer-generated projections of the triple-helical structure of hydrated crystalline curdlan (16), which is a (1→3)- β -D-glucan with no side chains. Each strand of the helix is shown in a different color. A cross-section through the triple-helix structure (**Figure 1b**) reveals the intrahelix hydrogen bonds that stabilize the triple helix. The triangles indicate hydrogen-bonded networks located at successive levels along the

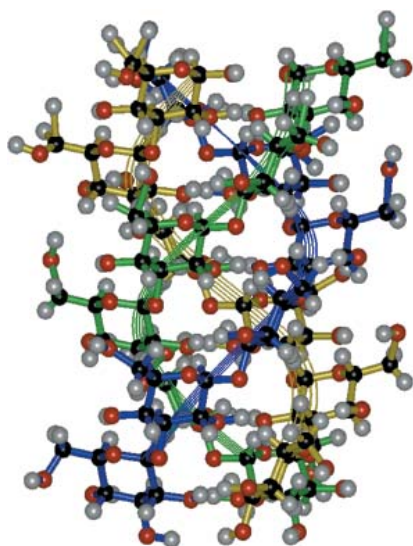


Figure 1a. The accepted triple-helical structure of hydrated crystalline curdlan.

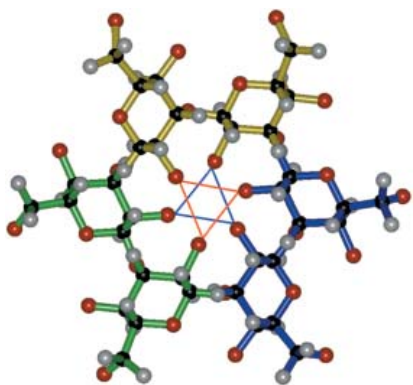


Figure 1b. A cross-section through the triple-helix structure reveals intrahelix hydrogen bonds.

(1→3)- β -D-glucan helix axis (16).

Materials and Methods

Crude scleroglucan powder (Actigum CS 6, Lot No. 26, CECA S.A., Velizy, France) was suspended in aqueous 0.02% sodium azide and stirred overnight at room temperature. The resultant turbid solution was centrifuged to remove cell debris and unsolubilized polymer. Aliquots of the clear supernatant were sonicated (20 kHz, 375 W, 0.5" tip) at 0 °C for 4 hours, centrifuged again, and filtered through 0.45- μ m filters. The resultant clear solution was then fractionally precipitated (17) with acetone to produce fractions of differing mean molar mass. Each fraction was redissolved separately in water, dialyzed exhaustively against distilled water, and recovered by freeze drying. Fractions chosen for study were redissolved in distilled water at 1 mg/mL.

For this study, one of the redissolved scleroglucan fractions was exposed to a denaturation-renaturation procedure:

Aliquots in sealed 10-mL micro-reaction vials were first heated to 160 °C for 15 minutes in a thermostat-controlled oil bath to dissociate the native triple helix (10, 18). The vials were then placed in a thermostat-controlled oil bath at 90°C, annealed at constant temperature for 4 hours, and quenched to room temperature in an ice bath (12, 15).

Results and Discussion

The denatured-renatured polysaccharide solution described was diluted with filtered distilled water to a final polymer concentration of about 10 μ g/mL. Aliquots (20 μ L) of the diluted solution were drop-deposited onto freshly cleaved mica and allowed to air dry. Approximately 200 μ L of n-propanol was placed onto the mica with the dried polysaccharide sample in the liquid cell of an Agilent atomic force microscope. The sample was imaged in MAC Mode with a small multi-purpose scanner. A piezoelectric scanner with a range up to 6.5 μ m was utilized for all images. The AFM tips used were MAClevers with a force constant of 2.0 N/m and resonant frequency ω_0 of approximately 80 kHz in air. Images were stored as 512 x 512-point arrays and analyzed using Agilent image-processing software.

Figure 2 shows a MAC Mode AFM image of a denatured-renatured scleroglucan sample (described below). Clearly, the image shows some aggregation of the molecules, which probably occurred during drying. The tiny spots in the background are due to impurities in the n-propanol. In air, the mean thickness of the renatured cyclic structures is 15–20% smaller than the mean thickness of the renatured linear structures. Both structures are independent of scleroglucan molecular weight and annealing temperature (15).

Conclusions

These studies have demonstrated that linear and circular polysaccharide molecules can be clearly visualized

in situ upon using MAC Mode AFM. With its high spatial resolution, MAC Mode offers a great opportunity to identify local structural differences at the surface of the linear and circular forms. Furthermore, easy setup makes MAC Mode AFM an excellent tool for optimizing imaging conditions for biological samples such as polysaccharides in buffered aqueous solutions. Future studies using MAC Mode AFM in conjunction with Agilent's temperature-controlled stage may reveal other interesting aspects of polysaccharide structure.

Acknowledgment

We thank Dr. Jennifer H.-Y. Liu for generation of the triple-helix structure of curdlan.

References

1. Binnig, G.; Quate, C. F.; Gerber, C. *Physical Review Letters* 1986, 56, 930-933.
2. Han, W.; Lindsay, S. M.; Jing, T. *Applied Physics Letters* 1996, 69, 4111-4114.
3. Hansma, P. K.; Cleveland, J. P.; Radmacher, M.; Walters, D. A.; Hillner, P. E.; Bezantilla, M.; Fritz, M.; Vie, D.; Hansma, H. G.; Prater, C. B.; Massie, J.; Fukunaga, L.; Gurley, J.; Elings, V. *Applied Physics Letters* 1994, 64, 1738-1740.
4. Han, W.; Lindsay, S. M.; Dlakic, M.; Harrington, R. E. *Nature* 1997, 386, 563.
5. *Industrial Gums: Polysaccharides and Their Derivatives*; 3 ed.; Whistler, R. L.; BeMiller, J. N., Eds.; Academic Press, Inc.: San Diego, CA, 1993, pp. 642.
6. Johnson, Jr., J.; Kirkwood, S.; Misaki, A.; Nelson, T. E.; Scaletti, J. V.; Smith, F. *Chemistry and Industry (London)* 1963, 820-822.
7. van der Valk, P.; Marchant, R.; Wessels, J. G. H. *Experimental Mycology* 1977, 1, 69-82.
8. Jong, S.-C.; Donovick, R. *Advances in Applied Microbiology* 1989, 34, 183-262.
9. Pretus, H. A.; Ensley, H. E.; McNamee, R. B.; Jones, E. L.; Browder, I. W.; Williams, D. L. *The Journal of Pharmacology and Experimental Therapeutics* 1991, 257, 500-510.
10. Kitamura, S.; Kuge, T. *Biopolymers* 1989, 28, 639-654.
11. Stokke, B. T.; Elgsaer, A.; Brant, D. A.; Kuge, T.; Kitamura, S. *Biopolymers* 1993, 33, 193-198.
12. McIntire, T. M.; Penner, R. M.; Brant, D. A. *Macromolecules* 1995, 28, 6375-6377.
13. Brant, D. A.; McIntire, T. M. In *Large Ring Molecules*; Semlyen, J. A., Ed.; Wiley: Chichester, UK, 1996; pp. 113-154.
14. McIntire, T. M.; Brant, D. A. *Biopolymers* 1997, 42, 133-146.
15. McIntire, T. M.; Brant, D. A. *Journal of the American Chemical Society* 1998, 120, 6909-6919.
16. Chuah, C. T.; Sarko, A.; Deslandes, Y.; Marchessault, R. H. *Macromolecules* 1983, 16, 1375-1382.
17. Kotera, A. In *Polymer Fractionation*; Cantow, M. J. R., Ed.; Academic Press: New York, NY, 1967; pp. 43-66.
18. Zentz, F.; Verchère, J.-F.; Muller, G. *Carbohydrate Polymers* 1992, 17, 289-297.

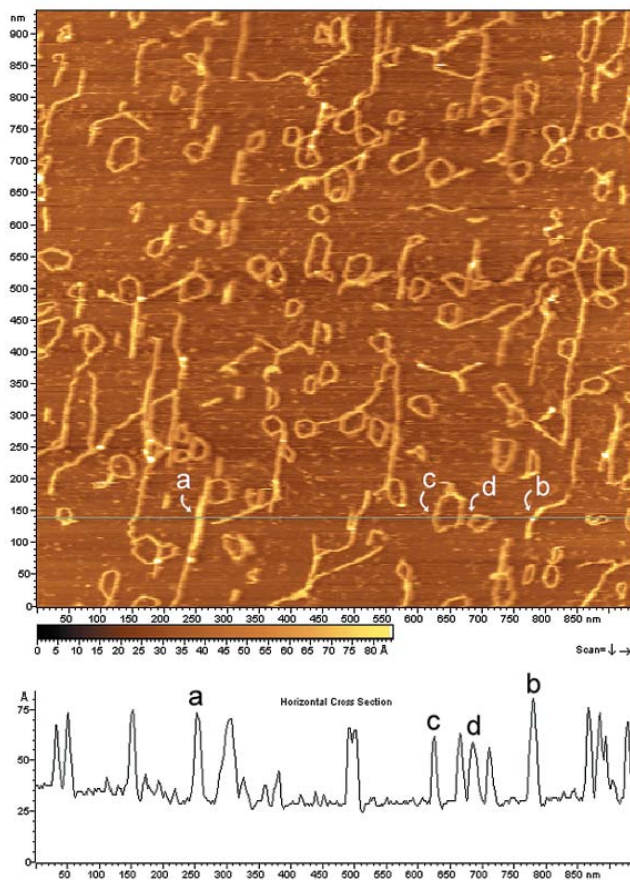


Figure 2. MAC Mode AFM image of renatured linear (a,b) and cyclic structures (c,d).

AFM Instrumentation from Agilent Technologies

Agilent Technologies offers high-precision, modular AFM solutions for research, industry, and education. Exceptional worldwide support is provided by experienced application scientists and technical service personnel. Agilent's leading-edge R&D laboratories are dedicated to the timely introduction and optimization of innovative and easy-to-use AFM technologies.

www.agilent.com/find/afm

www.agilent.com

For more information on Agilent Technologies' products, applications or services, please contact your local Agilent office. The complete list is available at:

www.agilent.com/find/contactus

Americas

Canada	(877) 894-4414
Latin America	305 269 7500
United States	(800) 829-4444

Asia Pacific

Australia	1 800 629 485
China	800 810 0189
Hong Kong	800 938 693
India	1 800 112 929
Japan	81 426 56 7832
Korea	080 769 0800
Malaysia	1 800 888 848
Singapore	1 800 375 8100
Taiwan	0800 047 866
Thailand	1 800 226 008

Europe

Austria	0820 87 44 11
Belgium	32 (0) 2 404 93 40
Denmark	45 70 13 15 15
Finland	358 (0) 10 855 2100
France	0825 010 700
Germany	01805 24 6333*
	*0.14€/minute
Ireland	1890 924 204
Italy	39 02 92 60 8484
Netherlands	31 (0) 20 547 2111
Spain	34 (91) 631 3300
Sweden	0200-88 22 55
Switzerland (French)	44 (21) 8113811(Opt 2)
Switzerland (German)	0800 80 53 53 (Opt 1)
United Kingdom	44 (0) 7004 666666

Other European Countries:

www.agilent.com/find/contactus

Revised: March 23, 2007

Product specifications and descriptions in this document subject to change without notice.

© Agilent Technologies, Inc. 2007

Printed in USA, April 27, 2007

5989-6610EN



Agilent Technologies