

Polystyrene Stars on Agilent PLgel 5 µm MIXED-C using Gel Permeation Chromatography

Application Note

Materials Testing and Research, Polymers

Author

Graham Cleaver Agilent Technologies, Inc.

Introduction

Recently, there has been increasing interest in the synthesis of star-branched polymers due to their unusual flow and viscosity properties compared to linear analogues. Star-branched polymers are constructed with several arms radiating from a central core, either by preparing the individual arms and attaching them to a central molecule, the arms first approach, or by growing the polymer arms from a central core, the core first approach. Many commercial polymers can be constructed with a star-branched morphology relatively easily, but their characterization is still a challenge to the analytical chemist. Gel permeation chromatography (GPC) employing a concentration detector (typically, a refractive index detector) combined with a viscometer can be used to measure not only the molecular weight of the materials, but also to investigate the star-branched structure.

Analysis of Polystyrene Stars

GPC/viscometry was used to analyze a series of star-branched polystyrenes that had been synthesized by a core first approach, giving theoretical 5-, 14- and 21-arm structures. Figure 1 shows a dual detection chromatogram of the 14-arm star-branched polystyrene.





Conditions

Columns 2 × Agilent PLgel 5 µm MIXED-C, 7.5 × 300 mm

(p/n PL1110-6500)

Eluent THF

Flow rate 1 mL/min

Temp 40 °C

Detector Agilent PL-GPC 220

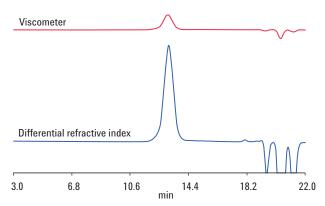


Figure 1 A 14-star-branched polystyrene detected by viscometry and refractive index.

The Universal Calibration approach was used to calculate the molecular weight averages for the star-branched polymers. The universal calibration curve was generated using linear polystyrene (PS) standards with narrow polydispersity (Figure 2).

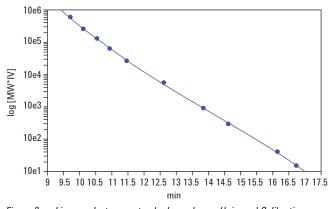


Figure 2 Linear polystyrene standards produce a Universal Calibration curve.

Based on this calibration, the molecular weight averages and weight average intrinsic viscosity (IVw) calculated for the PS star-branched polymers are given in Table 1.

Table 1. Molecular Weight Averages and Weight Average Intrinsic Viscosity for a 14-Star-Branched Polystyrene

Polystyrene

Mp

Molecular weight averages/gmol-					
Mn	Mw	Mz	Mz+1	Mv	ı

IVw

5-arm 56,120 10,460 64,856 98,594 134,877 46,292 6.20 0.28 14-arm 27,436 26,812 29,310 32,425 36,542 28,687 1.10 0.10 21-arm 149,752 111,377 157,884 201,225 256,977 141,293 1.42 0.21

Mark-Houwink plots of log intrinsic viscosity as a function of log molecular weight were calculated for the PS star-branched polymers and for a broad PS material that was known to contain no branching. Figure 3 shows an overlay of the Mark-Houwink plots obtained, indicating that increasing the number of arms on the star-branched PS results in a decrease in intrinsic viscosity at any given molecular weight when compared to the broad linear PS. However, for the 21-arm star-branched polymer, the change in intrinsic viscosity relative to the linear PS varied strongly with the molecular weight.

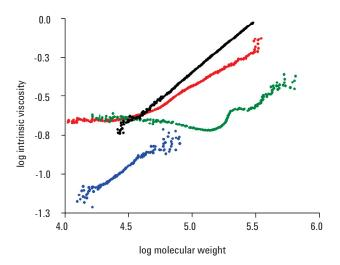


Figure 3 Overlaid Mark-Houwink plots for the linear PS (black) and the 5-arm (red), 14-arm (blue) and 21-arm (green) star-branched polymers.

Based on the linear regions in the Mark-Houwink plots for the stars, the intrinsic viscosity contraction factor for the stars, the intrinsic contraction factor g' was calculated as a function of molecular weight.

$$\frac{g'=[\eta] \text{ star}}{[\eta] \text{ linear}}$$

From the g' data, the radius of gyration contraction factor g was calculated.

$$g' = [a + (1a)g^p]g^b$$

Where a = 1.104, p = 7 and b = 0.906 (Weissmuller & Burchard (1997) Polymer Internat. 44, 380).

Figure 4 shows an overlay of Rg contraction g plots. Using the calculated g values, the functionality f for the stars (the theoretical number of arms) was calculated using a model based on an assumption that the arms were random, that is, polydisperse in molecular weight. For random stars with f number of polydisperse arms:

$$g = \frac{3f}{(f+1)^2}$$

(Burchard (1983) Adv. Polym. Sci. 48, 1; Burchard (1997) Macromolecules 10, 919)

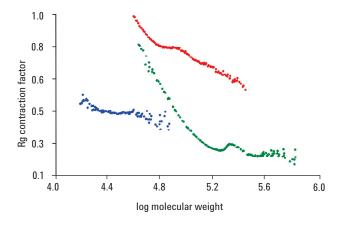


Figure 4. Rg contraction plots for the 5-arm (red), 14-arm (blue) and 21-arm (green) star-branched polymers.

Figure 5 shows an overlay of f for the stars as a function of the log molecular weight. The random model gave a prediction of the functionality f, the number of arms, which was in good agreement with the value expected from the synthesis. However, for all of the star-branched polymers, especially the 21-arm PS, the calculated value of f increased sharply with molecular weight, indicating that a considerable portion of the sample at low molecular weight contained components with fewer than expected arms.

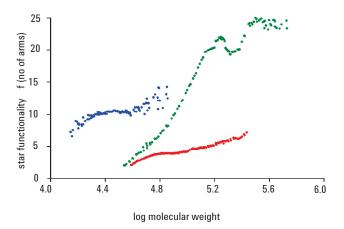


Figure 5. Functionality f plots for the 5-arm (red), 14-arm (blue) and 21-arm (green) star-branched polymers.

Conclusions

The results show that gel permeation chromatography employing refractive index and viscometry detectors can be used to investigate the structure of star-branched polymers. The variation in the functionality f (the number of arms on the polymers) with molecular weight gives valuable insight into the mechanism of the core first approach used to synthesize these materials.

For More Information

These data represent typical results. For more information on our products and services, visit our Web site at www.agilent.com/chem.

www.agilent.com/chem

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc., 2015 Printed in the USA April 30, 2015 5990-8487EN

