



Quantitative GPC Analysis of PVC and Plasticizers

Application Note

Author

Greg Saunders
Agilent Technologies, Inc.

Introduction

Poly(vinyl chloride) is a thermoplastic widely encountered in everyday life. It is light, non-flammable, robust and durable. PVC is permeable, does not deteriorate, is easy to maintain and its physical and mechanical characteristics make it ideal for many different uses. PVC application areas include the toiletry, food, water and car industries.

Unplasticized PVC has a high melt viscosity leading to some difficulties in processing. The finished product is also too brittle for many applications. In order to overcome these problems, it is routine to incorporate additives to the PVC. In addition to acting as impact modifiers, a number of polymeric additives may be considered as processing aids. Such materials are primarily included to ensure more uniform flow and hence improve surface finish. The properties of the final material are dependent on the molecular weight distribution of the PVC and the type and level of the added plasticizers. The analysis of the compounded material is, therefore, of primary importance, and GPC is the ideal analytical tool for its characterization.

With their linear resolving capability over a wide molecular weight range, PLgel MIXED columns provide resolution of both polymer and additives, particularly with the high efficiency 5 μm particle size columns.



Agilent Technologies

Conditions

Columns: PLgel 5 μ m MIXED-C, 600 x 7.5 mm (p/n: PL1110-8500)
Eluent: THF
Flow Rate: 1.0 mL/min
Detection: RI

Results and Discussion

Figure 1 shows the analysis of plasticized PVC. The high efficiency PLgel 5 μ m MIXED-C column provides sufficient scope to resolve the polymer distribution and the four phthalate plasticizers for quantitative analysis.

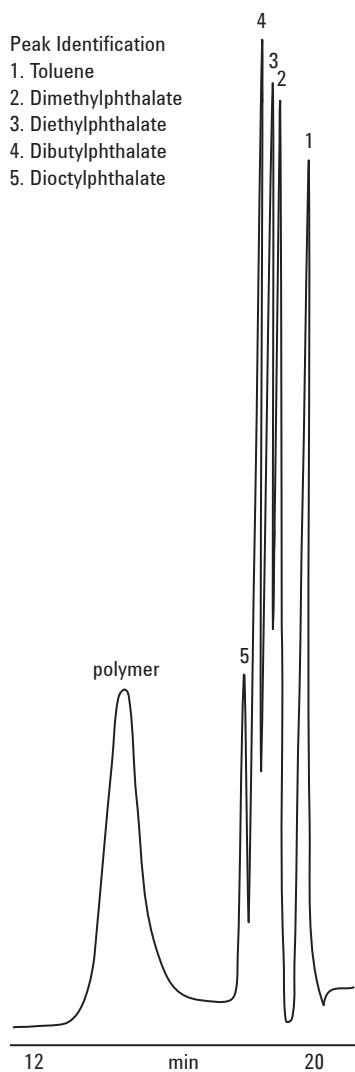


Figure 1. Plasticized PVC analysis.

Conclusion

GPC using PLgel MIXED-C columns permits the determination not only of the molecular weight distribution of PVC samples, but also the identification and quantification of their plasticizers.

PLgel 5 μ m MIXED-C columns are designed for rapid polymer analysis. With its linear calibration up to 2 million MW, this is the column of choice for highest resolution and accuracy in molecular weight distribution analyses. Rapid solvent change capability, excellent temperature stability and the high resolution of the PLgel 5 μ m MIXED-C also provide the versatility essential for the modern R&D laboratory.

www.agilent.com/chem

This information is subject to change without notice.

© Agilent Technologies, Inc. 2011

Published in UK, June 8, 2011

5990-8392EN



Agilent Technologies