

Polymer Degradation Mechanisms using Pyrolysis-GC/MS

Author:
Thomas Wampler

CDS Analytical, Inc.
465 Limestone Road
Oxford, PA 19363-0277

Introduction

The use of analytical pyrolysis¹ with gas chromatography/mass spectrometry to study the structure of polymeric material must be based on an understanding of how these large molecules behave at such elevated temperatures. The utility of the technique is based on the application of thermal energy to produce volatile fragments and products from a macromolecule – compounds capable of being analyzed using GC/MS. The production of these analytes must, of course, be reproducible and follow chemical principles which permit the interpretation of results applied to unknown polymers.

In general, the degradation mechanisms experienced by polymers are free-radical processes initiated by bond dissociation at the pyrolysis temperature. The specific pathway followed by a particular polymer is related to the relative strength of the polymer bonds, and the structure of the polymer chain. These mechanisms are generally grouped into three large categories – random scission, unzipping and side-group elimination.

This Field Application Report shows an example of each of these three general degradation processes. The analytical system consisted of a PerkinElmer Clarus 500 GC/MS interfaced with a CDS Analytical Pyroprobe 2500 Pyrolysis Autosampler*. Samples are rapidly pyrolyzed, automatically introduced into the GC carrier stream and transferred to the GC column for analysis. The specific polymers used are Polyethylene, Poly methylmethacrylate and Poly vinylchloride.

* Available for purchase directly from CDS Analytical Inc., 1-800-541-6593, www.cdsanalytical.com

Experimental

Table 1. Instrument Parameters

| <u>Clarus 500 GC</u> | | <u>Clarus 500 MS</u> | |
|-----------------------------|--|----------------------|-------------|
| Injector Temp | 300° C | Mass Range | 30 - 550 u |
| Oven Program | 40° for 2 min | Scan Time | 0.39 second |
| Rate | 6° C/min | InterScan Delay | 0.01 second |
| | 295° C for 10 min | Transfer Line | 250° C |
| Column | Elite-5* 30 m, 0.25 mm I.D., 1.0 µm film | Solvent Delay Time | 0 seconds |
| Carrier Gas | He (split ratio 50:1) | Source Temperature – | 280 ° C |
| | | Multiplier Voltage | 350 V |
| <u>Model 2500 Pyrolyzer</u> | | Trap Emission | 100 mA |
| Oven | 300° C | Threshold | xx |
| Transfer Line | 300° C | | |
| Pyrolysis Temperature | 750°C | | |
| Pyrolysis Time | 15 seconds | | |
| Heating rate ° C/ms | 10 | | |

* PerkinElmer part number (N9316077)

Samples of approximately 100 µg were analyzed using the conditions listed in Table 1. Sample preparation consisted of simply placing the sample into a quartz tube, which was introduced into the Pyroprobe automatically. No solvent is used, so there is no need for a solvent delay. In fact, some of the compounds found in the earliest eluting peaks may be important diagnostic products for the identification of a particular polymer. The GC column, carrier gas parameters, split ratio and so on are used in the same way as for samples introduced by any other means.

Results

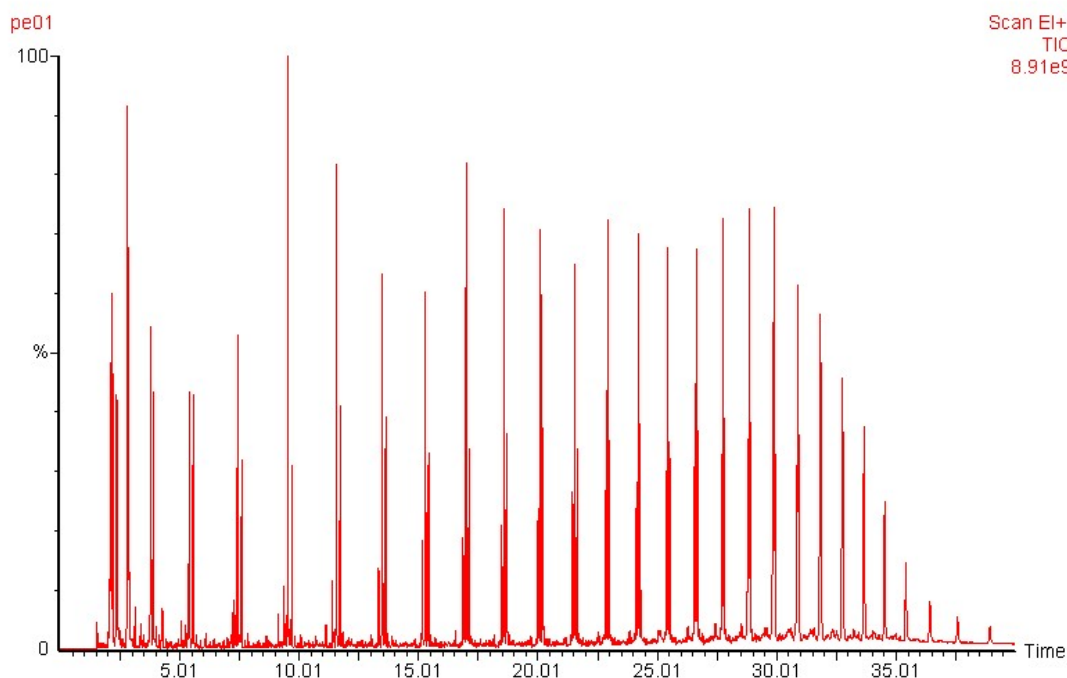


Figure 1. Total Ion Chromatogram resulting from pyrolyzing Polyethylene at 750°C for 15 seconds.

Random Scission – Polyethylene

The total ion chromatogram for the pyrolysis of polyethylene (the pyrogram) is shown in Figure 1. Polyethylene is very high molecular weight hydrocarbon, and the fragments produced via pyrolysis are just pieces of the original molecule small enough to go through the GC. This produces a pattern of peaks for successively longer oligomers, which is common for polyolefins. For polyethylene, the pattern consists of triplets of peaks, as shown in the expanded chromatogram in Figure 2. These are all normal hydrocarbons, and as seen in Figure 3, may be identified as the alkane, alkene and diene of increasing chain lengths. Other polyolefins, such as polypropylene, polybutylene and so on, behave in the same manner, producing characteristic patterns of oligomers specific for that polymer.

Unzipping – Poly methyl methacrylate

Some synthetic polymers, including the methacrylates, simply unzip at pyrolysis temperatures, generating mostly monomer. This is the case for poly methyl methacrylate, shown in Figure 3. Instead of a series of increasingly longer oligomers, the chromatogram shows a large peak for monomer, and little else. This behavior is seen in copolymers involving methyl methacrylate monomer as well as in the pure homopolymer, and acrylic copolymers in general produce monomer peaks for each of the monomers used in the formulation.

Side-group Elimination – Poly vinyl chloride

In the polymers discussed above, the weakest bonds are those holding the chain together, so pyrolysis produces either fragmentation to smaller oligomeric units, or unzipping all the way to monomer. For some polymers, however, the groups attached to the side of the chain are held by bonds which are weaker than the bonds connecting the chain. In such cases, the side groups are stripped off from the chain before it is broken into smaller pieces, so no monomer or higher oligomers are

seen. This is the case for some vinyl polymers, including poly vinyl chloride, poly vinylidene chloride and poly vinyl acetate. In PVC, the carbon-chlorine bond is the weakest, so a chlorine free radical is generated at relatively low temperatures. This chlorine takes a hydrogen from the polymer, making HCl gas and leaving behind a highly unsaturated chain. The chain then degrades to produce aromatics, so the pyrolysis products from PVC include HCl, benzene, toluene and other aromatics, even naphthalene.

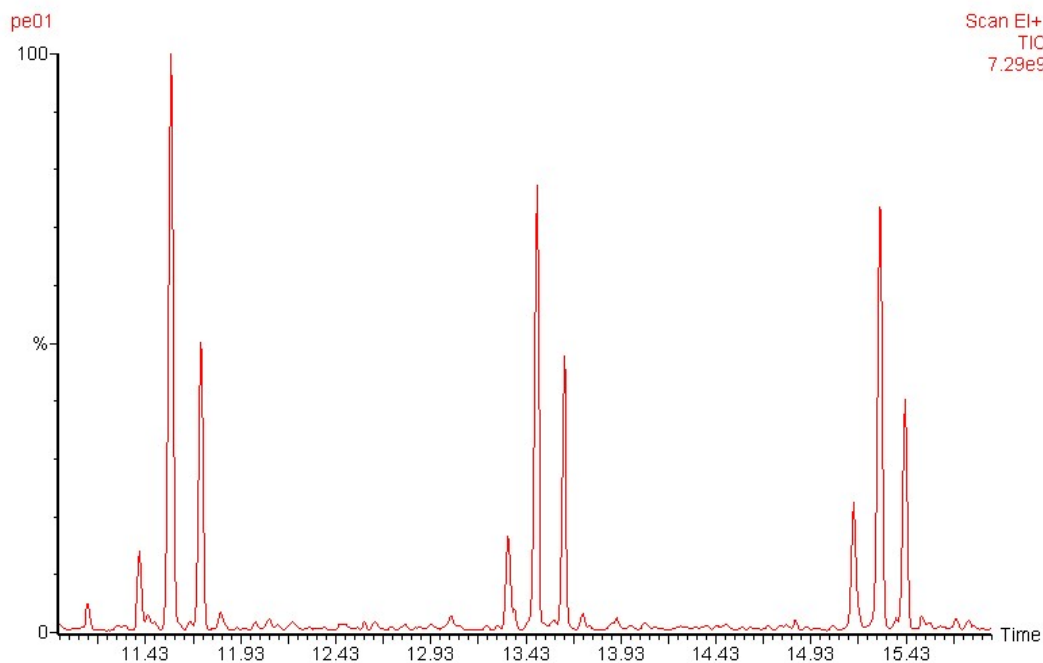


Figure 2. Expanded portion of the pyrogram shown in Figure 1.

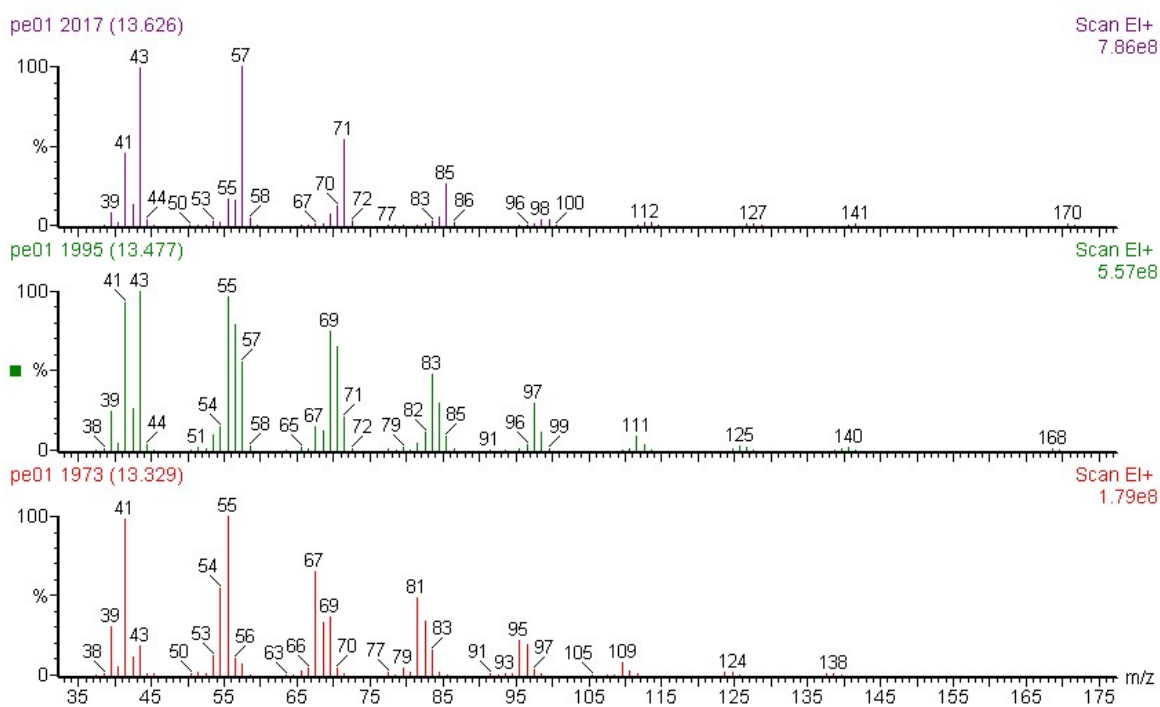


Figure 3. Mass spectrum from peaks at 13.3 to 13.6 minutes, identified using the Wiley 7 database as dodecadiene, dodecene and dodecane.

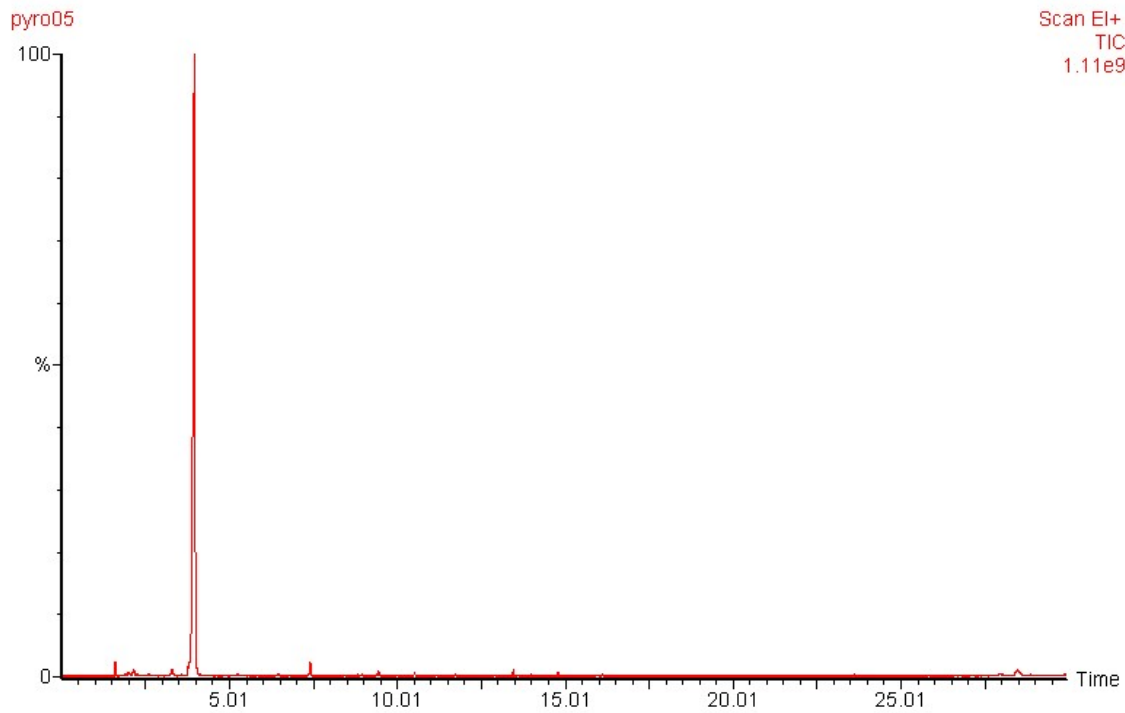


Figure 4. Pyrogram of poly methyl methacrylate.

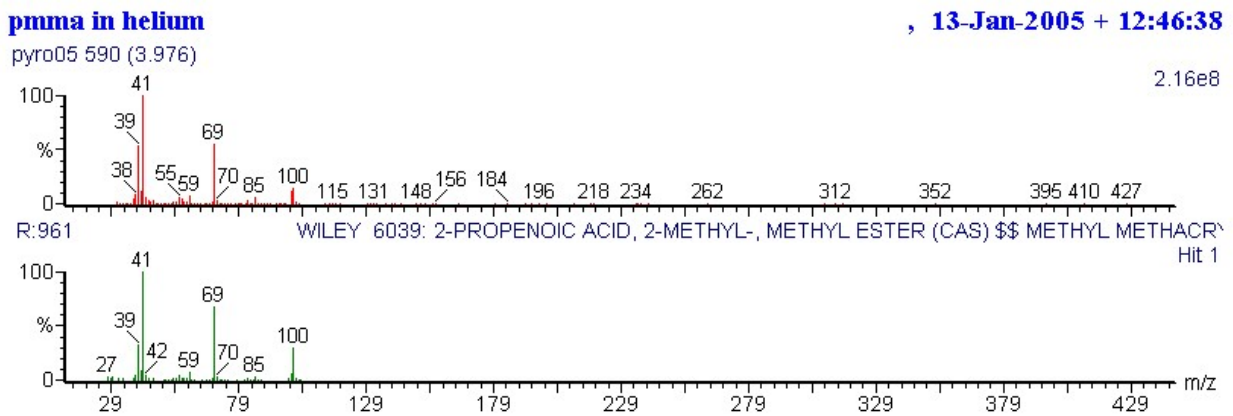


Figure 5. Spectrum of the peak at 4 minutes, identified as the monomer, methyl methacrylate.

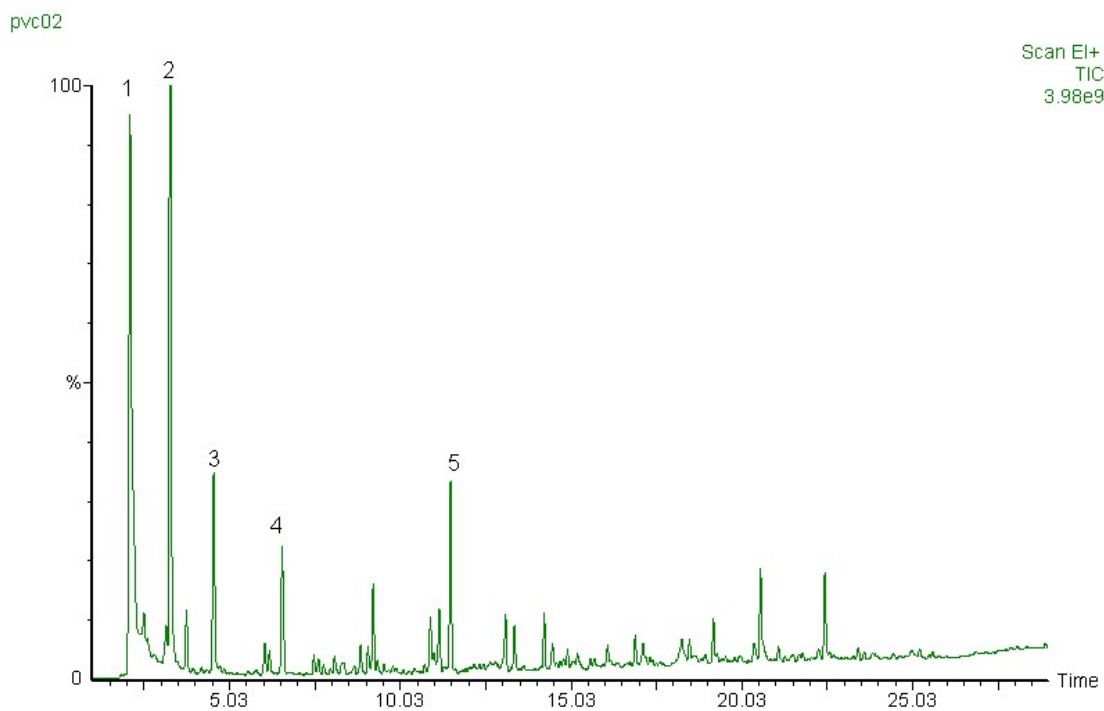


Figure 6. Pyrogram of poly vinyl chloride. Peak identification: 1)HCl, 2)Benzene, 3)Toluene, 4) Xylene 5) Naphthalene

Conclusions

Analytical pyrolysis is a simple technique which extends the use of GC and GC/MS to the analysis of polymeric materials. Interfacing a CDS Model 2500 Pyrolysis Autosampler to a PE Clarus 500 GC/MS produces a system capable of providing detailed analysis of up to 36 samples automatically. Frequently, the products generated by pyrolysis are simple molecules, such as the monomers of the original polymer, and may be identified readily by mass spectrometry. Sample preparation is minimal, and the reproducibility is excellent, with an RSD of just over 2%. The data provided may be used in product identification and reformulation, quality control assays of monomer content, forensic analysis, including the identification of paint fragments from crime scenes and polymer microstructure analysis.

References

1. Applied Pyrolysis Handbook, T. P. Wampler (Ed.) Marcel Dekker, New York (1995).
2. F. C.-Y. Wang, The Microstructure Exploration of Thermoplastic Copolymers by Pyrolysis-Gas Chromatography, J. Anal. Appl. Pyrolysis, 71 (2004) 83-106.

**PerkinElmer Life and
Analytical Sciences**
710 Bridgeport Avenue
Shelton, CT 06484-4794 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/lasoffices

©2004 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. (PKI product names) are trademarks and (PKI product names) are registered trademarks of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.

The data presented in this Field Application Report are not guaranteed. Actual performance and results are dependent upon the exact methodology used and laboratory conditions. This data should only be used to demonstrate the applicability of an instrument for a particular analysis and is not intended to serve as a guarantee of performance.

Enter Project ID number here.