

# Silane Oligomer Analysis with Agilent OligoPore and Gel Permeation Chromatography

## Application Note

Materials Testing and Research

### Authors

Greg Saunders and Ben MacCreath  
Agilent Technologies (UK) Ltd  
Essex Rd  
Church Stretton  
SY6 6AX  
UK

### Introduction

Alkyl silanes are a class of reactive molecules that can be hydrolyzed and polymerized by condensation reactions to form oligomeric and polymeric materials, including silica glasses and gels. This application note illustrates the use of Agilent OligoPore columns and gel permeation chromatography for the analysis of silane oligomers generated by the controlled hydrolysis of silanes under acidic conditions.

### Silane Oligomer Analysis

Dimethyldiethyl (DMDE) silane was mixed with a 0.1 N hydrochloric acid solution in 50:50 ratio and left to react for 12 hours. After the reaction period, the sample was made up at 0.2% (w/v based on the silane) in tetrahydrofuran and injected without further treatment. Figure 1 is the resulting chromatogram of dimethyldiethyl silane after the hydrolysis reactions.



**Agilent Technologies**

## Conditions

Samples	Hydrolyzed dimethyldiethyl silane (DMDES), 0.2% (w/v)
Columns	2 × Agilent OligoPore, 7.5 × 300 mm (p/n PL1113-6520)
Eluent	THF
Flow rate	1.0 mL/min
Inj vol	100 µL
Detector	RI
System	Agilent PL-GPC 50

## Conclusions

Agilent OligoPore columns resolve the individual low molecular weight DMDE silane oligomers after hydrolysis. Individual oligomers can be identified allowing a fingerprint of the silane oligomer distribution after hydrolysis for different times and under different conditions, a vital tool for quality and process control.

## 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](http://www.agilent.com/chem).

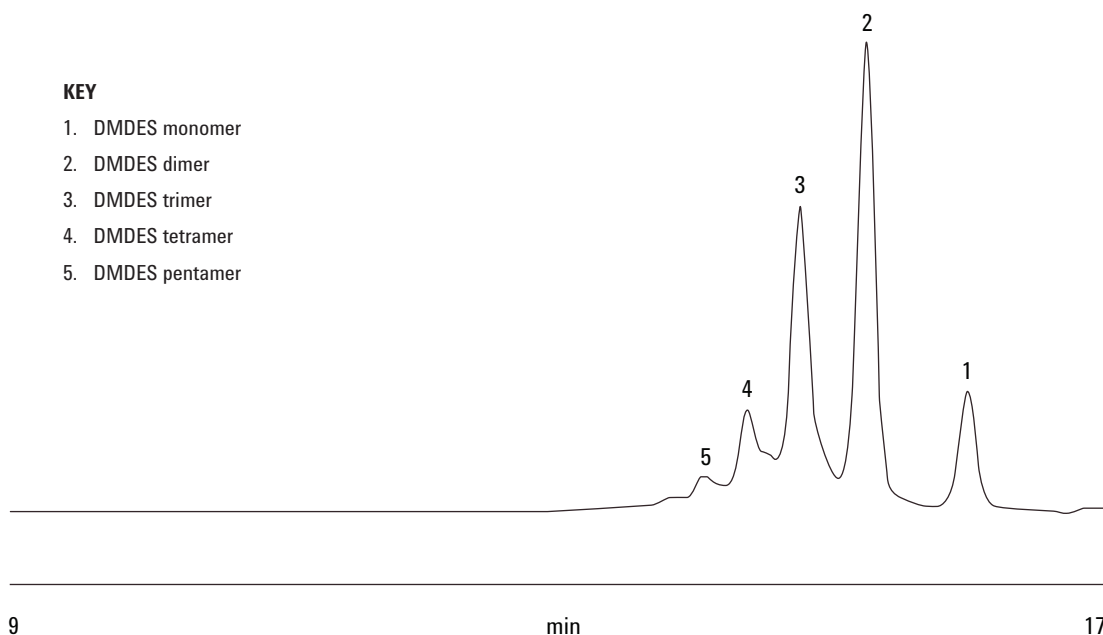


Figure 1. Fingerprinting silane oligomers with an Agilent OligoPore column.

[www.agilent.com/chem](http://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., 2011  
Printed in the USA  
July 12, 2011  
5990-8605EN