

Analysis of Triazine Pesticides in Wastewater by GC/MS

Anila I Khan, Rob Bunn, Tony Edge, Thermo Fisher Scientific, Runcorn, Cheshire, UK

Key Words

- US EPA Method 619
- Triazine
- pesticides
- waste water
- TG-5SiIMS
- 5% Phenyl

Introduction

US EPA 619 is an analytical GC/MS method used for determining certain triazine pesticides in municipal and industrial wastewater. This method was developed by the US Environmental Protection Agency to monitor industrial and municipal discharges under 40 CFR 136.1.

EPA method 619 was performed using a splitless injection mode on a Thermo Scientific TRACE GC coupled to a Thermo Scientific ion trap mass spectrometer.

The Thermo Scientific TraceGOLD TG-5SiIMS provides excellent performance for the analysis of triazine pesticides, in accordance with the EPA method 619.

Goal

To demonstrate the suitability and performance of TraceGOLD™ TG-5SiIMS for the analysis of EPA method 619; triazine pesticides in wastewater.

Experimental details

The triazine pesticides stated in the EPA method 619 were run on a TRACE™ GC fitted with a TriPlus autosampler. The ion trap mass spectrometer was used in a segmented mode to allow precise control of groups of ions for improved ion statistics and ion ratios. The column used for the analysis of a series of triazine compounds, was a low polarity silarylene phase with selectivity comparable to a 5% diphenyl/95% dimethyl polysiloxane phase. The data was acquired and processed using Thermo Scientific Xcalibur data handling software..

Sample preparation

A pre-mixed 1 ng/μL of each triazine pesticide standard solution prepared in acetone was used for the analysis.

Column	Part Number
TraceGOLD TG-5SiIMS, 30 m × 0.25 mm × 0.25 μm,	26096-1420
Guard Column 2 m × 0.32 mm	260RG497
Press-Fit Union	64000-001



Thermo Scientific TriPlus Autosampler

Sample volume	1 μL
---------------	------

TRACE GC Ultra

Oven Program	60 °C (5 min), 8 °C/min, 300 °C (10 min)
Equilibration Time	0.5 min
Injector	275 °C, Splitless (1 min)
Split Flow	30 mL/min
Column Flow	Helium, 1.5 mL/min (constant flow)
Transfer Line Temperature	300 °C

Thermo Scientific Ion Trap MS

MS Type	ITD 230 LT (250 L turbo pump)
MS Source Temperature	225 °C
MS Source Current	250 μA
Electron Energy	70 eV
Filament Delay	5 min
MS Acquisition Mode	El+, 45-450 amu Segmented Scan

Consumables

Consumables	Part Number
BTO 17 mm septa	31303211
3 mm ID Focus Liner, 105 mm long	45350032
Liner graphite seal	29033406
10 μL, 80 mm Syringe	36502019
Graphite ferrules to fit 0.32 mm id columns	29053487
Graphite/vespel 0.25 mm ID ferrules for GC/MS interface	29033496
2 mL clear vial and Si/PTFE seal	60180-599

Results

The requirements for the EPA method 619 were achieved with separation of all of the triazine analytes using the TraceGOLD TG-5SilMS on a GC/MS system. Triazine pesticides are generally difficult to analyse because any activity in the column results in peak tailing or reduced efficiency of the column. The stationary phase in the TraceGOLD TG-5SilMS provides excellent performance due to minimal interaction of active compounds with active sites on the column, therefore no significant peak tailing was observed. Figure 1 shows the TIC chromatogram for 1 ng/μL of triazine pesticides in acetone obtained using a TraceGOLD TG-5SilMS column. Table 1 shows the peak identification of triazine pesticides according to the retention times on the TraceGOLD TG-5SilMS column.

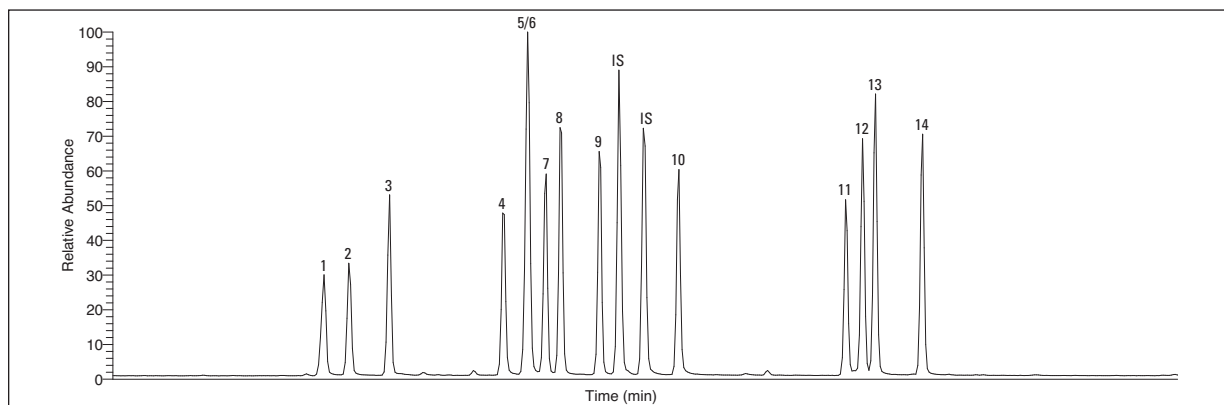


Figure 1: TIC chromatogram of 1 ng/μL of triazine pesticide mix separated on a TG-5SilMS column

Compound Name	Peak No.	Compound Name	Peak No.
Atrazine-desisopropyl	1	Terbutylazine	9
Atrazine-desethyl	2	IS	IS
Terbutylazine-desethyl	3	IS	IS
Atraton	4	Secbumeton	10
Simazine	5	Simetryn	11
Prometon	6	Ametryn	12
Atrazine	7	Prometryne	13
Propazine	8	Terbutryn	14

Table 1: List of triazine pesticides according to retention times

Conclusions

The TraceGOLD TG-5SilMS column demonstrated excellent performance for the analysis of triazine pesticides with no peak tailing of analytes. The chromatogram illustrates the superior performance of the TraceGOLD TG-5SilMS for the analysis of pesticides in accordance with the EPA method 619.

References

1. EPA method 619:
http://water.epa.gov/scitech/swguidance/methods/bioindicators/upload/2007_11_06_methods_method_619.pdf

Acknowledgement

Many thanks to the dedicated team at Cromlab S.L., Barcelona, Spain for all of their help providing the data for this application.

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

**North America
USA and Canada**
+1 800 332 3331

**Europe
France**
+33 (0)1 60 92 48 34

Germany
+49 (0) 2423 9431 -20
or -21

Switzerland
+41 56 618 41 11

United Kingdom
+44 1928 534110

**Asia
Japan**
+81 3 5826 1615

China
+86-21-68654588
or +86-10-84193588
800-810-5118

India
1800 22 8374 (toll-free)
+91 22 6716 2200

**Thermo Fisher
Scientific Australia
Pty Ltd**
1300 735 292 (free call
domestic)

**Thermo Fisher
Scientific New
Zealand Ltd**
0800 933 966 (free call
domestic)

All Other Enquiries
+44 (0) 1928 534 050

Technical Support

North America
800 332 3331

**Outside North
America**
+44 (0) 1928 534 440

www.thermoscientific.com/chromatography

©2011 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

ANCCSGPESTWWW 0311