

Generic HPLC-ELSD Method for Lipids

Dr. Eric Verette¹, Derek Hillbeck²,

¹SEDERE S.A.S., Alfortville Cedex, France, ²Thermo Fisher Scientific, Runcorn, Cheshire, UK

Abstract

This application note highlights the analysis of a mixture of twenty five fatty acids using a reversed-phase HPLC method with a 1.9 µm column and evaporative light scattering detection.

Introduction

Lipids are a diverse set of molecules, the analysis of which has come to the fore due to a greater interest in lipidomics and the drive to analyze the lipidome to provide further insight into disease states in clinical applications, as well as understanding normal physiological homeostasis.

Routine lipid analysis has historically been done by TLC. However, more recently faster technologies with better resolution have been developed including high throughput HPLC-MS and GC-MS systems. Establishing good separation, both analyte from analyte and analyte from matrix components is key to the success of these analyses. The Thermo Scientific Hypersil GOLD range of HPLC columns was developed to give reproducible and reliable chromatographic analysis with excellent peak shape. Greater resolution can be achieved through increased efficiency when using sub 2 µm particle size products. This application note demonstrates the effective separation of a number of typical compounds in the field of lipidomic analysis and their detection at low nanogram levels using a light scattering detector, the SEDEX LT-ELSD™.



Key Words

- Hypersil GOLD 1.9 µm
- Lipid
- Evaporative light scattering
- HPLC
- Fatty acids
- Glycerides

Chemicals and Reagents	Part Number
Fisher Scientific HPLC grade methanol	M/4056/17
Fisher Scientific HPLC grade acetonitrile	A/0626/17
Fisher Scientific HPLC grade formic acid	F/1900/PB15
Fisher Scientific HPLC grade acetone	A/0600/17
Fisher Scientific HPLC grade water	W/0106/17

Separation Conditions	Part Number
Column:	Hypersil GOLD® 1.9 µm, 200 x 2.1 mm

Mobile Phase													
Mobile phase:	A – 50:30:19.8:0.2 (v/v) methanol / acetonitrile / water / formic acid												
	B – 59.8:40:0.2 (v/v) methanol / acetone / formic acid												
Gradient:	<table border="1"> <thead> <tr> <th>Time (min)</th> <th>%A</th> <th>%B</th> </tr> </thead> <tbody> <tr> <td>0</td> <td>100</td> <td>0</td> </tr> <tr> <td>3</td> <td>100</td> <td>0</td> </tr> <tr> <td>43</td> <td>0</td> <td>100</td> </tr> </tbody> </table>	Time (min)	%A	%B	0	100	0	3	100	0	43	0	100
Time (min)	%A	%B											
0	100	0											
3	100	0											
43	0	100											
Flow rate:	0.3 mL/min												
Run time:	45 minutes												
Column temperature:	60 °C												
Injection volume:	2 µL												
Detector:	SEDEX 90LT evaporative light scattering, 28 °C, 3.5 bar pressure												

Experimental Details

Results

Good peak shape and resolution was achieved for the analysis of twenty five fatty acids, fatty alcohols, fat soluble vitamins, mono-, di- and tri-glycerides and related compounds. Detection was by evaporative light scattering detector. RSD values for retention time were less than 0.25 % and for detector response less than 5 % indicating good levels of reproducibility and assay stability, which is essential when comparing results across different data sets. Limits of detection (defined as signal/noise ratio >3) ranged from 0.5 ng on column to 5.7 ng on column with the exception of the semi-volatile component lauric acid where the limit of detection was 16.2 ng on column.

Thermo Fisher Scientific recognize the contribution of Dr Eric Verette (SEDERE, France), for providing the experimental data from which this application note is derived.

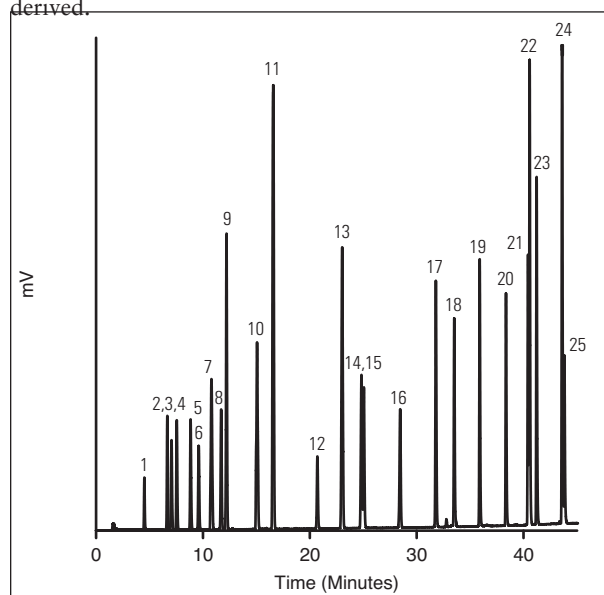


Figure 1: Chromatogram showing the analysis of twenty five lipid related compounds by HPLC-LT/ELSD. Compounds and peak numbering are provided in Table 1

Conclusions

This application note demonstrates that good chromatographic resolution can be obtained using a small particle size Hypersil GOLD 1.9 μm column. This coupled with the SEDEX LT-ELSD™ provides an effective way of analyzing and quantifying a broad range of lipid related compounds.

ID	Compound	t_r	% RSD (n=6)		LOD
		Minutes	t_r	Detector Response	ng (o.c.)
1	lauric acid	4.87	0.22	4.7	16.2
2	linolenic acid	7.17	0.21	3.3	4.1
3	myristic acid	7.58	0.21	2.1	1.6
4	retinol (vitamin A)	8.10	0.20	3.3	3.6
5	linoleic acid	9.43	0.20	2.1	5.1
6	monolein	10.21	0.14	3.3	4.8
7	palmitic acid	11.43	0.25	2.9	0.8
8	oleic acid	12.35	0.23	2.0	5.7
9	hexadecanol	12.88	0.12	4.6	2.1
10	stearic acid	15.77	0.16	2.2	0.5
11	octadecanol	17.32	0.11	2.6	0.5
12	eicosanol	21.63	0.06	3.1	0.7
13	cholesterol	23.80	0.17	2.8	1.3
14	docosanol	25.57	0.06	3.2	0.9
15	α -tocopherol (vitamin E)	25.80	0.05	2.9	3.8
16	vitamin K	29.80	0.02	3.6	3.8
17	squalene	32.54	0.12	2.0	2.4
18	diolein	34.13	0.05	2.8	2.3
19	trilaurin	36.50	0.10	3.1	2.1
20	trilinolenin	38.90	0.08	4.0	2.5
21	trimyristin	40.97	0.08	4.7	1.7
22	coenzyme Q10	41.09	0.03	2.7	1.8
23	trilinolein	41.73	0.06	3.6	1.9
24	tripalmitin	44.09	0.06	3.9	1.7
25	triolein	44.29	0.06	4.5	1.1

Table 1: List of compounds analyzed showing retention time, %RSD for retention time and response, and limits of detection. ng (o.c.): amount on column (ng)

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

United Kingdom
+44 1928 534110

**Asia
Japan**
+81 3 5826 1615

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

India
+91-22-6742 9494

**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. SEDEX LT-ELSD is a trademark of SEDER S.A. France. All other 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 rep for details.

ANCCSLCELSLDIIP 1211