

**Improving the Analysis
of Fatty Acid Methyl
Esters Using
Automated Sample
Preparation Techniques**

Rebecca Veeneman
Agilent Technologies

Pittcon 2011
March 16, 2011
1370-5

Outline

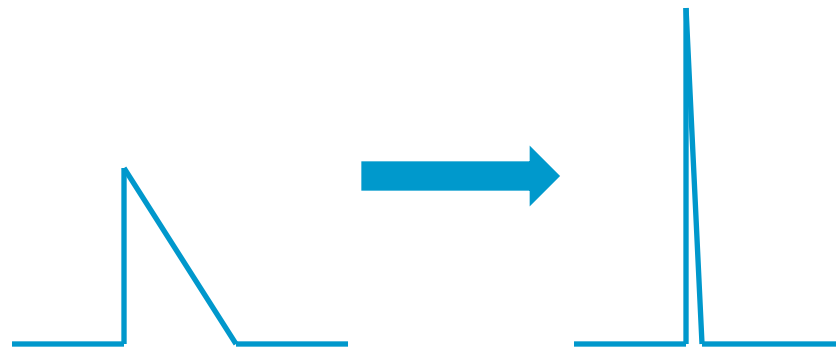
- Introduction
 - Fatty acids and the challenges associated with their analysis
 - Methods of derivatization
 - Automated solutions
 - 7696A Sample Prep WorkBench
- Development of an automated method for WorkBench
 - Calibration curves
 - Method validation
 - Oil sample analysis
- Conclusions

Analysis of Fatty Acids Is Common In Multiple Industries

- Fatty acids are found in many matrices
 - Food industry
 - Total lipid analysis (triglycerides, phospholipids, sterols): milk, eggs, meat, oils, seeds
 - Edible fat analysis (mainly fatty acids): oils
 - Biomedical applications
 - Fatty acid profiles as a diagnostic tool: blood, tissue
 - Chemical industry
 - Fatty acids found in cosmetics, surfactants, other household products

Fatty Acid Analysis Can Be Challenging

- Gas chromatography has been the predominant method of analysis
 - Underivatized fatty acids can be analyzed with polar columns but often have poor peak shape and long retention times
- Fatty acids are often derivatized to improve the peak shape and separation
 - More reproducible data



There Are Various Methods of Derivatization

- Acid catalyzed reactions forming methyl esters
 - Reagents include BF_3 , HCl , or H_2SO_4
 - HCl and H_2SO_4 often need long reaction times and high reaction temperatures
 - BF_3 can methylate fatty acids within 2 minutes
 - Works on free fatty acids, phosphoglycerides, and triglycerides
- Base catalyzed reactions forming methyl esters
 - Reagents include NaOH or KOH in methanol
 - Very fast
 - Does not work on free fatty acids
- Methyl esters produced with diazomethane
 - Toxic and explosive; can produce byproducts
- Silylation reactions to form trimethylsilyl esters
 - Sensitive to water

Preparation of Fatty Acid Methyl Esters with an Acid Catalyzed Reaction

Step 1: 100 mg sample in 20 mL test tube

Step 2: Add 2 mL 2N NaOH in methanol

Step 3: Heat 80°C for 1 hour

Step 4: Add 2 mL 25% BF₃ in methanol

Step 5: Heat 80°C for 1 hour

Step 6: Add 5 mL 1M NaCl in H₂O

Step 7: Add 5 mL Hexane

Step 8: Shake

Transfer supernatant to autosampler vial

F. David, P. Sandra, P. Wylie, "Improving the Analysis of Fatty Acid Methyl Esters Using Retention Time Locked Method and Retention Time Databases," Application Note 5990-4822EN, Agilent Technologies publication 5988-5871EN (2003).

AOAC Official Methods of Analysis (1990), method 969.33.



Preparation of Fatty Acid Methyl Esters with a Base Catalyzed Reaction

Step 1: 100 mg sample in 20 mL test tube

Step 2: Dissolve in 10 mL hexane

Step 3: Add 100 mL 2N KOH in methanol

Step 4: Mix 30s

Step 5: Centrifuge

Transfer supernatant to autosampler vial

F. David, P. Sandra, P. Wylie, "Improving the Analysis of Fatty Acid Methyl Esters Using Retention Time Locked Method and Retention Time Databases," Application Note 5990-4822EN, Agilent Technologies publication 5988-5871EN (2003).

Automated Procedures Can Simplify FAME Preparation

- Tecan Freedom Evo
- Gerstal MPS PrepStation
- Agilent 7696 Sample Prep WorkBench
 - 2 syringe modules
 - 150 vial rack
 - Single vial vortex mixer
 - Single vial heater (80°C)
 - Individually heated (80°C) and cooled (5°C) racks
 - Sample tracking



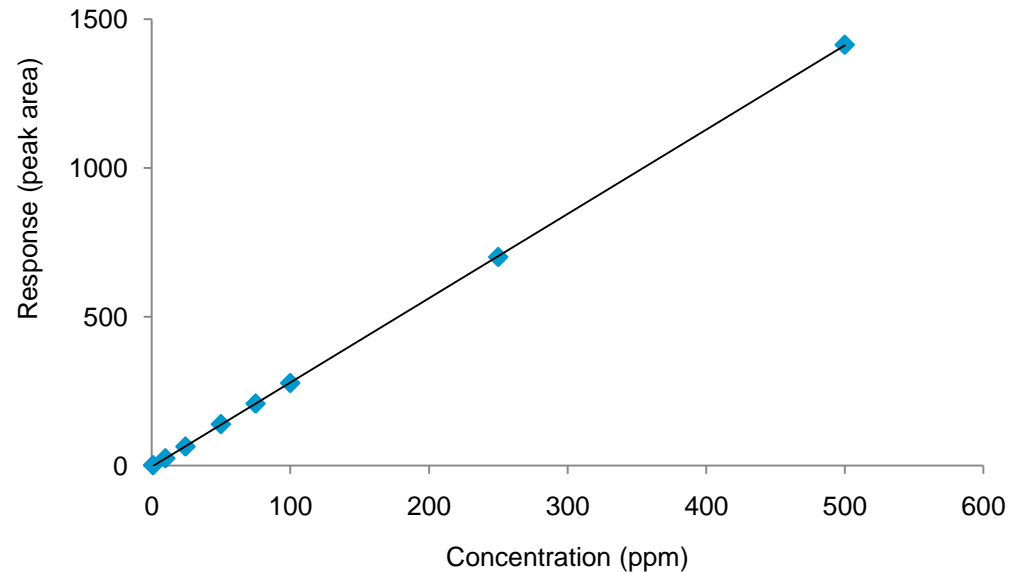
R. Perkins, K. Summerhill, and J. Angove, *Chromatography Today*, Sept/Oct, 17-19 (2008).
M. Athar Masood and N. Salem Jr., *Lipids*, **43**, 171-180 (2008).

Step-Wise Approach to Developing a Method for WorkBench

- Make calibration curve standards
- Validate the automated method
 - Translate manual AOAC method to WorkBench appropriate volumes
 - Verify WorkBench method gives the same (or better) results as a manual method with a fatty acid standard
- Acid reaction
 - Canola oil sample
- Base reaction
 - Canola oil sample

Calibration Curve Standard Preparation Is Fast and Yields Excellent Results

- Eight level calibration curve: 1 to 500 ppm
- Linear dilutions in ~100 μL of hexane
- Complete in 40 minutes
- Excellent linearity ($R^2=0.999$)



Calibration Curve Standard Preparation Is Fast and Yields Excellent Results

- Eight level calibration curve: 1 to 500 ppm
- Linear dilutions in ~100 μL of hexane
- Complete in 40 minutes
- Excellent linearity ($R^2=0.999$)

The screenshot displays the 'Setup Method' window for the Agilent 7696A Sample Prep Method. The interface includes an 'Actions' toolbar with icons for Add, Mix, Heat, Wait, and Flag as result. The 'Program' section shows a flowchart of 16 steps, with step 14 highlighted. The 'Steps' list on the right provides detailed instructions for each step, including the volume of Hexane or FAME 1 mg/mL added and the concentration at different towers. The 'Available Resources Tracked By Use' table is also visible.

Resource Name	Resource Type	Uses/Vial	Vial Range
New Vial	Empty Container	1	1-20

Resource Name	Resource Type	Usable Volume/Vial
BF3	Chemical Resource	1500
BSTFA	Chemical Resource	1500
Canola Oil	Chemical Resource	500
FAME 1 mg/mL	Chemical Resource	500
Hexane	Chemical Resource	1500

Translating the Acid Catalyzed FAME Preparation

- Step 1: Add 10 μL sample
- Step 2: Add 10 μL internal standard
- Step 3: Add 40 μL 2N NaOH in methanol
- Step 4: Mix 30s
- Step 5: Add 80 μL 14% BF_3 in methanol
- Step 6: Mix 30s
- Step 7: Heat at 65°C for 20 min
- Step 8: Cool 2 min
- Step 9: Add 100 μL 1M $\text{H}_2\text{O}/\text{NaCl}$
- Step 10: Add 100 μL hexane
- Step 11: Mix 20s
- Step 12: Transfer top 100 μL to a new vial

Translating the Acid Catalyzed FAME Preparation

Setup Method

Agilent 7696A Sample Prep Method | Agilent 7696A Configuration

Import | Export

Version 2.3.31.0

Actions

Add Mix Heat Wait Flag as result

Program

```
graph TD; S1[1. Mix] --> S2[2. Add]; S2 --> S3[3. Add]; S3 --> S4[4. Add]; S4 --> S5[5. Mix]; S5 --> S6[6. Add]; S6 --> S7[7. Mix]; S7 --> S8[8. Heat]; S8 --> S9[9. Wait]; S9 --> S10[10. Add]; S10 --> S11[11. Add]; S11 --> S12[12. Mix];
```

Steps

1. Mix Sample at 1000 RPM for 0 min 5 sec
2. Add 10 uL of Sample to oil 1 at Back Tower (washes, pumps)
3. Add 10 uL of ISTD to oil 1 at Back Tower (washes, pumps)
4. Add 40 uL of NaOH to oil 1 at Front Tower (washes, pumps)
5. Mix oil 1 at 1000 RPM for 0 min 30 sec
6. Add 80 uL of BF3 to oil 1 at Front Tower (washes, pumps)
7. Mix oil 1 at 1000 RPM for 0 min 30 sec
8. Heat oil 1 at 65 °C for 1 min 0 sec
9. Wait for 1 min 0 sec
10. Add 100 uL of H2O/NaCl to oil 1 at Front Tower (washes, pumps)
11. Add 100 uL of Hexane to oil 1 at Front Tower (washes, pumps)
12. Mix oil 1 at 1000 RPM for 0 min 20 sec

Available Resources Tracked By Use

Resource Name	Resource Type	Uses/Vial	Vial Range
New Vial	Empty Container	1	1-20

Available Resources Tracked By Volume

Resource Name	Resource Type	Usable Volume/Vial
BF3	Chemical Resource	1500
BSTFA	Chemical Resource	1500
Canola Oil	Chemical Resource	500
FAME 1 mg/mL	Chemical Resource	500
H2O/NaCl	Chemical Resource	1500

OK | Apply | Cancel | Help

WorkBench Method Validation with a Fatty Acid Standard

- WorkBench made 5 samples per day for 3 days
 - Determine day to day reproducibility
 - Sample was a 1 mg/mL fatty acid standard
 - 1 mg/mL alkane internal standard
- For any given day, the RSD for 5 samples was <2% and recovery was 103%
- For the 15 samples made across 3 days, the average RSD was 1.2% and recovery was 103%

Manually Prepared Samples Were Not as Reproducible

- 5 samples made on 3 days alongside the WorkBench made samples
- For a given day, the average RSD was $>4.5\%$ with an average recovery of $>110\%$
- For the 15 samples made over 3 days, the average RSD was 6.8% with an average recovery of 125%
- Conclusion: WorkBench made samples are as good as manually prepared samples

Dispensing Reagents with WorkBench Is Reproducible and Accurate

- Dispensing of the reagents was measured gravimetrically to determine the accuracy and precision of WorkBench as a liquid handling system

Dispensed Volume	RSD	Accuracy
10 μ L sample	0.84%	10%
40 μ L 2N NaOH in methanol	0.33	2.1%
100 μ L 2N NaOH in methanol	0.48%	1.1%
80 μ L 14% BF ₃ in methanol	0.30%	0.93%
100 μ L 1M H ₂ O/NaCl	0.55%	1.0%
100 μ L hexane	0.54%	1.9%
500 μ L hexane	0.27%	0.30%

WorkBench Prepared Samples – Acid Prep

Analyte	Amount (ppm)	RSD (%)	Recovery (%)
Methyl laurate	51	-	97
Methyl palmitate	1500	0.78	-
Methyl stearate	307	0.93	-
Methyl eicosanoate	227	1.1	-
Methyl behenate	112	0.86	-

- 11 samples prepared across 2 days
- Sample was mixture of the initial oil sample and a lauric acid standard
- RSD was calculated using methyl laurate as the internal standard
 - Average RSD=0.92%
 - Using an internal standard takes dilution inaccuracy into account
 - Using an external standard, the RSD was 4.0%

Base Catalyzed Preparation

Step 1: Add 10 μL sample

Step 2: Add 500 μL hexane

Step 3: Mix 30s

Step 5: Add 100 μL 2N NaOH in methanol

Step 6: Mix 60s

Step 7: Transfer top 100 μL to a new vial

Base Catalyzed Preparation

Setup Method

Agilent 7696A Sample Prep Method | Agilent 7696A Configuration

Import | Export

Version 2.3.31.0

Actions

Add | Mix | Heat | Wait | Flag as result

Program

```
graph LR; A[1. Mix] --> B[2. Add]; B --> C[3. Add]; C --> D[4. Mix]; D --> E[5. Add]; E --> F[6. Mix];
```

Steps

1. Mix Sample at 1000 RPM for 0 min 10 sec
2. Add 10 uL of Sample to oil 1 at Back Tower (washes, pumps)
3. Add 500 uL of Hexane to oil 1 at Front Tower (washes, pumps)
4. Mix oil 1 at 1000 RPM for 0 min 30 sec
5. Add 100 uL of NaOH to oil 1 at Front Tower (washes, pumps)
6. Mix oil 1 at 1000 RPM for 1 min 0 sec

Available Resources Tracked By Use

Resource Name	Resource Type	Uses/Vial	Vial Range
New Vial	Empty Container	1	1-20

Available Resources Tracked By Volume

Resource Name	Resource Type	Usable Volume/Vial
BF3	Chemical Resource	1500
BSTFA	Chemical Resource	1500
Canola Oil	Chemical Resource	500
FAME 1 mg/mL	Chemical Resource	500
U200 N-Cl	Chemical Resource	1500

OK | Apply | Cancel | Help

WorkBench Prepared Samples – Base Prep

Analyte	Amount (ppm)	RSD (%)	Recovery (%)
Hexadecane	9.7	-	99
Methyl palmitate	313	2.7	-
Methyl stearate	50	4.9	-
Methyl eicosanoate	41	2.2	-
Methyl behenate	18	2.8	-

- 10 samples prepared in 1 day
- Sample was mixture of the initial oil sample and an alkane standard
 - The base catalyzed reaction does not esterify free fatty acids
- RSD was calculated using hexadecane as the internal standard
 - Average RSD=3.2%
 - Using an external standard, the RSD was 4.5%

Conclusions

- FAME preparations can be easily translated to an automated system
 - Method development on WorkBench enabled a shorter reaction time
- Samples prepared with WorkBench are more reproducible than those prepared manually
- WorkBench can prepare FAMEs via acid or base catalyzed reactions
 - Acid catalyzed reactions result in an average RSD <1% with 97% recovery
 - Base catalyzed reactions result in an average RSD 3% with 99% recovery

Conclusions

- Preparing oil samples with WorkBench reduces the volume of reagents needed
 - Acid catalyzed reaction was reduced ~50-fold
 - Base catalyzed reaction was reduced ~10-fold
- Operator exposure is reduced
 - Reagents and samples are in capped vials
- Automated preps frees the operator to work on other tasks beyond sample preparation

THANK YOU