

Evaluation of SFC-MS Configurations for the Analysis of Lipids, Sterols, and PAHs

Introduction

Supercritical fluid chromatography (SFC) is a powerful separation technique, and especially for the separation of apolar compounds, it offers an interesting alternative to normal phase HPLC. Most SFC work has been performed using UV detection, although the coupling to mass spectroscopic detection can largely extend SFC applicability.

Initial experiments with the Agilent 1260 Infinity SFC/MS system showed that the hyphenation of SFC with MS results in high sensitivity and good mass spectral quality due to the fact that the major component of the mobile phase is CO₂, which evaporates upon entering the source, allowing for analysis at high flow rates. However, the carbon dioxide decompression and expansion after the backpressure regulator (BPR), used in all commercially available instruments, also causes cooling of the transfer line to the MS, which can result in solute deposition, baseline noise and irreproducible results.

For hyphenation of SFC with MS, different configurations can be used, including effluent splitting before or after the BPR [1]. Effluent splitting obviously reduces sensitivity and alternatives were studied. Best results in term of sensitivity, baseline stability, retention time and peak area repeatability were obtained using a configuration whereby a small (typically 0.2 mL/min) make-up flow is added before the BPR [2]. The total effluent is introduced in the MS (both ESI and APCI were tested). It was also found that heating of a part of the transfer line, just before the ionization source entrance, was needed to obtain optimal peak shape and sensitivity.

This SFC-MS configuration was used to analyze several different types of apolar solutes, including lipids, sterols, and polycyclic aromatic hydrocarbons (PAHs). These compounds are notoriously difficult to detect using LC-MS, and comparisons were made between the two techniques. In all cases, similar separation and MS conditions were used.

1 System Configuration and Initial Testing

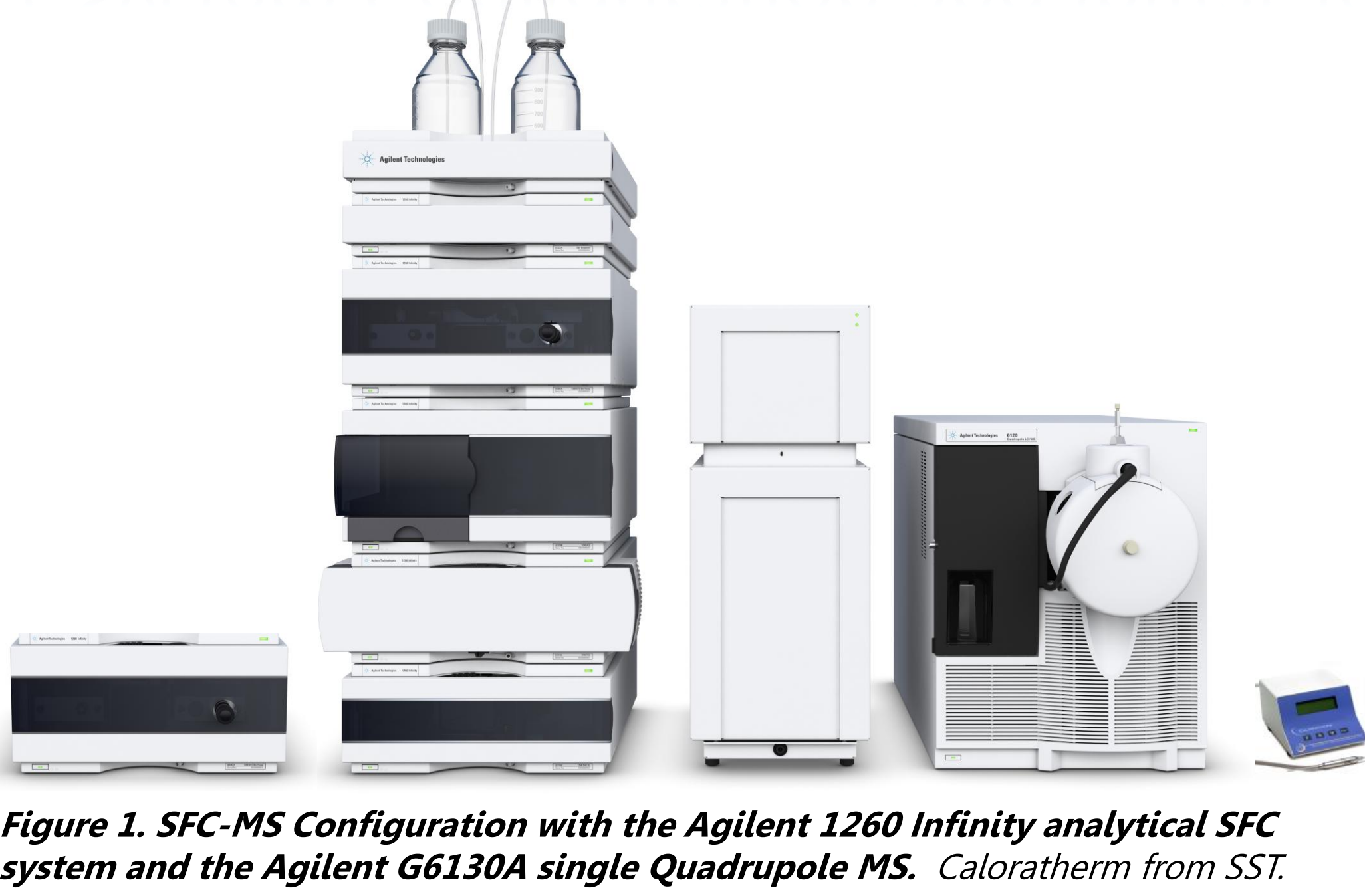
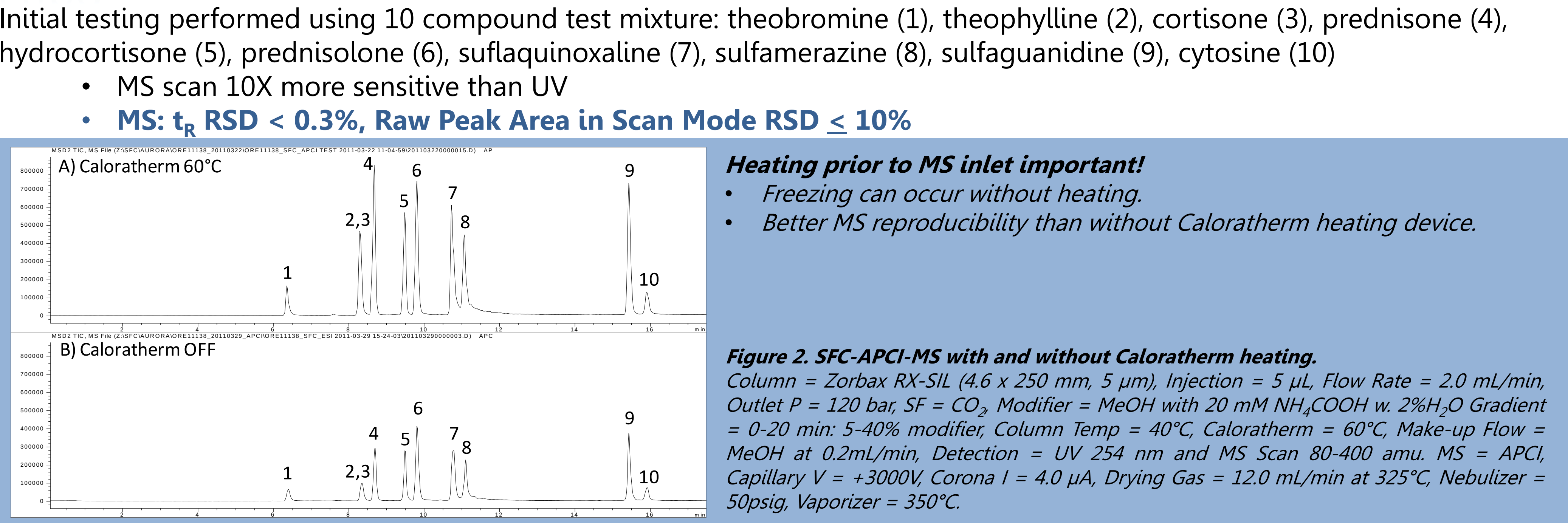
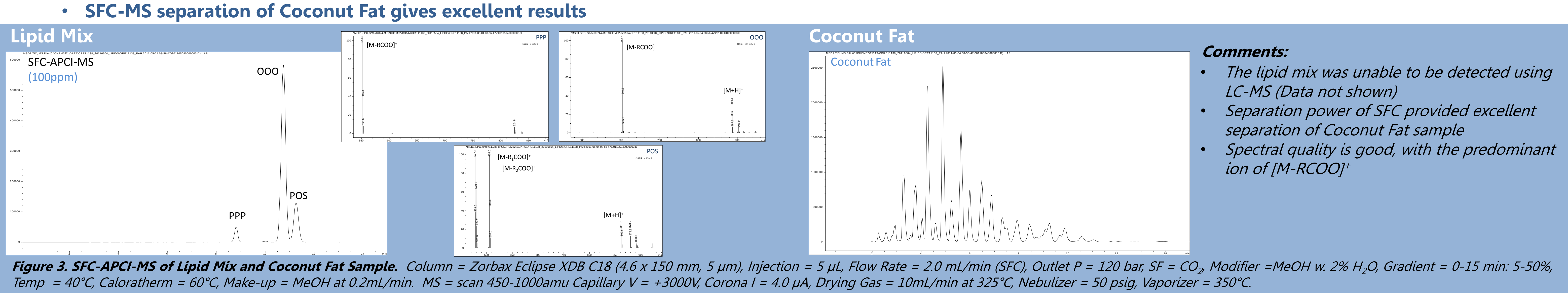


Figure 1. SFC-MS Configuration with the Agilent 1260 Infinity analytical SFC system and the Agilent G6130A single Quadrupole MS. Caloratherm from SST.



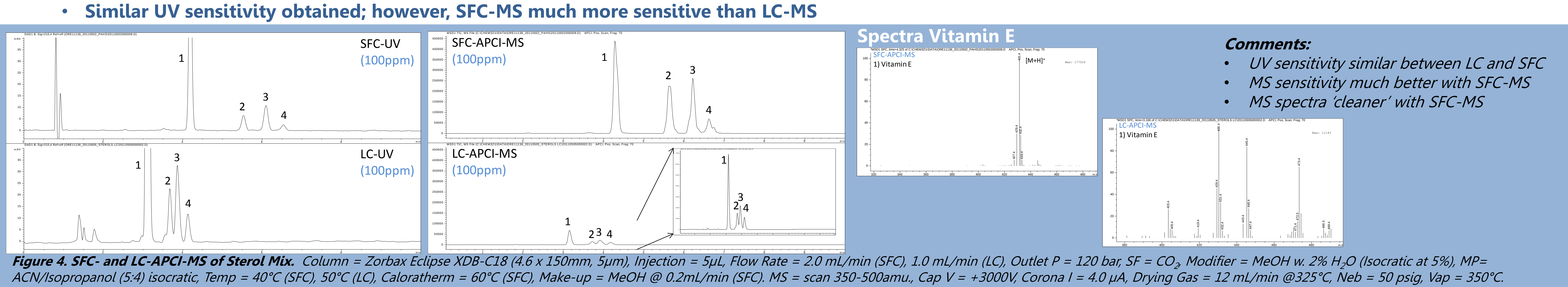
2 Lipids and Coconut Fat

A lipid mixture containing PPP, OOO, and POS was separated using LC-MS and SFC-MS. The same MS settings were used in both cases.



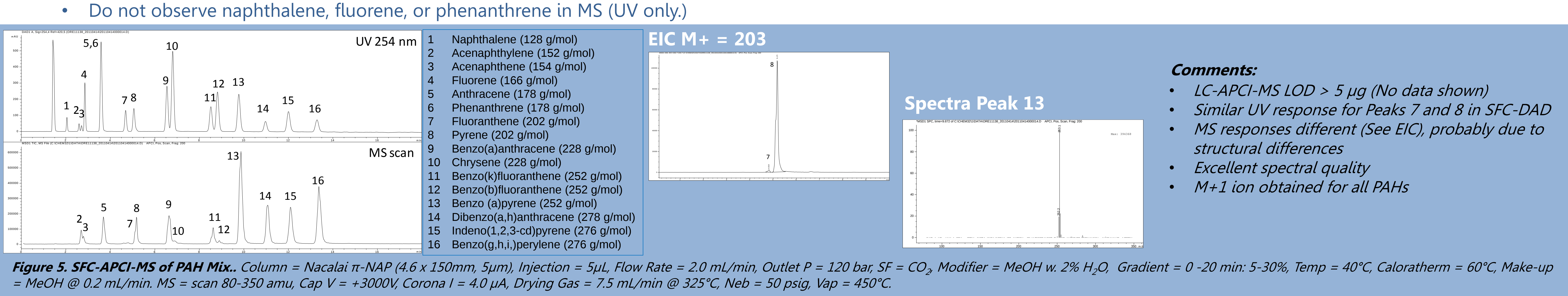
3 Sterols

A sterol mixture [vitamin E (1), cholesterol (2), stigmasterol (3), and β -sitosterol (4)] was separated using LC-APCI-MS and SFC-APCI-MS. The same MS settings were used in both cases.



4 PAHs

A PAH mixture containing 16 PAHs at 100ppm was separated using SFC-APCI-MS.



[1] Aurigemma and Farrell. Chromatogr A. 1217 (2010) 6110-6114
[2] Agilent Technical Note [5990-7972EN]

We acknowledge, Udo Huber, Hans-Gerog Härtl, and Markus Becker from Agilent for their help and support.