# Breath Biopsy: combining Thermal Desorption-Gas Chromatography with High Resolution Mass Spectrometry for improved sensitivity and selectivity in untargeted breath analysis

Lori Dolata<sup>1</sup>, Dominic Roberts<sup>1</sup>, Cristian Cojocariu<sup>1</sup>, Paul Silcock<sup>1</sup>, Jason Cole<sup>1</sup>, Max Allsworth<sup>2</sup>. <sup>1</sup>Thermo Fisher Scientific, Runcorn, UK. <sup>2</sup>Owlstone Medical Ltd., CB4 0GA UK.

## ABSTRACT

**Purpose:** The objective of this study was to employ a new analytical approach that allows reliable, sensitive and selective analysis of breath volatiles in untargeted metabolomics.

**Methods:** Breath samples were collected and pre-concentrated onto Breath Biopsy Cartridges using a ReCIVA® Breath Sampler. Samples were analysed using a thermal desorption (TD) autosampler interfaced with a Thermo Scientific<sup>™</sup> Exactive<sup>™</sup> GC Orbitrap<sup>™</sup> GC-MS system. Wide dynamic range (> 5 orders of magnitude) and ppt-level sensitivity of the Orbitrap mass spectrometer allow for the detection of both high-abundance volatile organic compounds (VOCs) (e.g. acetone) and tracelevel analytes within a single analysis. High mass resolution/accuracy and alternative ionisation methods were employed together with software-based features to improve identification of VOCs in high-density breath data.

**Results:** This study reports the first use of a thermal desorption-GC Orbitrap platform for preconcentrated breath sample analysis, proving it to be a powerful tool for metabolomics studies and biomarker discovery.

## INTRODUCTION

Exhaled breath contains a wide range of VOCs, products of metabolic activity and promising biomarkers for a range of diseases, making breath analysis a rapidly emerging technique employed in non-invasive metabolomics studies. The study of VOCs in breath has been challenging due to the complexity of the sample matrix and the wide concentration range of VOCs in breath. The ability to confidently identify and quantify low-abundance chemicals is crucial in untargeted approaches investigating small metabolic changes affecting a large number of biologically relevant VOCs. In the past, time-of-flight (TOF) analysers have been the mass analysers of choice in discovery studies as their 'full ions transmission all the time' detection mode results in good sensitivity over a wide mass range and their high resolution allows for accurate compound identification. Nonetheless, ion saturation is often observed at higher concentrations due the use of multichannel plate (MCP) detectors, potentially leading to inaccurate quantitation and identification. High-resolution, accurate mass (HRAM) analysers such as the Orbitrap address some of these issues with sub-ppm mass accuracy and high dynamic range by means of innovative technologies such as C-Trap's Automatic Gain Control (AGC). The instrument has extended compound identification capabilities as it is equipped with a variable-electron energy (VeV) electron ionization (EI) and easy to swap positive/negative chemical ionization (CI) source.

## MATERIALS AND METHODS

### **Breath collection**

Successful breath measurements require highly reproducible sampling and analysis techniques. The Breath Biopsy platform includes the ReCIVA Breath Sampler (Figure 1), which was designed to provide a standardized method to collect exhaled breath samples. Breath samples (1500 mL ontube) were collected using the ReCIVA Breath Sampler, which captures and pre-concentrates VOCs onto Breath Biopsy Cartridges (both Owlstone Medical Ltd.). Quality control samples consisted of a custom, 40-compound mixture prepared in methanol (1, 100 and 200 ppm median concentration).

Figure 1. The ReCIVA Breath Sampler (left and middle panel) enables reliable, reproducible collection of breath VOCs and pre-concentration for enhanced sensitivity. Pressure and CO2 sensors in ReCIVA provide real-time monitoring of the patient's breathing, allowing different breath fractions to be sampled in a single collection event if desired (right panel).



### Data Acquisition

Samples were dry purged to remove excess water, desorbed using a TD100-xr thermal desorption autosampler (Markes International) and transferred onto a 30m x 0.32 mm x 3.00 µm column (Quadrex) using splitless injection. Chromatographic separation was achieved via a programmed method on a Thermo Scientific<sup>™</sup> TRACE<sup>™</sup> 1310 GC oven. Mass spectral data were acquired using an Exactive GC Orbitrap mass spectrometer with both variable-EI and CI capabilities. The details of the TD-GC-Orbitrap setup used in this study as well as the experimental parameters are given in Figure 2 and Table 1.

Data Processing Data were processed using Thermo Scientific<sup>™</sup> Xcalibur<sup>™</sup> 4.0 software and Thermo Scientific<sup>™</sup> TraceFinder<sup>™</sup> 4.1 software, which allowed for both qualitative and quantitative analysis.

### **Table 1. Experimental parameters**

TD parameter name	TD parameter value	MS parameter name	MS parameter value
Pre-purge time (min)	0.1	Transfer line temperature (°C)	250
Trap in line (Y/N)	Yes	lonization type	El, PCI (methane)
Pre-purge trap flow (mL/min)	50	lon source temperature (°C)	280 (EI), 180 (PCI)
Tube desorb time (min)	10.0	Electron energy (eV)	35, 70
Tube desorb flow (mL/min)	50	C-Trap voltage (V)	0, 2
Tube desorb temperature (°C)	210	Acquisition mode	Full scan
Trap desorb time (min)	3.0	Mass range $(m/z)$	35 - 350
Trap purge flow (mL/min)	50	Mass resolution (FWHM at $m/z$ 200)	60,000
Trap high (°C)	250	Lock masses ( <i>m/z</i> )	207.03235; 218.05114; 355.06994
GC parameter name	GC parameter value		
Operation mode	Constant flow		
Carrier gas	Helium; 3.00 mL/min		
Temperature ramp (°C)	40 - 250		

## RESULTS



Peak deconvolution results in a features list containing hundreds of entries, which can be either true chromatographic peaks associated to chemicals in breath as well as false positives (i.e. noise). Reduction of the number of false positives is essential for generating high-quality data sets that can be used in untargeted metabolomics approaches. Together with typical spectral matching (e.g. against reference NIST spectra; SI score), TraceFinder calculates High Resolution Filtering (i.e. HRF) scores from high-accuracy EI data that allow for more precise compound identification. Similarly, retention index data via the use of e.g. alkanes results in an additional score (i.e.  $\Delta RI$ ) and leads to improved identification of unknowns (Figure 4).



### Figure 2. TD-GC-Orbitrap setup



### Mass Accuracy and Ion Ratios

The collection of HRAM data is crucial in metabolomics studies where low-concentration analytes are to be detected in a complex samples. Sub-1 ppm mass accuracy is achieved over the full chromatographic peaks (Figure 3). This significantly improves peak deconvolution and helps differentiation of analytes of interest from matrix ions. The high linear dynamic range of the Orbitrap analyser guarantees stable ion ratios even at high sample concentrations, which improves deconvolution, compound identification and the generation of custom libraries.

#### Figure 3. Sub-1 ppm mass accuracy across a chromatographic peak (100 ppm sample; left). Identical fragment ion ratios are observed even at high concentrations (200 ppm sample; right).

### **Deconvolution and Library Matching**

### Figure 4. TraceFinder data processing work flow with an example of true and false feature

Ionization and fragmentation The production of characteristic fragment ion techniques aiding compound identification (e.g. nspectra under EI conditions in GC-MS allows for Tetradecane). the identification of unknown analytes. For closely-related analytes with similar or identical fragmentation, EI data is often insufficient to lead to a conclusive identification. Alternative ionization methods like positive/negative CI can be explored to aid identification of unknowns as they lead to formation of higher m/z (e.g. [M+H+]). These can improve differentiation between compounds such as alkanes and terpenes. More complimentary data sets can be obtained by use of VeV ionisation. Using these methods EI-like fragment spectra are acquired and higher abundance is obtained for molecular ions [M +] (Figure 5).

### Smoking-related markers

A small set of 12 breath samples were analysed on the TD-GC-Orbitrap platform and above-mentioned tools were used for generating a list of reliable features for each sample. Samples were divided into three groups based on the subject's smoking history: never smoked (n=3), current smokers (n=4) and exsmokers (n=5). Based on smoking-related breath markers previously reported in literature, a custom 6compound database was created and used for quantitation of these markers in each of the breath samples. Fold changes relative to the non-smoker group for each marker and sample were calculated using extracted peak area responses.

Figure 6 shows low fold changes (blue Figure 6. Data (n=12) showing high fold changes for 2,5and white hue) for most of the targeted **Dimethylfuran and Toluene in samples from current** smoking-markers in all three groups, suggesting low correlation between smoking behaviour and the reported markers (e.g. ethylbenzene boxplot) High fold changes (red hue) are observed for 2,5-dimethylfuran and toluene in the current group of smokers, suggesting high correlation with smoking behaviour for these two markers (e.g. toluene boxplot). Nonetheless, one sample from the current smokers group gives low fold changes for these two markers.

## CONCLUSIONS

- compound identification and guantitation.
- generating dense, high-quality data sets.

## REFERENCES

- 8328-8335
- Chromatogr. 2018, 32 (4), 1-11.
- Perspect. 2002, 110 (7), 689-698.

## **TRADEMARKS/LICENSING**

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.

PO65784-EN 0422S



Figure 5. Comparison of various ionisation

smokers smoking	and beha	su viou	gge ır.	sts	cori	relat	ion	of	the	mai	'ker	s to
Compound	Never	Never	Never	Current	Current	Current	Current	Ex	Ex	Ex	Ex	Ex
Benzene	0.96	1.59	0.45	4.21	2.40	1.99	0.32	0.30	0.67	0.76	0.54	0.71
2,5-Dimethylfuran	0.55	1.61	0.84	83.74	59.50	42.66	0.74	0.78	0.50	0.32	0.80	1.31
Toluene	0.35	1.49	1.16	13.70	7.29	6.29	0.74	0.94	0.40	0.46	0.38	2.16
Ethylbenzene	0.53	0.91	1.57	2.17	2.28	2.38	2.07	0.71	1.68	1.46	1.09	1.14
m/p-Xylene	0.49	0.85	1.66	1.90	2.06	1.84	2.12	0.75	1.66	1.52	0.96	0.95
o-Xylene	0.49	0.91	1.60	1.84	2.15	1.98	2.08	0.72	1.69	1.51	0.97	1.01
2.50 2.00 1.50 0.50 0.50			NE	<ul> <li>Never</li> <li>Ex</li> <li>Curre</li> </ul>	, nt	Fold Change	90.00 80.00 70.00 50.00 40.00 30.00 20.00 10.00	2,5-D	IMETH		RAN	Never Ex Current

High mass accuracy, routine 60,000 resolving power and dynamic range improve peak deconvolution.

· Retention time alignment and High Resolution Filtering can be used to curate feature lists and

• Variable-EI at low/high energies together with chemical ionisation can be used to gain further insight into the chemical structure of VOCs, especially when studying compound classes showing similar EI data • Breath data shows high fold changes for two smoking-related markers, suggesting a high correlation between these markers and the smoking habits of the studied Subjects

1. N. W. Kwiecien et al. High-Resolution Filtering for Improved Small Molecule Identification via GC-MS. Anal. Chem. 2015, 87, 2. S. Capone et al. Chromatographic analysis of VOC patterns in exhaled breath from smokers and nonsmokers. Biomed. 3. S. M. Gordon et al. Volatile Organic Compounds as Breath Biomarkers for Active and Passive Smoking. Environ. Health 4. M. Castellanos et al. 2,5-dimethylfuran as validated biomarker of smoking status. Nicotine Tob. Res. 2018, Epub ahead of print.





