#### **Thermo Fisher** s c i e n t i f i c

## High throughput analysis of synthetic DNA using a compliant LC-MS based workflow

#### Andrew Williamson<sup>3</sup>

Angela Criscuolo<sup>1</sup>; Tabiwang N. Arrey<sup>2</sup>; Ken Cook<sup>3</sup>; Keeley Murphy<sup>4</sup>; Jennifer Sutton<sup>4</sup>; Cheryl Rhodick<sup>5</sup>; Melissa N. Sato<sup>6</sup>; Alexander Harder<sup>7</sup>

<sup>1</sup>Thermo Fisher Scientific GmbH, Dreieich, Germany; <sup>2</sup>Thermo Fisher Scientific, Bremen, Germany; <sup>3</sup>Thermo Fisher Scientific, Hemel Hempstead, United Kingdom; <sup>4</sup>Thermofisher Scientific, San Jose, CA; <sup>5</sup>Thermo Fisher Scientific, Paisley, United Kingdom; <sup>6</sup>Thermo Fisher Scientific, San Jose, CA; <sup>7</sup>Thermo Fisher Scientific (Bremen) GmbH, Bremen, Germany

The world leader in serving science



### Abstract

• The advances in high-speed chromatography, mass spectrometry, and software have been used in this study to develop high throughput workflows utilizing quantitation by mass spectrometry detection to match this growing requirement. In this abstract we present the analysis of synthetic oligonucleotides with respect to product identification, purity, and speed of analysis in a compliant environment.

# Introduction

#### **Primary Challenge**

- Throughput limitations due to LC method overhead times can account for up to 50% of the total analysis time
- Additional software is often required for data analysis

#### Novel approach

Rapid analysis of
oligonucleotides and
impurities using a dual UHPLC
system coupled to a HRMS
with automated sequence
annotation in a compliant
ready software.

#### **Overview content Summary**

An oligonucleotide method
confirming product identity in a
rapid-fire configuration, impurity
analysis and desalting using a
fast gradient in a tandem column
configuration, and high
throughput sequencing with
MS/MS enabled was developed.

## **Materials and Methods**

Sample prep

 5pmol of each oligonucleotide sample was analyzed using either a high throughput method or

chromatographically separated to perform desalting and impurity analysis, using a tandem column workflow.

#### LC- Parameters

The Oligonucleotide samples
were separated on a Thermo
Scientific<sup>™</sup> DNAPac RP
column using a Thermo
Scientific<sup>™</sup> Vanquish<sup>™</sup> Duo
UHPLC system. The columns
were maintained at 60° C,
with a flow rate of 400 µL/min.

MS and Data Analysis methods

•

The eluting analytes were measured on the Thermo Scientific<sup>™</sup> Orbitrap Exploris<sup>™</sup> 240 mass spectrometer in negative ion mode. Data analyses and system control were performed with Thermo Scientific<sup>™</sup> Chromeleon<sup>™</sup> 7.2.10 CDS, and any additional sequence data analysis was performed with Thermo Scientific<sup>™</sup> BioPharma Finder<sup>™</sup> 4.1 software.

# **Results: standardization and productivity**



- Rapid injections of 30 sec
- Short gradients possible or isocratic rapid injections
- Charger system available for long unattended sequences

### Maximizing productivity

• Dual tandem LC operation to re-equilibrate one set of columns while the other set is running the sample gradient



# **Results: standard methods and productivity**



- Stepped Normalized Collision Energy: 10-12-14 to 20-22-24 (18-20-22 for impurity analysis)
- High resolution MS and MS/MS

#### Maximizing productivity

- The MS only method was used for the rapid analysis of ٠ oligonucleotides with intact deconvolution.
- The ddMS2 method was used for sequence analysis of • oligonucleotides and was used on QE Plus, orbitrap Exploris 120 and 240
- Enables the identification and relative quantification of • oligonucleotides and their impurities, even those present at very low levels, in a single experiment.
- Easy to do with Chromeleon software for compliant intact • analysis and BioPharma Finder for automated annotation and identification with sequence analysis

### **Results: precision with accurate identification in BioPharma Finder software**

60

9.48

1.14

1.7

1.9

Quality of results at 120,000 resolution while maintaining sensitivity to see low abundant species 20200713 A1 21mer 0 1µL LowP A1 6288.088 CCACTCTGTTCTACTTAAATC 100 Matched Mass Error Modification Theoretical Mass (Da) Relative Abundance equence Name Monoisotopic Mass 80 • • • Relative Intensity = (NonBla... 🔻 🏹 🗛 = • A1 6288.088 6288.078 1.7 100.00

A1

A1

1xCnet(ps)

1xAdenine loss(A)

6341.115

6153.035

6341.104

6153.023



# **Results: precision with isotopic resolution in Chromeleon**

Quality of results at 120,000 resolution with UV and HRMS data collection in Chromeleon

TACTGGACCACCTGGCATCAAAGACAACTTTTCAGAGC 38mer



# **Results: sequence analysis**



### Reporting

.

- Target mass is compared with experimental mass to give a "pass/Fail" with Chromeleon in a compliant environment
- Can include fractional abundance •
- Component identification •
- UV and HRMS signal reported
- Oligonucleotide sequencing with automated annotation within . **BioPharma Finder software**

jectio	n Details						
).	Injection Name	Expected mass	Target Mass is most abundant component	TargetAccuracy	Fractional Abundance	Result Component Count	Calculated Mass
		Da			%		Da
	Sample D11	10802.0	Pass	5	73	5	10805.0
	Sample F04	13839.9	Pass	5	52	9	13843.9
	Sample H07	18394.9	Fail	5	28	19	18400.4
			1				1

### **Thermo Fisher**

### **Conclusions**

We developed high-throughput methods using the Vanquish Duo UHPLC system with DNAPac RP columns coupled with Orbitrap Exploris mass spectrometers. Rapid fire methods of 30 seconds are possible or tandem gradients for desalting and impurity separations. Using high resolution Orbitrap mass spectrometry methods, isotopic resolution can be achieved while maintaining the sensitivity required to confidently identify low abundant species.

MS/MS fragmentation methods allow sequence confirmation of oligonucleotides up to 100nt with automated annotation within BioPharma Finder software.

Rapid oligonucleotide primer QC testing is possible in a compliant environment using Chromeleon software for instrument control of the dual UHPLC and HRMS Orbitraps. Oligonucleotide deconvolution and "Pass/Fail" reporting within the same compliant software package.

#### **TRADEMARKS/LICENSING**

© 2021 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. PO66128