thermo scientific

High-throughput analysis of oligonucleotides using a single quadrupole mass spectrometer for quality control

Dennis Köhler, Mauro De Pra Thermo Fisher Scientific, Germering, Germany

ABSTRACT

Purpose: Show step-by-step the analysis, deconvolution, and reporting of oligonucleotide synthesis quality control with a single quadrupole mass detector.

Methods: Synthesized single length oligomers were analyzed without post synthesis purification. Thermo Scientific DNAPac RP (2.1x50mm, 4µm) was run on a Thermo Scientific[™] Vanguish[™] Flex Binary HPLC system with UV detection.

Results:

A step-by-step workflow including the analysis, deconvolution, and reporting of oligonucleotide synthesis quality control with a single quadrupole mass detector.

INTRODUCTION

Laboratories producing large arrays of customized DNA need to support this heightened throughput via increased automation and accuracy using intact mass determination for quality control. With this workflow from robotic DNA synthesis all the way through a confident pass/fail outcome for the expected sequence, Thermo Scientific[™] offers a complete package consisting of the Thermo Scientific[™] Vanquish[™] Flex UHPLC using the Thermo Scientific[™] DNAPac[™] RP column for the separation. Determination of the intact oligonucleotide mass uses the ISQ[™] EM Single Quadrupole Mass Spectrometer and is interpreted using the Thermo Scientific[™] Chromeleon[™] Chromatography Data System (CDS) with the inclusion of the Intact Protein Deconvolution (IPD) engine and oligonucleotide analysis capabilities. Minor method optimizations provide cost savings and the reduction of 1,1,1-3,3,3-hexa-fluoro-isopropanol (HFIP) and sodium adduct abundancy.

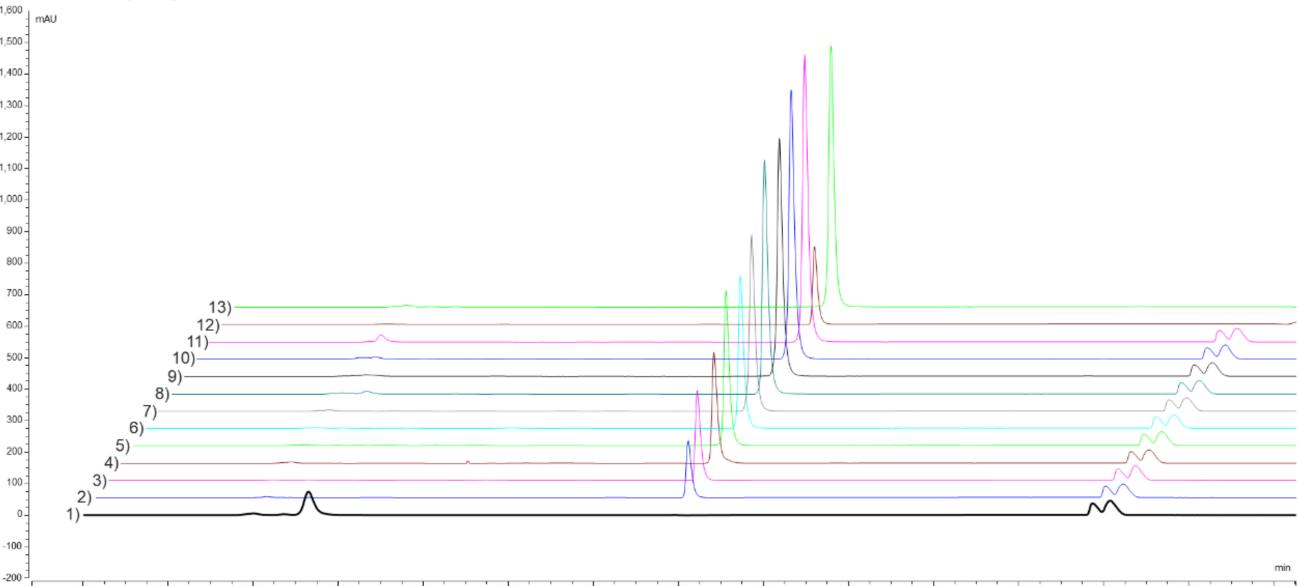
The vaporizer temperature, transfer tube temperature, sheath gas/auxiliary gas pressures, and spray voltage were optimized by maximizing the peak area associated with the most abundant charge state. The instrument source settings were optimized at the beginning of experiments using Custom Injection Variables in Chromeleon CDS in Figure 1. This order of optimization is represented in Table 4. It is important to note that the auxiliary gas pressure was always 10% that of the sheath gas pressure. Subsequently, the HFIP concentration was modified to improve the quality of the spectra. Finally, the source settings were optimized again at the new HFIP concentration.

Table 4. Variable source parameters in MS setting tuning

| | | <u>_</u> | |
|-------|----------------------------|---|----------------------|
| Order | Source Parameter | Optimization Range | Increments |
| 1 | Vaporizer temperature | 300 to 450 °C | 50 °C |
| 2 | Transfer tube temperature | 300 to 400 °C | 50 °C |
| 3 | Sheath gas (auxiliary gas) | 50 to 80 psig (5 to 8 psig; 10% of sheath gas) | 5 psig (0.5 psig) |
| 4 | Spray voltage | -1,000 to -5,000 V | 1,000 V |
| 4 | Spray voltage | -1,000 to -5,000 V | 1,000 V |

Using the optimal HFIP concentration of 0.1% and given in the LC method conditions presented in Table 2, the following chromatographic overlays represented in Figure 3. The results represented by the traces show the elution of the oligomers without the separation of impurities such as the N-1, N-2, N-3, etc. but removing all extraneous synthesizing reagents present during the oligomer synthesis. One can observe that a failed synthesis occurred, like the 10mer seen in chromatogram 1 (black), where the expected oligomer peak is absent.

Figure 3 UV chromatograms for the oligonucleotide array provided in Table 1 ranging from 10mer to 60mer.





MATERIALS AND METHODS

Sample Preparation

The samples were provided in a 96-well plate. They were collected directly from the DNA synthesizer and were injected neat.

Table 1. Oligomer sample array provided by GeneArt AG (part of Thermo Fisher Scientific), Regensburg, Germany. All oligomers are 10 mM in water and were not desalted.

| Oligo number | Sequence | Length (nt) | Theoretical average mass (Da) |
|-----------------|--|----------------|----------------------------------|
| 1 | AAGCCAGAGC | 10 | 3206.0 |
| 2 | CAATCTAAAGTATAT | 15 | 4559.0 |
| 3 | TCTCCCGGACGGAAACCGCC | 20 | 6047.9 |
| 4 | AGGTAATTTCGCCTCATTGGGGGCC | 25 | 7689.0 |
| 5 | CCGGCCTATGGCCCACAATGTAAAGAATTA | 30 | 9184.0 |
| 6 | GCCCGTGGTAAAGCAGTTCACGTGTACATAGTTGT | 35 | 10802.0 |
| 7 | GCCCATAATTGAGCCCCGCTGCCGACGAGCGGCTTTGTGC | 40 | 12249.9 |
| 8 | CCCTGAATTAAGGGGGGCAGCCCCTTAATGAATGCCCGGACTCGAA | 45 | 13839.9 |
| 9 | TAAACTGTTTATCGGGGGCTCAAATCTTAGGCCTAGGCAGGATCCCGTAAG | 50 | 15425.0 |
| 10 | ATAATCGAGAATTGGTATCGATTCGGGGGCCACCCACAAGTCCGGTACACCAACCG | 55 | 16897.9 |
| 11 | CACACCTCGAAGAGTATTCCGTCCCGGAGCTGGTTAGGTGACTAACACTGCAAATTCTCT | 60 | 18394.9 |
| 12 | GGGGCGCTCTATCTTCCATC | 20 | 6059.9 |
| 13 | CCCGAGCGGAGTTTTGCGATAGTACACCAACCGAGCATCTCGAATTAAAGGCCTG | 55 | 16928.9 |

Instrumentation

Thermo Scientific[™] Vanquish[™] Flex Binary HPLC System with Vanquish Variable Wavelength Detector F and ISQ EM single quadrupole mass detector.

Figure 1. Inserted custom variables are as follows: VaporizerTemp (orange), TransferTubeTemp (blue), SheathGas (purple), SprayVoltage (yellow).

| # | UV_VIS_1 | Name | Position | Volume [µl] | *VaporizerTemp [°C] | *TransferTubeTemp [°C] | *SheathGas [psig] | *SprayVoltage [V] | Instrument Method |
|----|----------|---|----------|-------------|---------------------|------------------------|-------------------|-------------------|-----------------------------|
| 1 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 300 | 300 | 75 | -3000 | HFIP Method v16 - ISQ Scout |
| 2 | l ma | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 350 | 300 | 75 | -3000 | HFIP Method v16 - ISQ Scout |
| 3 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 400 | 300 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 4 | l. | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 450 | 300 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 5 | | Sample G10 - 260nm with ISQ - 2μL injection - 55mer | G:F2 | 2.00 | 300 | 350 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 6 | I. | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 350 | 350 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 7 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 400 | 350 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 8 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 450 | 350 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 9 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 300 | 400 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 10 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 350 | 400 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 11 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 400 | 400 | 75 | -3000 | HFIP Method v16 - ISQ Scou |
| 12 | | Sample G10 - 260nm with ISQ - 2µL injection - 55mer | G:F2 | 2.00 | 450 | 400 | 75 | -3000 | HFIP Method v16 - ISQ Scou |

Chromatography Data System

Chromeleon 7.3 CDS was used for data acquisition and analysis.

The ISQ EM mass spectrometer is fully integrated into Chromeleon software, which was used for system operation, subsequent data analysis, and deconvolution using the integrated Intact Protein Deconvolution (IPD) feature. This feature is also intended for oligonucleotides specifically with the negative charge and peak model setting (Table 5). The obtained MS chromatograms were analyzed with the IPD settings shown in Table 5.

Table 5 Intact Protein Deconvolution settings

| | | Jottingo | |
|-----------------------|-------------------|---------------------------|-------|
| Parameter | Value | Parameter | Value |
| Peak retention window | 0.7-0.8 min | Low number adjacent | 3 |
| Algorithm | ReSpect™ | charges | |
| Output mass range | 2000-20000 Da | Intensity threshold scale | 0.01 |
| Deconvoluted spectra | Isotopic Profile | Min peak significance | 1 |
| display mode | | Negative charge | True |
| Model mass range | 2000-20000 Da | Noise compensation | True |
| Deconvoluted Mass | 100 ppm | Noise rejection | 95 |
| Tolerance | | Number of peak models | 1 |
| Peak model | Nucleotide | Peak model width scale | 1 |
| Resolution | Raw File Specific | Quality score threshold | 0 |
| Charge carrier | H+ | Relative abundance | 0 |
| Charge high | 30 | threshold | |
| Charge low | 1 | Target peak mass | 20000 |
| High number adjacent | 3 | Target peak shape left | 2 |
| charges | | Target peak shape right | 2 |

After the entire oligomer array was analyzed with the optimized HFIP concentration, LC method, and MS settings, data was analyzed using including the intact mass deconvolution, mass confirmation, and report. Using the deconvolution settings (Table 5) oligomer array spectra were analyzed for their respective intact masses (Table 1). The measured intact mass was then compared to the expected mass. This is performed with the Custom Injection Variables where the expected intact mass of the target oligomer and target mass accuracy is defined by the user within the injection sequence (Figure 5). The confirmation that the measured mass matched the expected mass within the specified target mass accuracy was automatically visualized as a pass/fail result in the sequence report (Figure 5).

Figure 4 Example of intact mass deconvolution using the 55-mer (sample 13). The identified charge states are overlaid to the original MS spectrum.

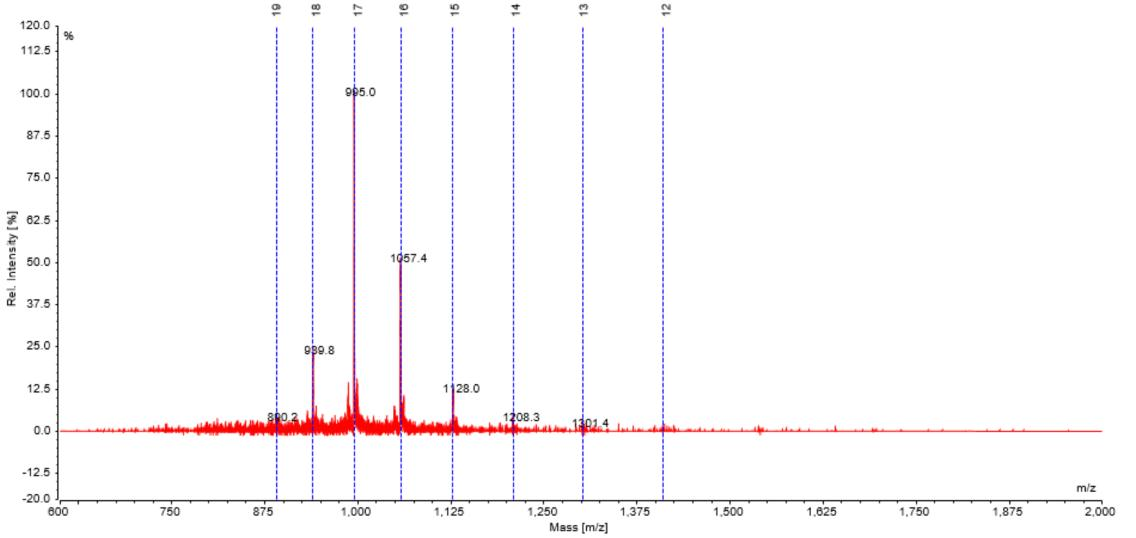


Table 2. Chromatographic conditions

| Column | Thermo Scientific DNAPac RP 2.1 × 50 mm, 4 µm (P/N 088924) | | | | | | | |
|-------------------------|---|----|-----|--|--|--|--|--|
| Flow rate: | 0.70 mL/min | | | | | | | |
| Mobile phase: | A: HFIP (0.01, 0.1, 0.5, 1.0, 2.0%), 0.1% TEA, in water | | | | | | | |
| | B: HFIP (0.01, 0.1, 0.5, 1.0, 2.0%), 0.1% TEA, in MeOH | | | | | | | |
| | Time (min) | %A | %B | | | | | |
| | 0.0 | 99 | 1 | | | | | |
| | 0.4 | 99 | 1 | | | | | |
| | 0.4 | 75 | 25 | | | | | |
| Gradient: | 1.0 | 75 | 25 | | | | | |
| | 1.0 | 0 | 100 | | | | | |
| | 1.6 | 0 | 100 | | | | | |
| | 1.6 | 99 | 1 | | | | | |
| | 4.0 | 99 | 1 | | | | | |
| | 70 °C, forced air mode | | | | | | | |
| Column temperature: | 70 °C, active pre-heater | | | | | | | |
| Injection volume: | 2 µL | | | | | | | |
| UV detector parameters: | λ=260 nm, 100 Hz | | | | | | | |

Table 3. MS Settings: Instrument and scan settings for the mass spectrometer used for the final sample analysis

| HESI Source Se | ttings | Scan S | ettings |
|-------------------------------|----------|--------------------|--------------|
| Vaporizer temperature | 350 °C | Mass range | 600-2000 m/z |
| Ion transfer tube temperature | 350 °C | Dwell/Scan Time | 0.5 s |
| Source voltage | -3000 V | Polarity | Negative |
| Sheath gas pressure | 75 psig | Spectrum Type | Profile |
| Aux gas pressure | 7.5 psig | Source CID voltage | 0 V |

RESULTS

Reversed-phase ion pairing chromatography was performed on the oligonucleotides. The method scope was to clean-up the sample from salt and other reagents and elute the target oligonucleotide and related impurities as single peak. Initial experiments focused on testing HFIP concentrations of 0.01, 0.1, 0.5, 1.0, and 2%. As seen in Figure 2, the HFIP concentration was incrementally increased from 0.01% to 2% to maximize oligo peak area and minimize HFIP adduction. The industry standard is 2%. For the ISQ EM, it was found that the adduct abundancy versus the maximum spectral intensity was the greatest at 0.1% HFIP which yielded the lowest HFIP adduct relative abundancy and the largest maximum charge state's intensity. This 20x reduction of HFIP usage has a notable cost-saving impact as well.

Figure 2 Impact of HFIP concentration on adduct abundance and signal intensity.

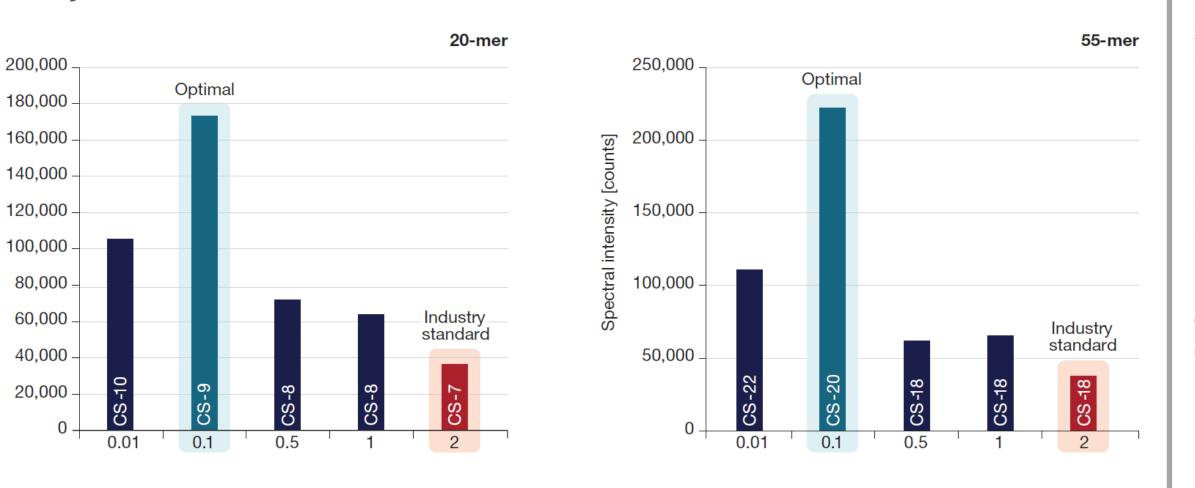


Figure 5 Expected Mass and Target Mass Accuracy with accompanying report. On the left using Custom Injection Variables in Chromeleon CDS allow the user to enter the expected target mass of the oligomer and define the target mass accuracy. This report template (right) confirms with an easy-to-read pass/fail result for the presence of the target mass. Red text "No Match": expected mass does not match any of the five most abundant deconvoluted masses. Green text "Yes, Most Abundant": expected mass matches the most abundant deconvoluted mass.

| | n Finished submit ▼ e ③ Studio ④ Print ▼ mb Up 3≈ Inse | rt Row 👻 🎛 Fill Dow | n 🔒 Lock 🛛 🍸 Filt | ering 🔚 Grouping $f_{\!X}$ | Custom Columns - | 🔻 🏭 Find Next 👻 | Inj. No. | Oligonucleotide Name | Position | TargetA ccuracy | ExpectedMass | Matches IPD Component? | Measured Mas |
|---|---|---------------------|---------------------|------------------------------|------------------------|--------------------------------------|----------|----------------------|----------|--------------------|--------------|------------------------|--------------|
| ‡ | UV_VIS_1 Name | *ExpectedMAss [Da] | *TargetAccuracy [Da |] Position Volum | [µl] Instrument Method | Processing Method | | | | Da | Da | | Da |
| ٠ | Sample A11 - 15mer | 4559.0 | 5 | G:A3 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 1 | 1 | G:A2 | 10.0 | 3206 | No Match | 12658.2 |
| | Sample A11 - 15mer | 4559.0 | 5 | G:A3 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 2 | 1 | G:A2 | 10.0 | 3206 | No Match | 12951.8 |
| | Sample A11 - 15mer | 4559.0 | 5 | G:A3 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 3 | 1 | G:A2 | 10.0 | 3206 | No Match | 8378.4 |
| | Sample B08 - 20mer | 6047.9 | 5 | G:A4 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 4 | 2 | G:A3 | 10.0 | 4559 | Yes, Most Abundant | 4559.8 |
| | Sample B08 - 20mer | 6047.9 | 5 | G:A4 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 5 | 2 | G:A3 | 10.0 | 4559 | Yes, Most Abundant | 4559.8 |
| | Sample B08 - 20mer | 6047.9 | 5 | G:A4 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 6 | 2 | G:A3 | 10.0 | 4559 | Yes, Most Abundant | 4560.0 |
| | Sample C08 - 25mer | 7689.0 | 5 | | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 7 | 3 | G:A4 | 10.0 | 6047.9 | Yes, Most Abundant | 6049.6 |
| | Sample C08 - 25mer | 7689.0 | | | | | 8 | 3 | G:A4 | 10.0 | 6047.9 | Yes, Most Abundant | 6049.2 |
| | × . | | | | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 9 | 3 | G:A4 | 10.0 | 6047.9 | Yes, Most Abundant | 6049.2 |
| | Sample C08 - 25mer | 7689.0 | 5 | G:A5 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 10 | 4 | G:A5 | 10.0 | 7689 | Yes, Most Abundant | 7690.6 |
| | Sample D06 - 30mer | 9184.0 | 5 | G:A6 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 11 | 4 | G:A5 | 10.0 | 7689 | Yes, Most Abundant | 7690.7 |
| | Sample D06 - 30mer | 9184.0 | 5 | G:A6 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 12 | 4 | G:A5 | 10.0 | 7689 | Yes, Most Abundant | 7690.7 |
| - | Sample D06 - 30mer | 9184.0 | 5 | G:A6 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 13 | 5 | G:A6 | 10.0 | 9184 | Yes, Most Abundant | 9186.2 |
| | Sample D11 - 35mer | 10802.0 | 5 | G:A7 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 14 | 5 | G:A6 | 10.0 | 9184 | Yes, Most Abundant | 9186.4 |
| | Sample D11 - 35mer | 10802.0 | 5 | | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 15 | 5 | G:A6 | 10.0 | 9184 | Yes, Most Abundant | 9186.2 |
| _ | × | | | | | | 16 | 6 | G:A7 | 10.0 | 10802 | Yes, Most Abundant | 10805.2 |
| | Sample D11 - 35mer | 10802.0 | 5 | G:A7 | .00 Oligo QC Method | Oligonucleotides ISQEM deconvolution | 17 | 6 | G:A7 | 10.0 | 10802 | Yes, Most Abundant | 10805.0 |

CONCLUSIONS

This work provides a complete workflow for the analysis of oligonucleotides via a high-throughput robust LC method, intact targeted mass confirmation, and a user-friendly report confirming that the expect oligonucleotide has been synthesized. The following features are included with this workflow:

• Optimal ISQ EM spectra quality is observed with 0.1% HFIP, much below the concentration typically found in the literature of 2% HFIP. Therefore, it reduces the consumption of HFIP by a factor of 20. In the case that 192 samples are run per day, a year's savings could amount to over \$3,500 in HFIP consumption.

Reduction of HFIP adducts and no sodium adducts are observed.

• Samples are collected directly from the DNA synthesizer and injected neat. No sample preparation is needed.

• The ISQ EM parameters have been optimized for oligomers in the range 10-60 chain lengths.

• Suggested deconvolution parameters provide for a reliable and automated recognition of the oligomer mass. For oligomers with mass outside the described range and/or different spectra quality, different parameters for the deconvolution method may be required

Quality control laboratories screening large arrays of synthesized oligonucleotides can now, with a high level of confidence, easily confirm the quality of their oligonucleotide syntheses.

Trademark/Licensing



© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific Inc. products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others

