

Poster Reprint

ASMS 2024
Poster number TP 556

A Novel and Highly Effective LC/MS Method for Phosphorodiamidate Morpholino Oligomer (PMO) Impurity and Sequencing Analysis

Guannan Li¹, Peter Rye¹ and Vincent Guerlavais²

¹Agilent Technologies, Inc., Santa Clara, CA

²Sarepta Therapeutics, Inc., Cambridge, MA

Introduction

Phosphorodiamidate morpholino oligomers (PMOs) are a type of Antisense Oligonucleotides (ASOs), a rapidly growing class of RNA modalities. Currently, four ASOs based on PMO chemistry are approved by the FDA.^{1,2} These PMOs possess a non-ionic backbone in which the ribose is replaced by a morpholine moiety and the phosphorodiester intersubunit bonds are replaced with phosphorodiamidate linkages (Figure 1). The PMO structure confers distinctive properties compared to other antisense strategies.^{3,4} These structural differences also make PMO characterization by LC/MS unique to other oligonucleotides. In this work, we describe the characterization of PMOs by LC/MS using a unique method. Purity analysis, impurity identification and PMO sequence confirmation are described.

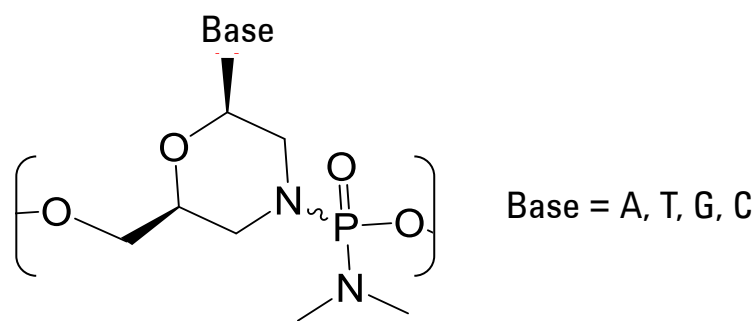


Figure 1. Chemical Structure of Phosphorodiamidate morpholino oligomers (PMOs)

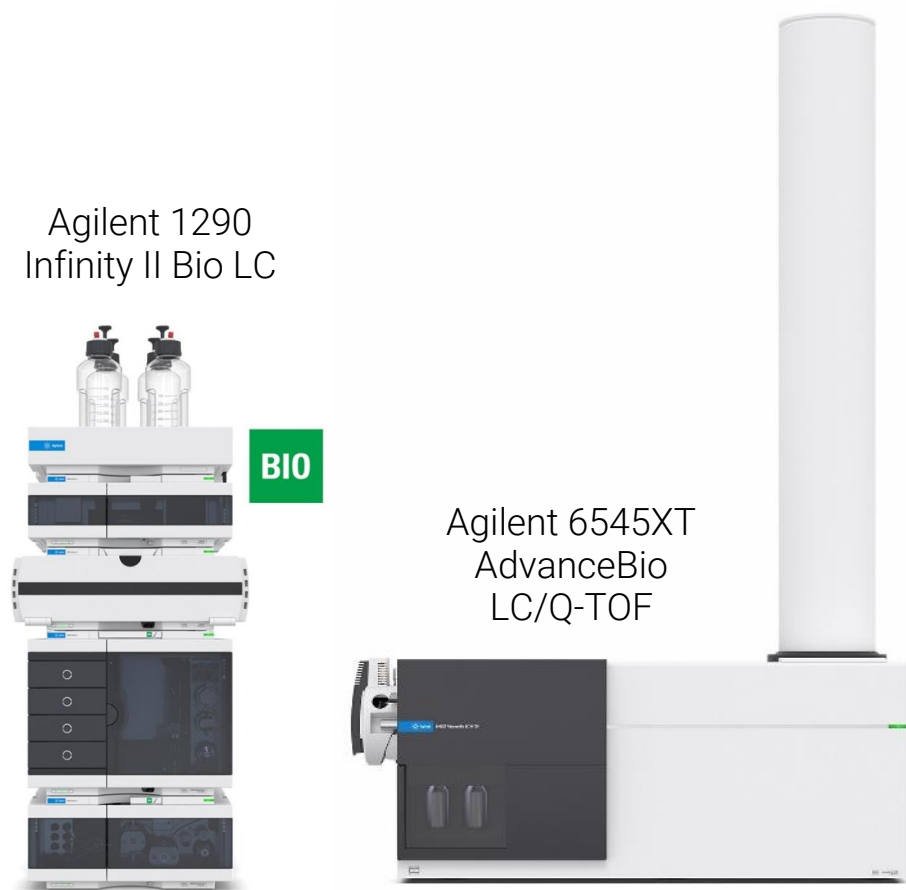


Figure 2. LC/MS Configuration

Experimental

Oligonucleotide Samples Analysis

LC/MS analyses were conducted on a 1290 Infinity II Bio LC system coupled with a 6545XT AdvanceBio LC/Q-TOF equipped with an Agilent Dual Jet Stream ESI source (Figure 2). LC separation was obtained with a ZORBAX RRHD 300Å StableBond C3 column, 1.8 µm, 2.1 x 50mm at 40 °C. For characterization and impurity analysis, data was acquired with a negative MS spectrum acquisition from 400 to 3200 *m/z*. For sequence confirmation analysis, data was acquired with targeted MS/MS mode on -3 charged ion. (Table 1). The resulting data were processed in Agilent MassHunter BioConfirm software 12.1 using both targeted and untargeted methods.

Table 1. LC/MS method used in the study

| Agilent 1290 Infinity II Bio LC Conditions | | |
|--|---|-------|
| Column | ZORBAX RRHD 300Å StableBond C3 column, 1.8 µm, 2.1 x 50 mm (pn: 857750-909) | |
| Injection volume | 1 or 2 µL | |
| Mobile phase | A = Water + 5 mM Ammonium Acetate B = Acetonitrile | |
| Flow rate | 0.4 mL/min | |
| Gradient program | Time (min) | B (%) |
| | 0.00 | 5 |
| | 4.00 | 90 |
| Stop time | 4.00 min | |
| Post time | 1.50 min | |

| 6545XT AdvanceBio LC/Q-TOF Source Conditions | |
|--|-------------------|
| Ion Polarity | Dual AJS Negative |
| Gas temperature | 325 °C |
| Drying gas flow | 12 L/min |
| Nebulizer gas | 35 psi |
| Sheath gas temperature | 350 °C |
| Sheath gas flow | 12 L/min |
| Capillary voltage | 3500 |
| Nozzle voltage | 2000V |
| Fragmentor | 175 V |

Experimental

Table 1. LC/MS method used in the study (continued)

| 6545XT AdvanceBio LC/Q-TOF Acquisition Conditions | |
|---|--------------------------|
| Characterization and Impurity Analysis | |
| Mass Range | 400 – 3200 <i>m/z</i> |
| Acquisition Rate | 1 spectra/sec |
| Sequence Confirmation | |
| MS Mass Range | 400 – 3200 <i>m/z</i> |
| MS Acquisition Rate | 4 spectra/sec |
| MS/MS Mass Range | 100 – 3200 <i>m/z</i> |
| MS/MS Acquisition Rate | 1 spectra/sec |
| Isotope Width | Medium (~ 4 <i>m/z</i>) |
| Collision Energies | 55 V, 60 V, 65 V or 70 V |
| Targeted Mass | 2068.7238 <i>m/z</i> |
| Retention Time | 2.16 min |
| Delta Retention Time | 1 min |

Results and Discussion

Characterization and Impurity Analysis

Figure 3 shows the raw spectrum of PMO (200 ng on column), and the -3 charged ion dominated.

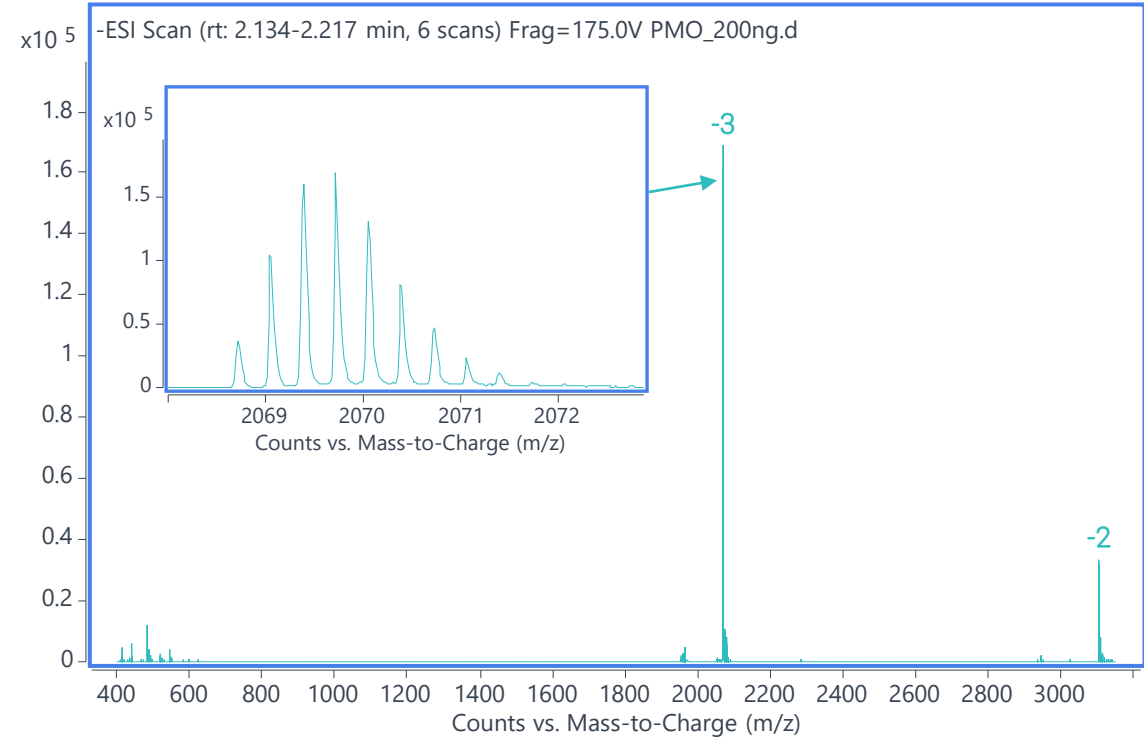


Figure 3. Raw spectrum for PMO.

Results and Discussion

Characterization and Impurity Analysis

Figure 4 shows the deconvolution result of PMO, and mass accuracy of PMO was 0.86 ppm. Multiple low-abundance impurities were identified: A deletion, C deletion and T deletion.

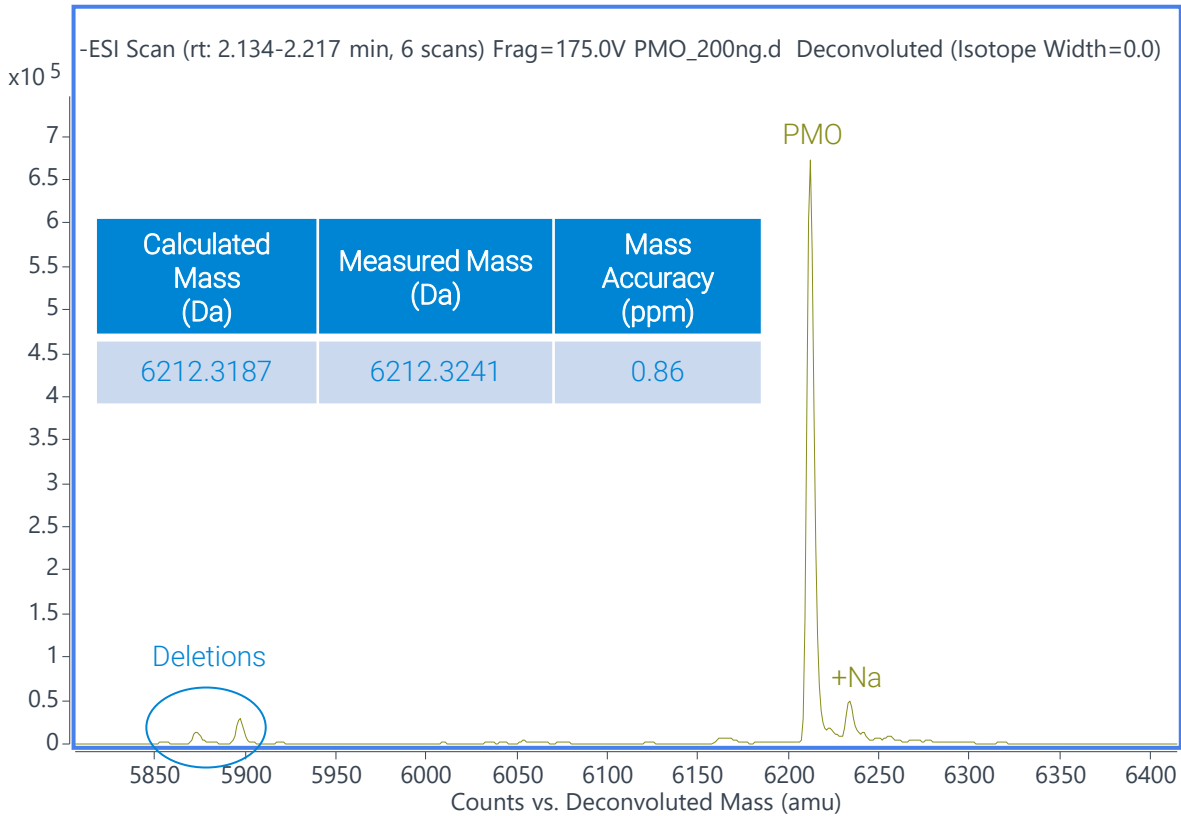


Figure 4. The deconvoluted spectrum for PMO.

Sequence Confirmation Analysis

For sequence confirmation analysis, targeted MS/MS was conducted on -3 charge state precursor (Figure 3) using 4 different collision energies.

Table 2. MassHunter BioConfirm method parameters used for PMO sequence confirmation analysis

| Agilent MassHunter BioConfirm 12.1 Parameters | |
|---|--|
| Workflow | Oligonucleotides |
| Experiment | Sequence confirmation |
| Match Tolerance | Tolerance: 5 ppm Theoretical profile relative abundance ≥ 12% |
| Absolute Height Threshold | 350 |
| Matching Criteria | Warn if score is < 90 Do not match if score is <85 |
| Extraction MS/MS | Group by collision energy 2 scans averaged |

Sequence Confirmation Analysis

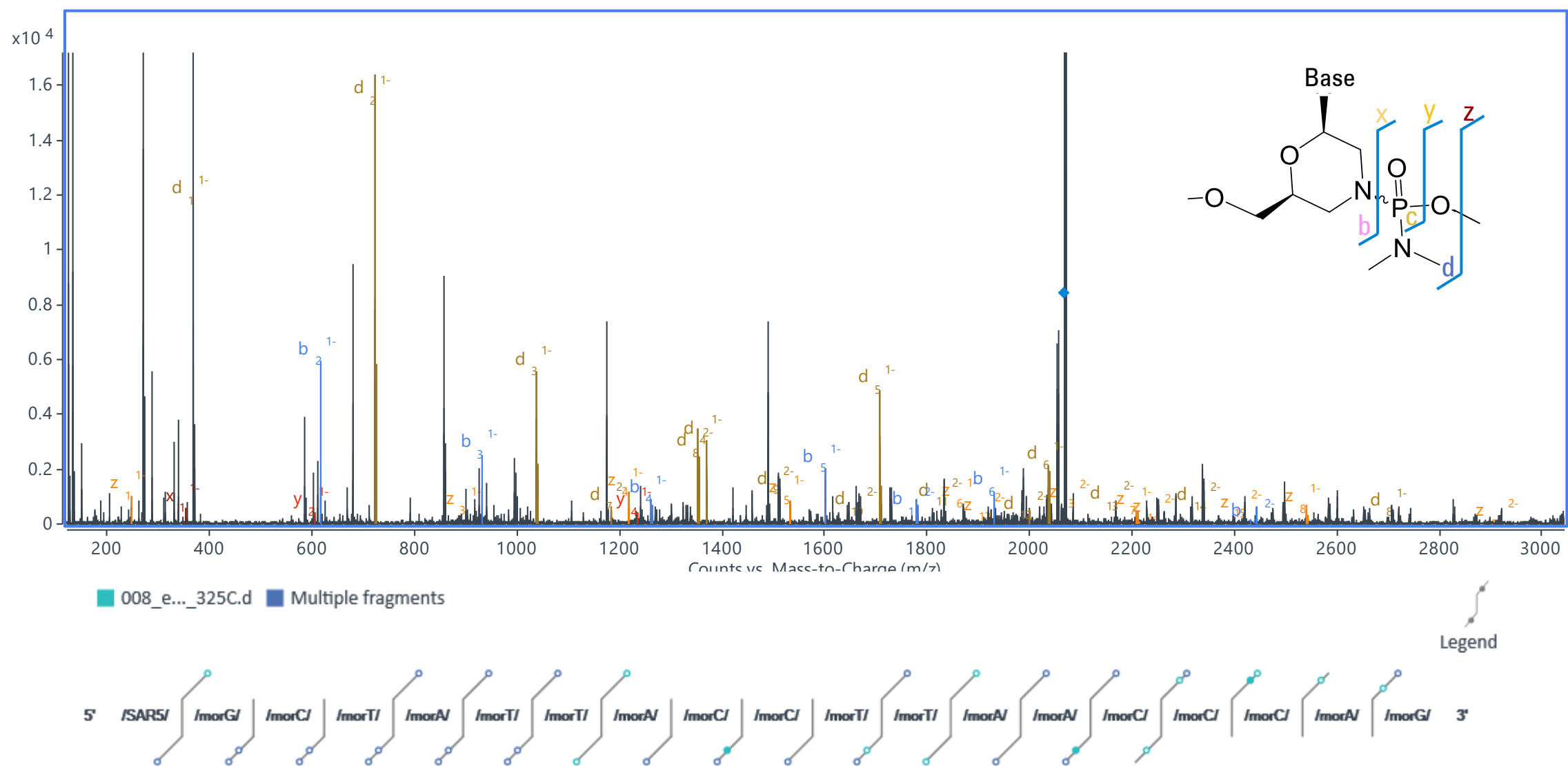


Figure 6. PMO sequence confirmation data. 100% sequence coverage achieved.

Complete sequence coverage was achieved in Agilent MassHunter BioConfirm software (Table 2) using the data from 2 μ L injection of 0.2 mg/mL PMO sample (Figure 6). From the fragment confirmation ladder in Figure 6, b, d and z ions dominated. Based on PMO structure, a and w ions should not be seen.

Conclusions

- This work describes an original LC/MS method for PMO impurity and sequencing analysis.
- Excellent mass accuracy for PMO was achieved.
- Multiple low-abundance impurities were identified.
- 100% sequence coverage was obtained with a single injection.

References

¹Oligonucleotide Therapeutics: From Discovery and Development to Patentability; Pharmaceuticals. 2022, 14 (2), 260,
²Next steps for the optimization of exon therapy for Duchenne muscular dystrophy, Expert Opin Biol Ther. 2023, 1-11,
³Pharmacokinetics and biodistribution of phosphorodiamidate morpholino antisense oligomers; Curr Opin Pharmacol. 2005, 5(5), 550-5.
⁴Invention and Early History of Morpholinos: From Pipe Dream to Practical Products, Methods Mol Biol. 2017, 1565, 1-15.

<https://www.agilent.com/en/promotions/asms>

This information is subject to change without notice.
DE15655762