

Poster Reprint

ASMS 2026
WP 589

An Integrated Software Ecosystem for End-to-End Oligonucleotides DataAcquisition, Analysis, and Reporting

Guannan Li, Stephen Sciuto, Jordy Hsiao, Linfeng Wu, David
Wong, Ruoji Luo

Agilent Technologies, Santa Clara, California

Introduction

Synthetic oligonucleotides, including small interfering RNA (siRNA), antisense oligonucleotides (ASO), guide RNA (gRNA) for CRISPR/Cas, and aptamers, have been emerging as therapeutics against a wide range of disease conditions.

Comprehensive characterization of oligonucleotide samples can be challenging and, in some cases, time-consuming, as it not only requires the profiling of the target oligonucleotide but also identifying and quantifying all related impurities.

This work focuses on developing a comprehensive software ecosystem that enables end-to-end analysis of siRNA for sequence confirmation and impurity profiling.

The results demonstrate the software’s ease of use and its capability to provide a complete solution for analyzing synthetic oligonucleotides.

Experimental

Two siRNA molecules, inclisiran and givosiran were used to evaluate the software’s performance (Table 1). They were dissolved in water without further purification. Ion-pair reversed-phase liquid chromatography coupled with time-of-flight mass spectrometry, along with novel inert column hardware, was employed to acquire data for both denatured and non-denatured forms. The software was implemented

Experimental

in a client–server architecture to enable end-to-end data acquisition, analysis, and reporting. The LC and Q-TOF conditions are presented in Table 2.

Table 1. siRNA sample sequences.

siRNA name	Length	Sequence
Inclisiran	21 S	mC*mU*mAmGmAmCfCmUfGmUdTmUmUmGmCmUmUmUmGmU/L96/
	23 AS	mA*fC*mAfAfAmGfCmAfAmAfAmCfAmGfGmUfCmUmAmG*mA*mA
Givosiran	21 S	mC*mA*mGmAmAmAfGmAfGmUfGmUfCmAmUmCmUmUmA/L96/
	23 AS	mU*mG*mGfUmCfUmUfUfCmUfCfAmCfAmGfAmGfUmAmGfA*fA*mU

Code	Description	Code	Description	Code	Description	Code	Description	Code	Description
*	Phosphorothioate bond	mA	2'-O-methyl A	mG	2'-O-methyl G	fA	2-fluoroadenosine	fG	2-fluoroguanadine
/L96/	C ₁₈ H ₁₃₉ N ₁₁ O ₃₁	mC	2'-O-methyl C	mU	2'-O-methyl U	fC	2-fluorocytidine	fU	2-fluorouridine

Table 2. LC and Q-TOF conditions.

	Target plus impurities		Sequence confirmation
	denatured	non-denatured	
Column	Agilent Altura Oligo HPH-C18, 2.1 × 50 mm, 2.7 μm		
Flow rate	0.4 mL/min		
Column temperature	65 °C	25 °C	65 °C
Injection volume	1 μL		
Mobile phase	A: 15 mM TEA and 50 mM HFIP in water B: Methanol		
Gradient	Time (min)	%B	Time (min) %B
	0	10	0 10
	7	40	1 10
	7.5	90	7 40
	8	90	7.5 90
Stop time	8 min		8 90
Post time	1.5 min		
UV detection	260 nm with 4 nm bandwidth		-
MS instrument	6545XT		
Drying gas temperature	350 °C	250/350 °C	350 °C
Drying gas flow	12 L/min		
Nebulizer	45 psi		
Sheath gas temperature	400 °C	250/400 °C	400 °C
Sheath gas flow	12 L/min		
Capillary voltage	3500 V		
Nozzle voltage	2000 V		
Fragmentor	175 V		
Skimmer	65 V		
Instrument mode	System tune		
Acquisition mode	MS		Targeted MS/MS
Mass range	m/z 400-3200		m/z 400-3200
Acquisition rate	1 spectra/s		4 spectra/s
MS/MS mass range	-		m/z 50-2000
MS/MS acquisition rate	-		2 spectra/s
Collision energy	-		12 V, 15 V, 18 V, 20 V
Targeted m/z	-		Calculated monoisotopic mass of the charged ion
Delta retention time	-		0.6 min
Isotope width	-		Medium (~4 m/z)

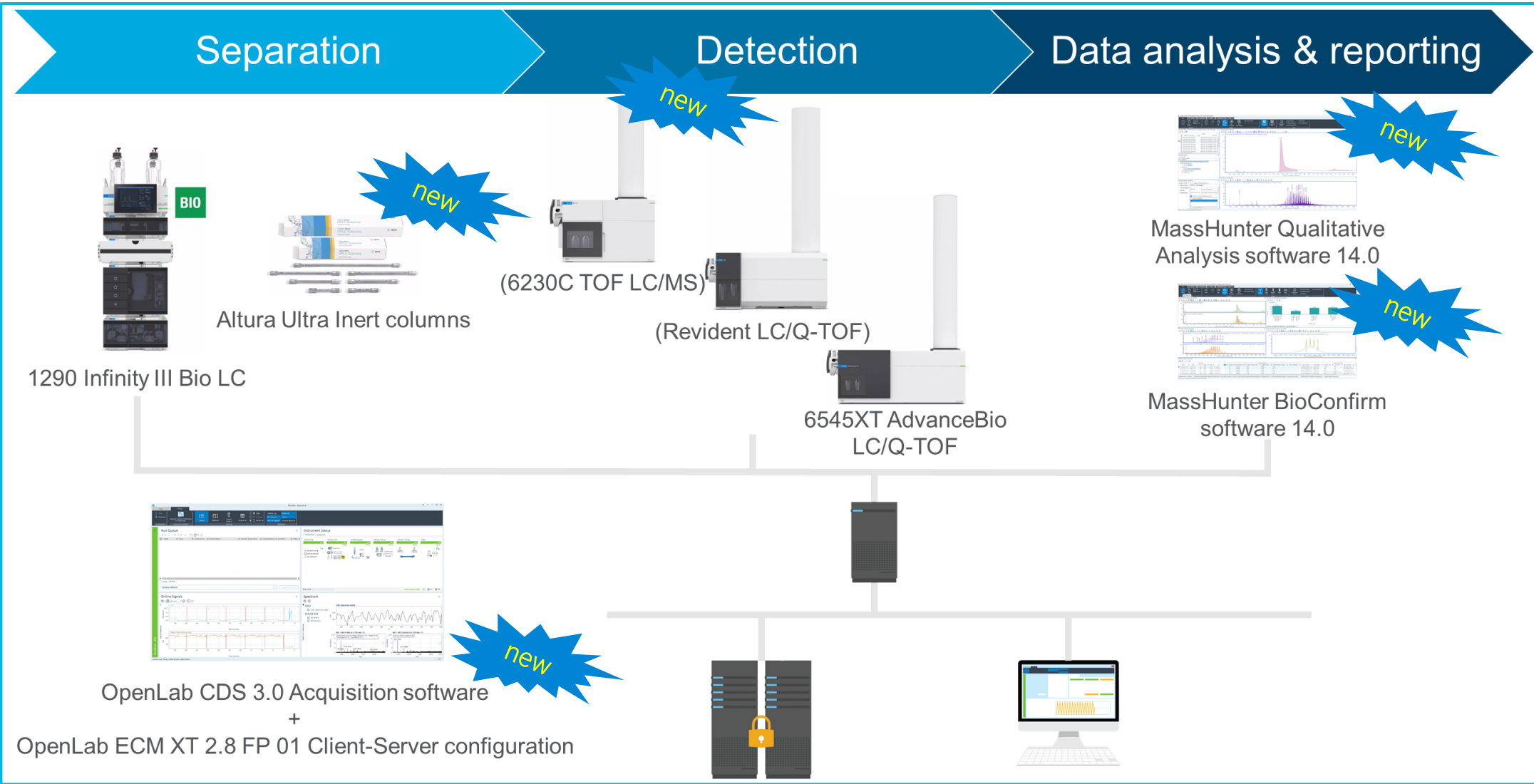


Figure 1. Solution overview.

Results and Discussion

Seamless sample processing: from data acquisition through data analysis to reporting

A combination of task actions, wait actions, and custom sample parameters enables the sample sequence to run without human intervention, potentially for several days (see Table 3). Unattended processing allows samples to be processed automatically immediately after data acquisition is complete, using predefined data analysis and reporting methods.

Table 3. A sequence for automated data acquisition, processing, reporting, and instrument control.

Action	Sample name	Inj/Vial	Acquisition method	Processing + Reporting method	Sample Custom Parameters	
					BioConfirm Sequence	BioConfirm Modification Profiles, and more...
Inject	Water	3	Oligo_TPI.amx	Oligo_TPI.bcpmx		
Inject	Inclisiran	3	Oligo_TPI.amx	Oligo_TPI.bcpmx	Inclisiran S, Inclisiran AS	
Inject	Inclisiran	5	Oligo_SC.amx	Oligo_SC.bcpmx	Inclisiran S	
Inject	Inclisiran	5	Oligo_SC.amx	Oligo_SC.bcpmx	Inclisiran AS	
Inject	Water	3	Oligo_TPI.amx	Oligo_TPI.bcpmx		
Inject	Givosiran	3	Oligo_TPI.amx	Oligo_TPI.bcpmx	Givosiran S, Givosiran AS	
Inject	Givosiran	5	Oligo_SC.amx	Oligo_SC.bcpmx	Givosiran S	
Inject	Givosiran	5	Oligo_SC.amx	Oligo_SC.bcpmx	Givosiran AS	
Inject	Water	3	Oligo_TPI.amx	Oligo_SC.bcpmx		
Task	Turn off the LC modules, switch MS to Eco Standby					
Wait	Wait until next morning, e.g. 8:00					
Task	Turn on all instruments, run mass calibration for MS					

Analysis of denatured and non-denatured siRNA

Both inclisiran and givosiran were analyzed in their denatured and non-denatured forms using ion-pairing HPLC coupled with DAD and Q-TOF detection. The results are shown in Figures 2 and 3.

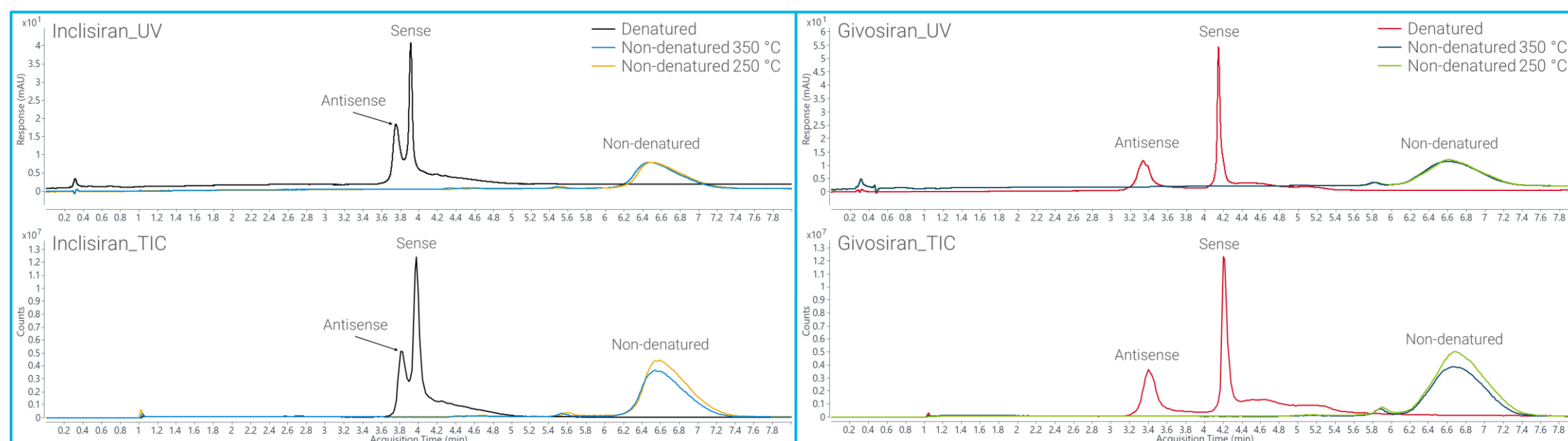


Figure 2. UV and total ion chromatograms of denatured and non-denatured inclisiran (left) and givosiran (right).

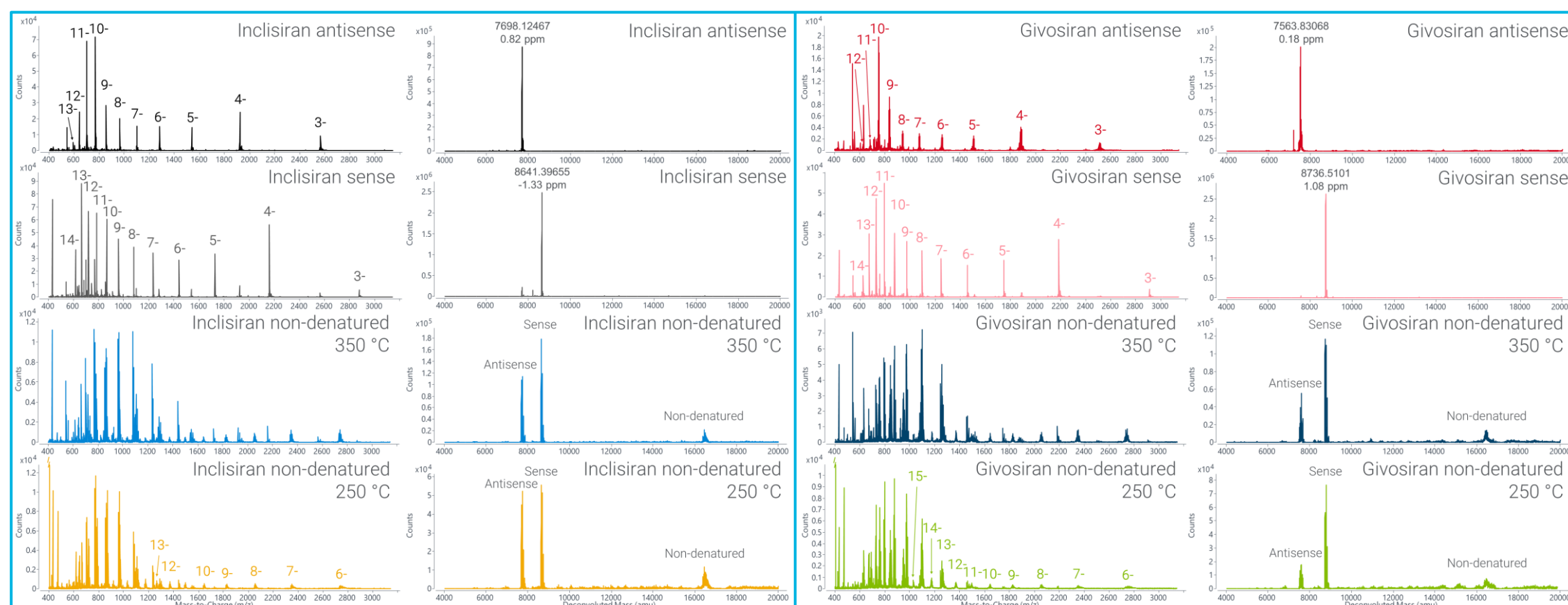


Figure 3. Raw and deconvoluted MS spectra of inclisiran (left) and givosiran (right).

Results and Discussion

Sequence confirmation of siRNA

As one example, inclisiran was analyzed using targeted MS/MS. 100% sequence coverage was achieved for both the antisense and sense strands of inclisiran, as shown in Figure 4.

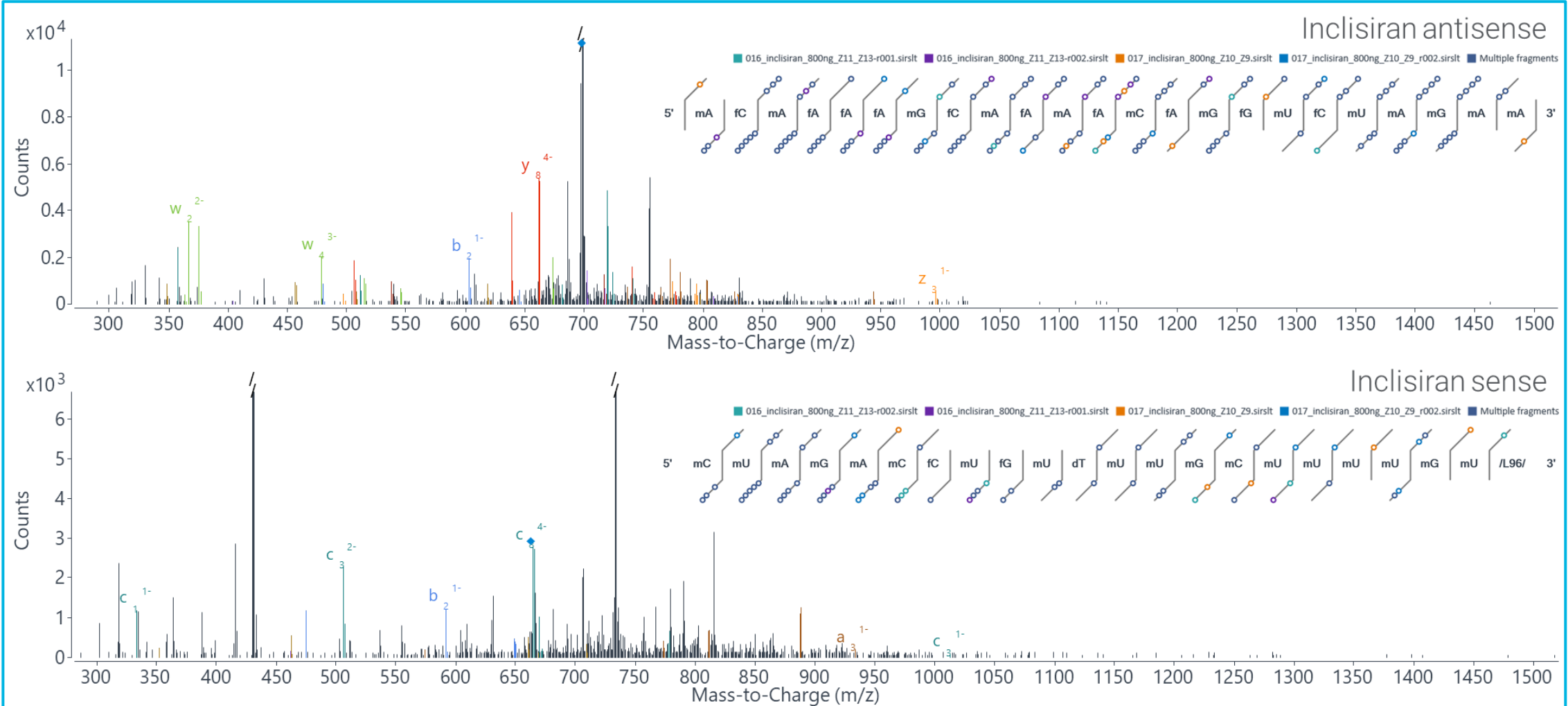


Figure 4. MS/MS analysis of inclisiran, measured by 6545XT. Sequence coverage was achieved by targeting the -10 and -11 charge states of the antisense strand, and the -9 and -13 charge states of the sense strand, in duplicate.

Reporting

Ready-to-use report templates are available for the analysis of oligonucleotides, including applications such as target and impurity analysis as well as sequence confirmation. A built-in report builder supports customized report design, enabling reports to be generated automatically and efficiently. An example of a sequence confirmation report for the antisense strand of inclisiran is shown in Figure 5 (right).

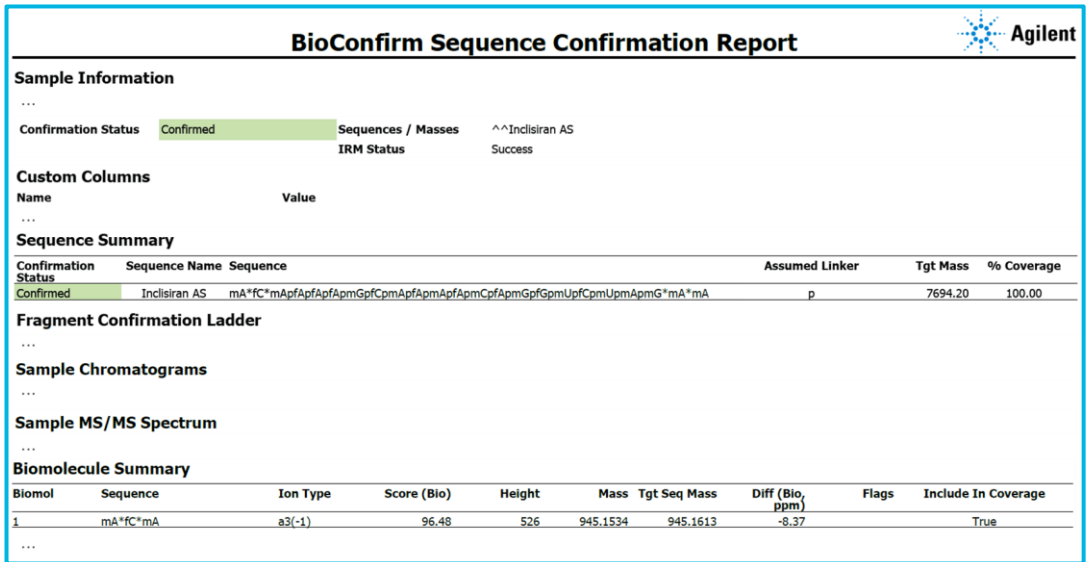


Figure 5. Report generated using the existing report templates in BioConfirm..

Conclusions

- An integrated software platform, together with new hardware and consumables, enables end-to-end data acquisition, analysis, and reporting for siRNA.
- Intelligent sample sequence management is supported through task actions, wait actions, sample custom parameters, and additional advanced features.

Table 4. Consumables for the solution.

Description	Part Number
Agilent Altura Oligo HPH-C18, 2.1 × 50 mm, 2.7 μm	227205-702
Agilent InfinityLab bio solvent bottle, PP, 1 L, with cap	9301-6028
Agilent InfinityLab bio solvent inlet filter, 10 μm, PEEK	5320-0070
Agilent InfinityLab Quick Connect LC fitting	5067-5965
Agilent InfinityLab Methanol for LC/MS	5191-5111
Agilent InfinityLab Water for LC/MS	5191-5121
Agilent Polypropylene Vials	5188-2788
Agilent Snap Top Caps	5182-3458

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

This information is subject to change without notice.

DE-014873