



## Sanger sequencing and fragment analysis

# SeqStudio series genetic analyzers – enabling simple, fast, flexible genomic analysis

### In this application guide, we demonstrate:

- Key features and connectivity of the Applied Biosystems SeqStudio™ and Applied Biosystems SeqStudio™ Flex Genetic Analyzers
- High data quality and comparability to other Applied Biosystems™ genetic analyzers
- A selection of the broad spectrum of genetic analysis applications that can be run on the SeqStudio family of instruments such as:
  - Cell line authentication
  - Microsatellite instability
  - dsDNA QC
  - Multiplex PCR analysis
  - Gene editing confirmation
  - SARS-CoV-2 sequencing
  - Minor allelic variant sequencing
  - Next-generation sequencing confirmation
  - Plasmid sequencing
  - Difficult template sequencing
  - Detection of KRAS mutations by Sanger sequencing
  - mRNA vaccine quality control using Sanger sequencing
  - Analyzing methylated DNA by Sanger sequencing
  - SNaPshot Multiplex System for SNP genotyping

### Introduction

The Applied Biosystems™ brand is a leader in capillary electrophoresis (CE) reagents, instrumentation, and software, and continues to provide innovative solutions that meet modern research needs. Capillary electrophoresis (CE) is a genomic analysis method that separates fluorescently labeled DNA fragments based on size by a single nucleotide, and is the foundation of Sanger sequencing and fragment analysis. Sanger sequencing is the gold-standard technology for accurate sequence determination with single-base resolution, ideal for targeted sequencing and confirmation analysis. Fragment analysis can provide sizing, relative quantitation, and genotyping information using fluorescently labeled DNA fragments produced by PCR using primers designed for a specific DNA target.

Applied Biosystems SeqStudio Family Genetic Analyzers are the trusted solutions in CE instrumentation. SeqStudio Family genetic analyzers are built upon reliability and trusted results. The SeqStudio Series Genetic Analyzers were purpose-built from the ground up to be easy to use, maintain, access and share data. Both the Applied Biosystems SeqStudio™ and Applied Biosystems SeqStudio™ Flex Genetic Analyzers are standalone instruments that offer an integrated touch screen computer, advanced connectivity features, fast run times and comprehensive primary and secondary analysis software. These flexible genetic analyzers are capable of performing both Sanger sequencing and fragment analysis.

The low-throughput, SeqStudio genetic analyzer is an easy-to-use system with the ability to run a standard 96-well plate. The SeqStudio Flex system, the latest addition to our CE portfolio, is our most advanced medium-throughput genetic analyzer with the ability to run 4-plates in either 96-well or 384-well varieties. The addition of features like continuous plate loading, voice command, streamlined one-button start up and remote start up and troubleshooting improve lab efficiency.

## Compare the instruments

### SeqStudio Genetic Analyzer

### SeqStudio Flex Series Genetic Analyzers



Number of capillaries	4	8 or 24
Capacity	96-well standard plate and standard 8-tube strips	4-plate capacity; 96-well plates, 384-well plates, 8-tube strips
Dimensions	49.5 x 64.8 x 44.2 cm	70 x 67.5 x 86.5 cm
Weight	53.6 kg	115 kg
Polymer type	POP-1, integrated into a simple click-in cartridge	POP-6, POP-7, POP-4
Continual plate loading	No	Yes
Sample reprioritization	No	Yes
Radio Frequency ID	Yes	Yes
Configuration	Integrated computer with touchscreen; optional desktop	Integrated computer with touchscreen; optional desktop
Integrated remote troubleshooting tools	No	Yes
Remote monitoring and data sharing	Yes	Yes
Voice command	No	Yes
Connectivity	USB, Ethernet ports, and Wi-Fi dongle	USB, Ethernet ports, and Wi-Fi dongle

\*Only the 24-capillary SeqStudio Flex instrument can use the 384-well plate

Customers looking for high-throughput capacity should consider the Applied Biosystems 3730 and 3730xL Genetic Analyzers. These systems are built for customers who require an economical, automated, and scalable instrument to process large numbers of samples with the option to process 96- or 384-well plates with plate stacking capabilities.

Learn more at [thermofisher.com/3730](https://thermofisher.com/3730)

## A connected experience

The integrated wireless connectivity make the SeqStudio series of genetic analyzers accessible via the onboard interface, a remote computer, or a mobile device app. (Figure 1).

Runs can be set up using either the onboard computer or by using PlateManager, the stand-alone software that operates within the cloud-based Connect platform or on a separate computer. By using web browser-based software, access to run setup, plate maps, run conditions, and analysis settings are all available from anywhere you have Internet access. Injection conditions, reinjections, and reordering of injections can all be modified during the run, maximizing the ability to collect quality data from each plate. After data collection, the web browser-based suite of applications makes analysis simple and accessible from any location, ideal for remote collaboration.

In this applications guide, we demonstrate how the SeqStudio family Genetic Analyzers integrate seamlessly into some of the most commonly used Sanger sequencing and fragment analysis applications. The ease of use and ease of data analysis provided give investigators flexibility and make either system an ideal addition to your genetic analysis toolbox.

Learn more at [thermofisher.com/connect](https://thermofisher.com/connect)



**Figure 1. SeqStudio Family Genetic Analyzers integrate seamlessly into Connect with cloud-enabled systems.** By registering your instrument and logging into your cloud account, remote features are accessible, including instrument monitoring, data analysis using cloud apps, and data sharing with your colleagues.

# Results

## Fragment analysis applications

Fragment analysis is a highly flexible method that involves the separation of different-sized and differentially labeled DNA fragments by CE. One of the most common methods of generating fragments for analysis is polymerase chain reaction (PCR); because of the flexibility afforded with the choice of PCR primers, a specifically sized fragment corresponding to a PCR target sequence is straightforward to generate. Along with the ability to label fragments with up to four different fluorophores, researchers have a large degree of flexibility in experimental design. Below we will review examples of how the SeqStudio and the SeqStudio Flex have been used to support various fragment analysis applications.

### Cell line authentication

Disease research relies on the analysis of samples obtained and manipulated *ex vivo*. These include human cell lines, induced pluripotent stem cells (iPSCs), and chimeric antigen receptor (CAR) T cells. Contamination and human error can occur and impact research results or put lives in danger.

For example, a study from 2019 found that that 32,755 articles reported results obtained with misidentified cells— these were in turn found to be cited in an estimated half a million other publications [2].

Finally, as research progresses into *ex vivo* cell therapies— cells that are removed from a donor, manipulated in a lab, and returned to a host—it is critical to confirm that donor cells have the expected genotype and match the intended recipient. Therefore, it is crucial to know the provenance of human cells and confirm that they have the desired identity.

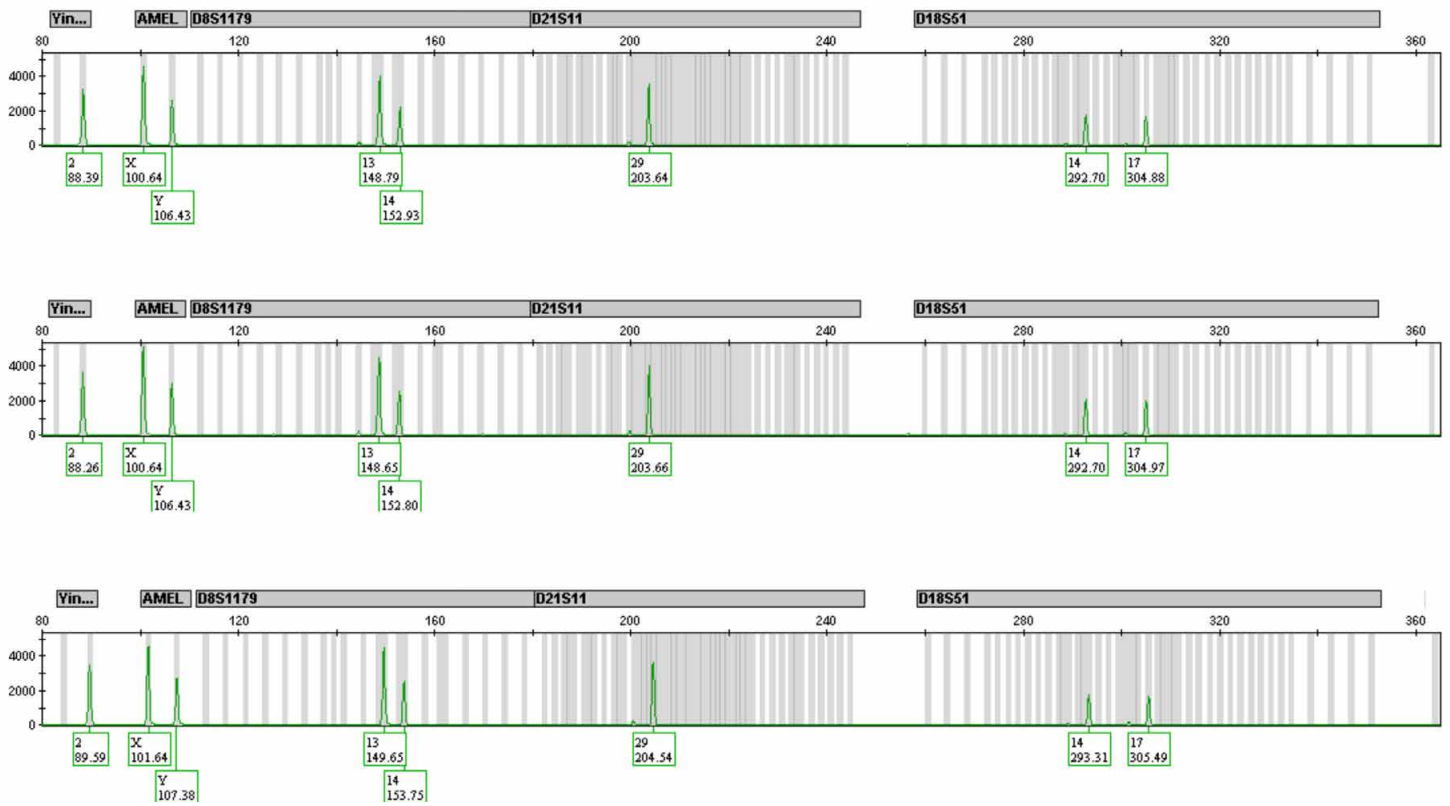
An array of Applied Biosystems™ solutions for identifying human samples are based on fragment analysis of short tandem repeats (STRs). STRs are microsatellite sequences that are highly variable and provide a unique molecular fingerprint for a human sample. The Applied Biosystems™ GlobalFiler™ cell line authentication (CLA) kit generates a molecular fingerprint for 24 different STR loci, while the Applied Biosystems™ Identifiler™ Plus CLA kit analyzes 16 STR loci. Both of these easy-to-use kits are optimized for use on the Applied Biosystems family of CE-based genetic analyzers. For more information, see reference 3.

To illustrate the performance of the SeqStudio Flex instrument relative to other Applied Biosystems genetic analyzers, DNA from six commonly used human cell lines (A549, Jurkat, U2OS, HEK293, M4A4GFP, and HeLa) was purified using Invitrogen™ RecoverAll™ kits. For STR analysis, 1 ng of genomic DNA was used with either the GlobalFiler CLA or Identifiler CLA kit according to the supplied protocols. Each sample was run in triplicate. Samples were run on the SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers. Allelic calls were made using the Applied Biosystems™ GeneMapper™ software. The authenticity of the cell lines was confirmed by checking the results against databases for STRs of common cell lines [4].

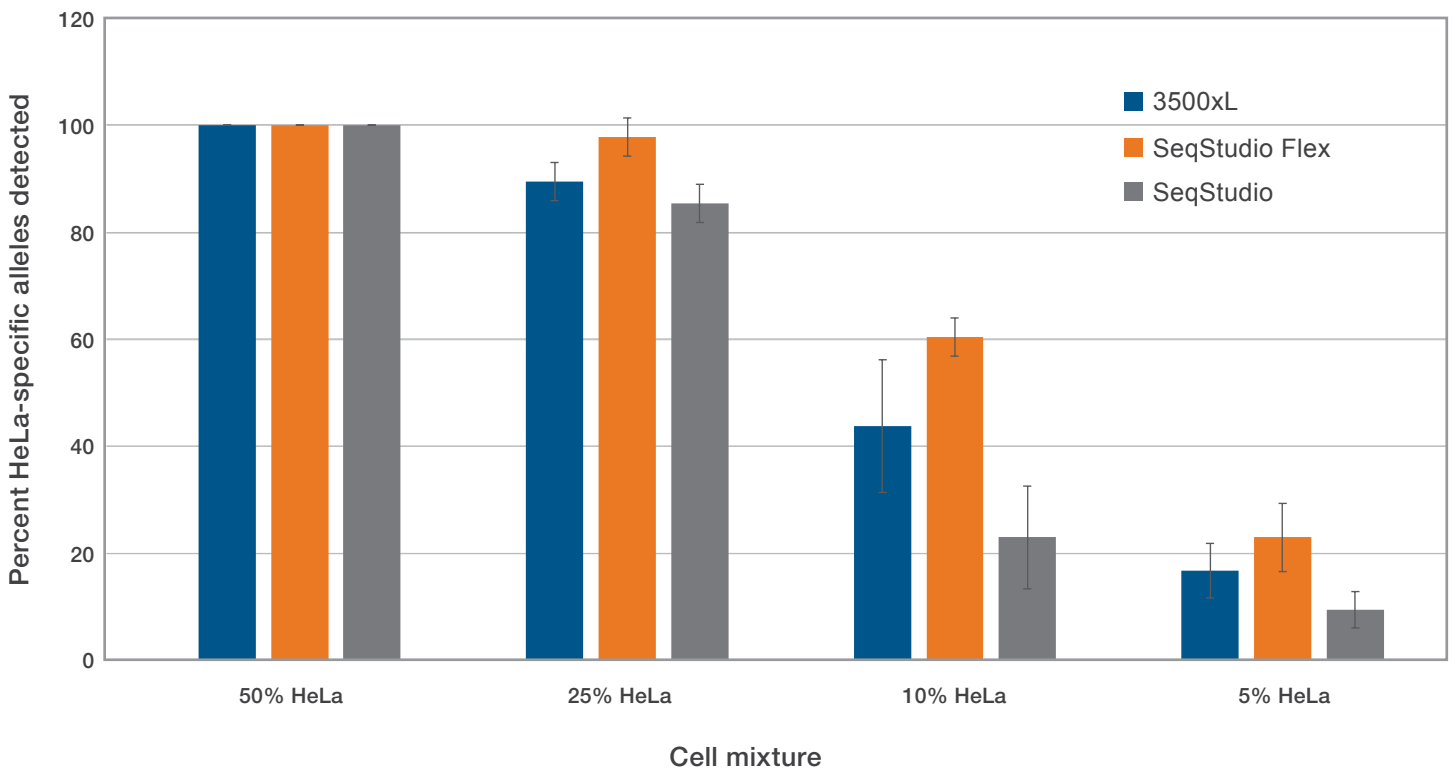
The STR peak profiles at all loci for all five cell lines were comparable across all instruments (Figure 2A). Using the GlobalFiler CLA kit, we found a 100% match to the expected STR allelic profile for all cells using the SeqStudio Flex and 3500xL genetic analyzers, and a 98.7% match with the SeqStudio genetic analyzer (Tables 1 and 2). Similarly, using the Identifiler CLA kit, we obtained a 100% match with the SeqStudio Flex system, a 99.6% match with the 3500xL system, and a 98.8% match with the SeqStudio system.

To examine the ability to detect cell line contamination, we mixed M4A4GFP and HeLa gDNA at 50%, 25%, 10%, and 5% HeLa DNA. The Identifiler CLA kit was used to analyze 1 ng of each gDNA mixture. HeLa cells have 16 alleles that are not found in the M4A4GFP profile. We found that even when the contaminating HeLa cells make up 5% of a gDNA sample, the genetic analyzers can detect between 9% and 23% of the unique HeLa alleles (Figure 2B). Together, these results demonstrate that the SeqStudio Flex instrument generates cell line authentication results of similar or better quality compared to the rest of the genetic analyzer family.

Read more about this application at [thermofisher.com/cla](https://thermofisher.com/cla)



**Figure 2A. CLA peak morphologies on three genetic analyzer platforms.** The SeqStudio Flex (top), 3500xL (middle), and SeqStudio (bottom) genetic analyzers produce similar peak morphologies and allele calls using the GlobalFiler CLA kit. A single channel is shown; similar results were seen in all five colored channels and with additional samples.



**Figure 2B. Detecting contaminating alleles at low concentrations.** Detecting contaminating alleles at low concentrations. Decreasing amounts of HeLa gDNA were added to M4A4GFP gDNA. The total amount of DNA was kept to 1 ng. The mixture was analyzed in triplicate using the Identifiler Plus CLA kit. The number of HeLa-specific alleles detected in each titration versus number expected was determined. Note that HeLa-specific alleles were detected in the 5% sample on all three instruments.

**Table 1. Allelic profiles were collected on SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers.** The example shown is from human A549 cells analyzed using the GlobalFiler CLA kit.

Locus	Channel	SeqStudio Flex		3500xL		SeqStudio	
AMEL	G	X	Y	X	Y	X	Y
CSF1PO	B	10	12	10	12	10	12
D10S1248	P	13	16	13	16	13	16
D12S391	P	18	18	18		18	
D13S317	R	11	11	11		11	
D16S539	B	11	12	11	12	11	12
D18S51	G	14	14	14		14	
D198433	Y	13	13	13		13	
D1S1656	P	17	18.3	17	18.3	17	18.3
D21S11	G	29	29	29		29	
D22S1045	R	15	15	15		15	
D2S1338	P	24	24	21		21	
D25441	Y	10	13	10	13	10	13
D3S1358	B	16	16	16		16	
D5S818	R	11	11	11		11	
D7S820	R	8	11	8	11	8	11
D8S1179	G	13	14	13	14	13	14
DYS391	G	10	10	10		10	
FGA	Y	23	23	23		23	
SE33	R	19	25.2	19	25.2	19	25.2
TH01	Y	8	8	8		8	
TPOX	B	8	11	8	11	8	11
VWA	B	14	14	14		14	
Yindel	G	2	2	2		2	

**Table 2. The total number of alleles that were correctly called using the GlobalFiler CLA kit was totaled across three replicates for six cell lines.** Each cell line has a slightly different combination of homozygous and heterozygous allele numbers, so the total number of alleles will vary between cell lines.

	SeqStudio Flex	3500xL	SeqStudio
A549	102/102	102/102	102/102
Jurkat	153/153	153/153	153/153
U2OS	105/105	105/105	105/105
HEK293	111/111	111/111	104/111*
M4A4GEP	96/96	96/96	96/96
HeLa	117/117	117/117	115/117*

\* Technical issues during one of three injections resulted in some anomalous calls in these samples.

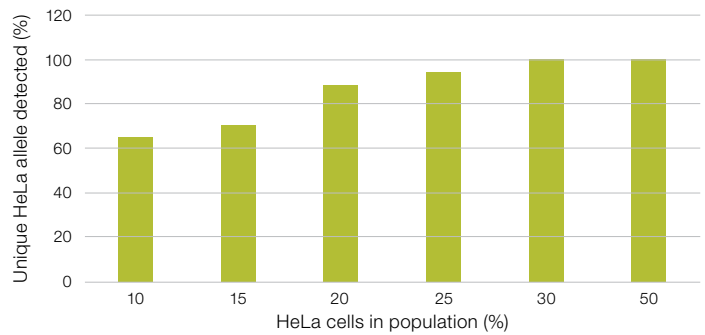
## Human cell line authentication

The study of development of human diseases relies heavily on the analysis of dissociated human cell lines grown in culture. However, an increasingly acknowledged problem is that cells grown in vitro can be misidentified or contaminated with other unrelated cell lines. The identity of cell lines can be verified by analysis of a highly specific genetic “fingerprint” of highly variable short tandem repeats (STRs). The SeqStudio platform integrates well with the Thermo Fisher Scientific cell line authentication solution.

The Applied Biosystems™ CLA Identifier™ Plus and CLA Identifier™ Direct kits can be used on purified and crude DNA preparations, respectively, for analyzing 16 highly variable human STR loci commonly used for verifying cell line authenticity. The Applied Biosystems™ GeneMapper™ Software and the cloud-based Microsatellite Analysis (MSA) software, used for analyzing alleles identified by CLA Identifier kits, are compatible with data produced by the SeqStudio instrument, and the results can be used to query ATCC or other STR databases to verify authenticity [16].

To demonstrate the utility of the SeqStudio instrument in a cell line authentication workflow, allelic information on STRs was obtained from five different, commonly used human cell lines. The identity of the cell lines was confirmed even with as little as 300 pg of gDNA. To show the ability to detect contaminating cells, a population of M4A4GFP cells was spiked with varying amounts of HeLa cells and analyzed using the CLA Identifier Direct kit. HeLa cell-specific alleles could be detected even if only 10% of the population had HeLa cells (Figure 3). Therefore, when coupled with the CLA Identifier kits, the SeqStudio instrument can be a central component for a cell line authentication solution.

Read more about this application at [thermofisher.com/cla](https://thermofisher.com/cla)



**Figure 3. Analysis of cell line contamination on the SeqStudio instrument.** HeLa cells and M4A4GFP cell suspensions were diluted to  $5 \times 10^5$  cells/mL, mixed in the indicated proportions, and spotted onto NUCLEIC-CARD™ Sample Collection Device. Contaminating HeLa cells can be detected with high confidence on the SeqStudio instrument if they make up approximately 20% of a population; however, some alleles unique to HeLa can be detected if they make up as little as 10% of a population.

## Microsatellite instability

Many types of cancer display deficiencies in DNA mismatch repair (MMR), producing an overall higher mutation rate across the genome [5]. A higher mutation rate often means a higher rate of neoantigen production, providing opportunities for immune therapy research [6].

There have been at least 11 different loci implicated in MMR [7]. Looking for an inactivating event in the sequence of all these loci can be complicated, time consuming, and expensive. Examining the outcome of perturbations in the MMR pathway by microsatellite instability (MSI) analysis provides a practical alternative. MSI analysis can be difficult on NGS systems due to the highly repetitive nature of microsatellite sequences. Thus, fragment analysis of microsatellite loci length using CE is a widely used method for detecting MSI.

The Applied Biosystems™ TrueMark™ MSI analysis system analyzes a panel of 13 microsatellite loci, including the Bethesda set of markers [8], as a measure of microsatellite instability. TrueMark MSI assays also provide STR information at two highly variable loci, allowing for sample identity confirmation. To ease the analysis of MSI raw data, we developed Applied Biosystems™ TrueMark™ MSI Analysis Software to simplify the calls at each locus (Figure 3A). Furthermore, we incorporated analysis algorithms that do not require side-by-side analysis of normal, non-tumor tissue, reducing the number of samples needed and the expense of an MSI analysis. For more information, see reference 9.

To illustrate the performance of the SeqStudio Flex instrument relative to other genetic analyzers, nine tumor/normal adjacent pairs and one tumor-only FFPE sample in blocks were obtained. Sections (10 µm) were cut from the FFPE blocks and gDNA was extracted from the sections, and 1–2 ng of genomic DNA was used in the TrueMark assay according to the published protocol. Fragments were analyzed on the SeqStudio Flex and 3500xL genetic analyzers, and results were interpreted using the TrueMark MSI Analysis Software (Figure 4A).

To compare the performance between genetic analyzers, we tallied the number of loci that could be called either stable, unstable, or no call (Figure 4B). The total number of loci that fell into each class was highly concordant. One sample (S07-001886-A5) had very little gDNA recovered, so using 2 ng of the recommended input amount was not possible. Nevertheless, the two instruments were able to make similar calls for the loci that were detectable.

Samples from the same FFPE blocks were used for standard immunostaining for MMR proteins, and the overall MSI determinations using results from the SeqStudio Flex and 3500xL systems were completely concordant with the immunostaining results. These results demonstrate that the SeqStudio Flex system generates MSI analysis data that are equivalent to those obtained with the 3500xL system.

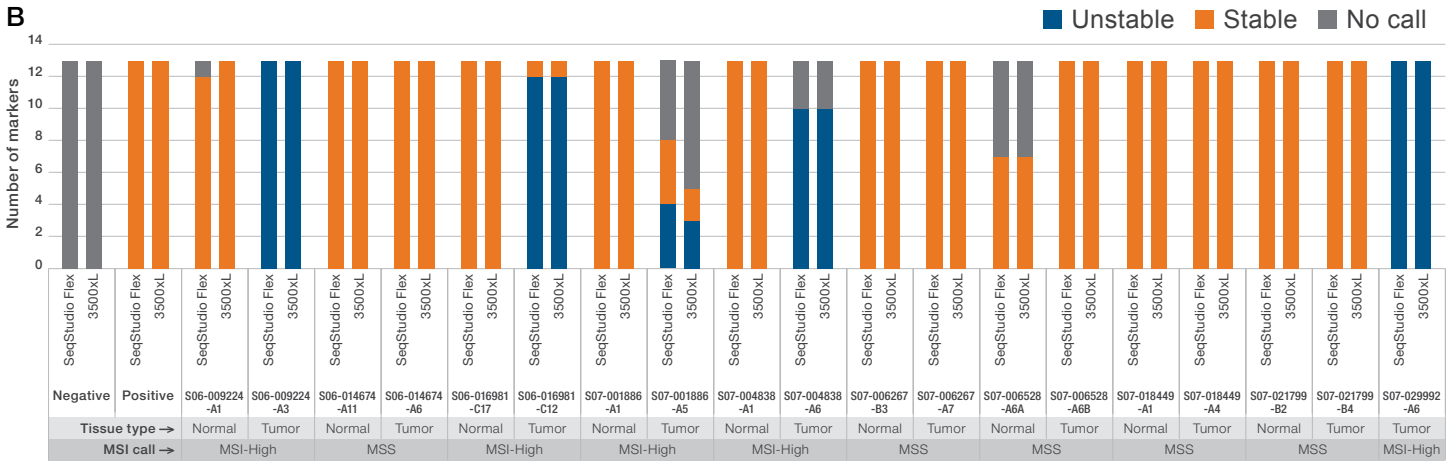
**Read more about this application at [thermofisher.com/msi](https://thermofisher.com/msi)**



A



B



**Figure 4. The SeqStudio Flex and 3500xL instruments produced similar data in MSI analysis.** (A) The TrueMark MSI assay analyzes 13 microsatellite loci for instability, including the widely used Bethesda standards. Loci that are determined to be unstable can be autocalled; the software will then use the totality of the calls to make an overall call for the sample. The assay includes two highly variable short tandem repeat (STR) sequences (TH01 and PentaD) that can be used to confirm sample identity. The proprietary algorithms used by the software do not require side-by-side analysis of normal, non-tumor tissue in order to make stable/unstable calls. (B) Nine tumor/normal adjacent pairs and one tumor-only sample were analyzed using the TrueMark MSI assay. The number of loci called by the software was highly similar using data from both instruments. Sample S07-001886-A5 had suboptimal gDNA recovered; not all of the loci amplified equivalently and thus produced slightly different results on the two instruments.

## dsDNA QC

Many protocols in genomic analyses start from a pool of DNA. Frequently the size and amount of DNA in that pool are confirmed prior to initiating further studies. For example, extracting DNA from FFPE-preserved tissue can result in fragmented DNA that may be too small to provide useful information. Thus, it is beneficial to know the overall size distribution before moving experiments forward.

In next-generation sequencing (NGS) experiments, libraries of pooled fragments are often generated before loading onto the sequence-reading chip. Knowing the size and abundance of DNA fragments in that library is important for predicting the success of the NGS result.

Finally, it is beneficial to confirm whether a specific amplicon was generated from a PCR reaction. Agarose gels are often useful for this, but while they are inexpensive, they are not extremely precise or high in throughput.

The size and relative abundance of double-stranded DNA fragments can be analyzed on Applied Biosystems genetic analyzers. For this application, Invitrogen™ TOTO™-1 DNA-intercalating fluorescent dye is incubated with a sample. The sample is subsequently run using Applied Biosystems™ POP-7™ polymer with the capillary heater turned off. This helps maintain the DNA in a mostly double-stranded, nondenatured state. By comparing the resulting migration of a test fragment with the migration of known dsDNA standards (e.g., phage lambda DNA digested with HindIII, phage phiX174 DNA digested with BsuRI), the size of the test fragment(s) can be determined.

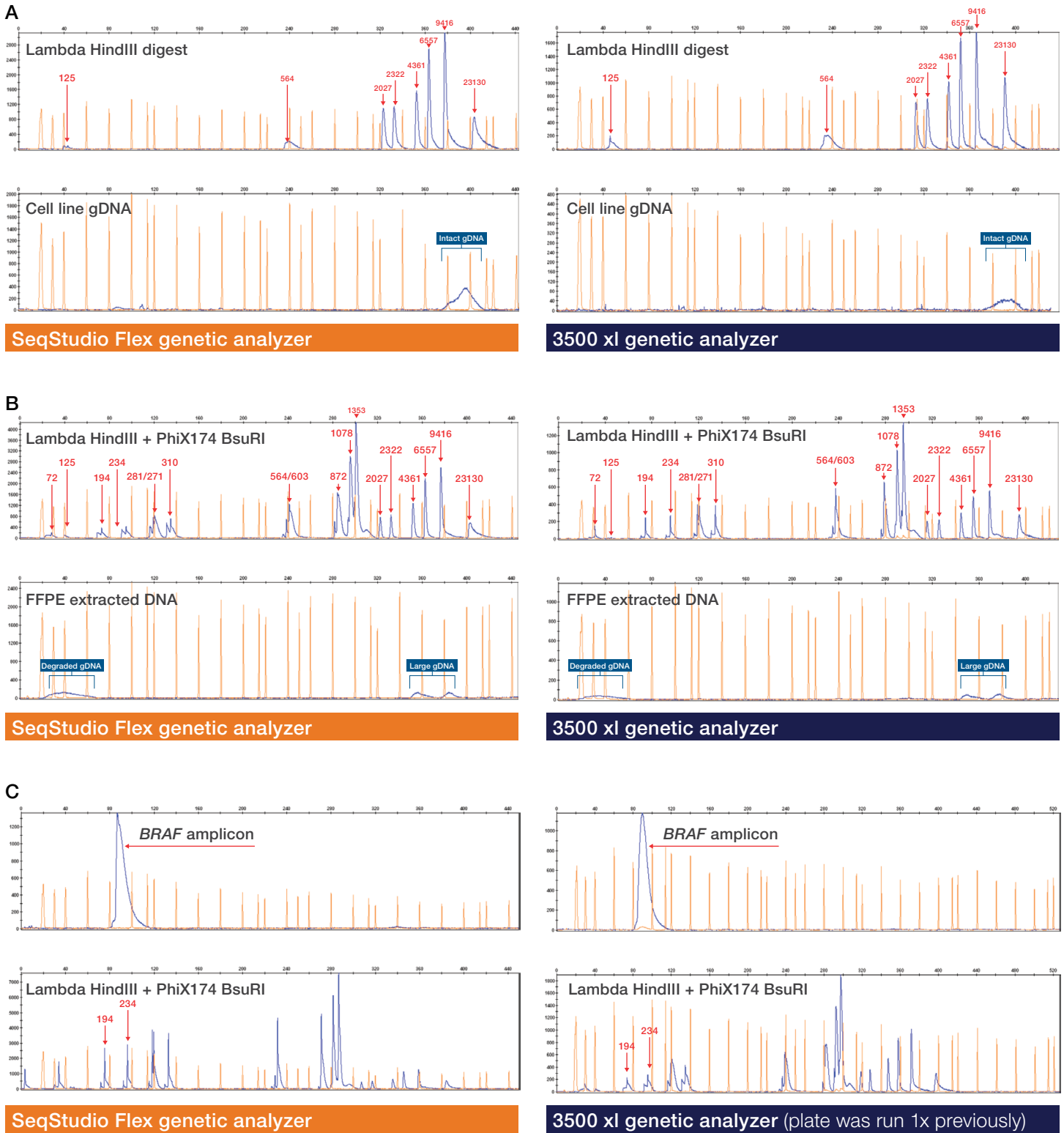
To assess the performance of the SeqStudio Flex system relative to the 3500xL system for analyzing dsDNA, we obtained samples of extracted genomic DNA from cell lines, extracted genomic DNA from FFPE slides, a BRAF amplicon for cycle sequencing, or lambda HindIII and phiX174 BsuRI digested DNA. dsDNA samples were mixed with a 50 nM solution of TOTO-1 dye and single-stranded (ss) Applied Biosystems™ LIZ™ size standards in water. Samples were incubated at 37°C for 10 minutes and separated on the genetic analyzer calibrated with the E5 dye set and capillary heating off. Fragments were analyzed using GeneMapper 6.1 software.

When lambda HindIII standards are analyzed using this method, the well-known pattern of eight fragments, ranging from 23,130 bp to 125 bp, can be seen (Figure 5A). In the same injection but in a different capillary, we analyzed 50 ng of genomic DNA purified from a cell line. Comparing the migration of the gDNA to the lambda standard shows that the majority of the intact DNA ranges between approximately 9 kb and 20 kb.

Similarly, we ran a mixture of lambda HindIII and phiX174 BsuRI standards in another capillary. In addition to the eight lambda fragments, 10 more fragments that range from 1,353 to 72 bp can be seen (Figure 5B). DNA extracted from an FFPE slide runs from about less than 70 bp to about 200 bp, plus some large DNA fragments in the 4 kb–10 kb range.

Finally, we checked the quality of a BRAF amplicon before cycle sequencing (Figure 5C). According to the lambda and phiX174 standards, the amplicon migrates in the range of around 200–230 bp. The expected size of this amplicon is 208 bp, matching well with the observed size. Note that in each example, the profiles obtained on the SeqStudio Flex and 3500xL analyzers were very similar, demonstrating that the SeqStudio Flex and 3500xL systems generate equivalent dsDNA fragment data.

This method provides a path to fast and simple QC of dsDNA samples. There are several enhancements that would increase the information gained from these runs. For example, the ss LIZ internal standards can be recalibrated to the lambda/phiX marker set. Once they have been redefined, the LIZ standards can act as an internal size standard when run in the same capillary as the unknown. Thus, it may be possible to quantify the amount in a peak, provided a standard curve is run at the same time. Finally, we have found that this application also works on the 4-capillary SeqStudio instrument (data not shown).



**Figure 5. Analysis of double-stranded DNA by CE.** (A) Different DNA preparations were electrophoresed through capillaries containing POP-7 polymer at reduced temperature. Single-stranded LIZ-labeled fragments were included in the capillary. Lambda HindIII digest fragments and genomic DNA from A549 cells were analyzed in different capillaries. Note the typical pattern of lambda fragments (top panels). These can be used to estimate the size of the gDNA population of about 9–23 kb (bottom panels). Equivalent performance was seen on both instruments. (B) A mixture of lambda HindIII and phiX174 BsuRI digests and DNA extracted from an FFPE tumor sample were analyzed in different capillaries. Note the typical pattern of phage fragments (top panels); these can be used to estimate the size of the fragmented DNA (75–190 bp) and the large genomic DNA (bottom panels). Equivalent performance was seen on both instruments. (C) An amplicon containing the BRAF V600E region was analyzed before cycle sequencing, along with a mixture of lambda HindIII and phiX174 BsuRI digests in different capillaries. Note the typical pattern of phage fragments (top panels); these can be used to estimate the size of the BRAF amplicon (194–230 bp). This is very similar to the expected size of 208 bp. Equivalent results were seen on both instruments.

## Multiplex PCR analysis

Amplification of DNA sequences by PCR remains a workhorse of almost all aspects of molecular biology and genomic research. As the depth of knowledge and scientific questions become more complex, there is an increasing need to analyze many targets in a single sample. However, singleplex PCR can be tedious, involving setting up individual reactions for each of the desired targets.

Methods that can analyze multiple PCR amplicons from a single sample would streamline workflows and conserve precious samples. One method to address this is fragment analysis by CE, using multiple fluorescent dyes for multiplex capabilities. Here, a very large number of targets can be analyzed from a single sample, since it can separate unique and specific amplicons based on both size and fluorescent dye in a single capillary. In fact, the CLA and MSI analysis kits described above are examples of multiplex PCR analysis of fragments.

We demonstrated how fragment analysis by multiplex PCR works using a custom panel for detecting respiratory viral pathogens. We designed a set of PCR primers that could detect 12 different respiratory RNA viruses. The panel was designed such that each organism generated a different sized amplicon. Oligos contained 6-FAM™ labeling at the 5' end, and the set of primers was pooled so the panel could be mixed with sample in a single step. The resulting amplicons could be separated and examined using any of the CE genetic analyzers.

To illustrate the performance of the SeqStudio Flex instrument compared to the other genetic analyzers, we obtained Invitrogen™ GeneArt™ synthetic DNA targets or purified genomic RNA targets (ATCC) for common respiratory viruses. These targets, either single or mixed, were amplified using Applied Biosystems™ TaqMan® Multiplex Master Mix and the custom panel of pooled primers described above. Following PCR, amplicons were mixed with Applied Biosystems™ Hi-Di™ formamide and LIZ standards and separated on SeqStudio Flex and 3500xL genetic analyzers. The resulting data were analyzed using GeneMapper v6.1 software.

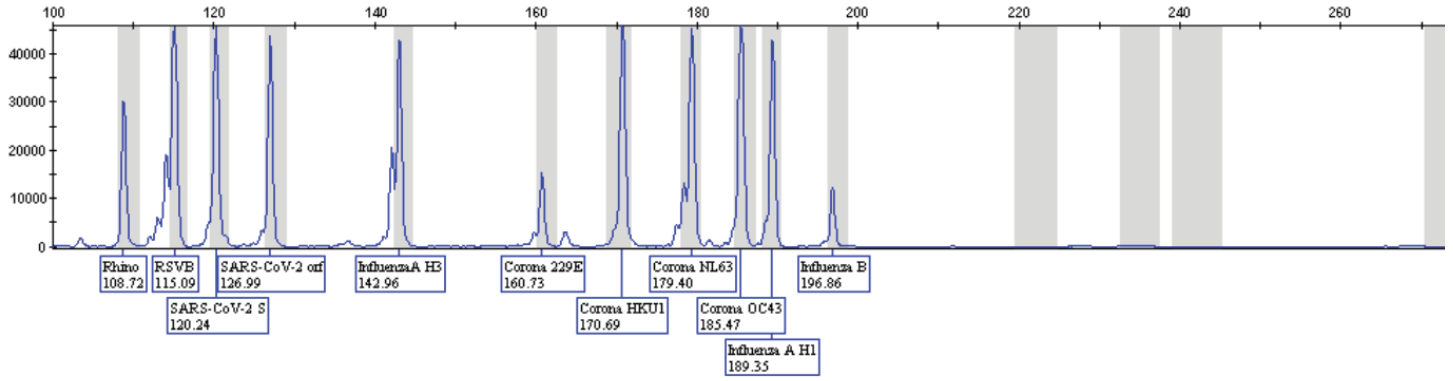
The functionality of the primers was confirmed by testing each primer pair individually against its cognate target (data not shown), and against the mixture of all DNA targets with the pooled set of primers (Figure 6A). Each of the primer pairs was able to recognize its target if present, in either the singleplex or multiplex assay. We confirmed that the approach could work with RNA genomes by first synthesizing cDNA from either single viral genomes or a mixture of all viral genomes, then PCR amplifying with either individual primers or the pooled primer set. Each of the organisms was detected if the appropriate target was in the sample (Figure 6B).

Finally, we compared the results obtained with the SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers. Each of the instruments produced comparable results (Figure 6C). These results therefore confirm that the SeqStudio Flex system generates multiplex PCR data that is equivalent to that obtained with the other members of our CE family of instruments.

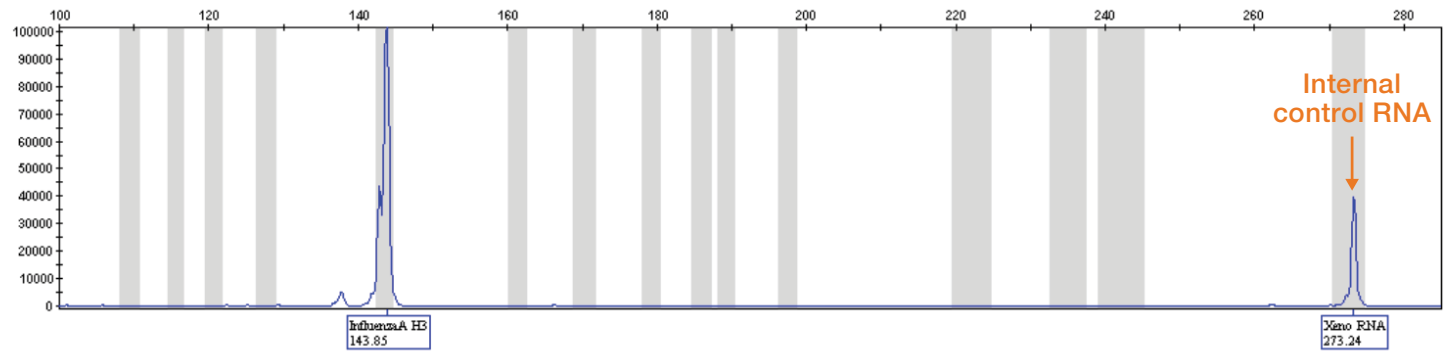
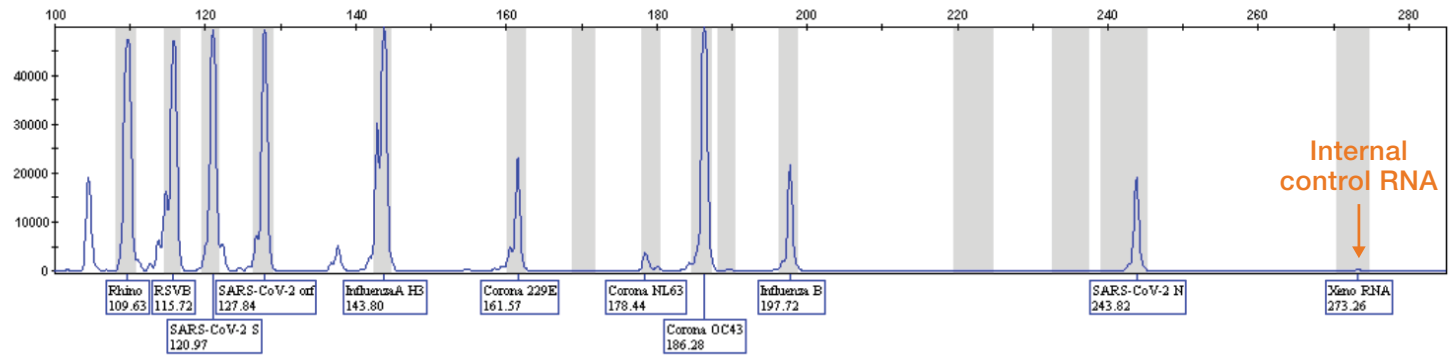
Multiplex PCR coupled with fragment analysis presents a tremendous opportunity to perform complex analyses with minimal effort. Designs of a panel are limited only by the ability to design primer sets against the targets of interest. It can therefore be adapted and utilized in nearly any scenario where detection of many target sequences in a single sample is desired.

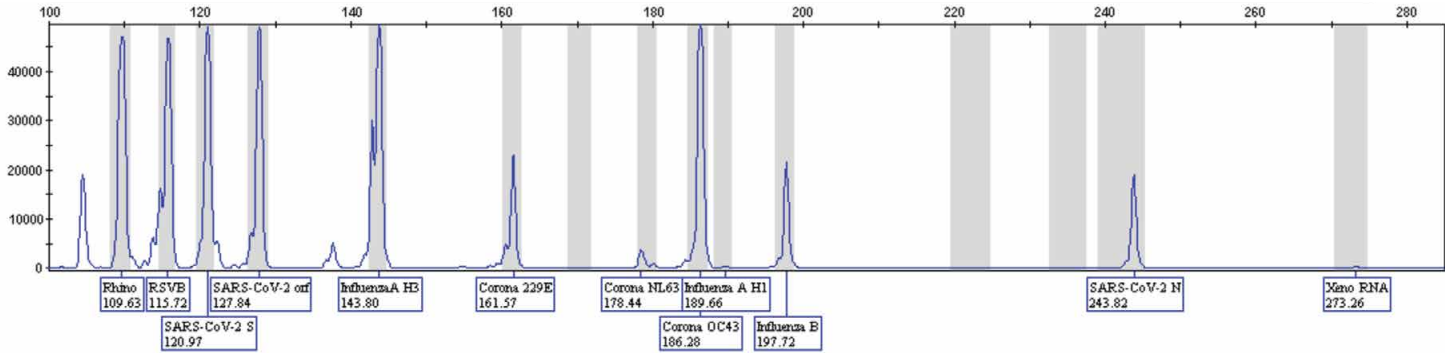
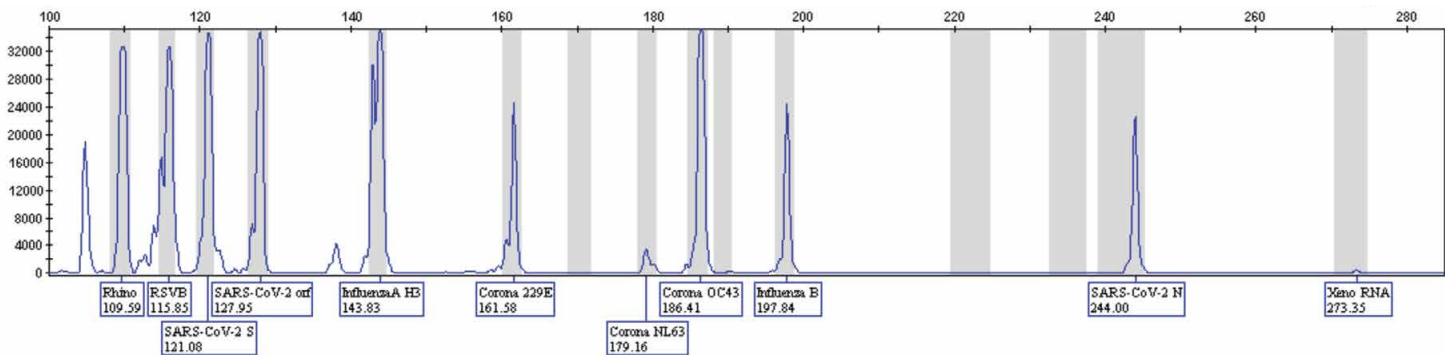
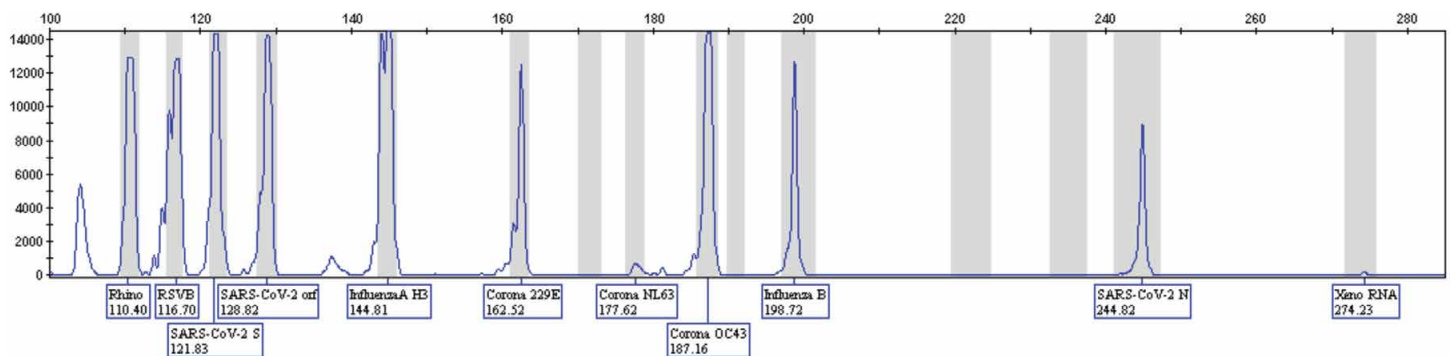
**Read more about this application at**  
[thermofisher.com/ce-clinical](https://www.thermofisher.com/ce-clinical)

A



B



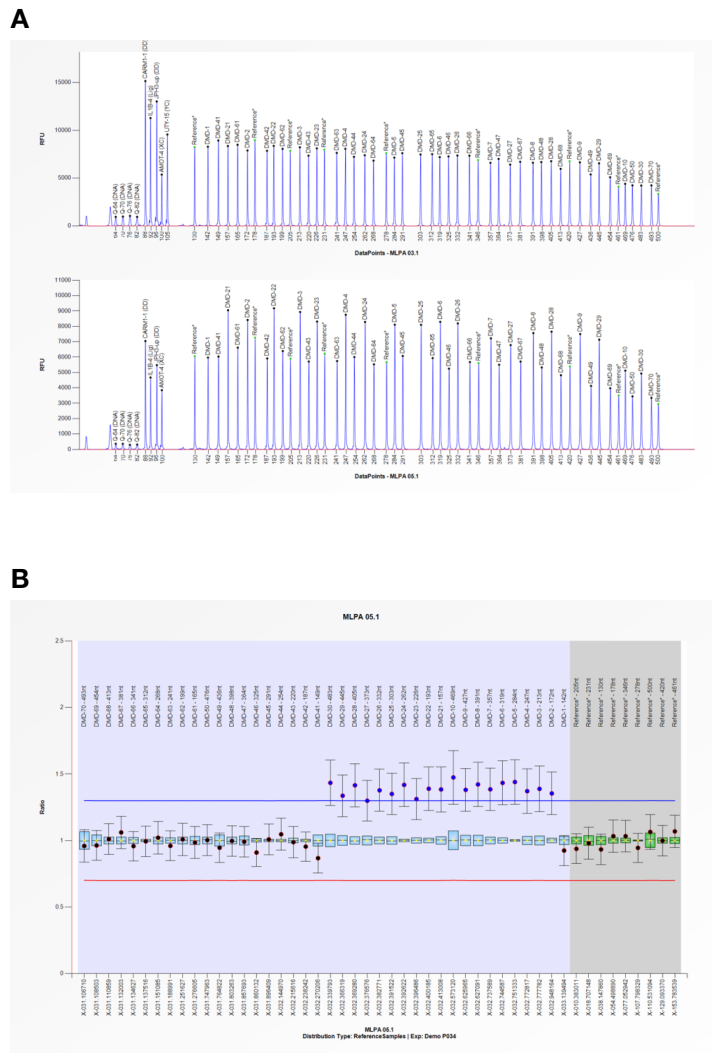
**C****SeqStudio Flex genetic analyzer****3500xL genetic analyzer****SeqStudio genetic analyzer**

**Figure 6. Analysis of multiple targets by multiplex PCR and fragment analysis. (A)** DNA targets of the RNA viruses shown above were amplified in a single reaction using a panel of 6-FAM-labeled PCR primers. Each primer pair will produce a virus-specific amplicon size that can be separated and defined by fragment analysis. The grey bins were defined to detect the targets of this particular panel. **(B)** Purified RNA genomes of the viruses shown above were either pooled (top) or left single (bottom). cDNA synthesis followed by amplification using the pooled primers confirmed that multiple targets or a single target can be picked up in a single reaction. As an internal control for cDNA synthesis and amplification, an exogenous RNA (Applied Biosystems™ Xeno™ RNA) was spiked into the RNA pool and the primer pool contained primers for the exogenous RNA. **(C)** Equivalent ability to detect the multiplex amplicons was seen on all instruments; an example from the RNA genomic pool is shown here.

## Multiplex ligation-dependent probe amplification (MLPA)

One widely used method for studying inherited human diseases arising from variations in copy number of a locus is multiplex ligation-dependent probe amplification [18]. This method, developed and commercialized by MRC Holland, can analyze up to 50 multiplexed pairs of adjacently located probes hybridizing to the loci of interest. The high dynamic range, sizing precision, and peak-height fidelity necessary for analyzing MLPA probe amplicons make the SeqStudio system an ideal platform for performing MLPA analyses. Results obtained on the SeqStudio instrument are compatible with MRC Holland's Coffalyzer.Net software for analyzing MLPA data.

MLPA on the SeqStudio instrument was used to analyze a DNA sample from a probe that is known to carry a duplication of exons 2–30 in the Duchenne muscular dystrophy (DMD) gene and a normal sample using the P034 DMD assay set from MRC Holland. The peak heights and relative sizes of these samples can readily be translated into an accurate detection of the region containing the duplication (Figure 7). Similar results were obtained using probes for large and small deletions. Therefore, the SeqStudio instrument can be an integrated tool for MLPA investigations of regions containing CNVs.



**Figure 7. Analysis of CNV on the SeqStudio instrument using MLPA.** (A) Fragment profile of a sample analyzed for CNVs at the DMD gene. Note that peak heights of some of the fragments are larger on the bottom sample versus the top, indicating an overrepresentation of DNA in that region. (B) The corresponding ratio chart, aligning the probes on the DMD locus, and clearly displaying increased ratios for DMD probes in exons 2–30. Probes that cross the blue threshold are indicative of a gain in copy number from 2 to 3.

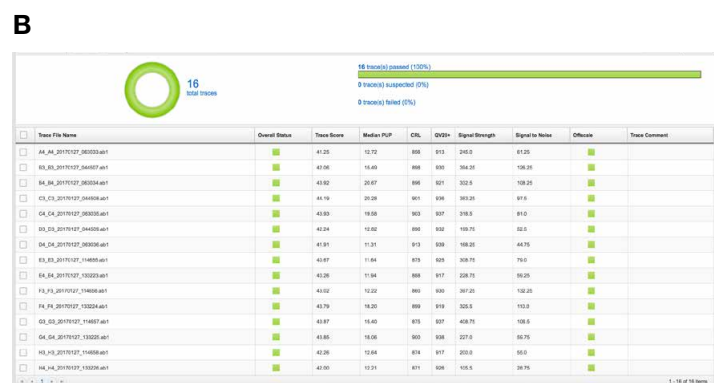
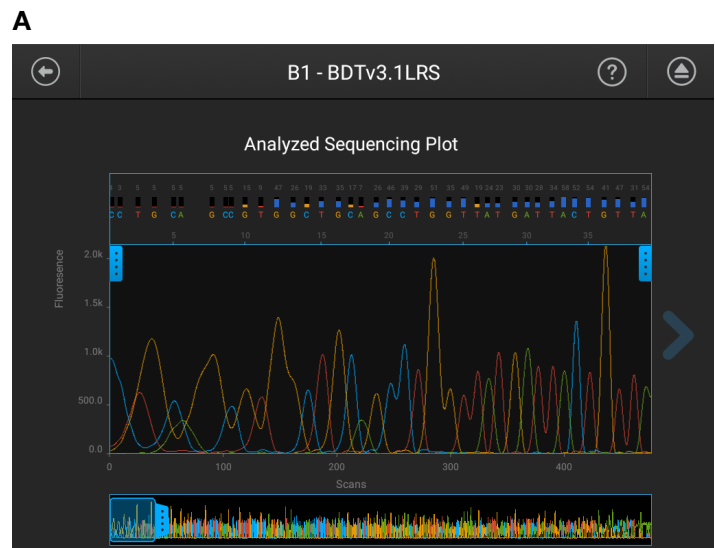
## Sequencing applications

Sanger sequencing is the trusted standard for obtaining DNA sequence information. It powered the Human Genome Project and investigators continue to rely on this method to generate highly accurate, targeted sequencing results. Applied Biosystems™ products support fast and straightforward Sanger sequencing workflows that enable a high degree of accuracy, long-read capabilities, and simple data analysis. An entire workflow can be completed in less than one workday, from sample to answer, providing the flexibility to support a diverse range of applications in many research areas. We describe below several popular applications that commonly use Sanger sequencing methods enabled by the SeqStudio and the SeqStudio Flex.

### Plasmid sequencing

One of the most common applications of Sanger sequencing is the analysis of inserts subcloned into plasmids. Applied Biosystems™ BigDye™ chemistries are widely used for Sanger sequencing and an integral part of plasmid sequencing workflows. Several of the features on the SeqStudio platform offer benefits to researchers performing basic plasmid sequencing methods. The instrument is preloaded with sequencing modules optimized for short (<300 bp), medium (500 bp), and long (>600 bp) read lengths, and can also be customized on the instrument to meet specific needs. The swappable cartridges can be associated with individual projects and users. The cloud-based Sanger sequencing Quality Check application provides an intuitive set of tools to analyze sequencing traces. Finally, the cloud connectivity for remote monitoring, accessing, and sharing sequencing information can help collaborators rapidly analyze the same data sets.

The performance of the SeqStudio instrument for plasmid sequencing was determined by sequencing the pGEM7zf+ plasmid with M13 primers and Applied Biosystems™ BigDye™ Terminator v3.1 chemistry. Results were obtained by analyzing the sequencing traces using the Sanger Quality Check module on Connect (Figure 8). In the example shown, the same plasmid was sequenced in 16 wells and analyzed on the SeqStudio Genetic Analyzer in four different injections. Note that the trace score, peak under peak (PUP) values, contiguous read length (CRL), and QV20+ (length with quality values >20) are similar for each sample. Similar results were obtained in traces on the other strand, and in other experiments by using Applied Biosystems™ BigDye™ Terminator v1.1 chemistry. These data demonstrate that the SeqStudio platform can generate plasmid sequencing results of very high quality.



**Figure 8. Analysis of sequencing quality using the Sanger sequencing Quality Check Connect app.** (A) Once a run is completed, the SeqStudio instrument displays the resulting sequence file as well as the quality scores for each base. (B) Sixteen separate pGEM7zf+ sequencing reactions were run on the SeqStudio instrument, and the .ab1 files were uploaded to the cloud and analyzed. Note that the sequencing metrics were very similar in the sixteen different reactions. CRL = contiguous read length, QV20+ = number of nucleotides with a quality value >20.



## Plasmid sequencing (long read length)

Modern genetic analyses often require extensive manipulation of DNA sequences. Pieces of DNA are cut, amplified, joined, propagated, and purified in tasks ranging from subcloning inserts into plasmids, through generating libraries, to constructing viral vectors for gene therapy.

Because many of these steps involve DNA polymerases, misincorporation of bases and genetic drift can lead to undesired mutations in the target sequences. Confirmation of manipulated DNA sequence is therefore necessary to help ensure the intended sequences in the final product are correct.

Sequence confirmation is also advisable for any genetic engineering experiment, such as AAV viral vector construction or gene synthesis experiments. The ability to have long Sanger sequencing reads simplifies the chemistry and analysis workflow needs.

Applied Biosystems™ BigDye™ Terminator v1.1 and v3.1 cycle sequencing chemistries are widely trusted for Sanger sequencing by CE. After cycle sequencing, various options exist for clean-up before electrophoresis, including Centri-Sep™ purification columns and plates and Applied Biosystems™ ExoSAP-IT™ enzyme mix. Finally, the genetic analyzers have running modules that have been optimized for long reads.

To illustrate the performance of the SeqStudio Flex system against our other genetic analyzers, we performed a standard sequencing experiment. Plasmid pGEM™-3Zf(+) was sequenced using both v1.1 and v3.1 BigDye Terminator chemistries in forward and reverse directions. Sequencing reactions for each direction were performed with 8 replicate reactions in each direction. Reactions were purified using Centri-Sep plates, dried, and resuspended in Hi-Di formamide.

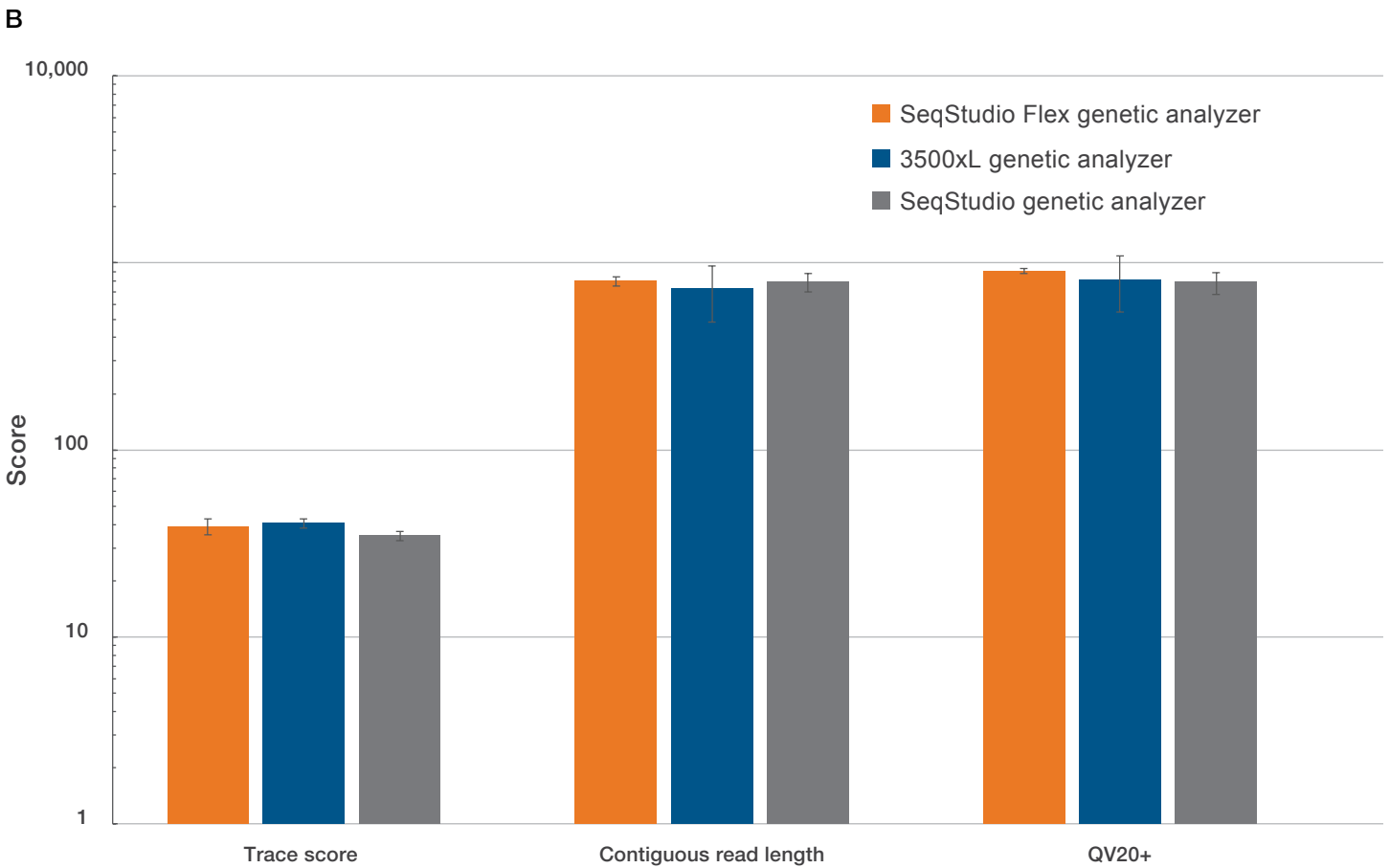
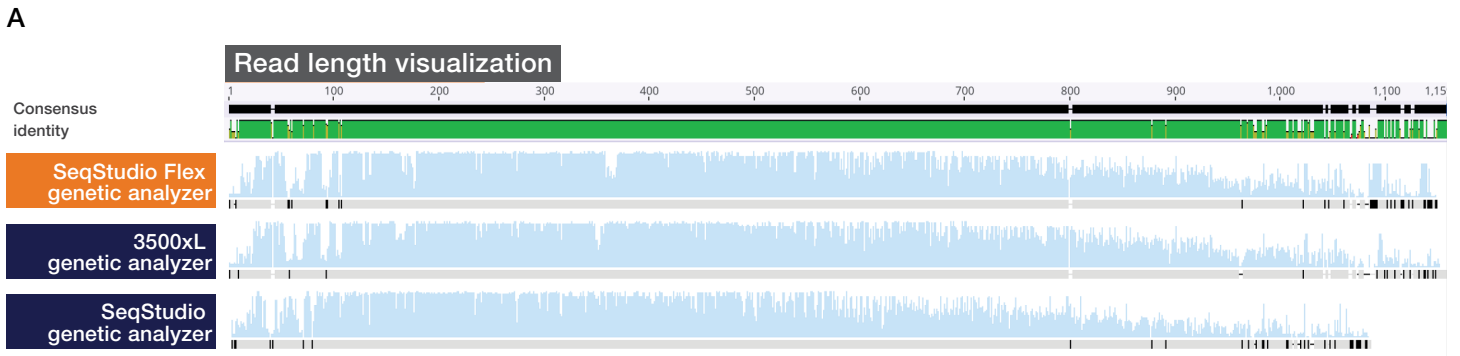
The resulting reactions were run on the SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers. Sequencing traces were analyzed using Applied Biosystems™ Sequence Scanner and Biomatters Geneious™ software. Using BigDye Terminator v1.1 chemistry, we were able to obtain over 1,000 bp of high-quality sequence on the SeqStudio Flex and 3500xL systems, and slightly less on the SeqStudio system (Figure 9A).

There are several quality control metrics that are generated by the Sequence Scanner Software. One measurement is the basecalling quality value (QV); it represents the shape and signal-to-noise ratio of a peak in the chromatogram. High-quality peaks typically have a QV greater than 20. The QV provides the most objective metric to evaluate the confidence of the base call and is used to calculate the trace score.

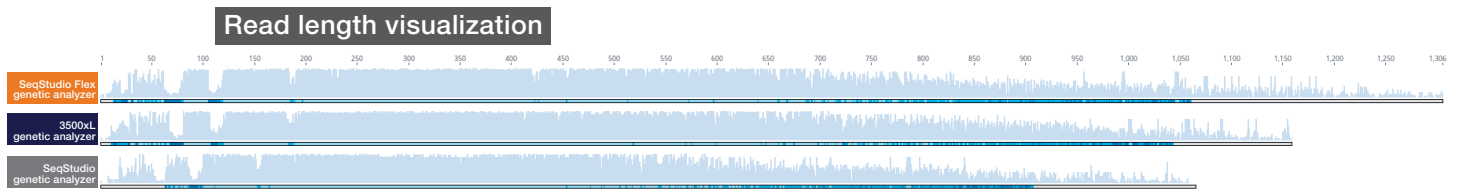
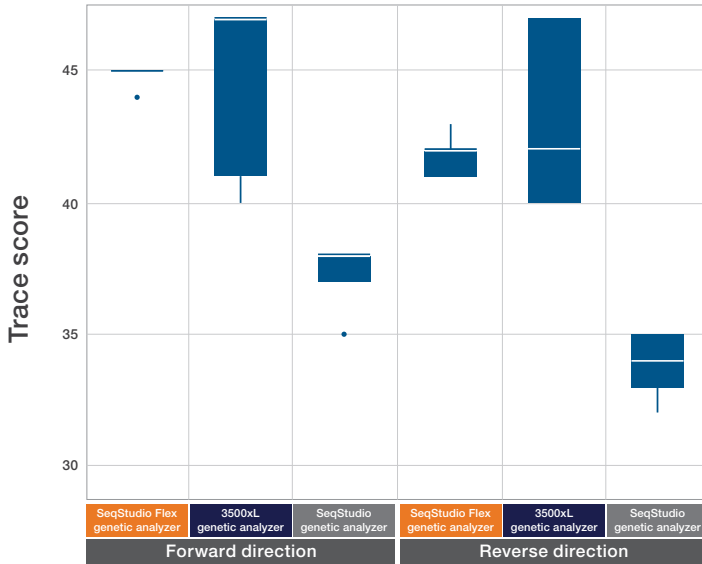
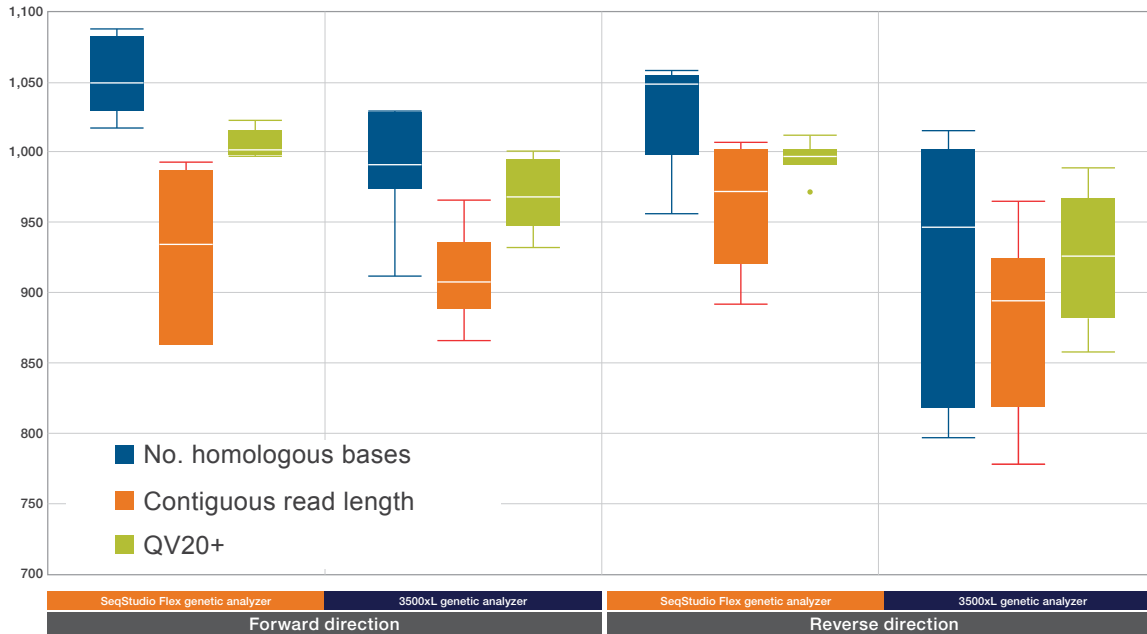
The trace score is the average of basecalling QVs for bases in the legible range of the chromatogram. The contiguous read length (CRL) is the longest read generated without poor-quality breaks in the chromatogram. The QV20+ value is the number of bases that had a quality value greater than 20 (high-quality reads) in the entire chromatogram. We used these metrics to compare the performance (Figure 9B).

The values for each of these QC metrics was nearly identical across the three instruments. Similarly, we sequenced the pGEM-3Zf(+) cloning vector using BigDye Terminator v3.1 chemistry, and obtained very similar read lengths on the SeqStudio Flex and 3500xL systems (Figure 10A). Furthermore, the trace score (Figure 10B), CRL, and QV20+ (Figure 10C) values were very similar.

Finally, we aligned the total read, which may include some regions that were homologous but not contiguous, to the pGEM-3Zf(+) reference sequence. We found an excellent degree of homology for the entire long read obtained on all instruments (Figure 10C). These results confirm that the SeqStudio Flex system generates high-quality, long reads that are equivalent to the quality obtained with our other CE instruments.



**Figure 9. Analysis of plasmid sequencing results of long reads using BigDye Terminator v1.1 chemistry. (A)** The read lengths are depicted in the Manhattan plot; the alignment with the consensus pGEM-3Zf(+) sequence is shown on the green bar on the top. The light blue bars indicate the quality of the read at that base; taller bars are better-quality reads. The SeqStudio Flex and 3500xL genetic analyzers produced almost identical results and were able to read more than 1,000 bp; the SeqStudio instrument had a slightly lower read length (~1,000 bp). **(B)** Quality metrics for reads were obtained on all three instruments. The sequence quality metrics produced on all three were nearly identical.

**A****B****C**

**Figure 10. Analysis of plasmid sequencing results of long reads using BigDye Terminator v3.1 chemistry. (A)** The read lengths are depicted in the Manhattan plot. The light blue bars indicate the quality of the read at each base; taller bars are better-quality reads. The SeqStudio Flex and 3500xL genetic analyzers produced almost identical results and were able to read more than 1,100 bp; the SeqStudio instrument had a slightly lower read length (~950 bp). Trace score **(B)**, contiguous read length, and QV20+ values **(C)** were obtained on all three instruments. In addition, the number of homologous bases in the reads was also obtained. The SeqStudio Flex and 3500xL systems gave similar results, and the SeqStudio system had slightly lower values (data not shown).

## SARS-CoV-2 sequencing (medium read-length)

Some investigations require focused sequence information for a specific region of a genome. While genome-wide discovery tools such as NGS provide vast amounts of data over very large regions, such complex approaches are not always needed in focused-based research. Sanger sequencing provides a simple, inexpensive, and easy-to-interpret solution when focusing on a region.

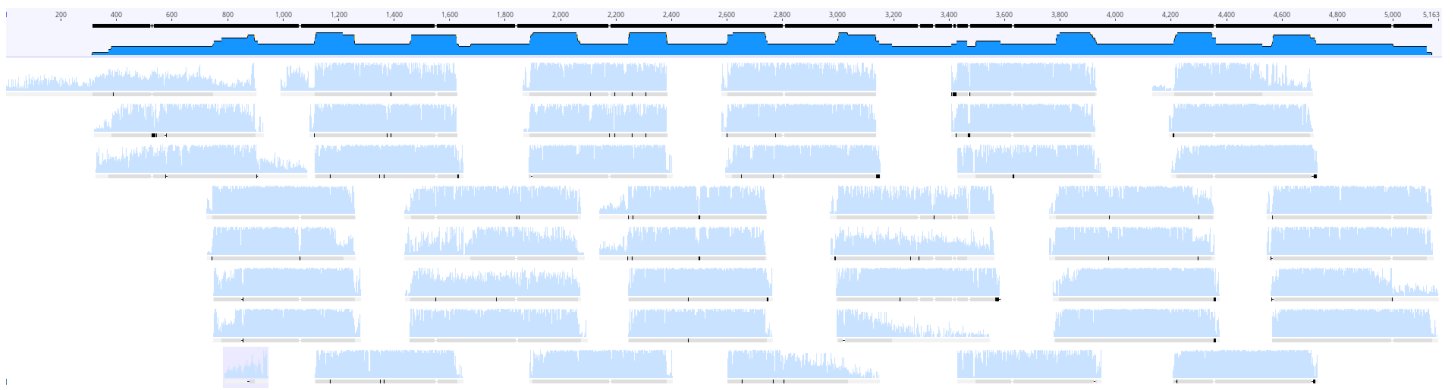
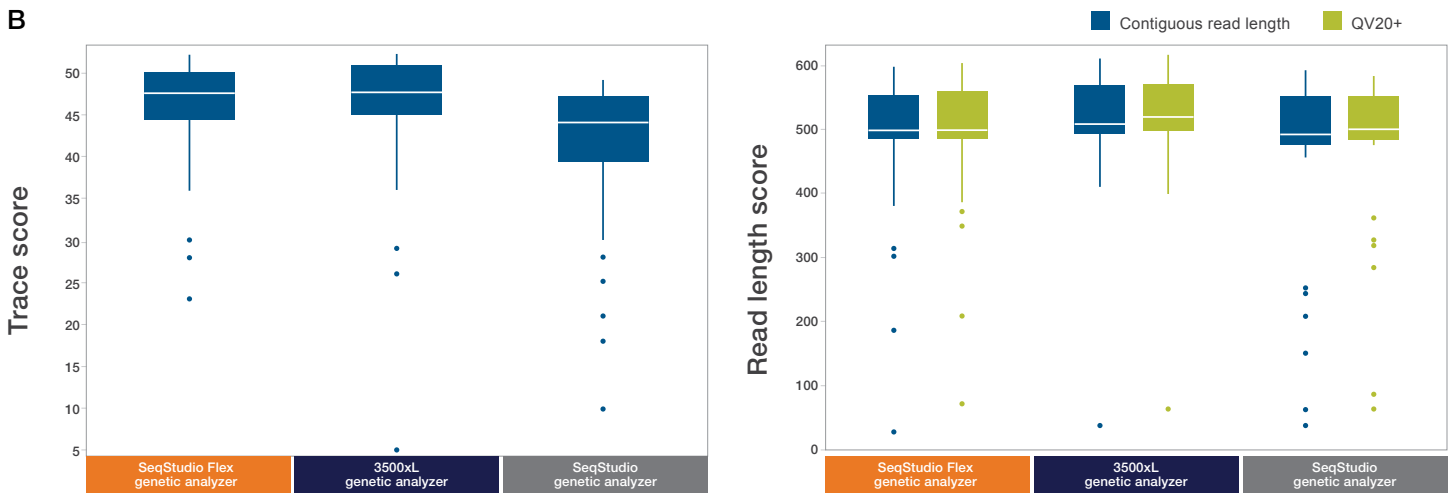
The Applied Biosystems™ BigDye™ Terminator v1.1 and v3.1 chemistries are also widely used for medium read-length sequencing. Another option, Applied Biosystems™ BigDye™ Direct chemistry, makes use of universal M13 primers to further simplify the cycle sequencing workflow.

When coupled with clean-up using the Applied Biosystems™ BigDye XTerminator™ Purification Kit, the entire cycle sequencing workflow can be performed in a single well of a 96-well plate. The genetic analyzer family has sequencing modules that have been optimized for BigDye Direct and BigDye Terminator sequencing reactions.

To compare targeted sequencing results, we sequenced the region of the SARS-CoV-2 genome that encompasses the Spike gene and adjacent sequences. Overlapping M13-tailed sequencing primers (12 pairs) for the entirety of the SARS-CoV-2 Spike gene have been previously described [10]. One thousand copies of SARS-CoV-2 genomic RNA were reverse-transcribed using Invitrogen™ SuperScript™ IV VIL0™ Master Mix. Amplicons (500–600 bp) were generated using 1 µL of cDNA and amplicon pairs, cycle sequenced using the BigDye Direct kit, and purified using the BigDye XTerminator kit. The resulting reactions were run on SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers. Chromatograms were analyzed using Sequence Scanner and Geneious software.

The Spike gene region of SARS-CoV-2 is covered by the 12 overlapping amplicons spanning over 5 kb. High-quality Sanger sequencing results were obtained from nearly all 12 amplicons in both directions (Figure 11A). These results could be assembled into a single contiguous read (contig) that was confirmed to be the SARS-CoV-2 Spike gene by BLAST alignment. The read quality metrics were nearly identical on all three genetic analyzers (Figure 11B). Consistent basecalling quality, as demonstrated by the data obtained on the three instruments, facilitates using Sanger sequencing to assemble the sequence of focused regions using overlapping mid-length amplicons, generating contigs that cover a locus of interest.

**Read more about this application at**  
[thermofisher.com/cecoronavirus](https://thermofisher.com/cecoronavirus)

**A****B**

**Figure 11. Sequencing the SARS-CoV-2 Spike gene by overlapping amplicons.** (A) Twelve amplicons were used to sequence SARS-CoV-2 RNA in duplicate. Reads produced by the amplicons were mapped to each other to recapitulate the Spike gene sequence. The read quality in each amplicon is indicated by the height of the light blue bars. Data were collected on the SeqStudio Flex system; similar alignments were performed with data from the 3500xL and SeqStudio systems, with highly similar results. Note that some of the amplicons had some quality drop-off; this was also observed with those runs on the other instruments. (B) Trace score (left) and contiguous read length and QV20+ values (right) were compared on all three instruments. Sequencing quality results were almost identical.

## Minor allelic variant sequencing and NGS confirmation

Discovery-based genomic research, such as next-generation sequencing (NGS), often uncovers novel or unexpected variants or other sequence anomalies. Investigators look for ways to verify these new discoveries using orthogonal methods. Sanger sequencing is the method of choice for confirming NGS results studies because of its simple workflow and unambiguous results.

For these confirmatory studies, short amplicons, usually covering only the region to be confirmed, need to be sequenced. Moreover, minor allelic variants, present in a heterogeneous sample, can be identified by Sanger sequencing and software processing from an electropherogram that contains mixed base calls.

For these types of NGS confirmation studies, BigDye Terminator or BigDye Direct cycle sequencing chemistries and the BigDye XTerminator purification system are frequently used. The genetic analyzer instruments have run modules optimized for short amplicons, so that confirmation data can be rapidly obtained. Finally, we have developed the specialized Applied Biosystems™ Minor Variant Finder software that can detect allelic variants that are present down to 5% in Sanger sequencing reactions.

To illustrate the performance of the SeqStudio Flex instrument for minor allele confirmation experiments, we focused on the BRAF gene. Standardized samples with *BRAF V600E* variants at 50%, 10%, 5%, and 2.5% minor allele frequencies were used as templates. Primers covering 172 nt surrounding the *BRAF V600E* mutation were used to generate sequencing amplicons

from these samples in duplicate. The resulting amplicons were processed using BigDye Direct cycle sequencing and BigDye XTerminator purification chemistries. The purified reactions were run on SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers, and analyzed using Sequence Scanner and Minor Variant Finder software.

We successfully detected the *BRAF V600E* variants at each of the concentrations tested (Figure 12A). In the results view of Minor Variant Finder, the frequency of the minor variant calculated by the software can be seen in each of the samples. In the chromatogram view of Minor Variant Finder, the forward and reverse traces for the control and test sample, as well as the software-processed traces, can be used to verify the software call of a minor allele.

Each of the genetic analyzers produced very similar results in the frequency of the minor allele calculated (Table 3), even below 5% minor allele frequency. We also compared the sequencing quality metrics from these reactions and found that the three instruments produced highly similar quality run data (Figure 12B). These results demonstrate that the SeqStudio Flex instrument can confirm the presence of minor allelic variants by Sanger sequencing with a performance equivalent to 3500xL and SeqStudio genetic analyzers.

**Table 3. Quantification of minor allele determinations on the three genetic analyzer platforms.** Note that although the measured frequency was higher than expected, it was equivalently high on all instruments, suggesting an anomaly of the sample rather than the instrument.

Sample	Expected	Measured: SeqStudio Flex instrument	Measured: 3500xL instrument	Measured: SeqStudio instrument
50% <i>BRAF V600E</i>	50	62.4	62.6	62.8
10% <i>BRAF V600E</i>	10	14.8	13.9	15.1
5% <i>BRAF V600E</i>	5	7.8	6.9	7.0
2.5% <i>BRAF V600E</i>	2.5	4.3*	4.8*	4.2*

\* Although a 2.5% concentration was tested and detected in this experiment, the ability to call at this low frequency may vary in other samples and with other alleles.

## A Results view

SeqStudio Flex genetic analyzer

Amplicon	Test Specimen	Chrom...	Position	Variant	Edi...	Review status	Origin	Review Indicator	For...	Rev...	Type	Variant Comments
BRAF	V600E-50	BRAF	59	A->T	N	Accepted	MVP	54.4	69.8	SNV	<MVP_UbernA	
BRAF	V600E-50	BRAF	59	A->T	N	Accepted	MVP	55.4	70.1	SNV	<MVP_UbernA	
BRAF	V600E-10	BRAF	59	A->T	N	Accepted	MVP	11.2	18.2	SNV	<MVP_UbernB	
BRAF	V600E-10	BRAF	59	A->T	N	Accepted	MVP	10.1	19.7	SNV	<MVP_UbernB	
BRAF	V600E-05	BRAF	59	A->T	N	Accepted	MVP	5.6	13.2	SNV	<MVP_UbernC	
BRAF	V600E-05	BRAF	59	A->T	N	Accepted	MVP	4.3	7.8	SNV	<MVP_UbernC	
BRAF	V600E-2.5	BRAF	59	A->T	N	Accepted	MVP	3.3	5.2	SNV	<MVP_UbernC	

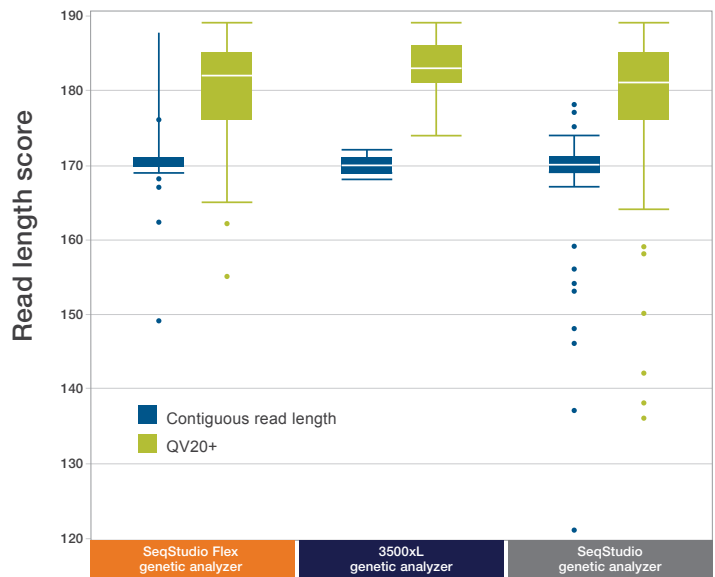
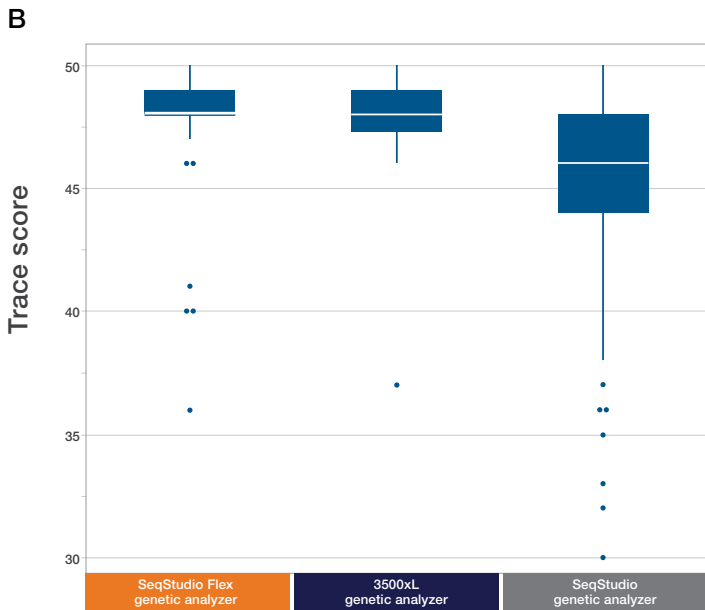
3500xL genetic analyzer

Amplicon	Test Specimen	Chrom...	Position	Variant	Edi...	Review status	Origin	Review Indicator	For...	Rev...	Type	Variant Comments
BRAF	V600E-50	BRAF	59	A->T	N	None	MVP	57.2	68.8	SNV	<MVP_UbernA	
BRAF	V600E-50	BRAF	59	A->T	N	None	MVP	57.6	65.5	SNV	<MVP_UbernA	
BRAF	V600E-10	BRAF	59	A->T	N	None	MVP	11.2	15.6	SNV	<MVP_UbernB	
BRAF	V600E-10	BRAF	59	A->T	N	None	MVP	12.6	16.1	SNV	<MVP_UbernB	
BRAF	V600E-05	BRAF	59	A->T	N	Accepted	MVP	4.3	9.2	SNV	<MVP_UbernC	
BRAF	V600E-05	BRAF	59	A->T	N	None	MVP	5.2	9.0	SNV	<MVP_UbernC	
BRAF	V600E-2.5	BRAF	59	A->T	N	None	MVP	3.6	6.2	SNV	<MVP_UbernC	

SeqStudio genetic analyzer

Amplicon	Test Specimen	Chrom...	Position	Variant	Edi...	Review status	Origin	Review Indicator	For...	Rev...	Type	Variant Comments
BRAF	V600E-50	BRAF	59	A->T	N	None	MVP	56.8	69.9	SNV	<MVP_UbernA	
BRAF	V600E-50	BRAF	59	A->T	N	None	MVP	56.5	68.2	SNV	<MVP_UbernA	
BRAF	V600E-10	BRAF	59	A->T	N	None	MVP	13.9	17.1	SNV	<MVP_UbernB	
BRAF	V600E-10	BRAF	59	A->T	N	None	MVP	10.9	16.6	SNV	<MVP_UbernB	
BRAF	V600E-05	BRAF	59	A->T	N	Accepted	MVP	5.6	8.4	SNV	<MVP_UbernC	
BRAF	V600E-05	BRAF	59	A->T	N	None	MVP	8.2	8.9	SNV	<MVP_UbernC	
BRAF	V600E-2.5	BRAF	59	A->T	N	None	MVP	3.1	6.2	SNV	<MVP_UbernC	

## Chromatogram view



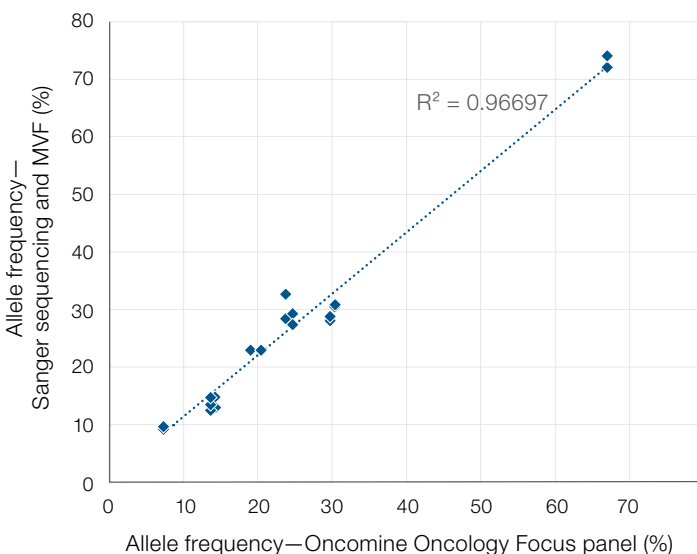
**Figure 12. Detecting minor alleles by short amplicon sequencing.** (A) Minor Variant Finder software can analyze Sanger sequencing traces and detect sequence variants that are present at 5% of the total. Templates containing *BRAF V600E* variants at 50%, 10%, 5%, and 2.5% were analyzed on the SeqStudio Flex, 3500xL, and SeqStudio instruments. In each case, the software was able to detect the 5% variant reproducibly (yellow highlights, left). The Sanger traces of control, test, and software-processed samples can be viewed in the chromatogram view (right). The traces looked identical on the three platforms; the allele quantity determined by the software in each set is indicated with the yellow highlight. (B) Trace scores (left) and contiguous read length and QV20+ values (right) were compared on all three instruments. Sequencing quality results for these small amplicons were almost identical.

## Oncology research, next-generation confirmation

The SeqStudio Genetic Analyzer can be used by clinical researchers to maintain high quality for detecting and verifying the presence of mutant alleles in tumor tissue. The SeqStudio system integrates with the following tools to simplify Sanger sequencing workflows:

- The SeqStudio Genetic Analyzer comes preloaded with running modules optimized for fragmented DNA extracted from formalin-fixed, paraffin-embedded tissue.
- The cloud-based NGC module allows investigators to rapidly verify variants identified in next-generation sequencing (NGS) .vcf files using Sanger sequencing traces.
- Allelic variants at frequencies down to 5% can be detected using the Applied Biosystems™ Minor Variant Finder (MVF) Software and Sanger traces generated by the SeqStudio instrument.
- Applied Biosystems™ BigDye™ Direct and BigDye XTerminator™ chemistries simplify the Sanger sequencing workflow by providing one-tube sequencing and cleanup.

The performance of the SeqStudio Genetic Analyzer for detecting mutant alleles in tumor samples was determined by analyzing genomic DNA extracted from 10 different formalin-fixed, paraffin-embedded (FFPE) tumor samples, and determining variant frequencies at four different hotspot regions. The frequency of mutant alleles was determined by NGS using the Ion Torrent™ Oncomine™ Oncology Focus Panel, and Sanger sequencing using BigDye Direct/BigDye XTerminator chemistries and MVF Software. The correlation between the frequencies measured by the SeqStudio Genetic Analyzer was excellent when compared to NGS at allele frequencies—from about 9% to about 70% (Figure 13).



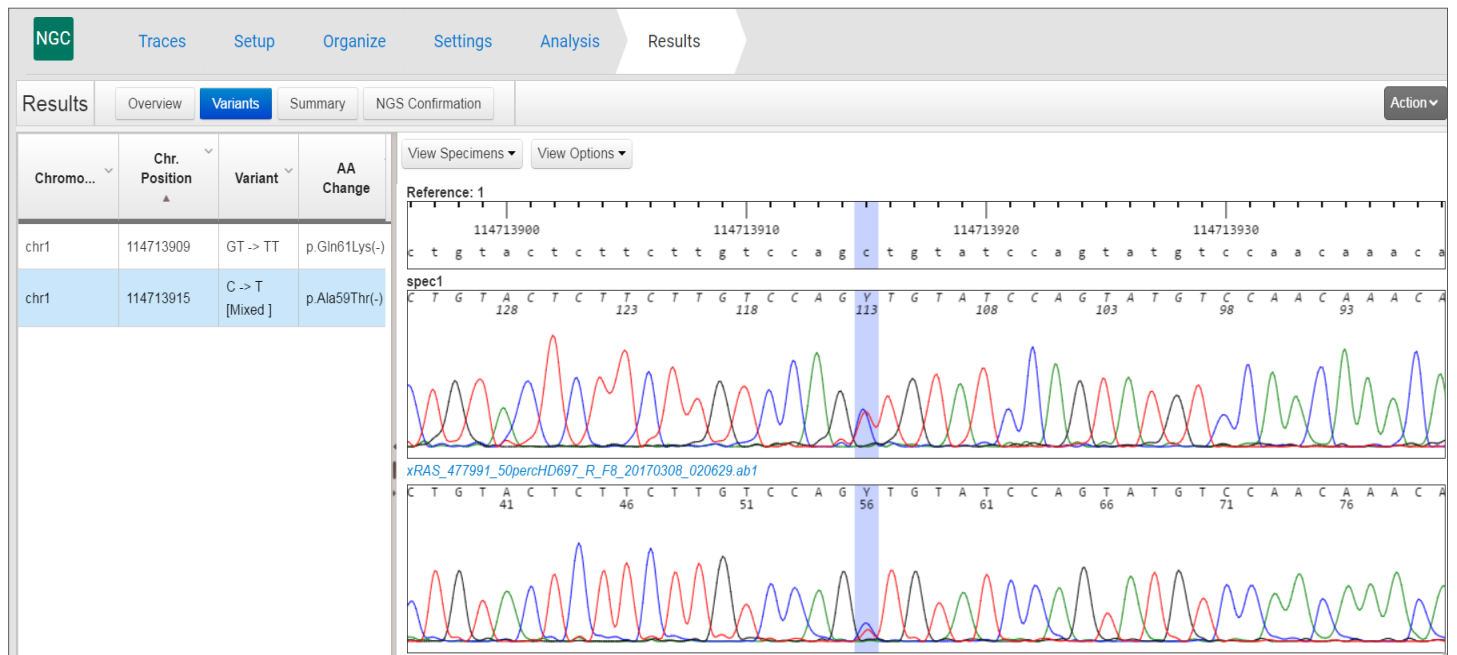
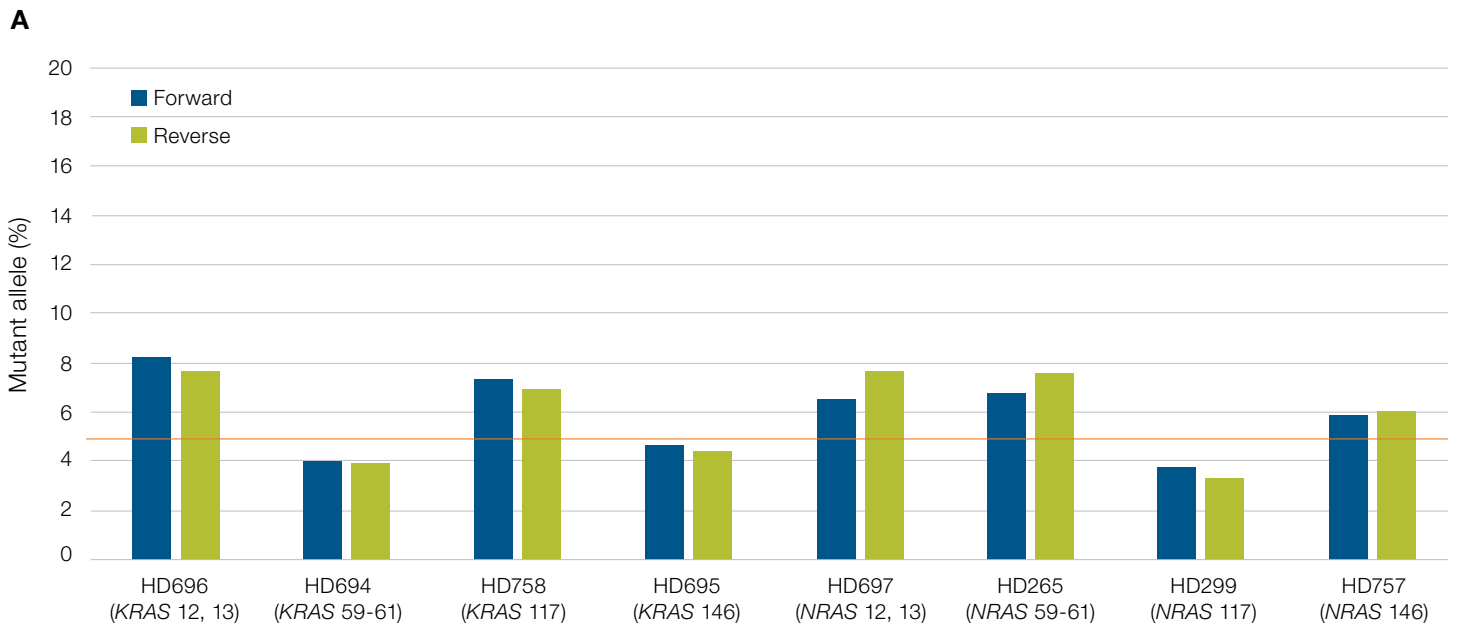
The ability of the SeqStudio Genetic Analyzer to analyze variant frequencies was also determined using a 96-well plate containing Sanger sequencing primers that query the most common tumorigenic mutations in KRAS and NRAS. The minor allele frequency analysis of SeqStudio instrument traces accurately measured the allele frequencies in 1 ng of diluted FFPE-extracted DNA (Figure 14A). Therefore, researchers needing to detect rare alleles can be confident that the SeqStudio Genetic Analyzer will produce accurate results on FFPE tissues. For more details, see reference 11.

Finally, the cloud-based NGC application simplifies the confirmation of variants identified by NGS by organizing Sanger sequencing traces by amplicons and specimens, and aligning them in the proper orientation to the candidate variant sequences in a .vcf file. To show the utility of the NGC app in an oncology workflow, we confirmed the presence of an NRAS mutation identified using the Oncomine Oncology Focus panel by Sanger sequencing (Figure 14B). The SeqStudio results verified that the mutation in NRAS (p.Ala59Thr) was present. Therefore, focused and rapid examination of the most meaningful portions of sequencing traces by the NGC app facilitates NGS variant confirmation.

**Read more about next-generation sequencing confirmation with Sanger sequencing at [thermofisher.com/ngsconfirmation](https://thermofisher.com/ngsconfirmation)**

**Figure 13. Analysis of FFPE samples using the SeqStudio instrument at allele frequency of 9–70%.** Ten different tumor samples were analyzed for mutations at four different hotspots by Ion Torrent™ sequencing and Sanger sequencing on the SeqStudio instrument. The mutant allele frequency correlated extremely well between the two methods across a wide range of allele frequencies.





**Figure 14. Analysis and confirmation of variants by SeqStudio Genetic Analyzer and the NGC application, respectively. (A)** Eight different FFPE samples with mutations at known RAS hotspots were diluted to 5% allele frequency, then analyzed using a 96-well plate containing Sanger sequencing primers that query the most common tumorigenic mutations in KRAS and NRAS, and using the SeqStudio Genetic Analyzer. Each of the allele queries accurately measured the allele frequencies; deviations from 5% reflected slight inconsistencies in starting concentration of the samples. Orange line is 5% frequency. Similar results were seen with 10% and 50% dilutions. **(B)** Confirmation of variants identified by NGS. From a .vcf file generated using Ion Reporter™ Software, Sanger sequencing primers targeting loci of interest were ordered from Primer Designer, samples were sequenced on the SeqStudio instrument, and variants common to the .vcf file and the Sanger sequencing traces were highlighted using the NGC cloud app.

## Microbial Species identification

The ready availability of genomic data opens the opportunity to identify species in an unknown sample by sequencing DNA of “fingerprint” loci. In the Applied Biosystems™ family of kits, for example, MicroSEQ™ kits have simplified the identification of prokaryotes and fungi by Sanger sequencing ribosomal DNA (rDNA) sequences [12]. Similarly, eukaryotic organisms can be identified using the mitochondrial-specific loci as the identifying locus. This strategy has been exploited in the Barcode of Life project (barcodeoflife.org, for review see reference 13), providing a means for rapidly establishing the identity of unknown eukaryotic samples.

To illustrate the performance of the SeqStudio Genetic Analyzer for microbial identification, we obtained genomic DNA samples from ATCC for a variety of microorganisms, and sequenced them using the Applied Biosystems™ MicroSEQ™ 500 PCR kit and the SeqStudio instrument. The resulting sequences were queried against the BLAST database. For each sequencing reaction, the correct organism was identified with the highest BLAST confidence. Similarly, using primers for fish mitochondrial sequences (CO1 gene) and fish samples, the fish species was correctly identified as the top BLAST hit. The accurate identification of the species queried with BLAST illustrates how well the SeqStudio platform can be used for species identification.

Learn more at [thermofisher.com/microseq](http://thermofisher.com/microseq)

**Table 4. Analysis of species ID using the SeqStudio Genetic Analyzer.** Samples of microorganism DNA or genomic DNA extracted from fish were sequenced using primers for 16S rDNA and the MicroSEQ kit (BigDye Terminator v1.1 chemistry), or using primers for fish mitochondrial CO1 sequences and BigDye Terminator v3.1 chemistry.

	Number of organisms	Number of queries	Percent correct
Microorganisms	24	48	100
Piscine organisms	12	24	100

## Sequencing of difficult templates

Some organisms and regions of the human genome are difficult to sequence due to a high GC content. These regions cause polymerases to stutter, stop, or disengage entirely from a sequencing template, precluding high-quality results. Furthermore, some regions of genomes contain a large amount of highly repetitive sequences. NGS platforms can have difficulty aligning such sequences to a reference genome. Because of its robust chemistry and unambiguous results without the need of a reference genome, Sanger sequencing can overcome many of these difficulties.

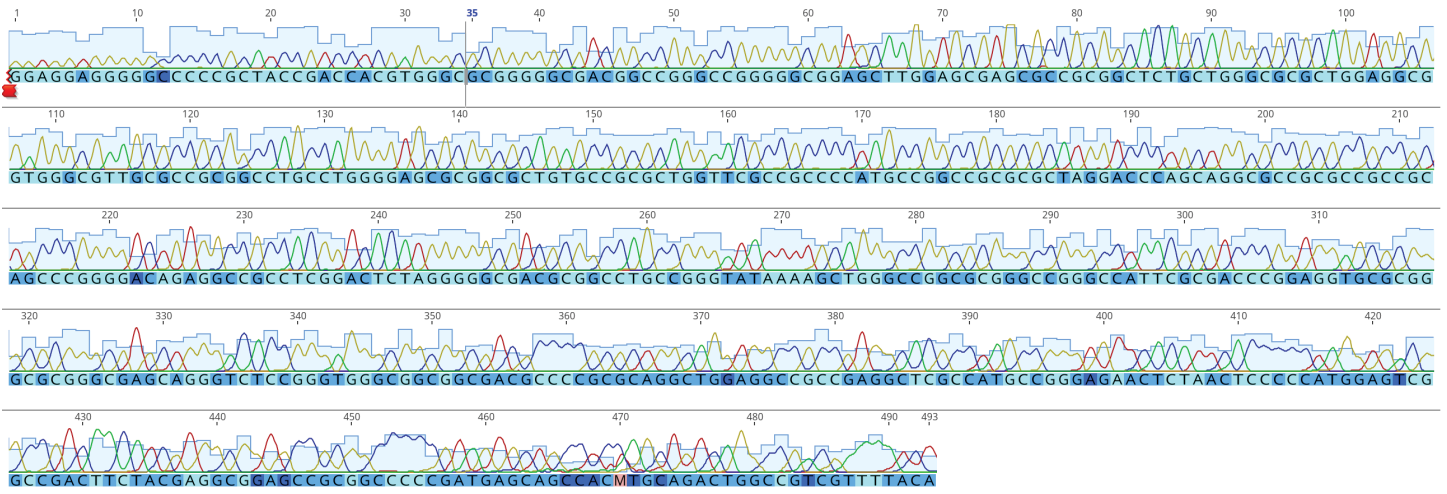
To assist with the PCR amplification of GC-rich sequences in a genome, we developed a GC enhancer reagent. This add-in reagent helps DNA polymerases generate templates from GC-rich sequences that can be used for cycle sequencing. BigDye Terminator or BigDye Direct sequencing chemistries are subsequently used, followed by purification and capillary electrophoresis on a genetic analyzer. To analyze GC-rich or highly repetitive sequences, the genetic analyzers require no special run modules.

To determine the quality of sequence obtained from difficult templates, we analyzed the human CEBPA gene, a gene with regions on average 66% GC. Primers were chosen to amplify eight different regions in the human CEBPA gene. These regions were amplified using Applied Biosystems™ AmpliTaq Gold™ polymerase and the GC enhancer reagent. Amplicons were sequenced using BigDye Direct cycle sequencing and BigDye XTerminator purification chemistries. The resulting reactions were run on SeqStudio Flex, 3500xL, and SeqStudio genetic analyzers.

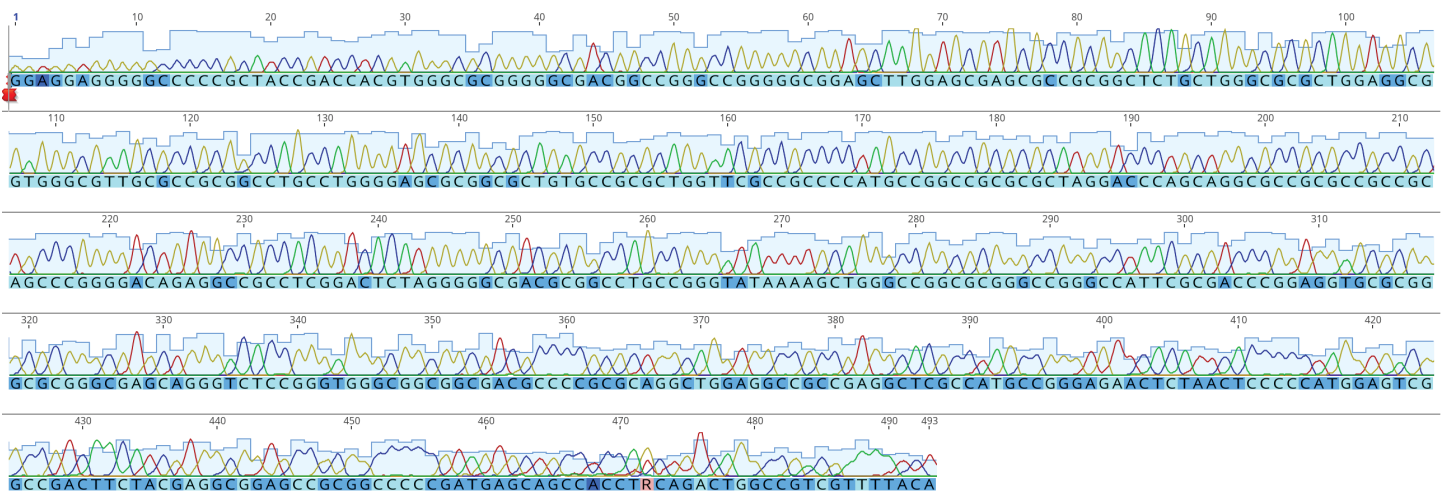
All of the instruments were able to produce high-quality results from these GC-rich amplicons. Comparing the electropherograms obtained on SeqStudio Flex and 3500xL instruments revealed nearly identical traces (Figure 15A). Although only one representative sequencing trace is shown, similar results were seen with other amplicons and in both sequencing directions. In support of these results, the QC analysis revealed nearly identical values for each of the metrics (Figure 15B). Together, these results show that even with difficult templates, the SeqStudio Flex instrument produces data equivalent to that generated on other Applied Biosystems genetic analyzers.

A

SeqStudio Flex genetic analyzer



3500xL genetic analyzer



B



**Figure 15. Sequencing through difficult regions.** Portions of the human CEBPA gene, which averages 66% GC content, were amplified with the GC enhancer reagent before sequencing. **(A)** Chromatograms from one amplicon, obtained on the SeqStudio Flex and 3500xL genetic analyzers, are shown. This particular region is 77.8% GC. Note that the chromatograms are nearly identical. **(B)** Trace scores, contiguous read length, and QV20+ values were compared for all amplicons on all three instruments. Sequencing quality results for these difficult amplicons were almost identical.

## Sanger sequencing for genome edit efficiency analysis

Genome editing has revolutionized the way biological research is performed. Because of its simplicity and efficiency, an investigator can make any type of a change to a genome, from engineering a specific SNP, to knocking out the function of a gene, to making large-scale genomic rearrangements.

Most researchers describe a continuous genome editing process that starts with design of a guide RNA and ends with a stable cell line containing the desired edit. However, throughout the process, there are steps where knowing the efficiency of the modification is necessary. Often, this involves monitoring a cell culture for the presence and frequency of a desired edit. In many cases, the efficiency challenge is linked to the efficiency of gRNA and CRISPR-related enzymes. Researchers therefore need unambiguous and straightforward methods for analyzing the efficiency of their genome editing reactions.

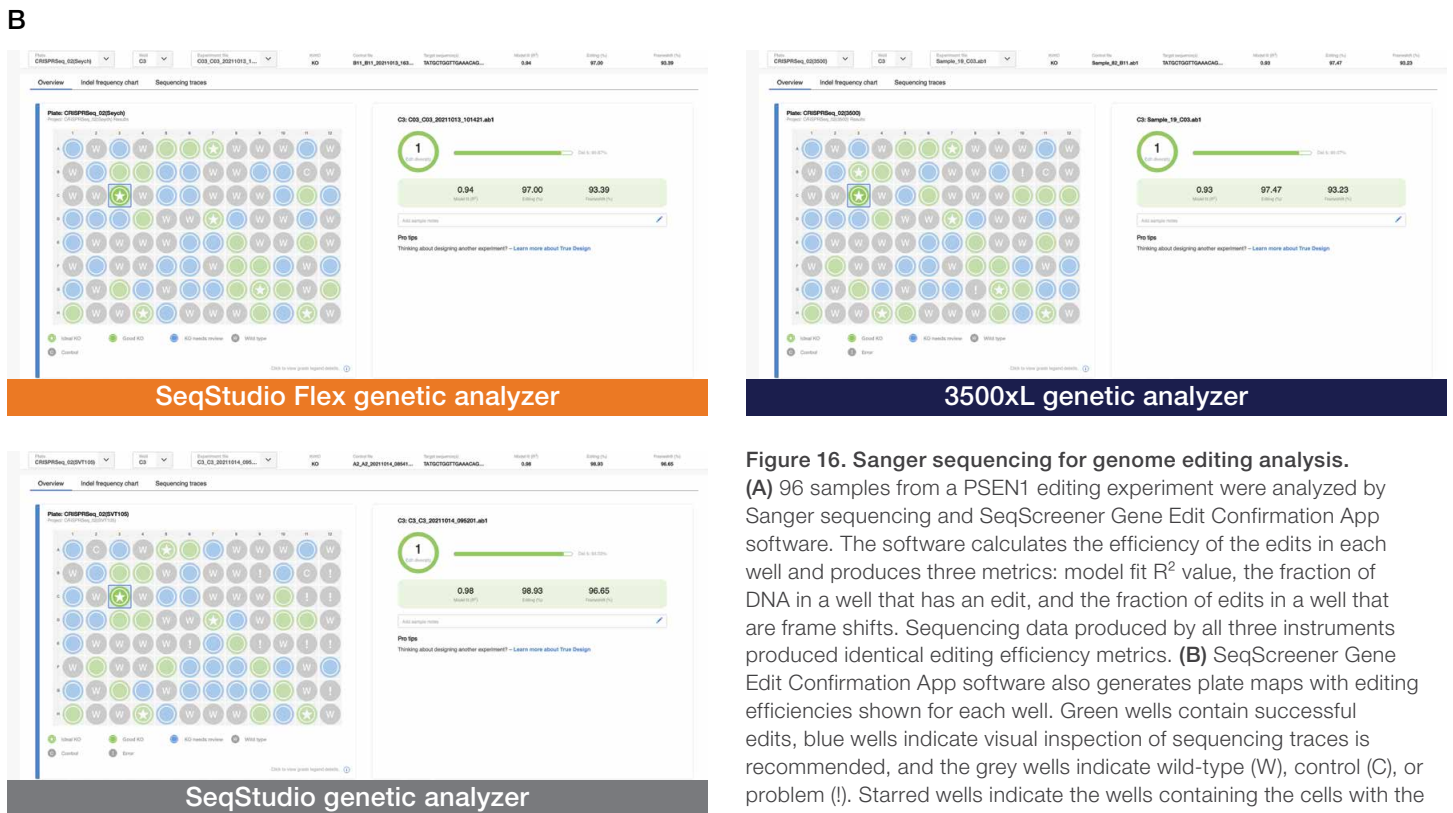
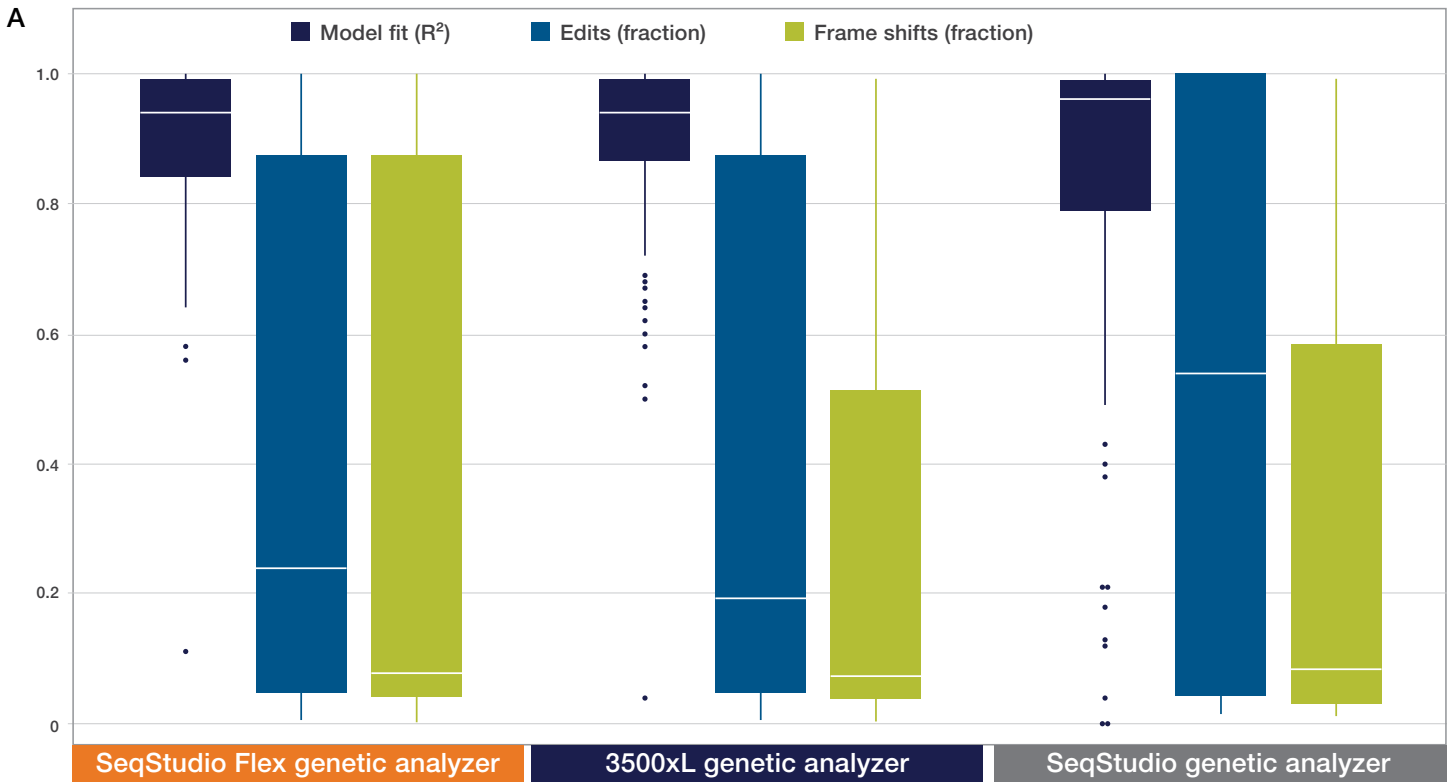
Analysis of the steps in a genome editing experiment often make use of cycle sequencing using BigDye Terminator or BigDye Direct chemistries, coupled with purification and electrophoresis on genetic analyzers. To facilitate the analysis of genome editing experiments, we developed the Applied Biosystems™ SeqScreener Gene Edit Confirmation App. This groundbreaking software is used to analyze the efficiency of edits and visualize subcultures that contain the optimal edit.

To characterize the genome editing analysis on the SeqStudio Flex instrument, we generated 96 CRISPR-directed edits in the human presenilin gene (PSEN1). For this experiment, we used different guide RNAs, transfection methods, and homology-directed repair templates, resulting in a diversity of editing efficiencies. gDNA from transfected cultures was isolated and sequenced using PSEN1 primers, BigDye Direct cycle sequencing, BigDye XTerminator purification, and electrophoresis on SeqStudio Flex, 3500xL, and SeqStudio instruments. The resulting sequencing traces were analyzed using Sequence Scanner and SeqScreener Gene Edit Confirmation App software.

The SeqScreener Gene Edit Confirmation software computes and displays results in several different ways. One way is by generating numerical metrics for the results (Figure 16A). The Model Fit ( $R^2$ ) is the overall confidence in the results. Typically, good-quality results have  $R^2$  values of 0.8 or greater; however, the threshold can vary for each sample. The Edits value is the cumulative fraction of all successful edits in the measured sample. The wide range of values indicates the wide range of efficiencies across all the different editing conditions tested. The Frame Shift value is the fraction of gene edits that are predicted to eliminate the expression of the encoded protein. This information is useful when screening for loss-of-function gene knockouts. The three genetic analyzers produced very similar genome editing efficiency results.

Another way SeqScreener software displays results is in an infographic showing the status of analysis in wells of a 96-well plate, modeling the results contained within the test plate (Figure 16B). In this representation, green wells contain successful edits, blue wells indicate visual inspection of sequencing traces is recommended, and the grey wells indicate wild-type (W), control (C), or problem (!). Starred wells indicate the wells containing the cells with the highest-efficiency, idealized edit. Clicking on a well brings up more information about the success in that well. The plate infographic produced by the three instruments was very similar. There were a few differences, mostly in the visual inspection vs. problem categories in the SeqStudio system compared to the SeqStudio Flex and 3500xL systems. This was likely due to the degradation in sequence read quality, since the plate was used multiple times. Nevertheless, these results demonstrate that the performance of the SeqStudio Flex Genetic Analyzer is equivalent to the 3500xL and SeqStudio systems for monitoring genome editing efficiencies.

**Read more about this application at**  
[thermofisher.com/ce-geneedit](https://www.thermofisher.com/ce-geneedit)

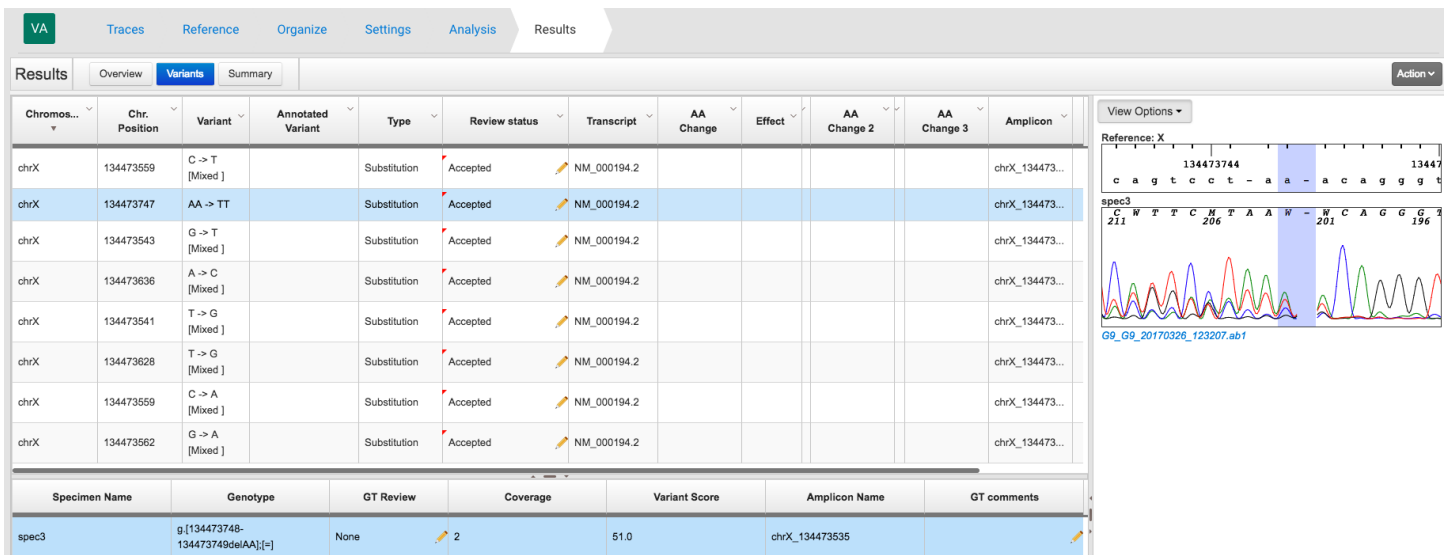


**Figure 16. Sanger sequencing for genome editing analysis.** (A) 96 samples from a PSEN1 editing experiment were analyzed by Sanger sequencing and SeqScreener Gene Edit Confirmation App software. The software calculates the efficiency of the edits in each well and produces three metrics: model fit  $R^2$  value, the fraction of DNA in a well that has an edit, and the fraction of edits in a well that are frame shifts. Sequencing data produced by all three instruments produced identical editing efficiency metrics. (B) SeqScreener Gene Edit Confirmation App software also generates plate maps with editing efficiencies shown for each well. Green wells contain successful edits, blue wells indicate visual inspection of sequencing traces is recommended, and the grey wells indicate wild-type (W), control (C), or problem (I). Starred wells indicate the wells containing the cells with the highest-efficiency, idealized edit. Clicking on a well will bring up more information (to the right of the plate map) about the success in that well. The plate infographics produced by the three instruments were very similar; most of the differences related to the need for visual inspection rather than outright discrepancies.

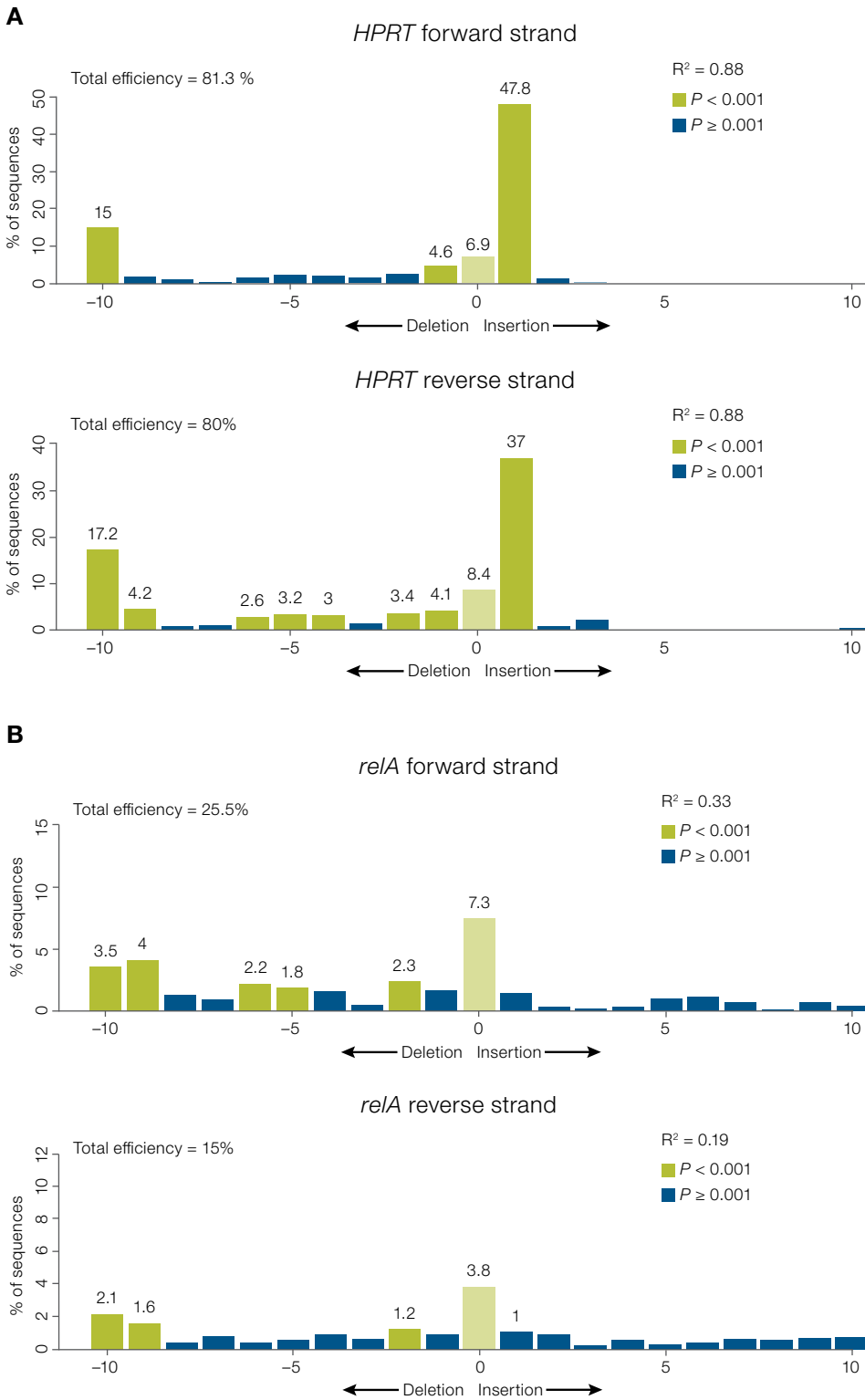
## CRISPR-Cas9 genome editing analysis

Genome editing technologies, including CRISPR-Cas9-mediated editing events, are rapidly becoming accessible to a majority of life science researchers and are poised to revolutionize all fields of biology and health care. Thermo Fisher Scientific offers all the tools necessary for a genome editing project. As an integral part of such a project, the features of the SeqStudio Genetic Analyzer facilitate Sanger sequencing analyses and fit well within a genome editing workflow. In particular, the data generated are compatible with Tracking of Indels by Decomposition (TIDE) software [14], a widely available tool for analyzing the efficiency of genome editing events.

The utility of the SeqStudio Genetic Analyzer in a genome editing project was shown by obtaining whole-cell lysates from HEK293 cells that were edited to introduce random deletions around a targeted site in the HPRT or the relA locus. To confirm the position of the edit, the Sanger sequencing traces were uploaded to the cloud and analyzed using the Sanger Variant Analysis module (Figure 17). Note that the position of the edit is clearly indicated and can be visualized by the abundant mixed-base peaks downstream of the break. The efficiency of the edits in this mixed primary culture was determined by analyzing these trace files using the TIDE software. In each case, the spectrum and frequencies of deletions at each locus was nearly identical using the data generated in the forward and reverse directions (Figure 18). These frequencies confirm results obtained using Invitrogen™ TOPO™ cloning and followed by Sanger sequencing results of the same edited cell populations [15].



**Figure 17. Analysis of genome-edited samples using the SeqStudio instrument.** A mixed population of cells with a genome editing event at the human HPRT locus was analyzed using the cloud-enabled Sanger Variant Analysis app. Note that this app finds single-nucleotide variants common to both forward and reverse strands, but is also able to detect where the genome cleavage event occurred, producing a population of mixed sequences downstream (to the right in this example) of the breakpoint.



**Figure 18. Analysis of two different genome editing events at the *HPRT* and *relA* loci using TIDE software and mixed population sequencing traces generated by the SeqStudio instrument.** The bars show the proportion of the population having the indicated number of nucleotides deleted or inserted. For (A) *HPRT*, the overall efficiency of the edit was around 80%, whereas the overall efficiency at the (B) *relA* locus was around 20%.



## Detection of KRAS mutations by Sanger sequencing

Kirsten rat sarcoma viral oncogene homologue (KRAS) is a hyper mutated oncogene across all cancer types. Ninety-eight percent (98%) of KRAS mutations occur within codons 12 or 13 (exon 2) and codon 61 (exon 3). A variety of mutations within these codons exist and their frequency varies by cancer type. Knowledge of specific KRAS mutations can inform cancer treatment and prognosis. Sanger sequencing is useful for the detection of known and unknown mutations within this region.

Using pre-designed primers and a simple workflow on the SeqStudio Genetic Analyzer, results can be obtained in as little as 6 hours. To test this assay reference DNA standards were purchased containing common KRAS mutations as well as wild type controls. Four primers were used for amplification, 2 forward and 2 reverse for exons 2 and 3. Samples were amplified then cycle sequenced using BigDye Direct Sequencing Master Mix. Reactions were cleaned up and run on the SeqStudio Flex genetic analyzer. Results were analyzed using SeqScape v4.1 software. All expected mutations were accurately detected. See sample results below (figure 19).

Read more about this application at [thermofisher.com/kras](https://thermofisher.com/kras)

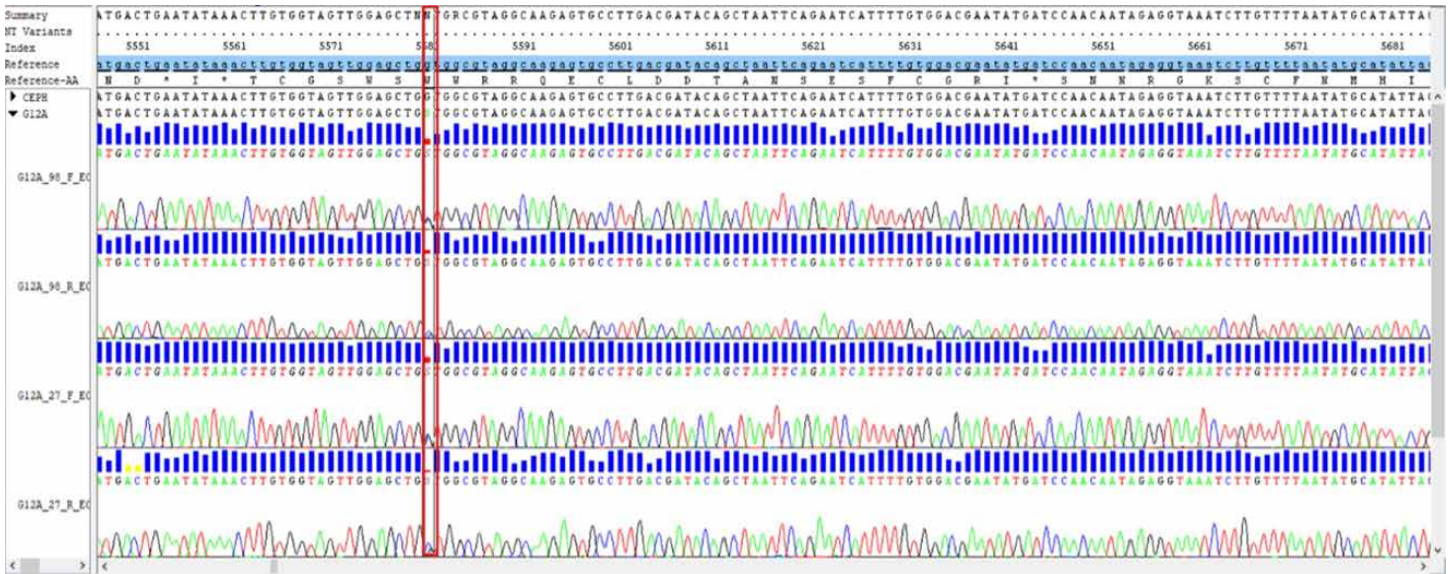


Figure 19. Detection of KRAS G12A mutation using two different PCR primer assays and forward and reverse sequencing primers.

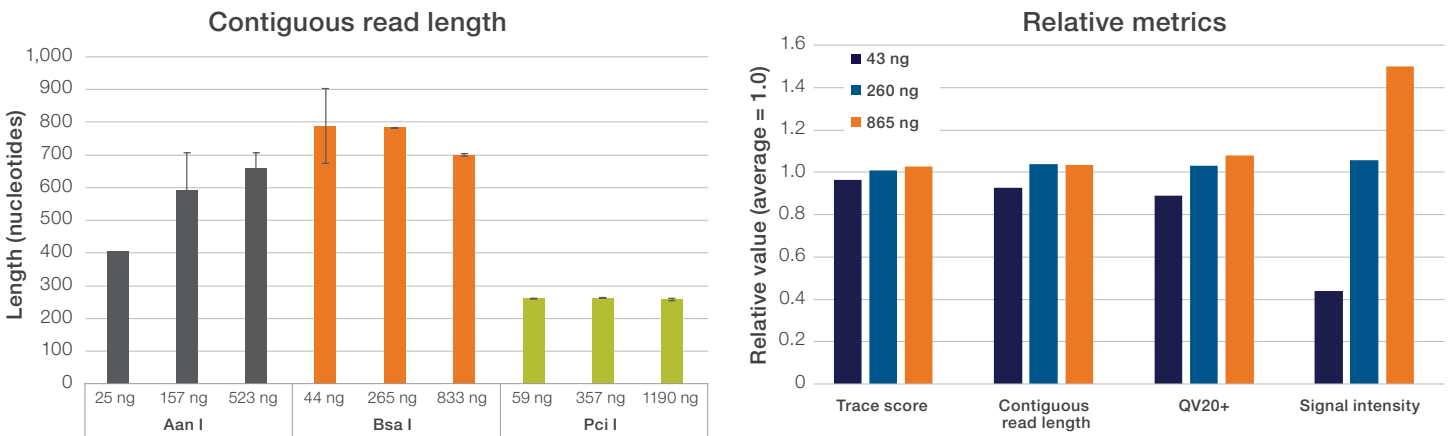
## mRNA vaccine quality control using Sanger sequencing

The use of mRNA for vaccines and therapeutics is an emerging area driven by the rapid and successful development of the COVID-19 vaccine. Sequence confirmation is a critical activity during mRNA therapeutic and vaccine manufacturing, both during plasmid production and mRNA identity testing. Sanger sequencing enables accuracy, workflow simplicity, scalability, and reproducibility to help ensure testing consistency and final product safety.

There are several methods that can be used in the QC process (20). One convenient method to QC the final mRNA product is sequencing directly from cDNA. The cDNA is generated from mRNA using Invitrogen Super Script IV VILO Master Mix. After removal of primers with ExoSAP-IT PCR Product Cleanup reagent, the reaction product can go into the Big Dye Terminator cycle sequencing reaction. It is then cleaned up and run on a SeqStudio Genetic Analyzer. Data can then be analyzed using Sequence Scanner Software 2.0. The entire protocol takes 4-6 hours.

To test this protocol, transcripts were generated from restriction digested pGEM plasmid. Different amounts of RNA were used to determine minimal inputs and the quality of data obtained. Good quality reads were generated with as little as 44 ng of RNA. While most metrics were constituent, signal intensity increased with increasing amounts of input mRNA.

Read more about this application at [thermofisher.com/ce-mrna](https://thermofisher.com/ce-mrna)



**Figure 20.** Left. Read length of indicated quantities of starting mRNA from transcripts generated from different restriction digest products from a pGEM vector. Right. Relative metrics from indicated quantities of mRNA transcript of BsaI restriction digested pGEM vector.

## Analyzing methylated DNA by Sanger sequencing

Changes in DNA methylation patterns on specific sequences are useful biomarkers for many pathological states including cancer (21). In eukaryotes, DNA methylation is usually found as 5-methylcytosine (5-meC) in CpG sites that are present in promoters and enhancers of genes. These methylated sites can affect the binding of DNA recognition proteins, and when DNA methylation status changes at these sites, it often results in changes in transcription of the associated gene.

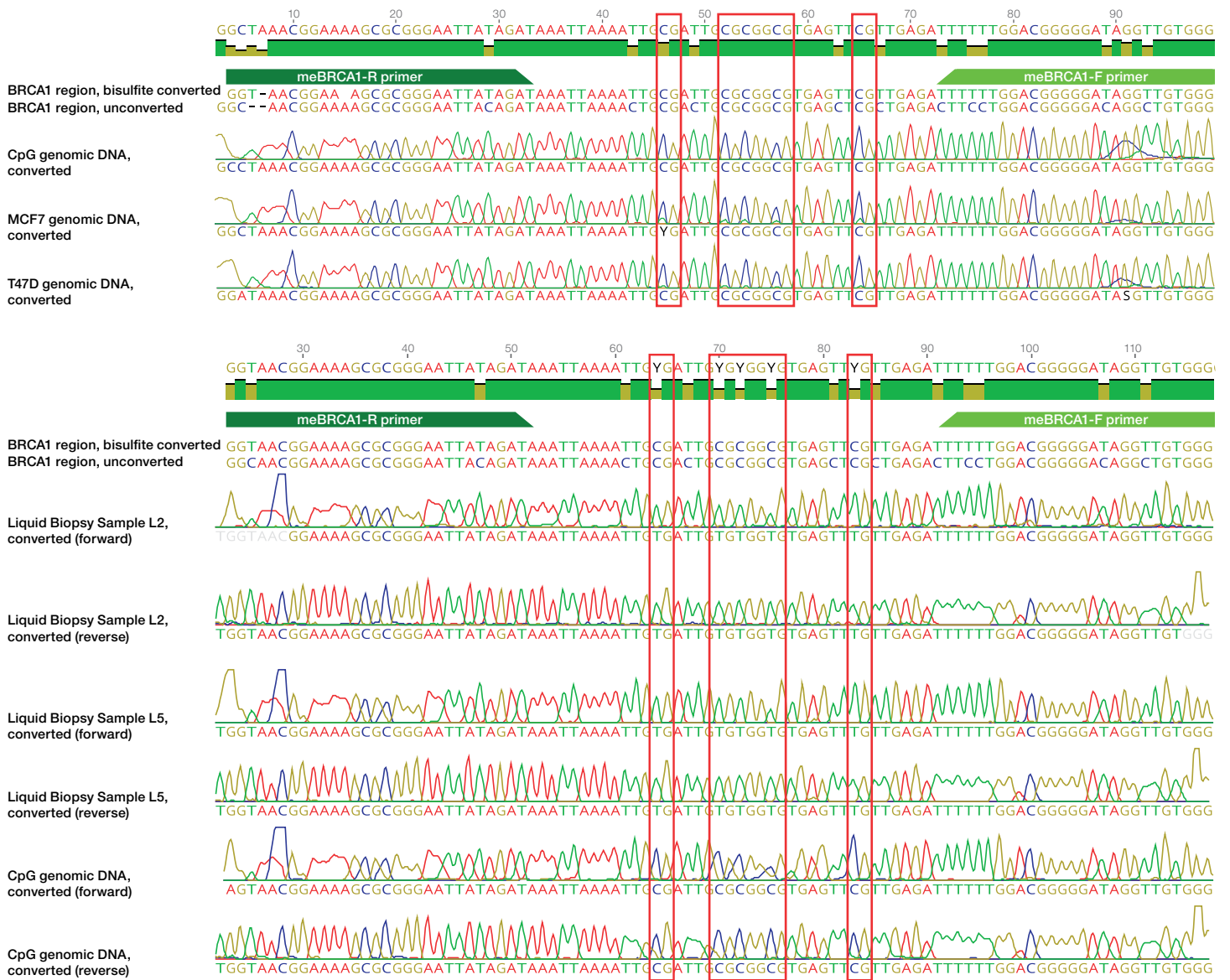
Sanger sequencing is a useful tool for the detection of DNA methylation from many sample types including circulating free DNA (cfDNA) and formalin-fixed, paraffin-embedded (FFPE). Sanger offers the benefit of a relatively simple, fast, and low cost workflow with straightforward data analysis. Additionally, a single reaction can give information about all positions in the amplicon, facilitating the detection of partially modified sequences.

Sequencing methods cannot detect the difference between 5-methylcytosine and normal cytosine – both pair equally well with guanine. However, unmethylated cytosine can be deaminated by treating with bisulfite, converting it to uracil that subsequently behaves as thymidine. Therefore, bisulfite conversion is commonly used for the detection of methylated DNA. Bisulfite conversion does not affect 5-methylcytosine, and it therefore continues to basepair as cytosine. Thus, after

bisulfite conversion, any unmethylated cytosines change from CG base pairs to TA base pairs in subsequent assays, while any methylated cytosines maintain the CG basepairing. Bisulfite conversion completely changes a sequence, resulting in two new non-complementary strands.

To show this application in an example, genomic DNA (gDNA) was isolated from cell lines and cfDNA. Isolated DNA was treated with bisulfite to convert unmethylated Cs to Us. Next, a region of the BRCA1 promotor was amplified using PCR primers that have been designed for the unique bisulfite-converted sequence. Note that DNA obtained from liquid biopsy (or FFPE samples) is fragmented, and the bisulfite conversion process fragments it further. Therefore, it is best to use primers that generate amplicons in the 80-120bp range. After PCR, the reaction was cleaned up by removing free primers using the ExoSapIt kit. A portion of the PCR reaction was then used in the cycle sequencing reaction, followed by cleanup with BigDye XTerminator. Finally, the reactions were run on the SeqStudio Flex Genetic Analyzer and peaks analyzed using standard software. In the data below, changes in BRCA1 methylation were detected in MCF7 cells but not T47D cells. Both liquid biopsy samples were unmethylated compared to controls.

**Find more on DNA methylation analysis for liquid biopsy and FFPE samples at [thermofisher.com/ce-methylation](https://www.thermofisher.com/ce-methylation)**



**Figure 21. Analysis of methylation sites in BRCA1 promoter.** In both panels, the predicted bisulfite converted and unconverted sequences are shown at the top, as well as the position of the primers used for PCR and cycle sequencing. Note that the position numbers are relative to the input sequence, and not the genomic coordinates. The electropherograms of each sample are shown below the reference sequences. **A.** Analysis in common breast cancer cell lines. In both cell lines, the CpGs highlighted by red box contained methylcytosine. In the MCF7 cells, the C at position 46 is partially methylated (indicated by a Y call). **B** Analysis in liquid biopsy samples. In both L2 and L5 liquid biopsy samples, this region is unmethylated. The same sequence was observed when sequenced either direction. For comparison, the CpG fully methylated control DNA sample is shown.

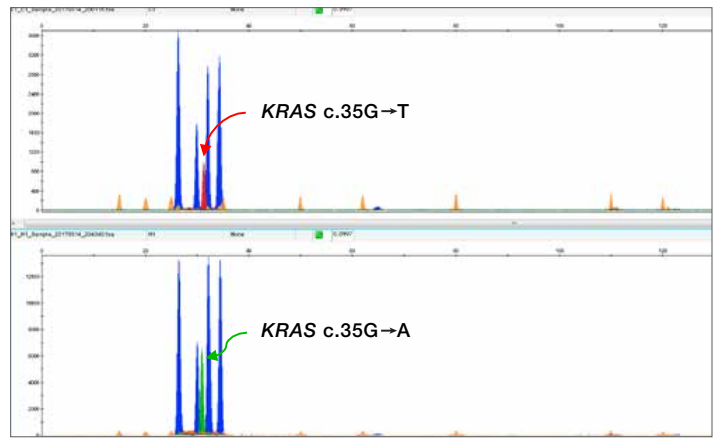
## SNaPshot Multiplex System for SNP genotyping

The ability to detect single-nucleotide polymorphisms (SNPs) plays a critical role in understanding how the genome influences biological phenotypes. To analyze SNP variants, the Applied Biosystems™ SNaPshot™ Multiplex System was developed [17]. This kit is customized for your target, offers multiplexing capability (up to 10-plex) and is compatible with all CE instruments.

Customizable, color-coded fragments of differing sizes, corresponding to specific alleles, are analyzed by fragment analysis. The SeqStudio system includes features that facilitate SNaPshot analysis, including built-in reporting of fragment analysis results of size and peak area. Additionally, the ability to mix fragment analysis and sequencing reactions on one plate enables investigators to perform SNP profiling and Sanger sequencing on a single run.

To illustrate the functional utility of the SeqStudio instrument in SNaPshot workflows, genomic DNA from FFPE-preserved tumor slices was collected and analyzed using probes targeting KRAS G12X and G13X alleles using the SNaPshot multiplex reagent kit. The SeqStudio instrument produced results that clearly showed the presence and accurate calls of the different alleles at this position (Figure 22). Note that although the detection of the alleles was accurate on SeqStudio instrument, the absolute migration of all peaks will differ slightly when compared to that in other platforms due to the different chemical nature of the different polymers. Therefore, to associate a peak with an allele without ambiguity, a calibration with known alleles should be performed before undertaking a large-scale analysis.

Read more about this application at [thermofisher.com/SNP](https://thermofisher.com/SNP)



**Figure 22. Analysis of SNPs using kit on the SeqStudio instrument.** One nanogram of FFPE-extracted DNA from two different tumors was processed using the SNaPshot multiplex kit and KRAS-specific primers, followed by fragment analysis on the SeqStudio system. The SNaPshot multiplex kit produces fragments with allele-specific colors and lengths. The four blue peaks represent wild type alleles at KRAS c.34, c.35, c.37, and c.38. The red peak (top graph) indicates that the sample also had a KRAS c.35G → T mutation present, whereas the green peak (bottom graph) results from a different allele at the same position in a different sample.

## Conclusions

Capillary electrophoresis is a powerful method for characterizing genetic information. Sanger sequencing and fragment analysis remain the gold-standard methods for sequencing and DNA analysis.

The SeqStudio family of instruments are intuitive, purpose-built systems that generate high-quality data for a wide variety of applications.

The SeqStudio genetic analyzer is an excellent low-throughput instrument for plasmid sequencing, NGS confirmation, genotyping, CRISPR-Cas9 genome editing confirmation, cell line authentication, and detection of single nucleotide polymorphisms (SNPs).

The SeqStudio Flex instrument is the most advanced Applied Biosystems genetic analyzer, featuring the latest design improvements and technological advances for increased flexibility, easier operation, enhanced connectivity, and remote serviceability. This medium-throughput system excels at the same applications as the SeqStudio genetic analyzer with the addition of multiplexed PCR, microbe or virus identification, microsatellite instability (MSI) analysis and dsDNA quality control.

## References

1. <https://iclac.org/databases/cross-contaminations/>
2. Horbach SPJM, Halffman W (2017) The ghosts of HeLa: How cell line misidentification contaminates the scientific literature. *PLoS ONE* 12(10): e0186281. <https://doi.org/10.1371/journal.pone.0186281>
3. Authenticating human cell lines using CLA Identifiler and CLA GlobalFiler kits on capillary electrophoresis platforms. <http://assets.thermofisher.com/TFS-Assets/GSD/Application-Notes/authenticating-human-cell-lines-str-kits-capillaryelectrophoresis-application-note.pdf>
4. ATCC STR database. <https://www.atcc.org/en/search-str-database>
5. Gologan A, Sepulveda AR (2005) Microsatellite instability and DNA mismatch repair deficiency testing in hereditary and sporadic gastrointestinal cancers. *Clin Lab Med* 25(1):179–196. doi: 10.1016/j.cll.2004.12.001. PMID: 15749237.
6. Zhao P, Li L, Jiang X et al. (2019) Mismatch repair deficiency/microsatellite instability-high as a predictor for anti-PD-1/PD-L1 immunotherapy efficacy. *J Hematol Oncol* 12:54. <https://doi.org/10.1186/s13045-019-0738-1>
7. Reyes GX et al. (2015) New insights into the mechanism of DNA mismatch repair. *Chromosoma* 124:443–462.
8. Boland CR et al. (1998) A National Cancer Institute workshop on microsatellite instability for cancer detection and familial predisposition: development of international criteria for the determination of microsatellite instability in colorectal cancer. *Cancer Res* 58:5248–5257.
9. TrueMark MSI Assay For Microsatellite Instability Testing User Guide. [https://assets.thermofisher.com/TFS-Assets%2FSLG%2Fmanuals%2FMAN0018868\\_MSI%20Assay\\_UG.pdf](https://assets.thermofisher.com/TFS-Assets%2FSLG%2Fmanuals%2FMAN0018868_MSI%20Assay_UG.pdf)
10. Protocol for Sanger sequencing of the SARS-CoV-2 spike (S) gene. <https://assets.thermofisher.com/TFS-Assets/GSD/brochures/sequencingsars-cov-2-spike-gene-protocol.pdf>
11. Low-level somatic variant detection in tumor FFPE samples by Sanger sequencing COL31176 0616.
12. Technical note: MicroSEQ Rapid Microbial Identification System.
13. Borisenko AV et al. (2009) The front-end logistics of DNA barcoding: challenges and prospects. *Mol Ecol Resour* 9:27–34.
14. Brinkman EK et al. (2014) Easy quantitative assessment of genome editing by sequence trace decomposition. *Nucleic Acids Res* 42:e168.
15. Application note: Using Sanger sequencing to facilitate CRISPR- and TALEN-mediated genome editing workflows.
16. [https://www.atcc.org/STR\\_Database.aspx](https://www.atcc.org/STR_Database.aspx).
17. Product bulletin: SNaPshot Multiplex System for SNP genotyping.
18. Schouten JP et al. (2002) Relative quantification of 40 nucleic acid sequences by multiplex ligation-dependent probe amplification. *Nucleic Acids Res* 30:e57.
19. Sanger sequencing Protocol guide: KRAS variant identification.
20. Sanger sequencing Protocol guide: mRNA vaccine quality control using Sanger sequencing.
21. Papanicolaou-Sengos A, Aldape K. DNA Methylation Profiling: An Emerging Paradigm for Cancer Diagnosis. *Annu Rev Pathol.* 2022 Jan 24;17:295-321. doi: 10.1146/annurev-pathol-042220-022304.

## Ordering information

Product	Quantity	Cat. No.
<b>Instruments</b>		
<b>SeqStudio 8 Flex System</b> Includes: SeqStudio 8 Flex Genetic Analyzer with data collection software, 1-year warranty, 1-day onsite SmartStart orientation, and DNA sequencing and fragment analysis reagent kits for system performance check		A53627
<b>SeqStudio 8 Flex Genetic Analyzer System plus 1-year extended warranty</b> Includes all items from A53627 plus additional 1-year warranty		A53789
<b>SeqStudio 24 Flex System</b> Includes: SeqStudio 24 Flex Genetic Analyzer with data collection software, 1-year warranty, 1-day onsite SmartStart orientation, and DNA sequencing and fragment analysis reagent kits for system performance check		A53630
<b>SeqStudio 24 Flex Genetic Analyzer System plus 1-year extended warranty</b> Includes all items from A53630 plus additional 1-year warranty		A53792
<b>SeqStudio 8 Flex for Pharma Analytics</b> Includes: SAE v2.1 and XE3 Computer		A57829
<b>SeqStudio 24 Flex for Pharma Analytics</b> Includes: SAE v2.1 and XE3 Computer		A57830
<b>SeqStudio™ Genetic Analyzer System with SmartStart orientation</b>		A35644
<b>SeqStudio™ Genetic Analyzer for HID, laptop, with training</b>		A46228
<b>SeqStudio™ Genetic Analyzer QST with SAE</b>		A49988
<b>Reagents</b>		
CLA Identifiler Plus PCR Amplification Kit	200 reactions	A65672
CLA Identifiler Direct PCR Amplification Kit	200 reactions	A65908
CLA GlobalFiler PCR Amplification Kit	200 reactions	A65909
TrueMark MSI Assay	100 reactions	A45295
BigDye Terminator v1.1 Cycle Sequencing Kit	100 reactions	4337450
BigDye Terminator v3.1 Cycle Sequencing Kit	100 reactions	4337455
BigDye Direct Cycle Sequencing Kit	100 reactions	4458687
BigDye XTerminator Purification Kit	100 preps	4376486
ExoSAP-IT PCR Product Cleanup Reagent	100 reactions	78200.200.UL
RecoverAll Total Nucleic Acid Isolation Kit for FFPE	40 reactions	AM1975
MicroSEQ 500 16S rDNA PCR Kit	1 kit	4348228
MicroSEQ 500 16S rDNA Sequencing Kit	1 kit	4346480
EpiJet Bisulfite conversion kit	50 reactions	K1461
SNaPshot Multiplex kit	200 reactions	4323151

 Learn more about CE applications at [thermofisher.com/ce](https://thermofisher.com/ce)

**applied biosystems**

**For Research Use Only. Not for use in diagnostic procedures.** © 2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TaqMan is a registered trademark of Roche Molecular Systems, Inc., used under permission and license. Geneious is a trademark of Biomatters Limited. pGEM is a trademark of Promega Corporation. MLPA is a trademark of MRC-Holland. NUCLEIC- CARD is a trademark of Copan Flock Technologies. Thermo Fisher Scientific's GlobalFiler and Identifiler kits are supplied for forensic and paternity use only in the following countries: Japan, China, and Hong Kong. **EXT6867 0624**