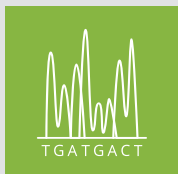


Laboratory Data Management Solutions

Genomics Workflows

Thermo Fisher™ Platform for Science™ Software

Applications provide functionality for many different pieces of the sequencing workflow



Sanger Sequencing



Nucleic Acid Purification



NGS Reagent Management



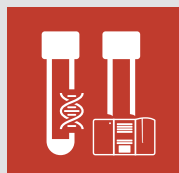
NGS Lab Operations



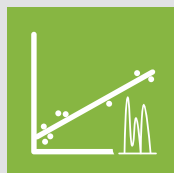
Sequencing Request Portal



Illumina DNA Library Prep



Axiom Genotyping Pre Lab



Sanger Lab QC



NGS Sample Accessioning



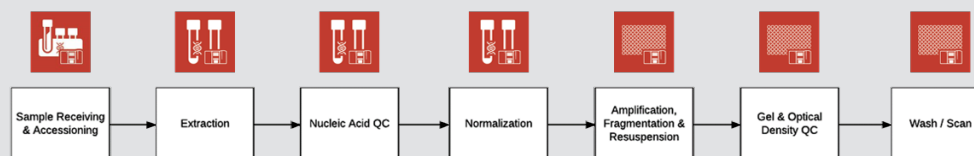
Reverse Transcription PCR

Axiom

Instrument supported



Axiom workflow



For use with

- 96 and 384 Well Plate Assay Formats
- Nimbus™ Target Prep Instrument
- Off the Shelf or Custom Arrays
- Gene Titan™ Multi-Channel Instrument


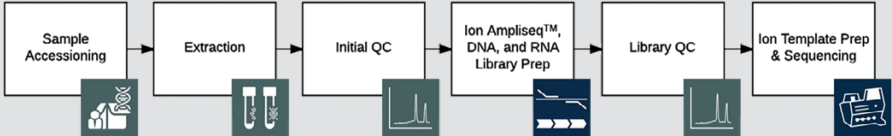
Key benefits


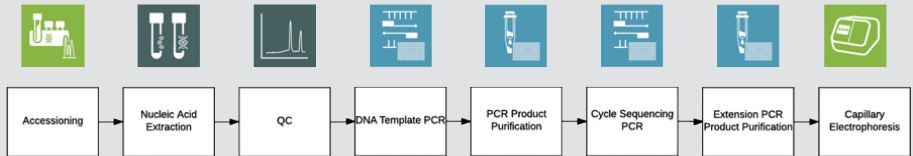
- Manage workflows and track all metadata generated.
- Easily navigate the system and track status of samples with dashboards.
- Automate capture of quantitation values after PicoGreen™ assay.
- Compare results captured from different instruments in the lab.
- Gain visibility into processing bottlenecks, sample failures, and reagent performance.
- Track projects, samples, users, instruments, consumables, and shipping requests.

Laboratory Data Management Solutions

Genomics Workflows

Thermo Fisher™ Platform for Science™ Software

Ion Torrent	
Instrument supported 	Ion Torrent workflows 
For use with <ul style="list-style-type: none"> • Ion Chef™ System • Ion OneTouch™ System • Ion GeneStudio™ Systems • Ion PGM™ System • Ion Proton System™ 	Key benefits <ul style="list-style-type: none"> • All data lives in the same platform architecture, streamlining data review. • Metadata generated from the creation of Ion libraries is captured centrally, including reagent kit, lots consumed, and equipment and instruments utilized. • Easily add capabilities over time through apps, solution sets, and configuration – with no custom code. • Process extracted DNA, RNA or FFPE samples through library prep and pooling with the Ion Chef System. • Execute sample enrichment and chip loading with the Ion OneTouch 2 and Ion Chef systems. Scientists can track reagents lots and plan runs for the Ion PGM, Ion Proton™ and Ion S5 sequencing instruments.

Capillary Electrophoresis and Sanger Sequencing	
Instrument supported 	Capillary Electrophoresis and Sanger Sequencing 
For use with <ul style="list-style-type: none"> • SeqStudio™ Genetic Analyzer • Applied Biosystems 3500 and 3700 Series Genetic Analyzer 	Key benefits <ul style="list-style-type: none"> • Provides scientists with a dashboard to manage samples ready to be run on the SeqStudio Genetic Analyzer. • Track run metadata such as instrument, SeqStudio cartridge, and application type • Automated generation of Plate Template File for import into the SeqStudio Plate Manager. • Connect other instruments and software systems by utilizing the OData Application Programming Interface.

Find out more at coreinformatics.com

For Research Use Only. Not for use in diagnostic procedures. © 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. **PSB80122-EN0618**

ThermoFisher
SCIENTIFIC