

Thermo Scientific Momentum—laboratory automation software that seamlessly connects to your software ecosystem

An evolving NEED...

- Labs need deep dynamic digital connectivity to software ecosystems
- Harmonized data must be stored with metadata (contextualized)

An evolved SOLUTION!

- Thermo Scientific™ Momentum™ scheduler dynamically controls automated workflows
- Momentum's Unite module built-in data tools connect to your lab's software ecosystem

Momentum Data Tools

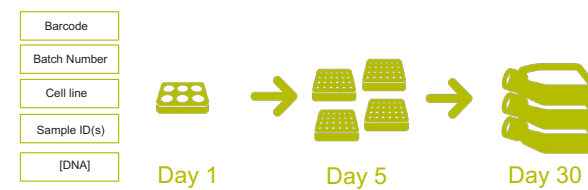
UNITE Module



- Real-time bi-directional communication
- Exchange data using Transform tools
- Compatible with your data interfaces:



Persistent Data

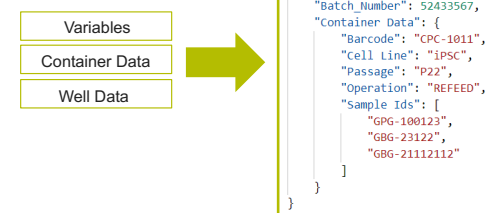


- Enables user defined plate/well data
- Finds & uses specific containers (machine logic)
- Shares containers across experiments

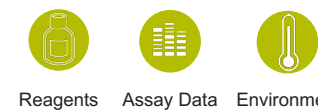


Data Builder

- Builds JSON or delimited formatted data

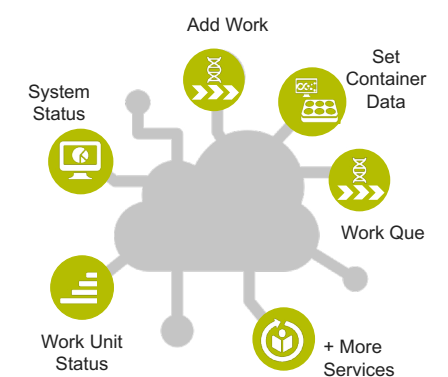


Metadata Capture



- Adds context to data for deeper insight & analysis

Momentum RESTful API



- Automates entire process
- Enables remote task & data management

Case Study—Momentum Unite & DataBuilder actions with gathered metadata: Thermo Fisher Scientific Amplitude™ COVID-19 Module 2 Testing Solution

Connected Amplitude Systems

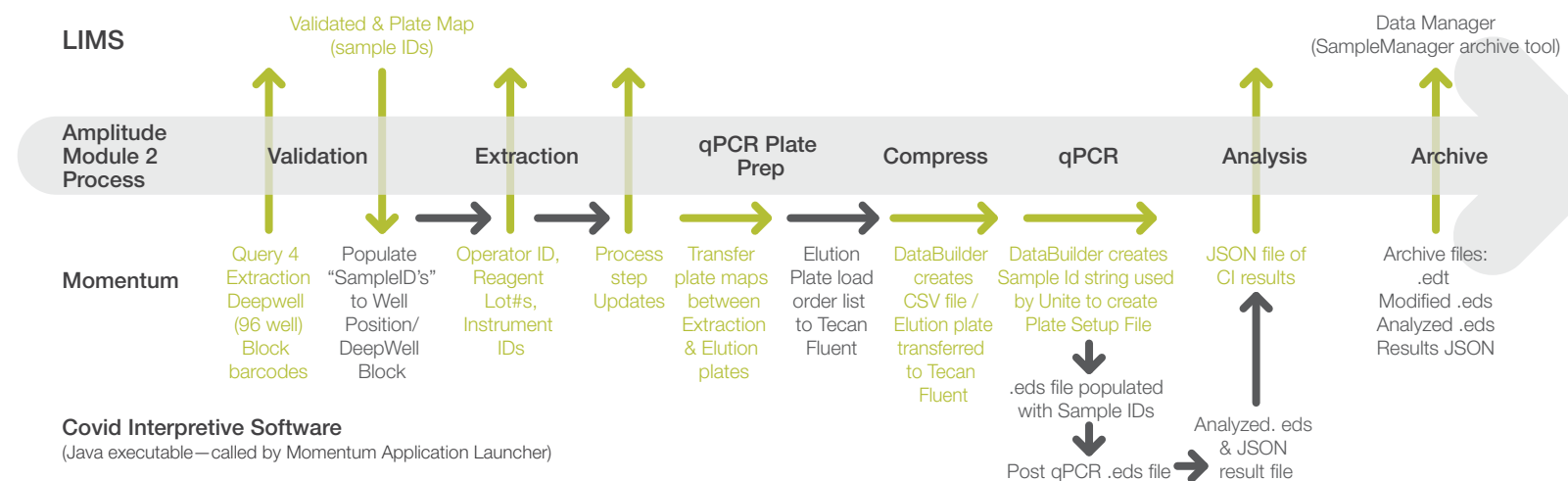


Module 1—Array samples

Input: 378 tubes
Output: Four 96 well DWP

Module 2—Extraction to result

Input: Four 96 well DWP
Output: Results from 384 qPCR plate



An evolved PRODUCT

Dynamic data driven workflows powered by digital connectivity

- Real-time bi-directional communication between Momentum and your lab's software ecosystem
- Actionable JSON or delimited formatted data
- Metadata available for use by ML/AI tools

Learn more at thermofisher.com/momentum

thermo scientific