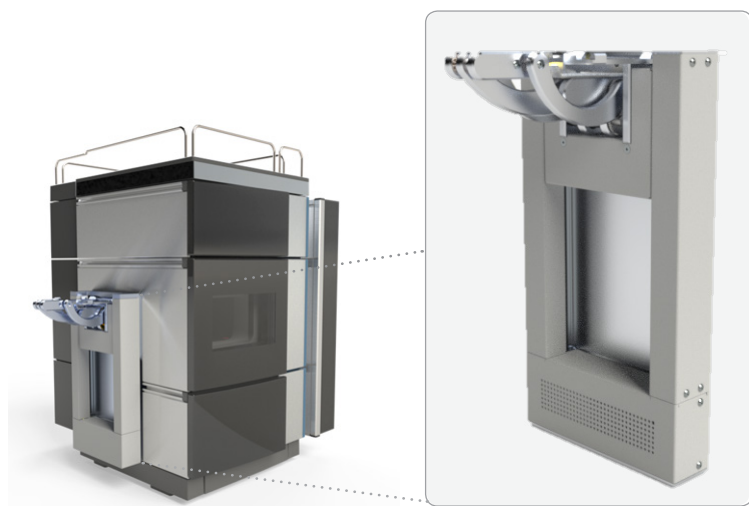


End-to-end analytical workflow solution—Thermo Scientific Vanquish UHPLC Loader

Introducing Thermo Fisher Scientific's new loader technology



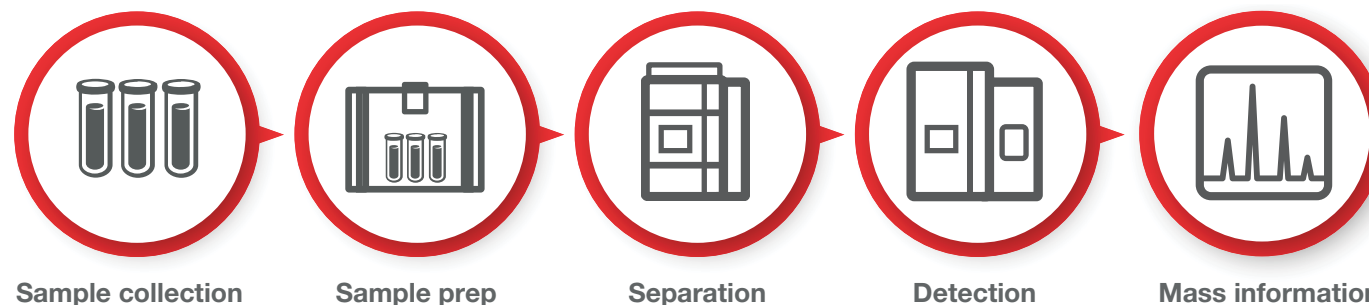
The field of Analytical Science is evolving with advancements in instrumentation and software. Accelerated by the recent pandemic, strategic discussions on how to embrace automation and digital transformation have consumed the industry and highlighted the need for diverse and connected solutions. It gave rise to new improvements in the field to ensure business continuity without compromising on productivity and data reproducibility.

Take your analytical science one step further with the Thermo Scientific™ Vanquish™ UHPLC Loader. Automate your entire workflow creating an end-to-end solution with little to no human intervention. Achieve seamless integration between upstream sample preparation and downstream analysis.

- All the steps in the Analytical workflow can be integrated using Thermo Fisher Scientific's Laboratory Automation products
- Prepped Samples can be transferred to Vanquish UHPLC by using one of our robotic movers and loaded through the Vanquish Loader
- A complete workflow automation solution powered by Thermo Scientific™ Momentum™ Workflow Scheduling Software along with Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS)

Why is this necessary?

CURRENT STATE



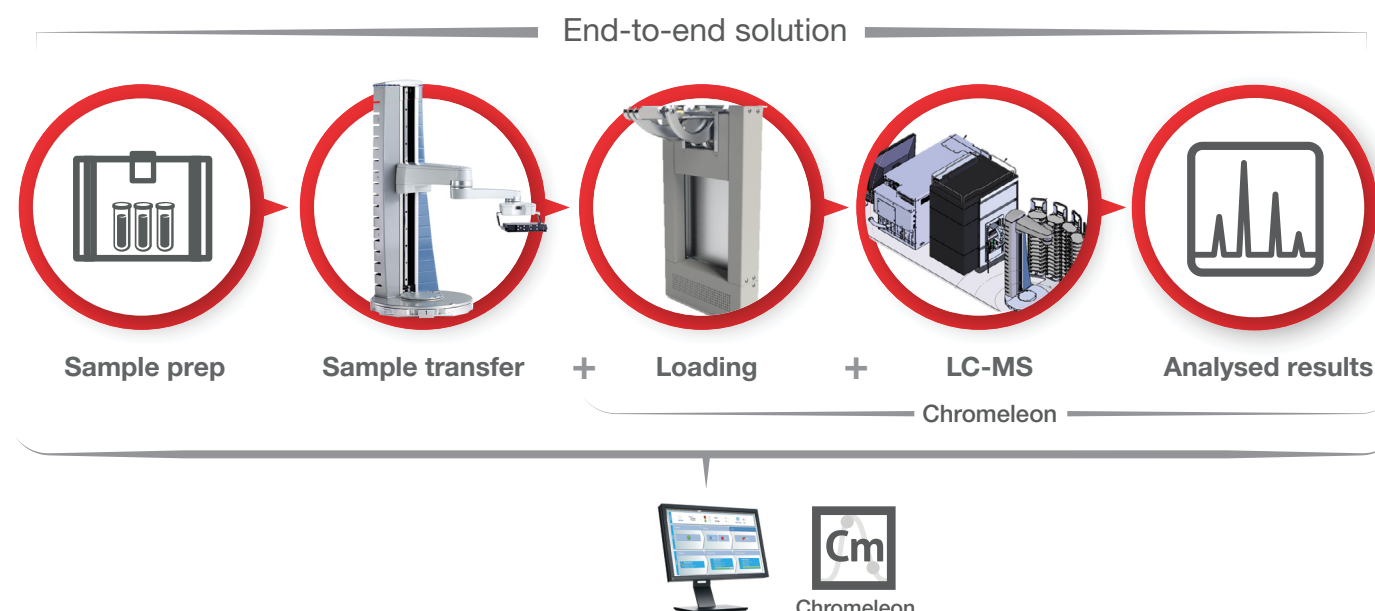
- Eliminate the need for human intervention
 - Stop worrying about time-consuming, labor intensive and error prone experimental workflows
- 
- Your downstream analytical workflow need not be disconnected from the upstream sample prep workflow

Technology benefits

-  Productivity
-  Reproducibility
-  Data tracking
-  **Shift the focus to data—not sample handling**
One step closer to an end-to-end workflow and a fully connected laboratory





How does workflow automation solve the problem?

FUTURE STATE



Thermo Scientific Momentum workflow scheduling software

Applications

-  Monoclonal antibodies
-  Biologics
-  Drug screening (small molecule)
-  Environmental

Learn more at thermofisher.com/labautomation