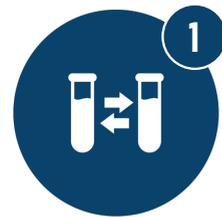


How the AQUIOS CL Load & Go flow cytometer helps to address the **8 WASTES OF LEAN PROCESSING** in a clinical flow cytometry laboratory

DEFECTS caused by the addition of a wrong reagent to your sample, sample mix-up during preparation or transcription errors can be largely eliminated from your process as the AQUIOS CL automates and traces the entire process from specimen input to sample ready to analyze.



OVERPRODUCTION such as time-consuming training of your staff to ensure process standardization will be mitigated as the AQUIOS CL simultaneously prepares and analyzes samples leveraging pre-programmed and standardized methods for Immunophenotyping and stem cell enumeration



WAITING times in quality control can be eliminated with IQAP. The AQUIOS CL's continuous loading and parallel processing, combined with IQAP's streamlined workflow and peer review capabilities, optimize time-to-result and ensure efficient outcomes for patients.



NON-UTILIZED TALENT is a real threat as skilled techs are hard to find. Full automation of low complexity assays including a comprehensive audit trail will allow your highly qualified staff to concentrate on more value adding tasks like the development of new assays or data analysis.



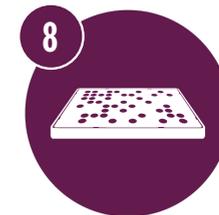
TRANSPORTATION of liquid reagents can be reduced by the AQUIOS CL capacity to store reagents and support inventory management.



INVENTORY management can be time consuming and complex. AQUIOS CL uses barcoded reagents and provides a full reagent audit trail to help address this.



MOTION like transcription of data, manual pipetting, opening and closing containers and refrigerators can be significantly reduced via the AQUIOS CL capacity to store reagents, pierce reagent caps as well as the option to integrate with your LIS.



EXTRA-PROCESSING like repeated testing performed to mitigate process variability or defects can be minimized as automation of the full sample preparation process will help standardization and reduce the risk of manual error-prone steps in your workflow.