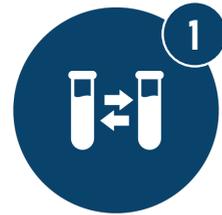


# How CellMek SPS Sample Preparation System helps to address the **8 WASTES OF LEAN PROCESSING**

**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 CellMek SPS **automates and traces the entire process** from specimen input to sample ready to analyze.



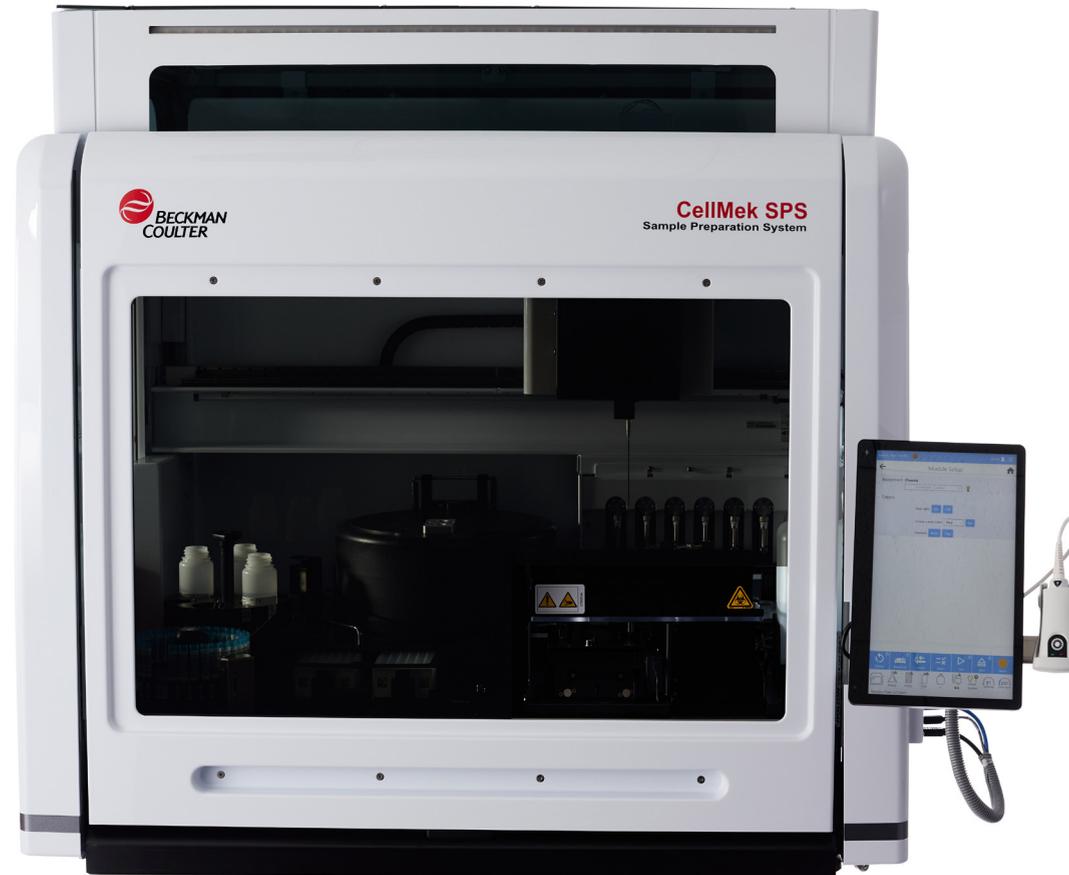
**OVERPRODUCTION** like a surplus of pre-mixed antibody cocktails or time consuming training of your staff to ensure process standardization will be mitigated as the CellMek SPS simultaneously prepares a **wide variety of samples** leveraging **pre-programmed and standardized methods** for your LDT workflows.



**WAITING** based on batch processing bottlenecks, centrifugation and incubation times can be eliminated by the CellMek SPS's **continuous loading and unloading capabilities, onboard cell wash** and **parallel processing** that empower the optimization of individual time to result for your patients.



**NON-UTILIZED TALENT** is a real threat as skilled techs are hard to find. **Fully automated sample preparation** including a **comprehensive audit trail** will allow your staff to concentrate on more value adding tasks like the development of new assays or data analysis.



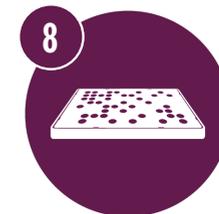
**TRANSPORTATION** of multiple refrigerated deliveries of liquid reagents can be addressed by the system's **ability to process DURACartridges**, an innovative, new format of customized dry reagent panels.



**INVENTORY** management can be time consuming and complex. CellMek SPS provides **on-board refrigerated reagent storage** as well as 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 CellMek SPS system's capacity to **store reagents, pierce reagent caps, wash samples and specimens**, and 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.