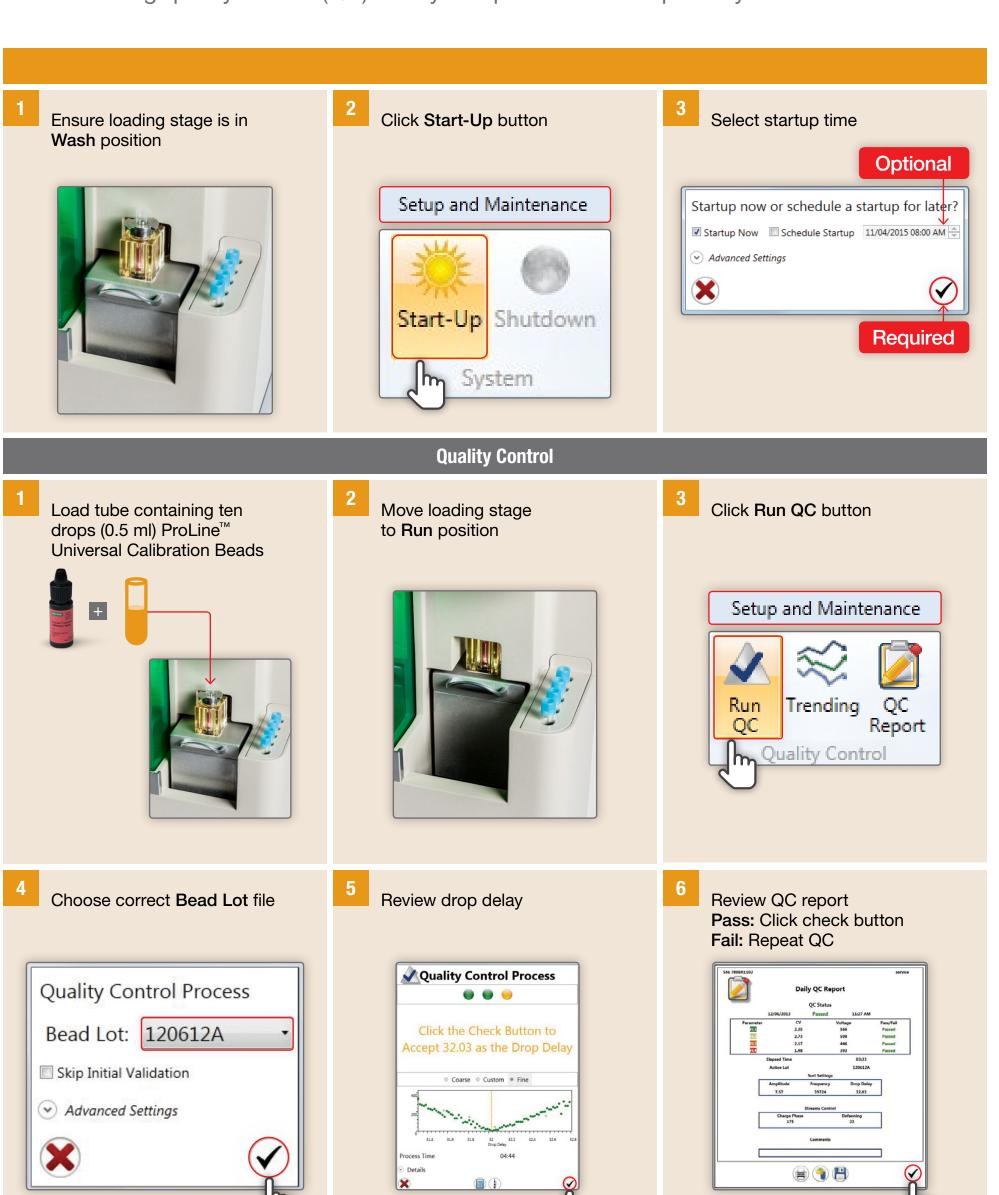
Daily Startup and Quality Control

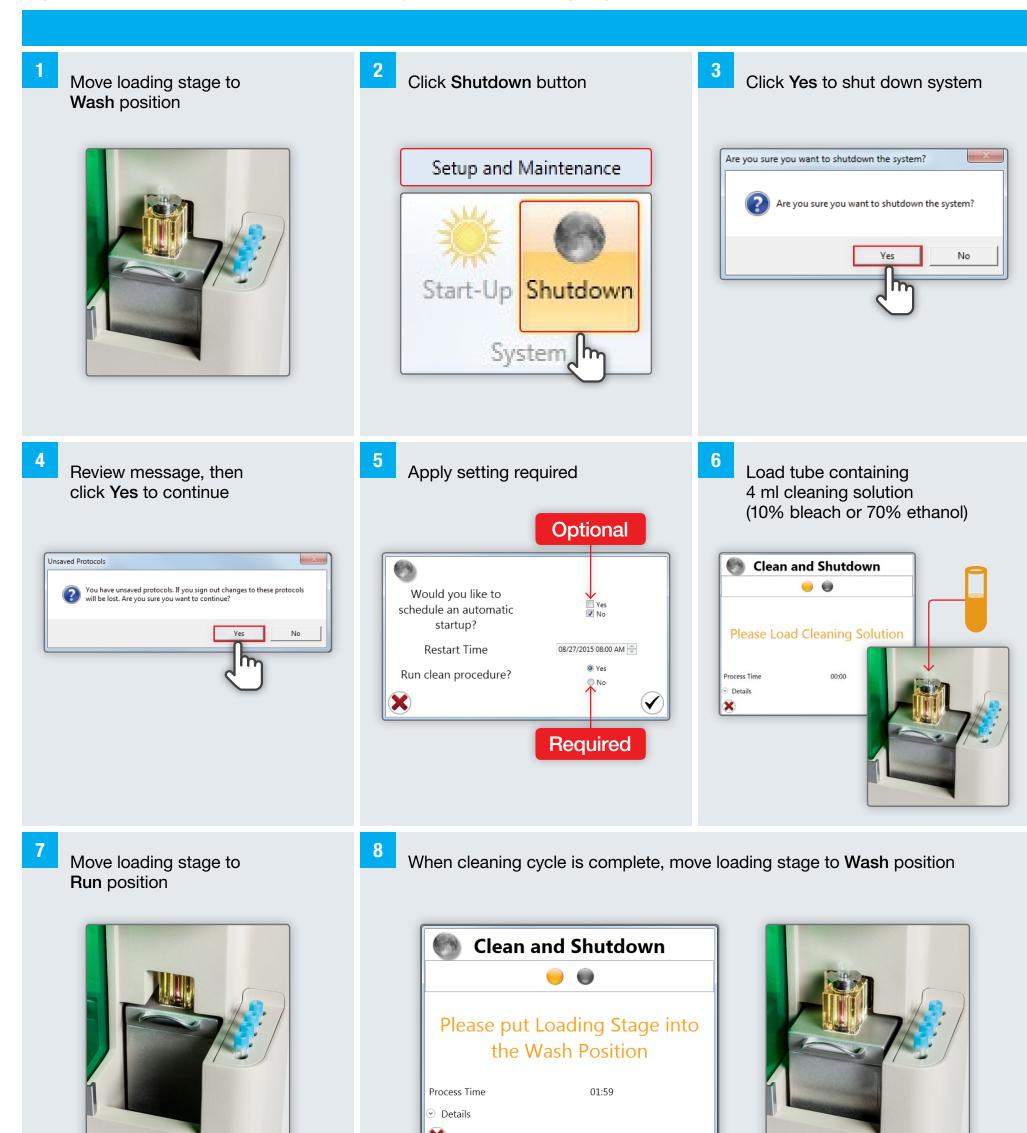
On a daily basis, the S3e Cell Sorter requires approximately 20 minutes of warm-up time before running quality control (QC) or any samples. To warm up the system:



Daily Shutdown Procedure

These steps should be performed after the last sort of the day:

Important! Prior to shutdown, make sure that you have saved all open protocols.



High Pressure Clean

For a more stringent cleaning process, these steps should be performed after the last sort of the day:

