



Cell Culture Automation in Biopharma

Cell Line Optimization and
Monoclonal Antibody Production

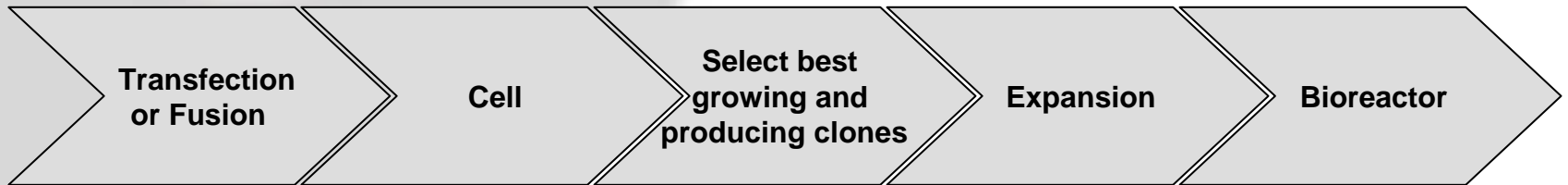
Dr. Christoph Krüll
Product Manager Automated Solutions
Beckman Coulter Europe

Agenda

- Cell Line Optimization and Expansion
- Monoclonal Antibody Production
- Software Control

Biopharmaceuticals

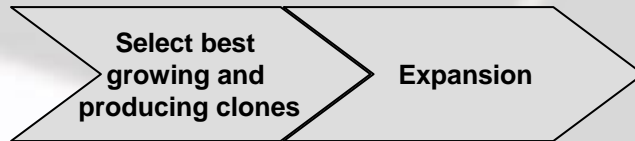
Basic Process



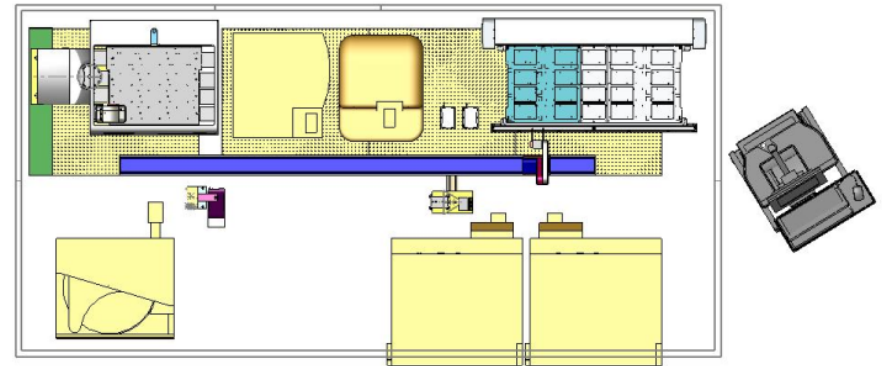
Monoclonal Antibodies

Cell Line Optimization and Expansion

Cell Line Optimization and Expansion System

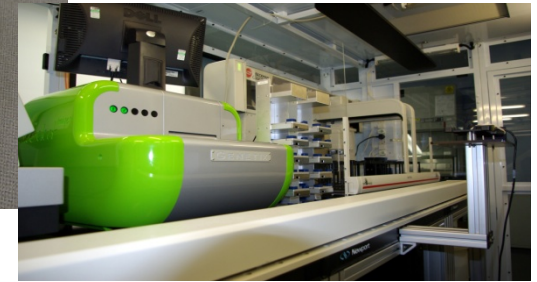
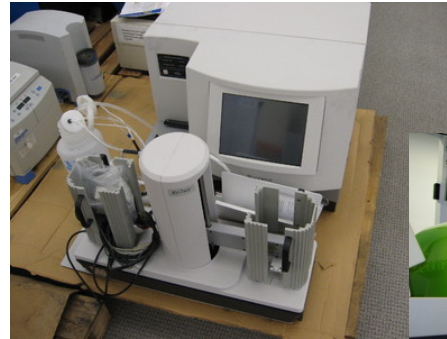


- Requirements :
 - On-line incubation
 - Cell Line passage
 - Media exchange
 - Cell imaging for confluency and monoclonality
 - Specific protein analysis (titre)
 - Hitpicking and Expansion (96 / 24 / 6 / 1 well plates and Autoflasks)
 - Cell counting/viability assay
 - Multiple experiments occur simultaneously
 - 2 months per batch
 - 5 batches in parallel

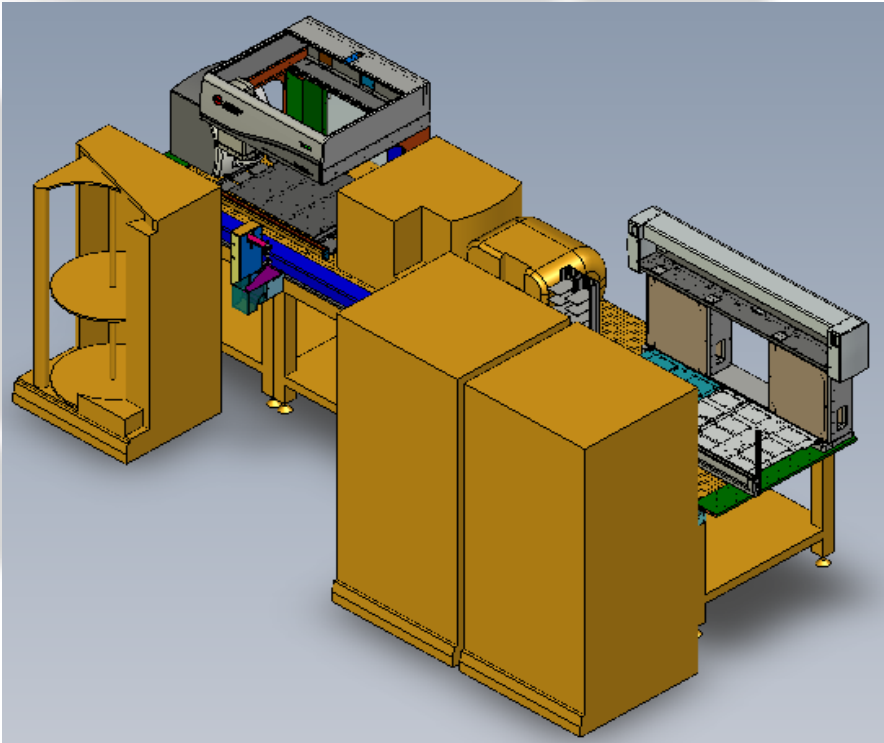


Cell Line Optimization and Expansion: Decision Making on the Fly!

- **96 well to 24 well Expansion**
 - combined M384 and Clone Select data used to generate pick list.
- **24 well to 6 well Expansion**
 - CloneSelect imager data generates pick list
- **Amplification in 96 well plates**
 - Vi-CELL[®] XR data used to plate cells back into 96 well plates at correct seeding density

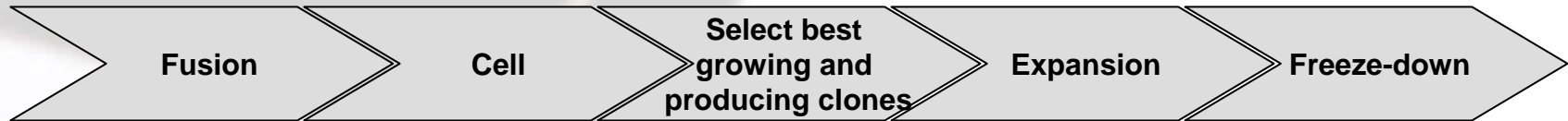


Cell Line Optimization and Expansion System



Monoclonal Antibody Production

Monoclonal antibody production – The Automation Mission



- Mission : Automate all non-animal tasks within the MAP
 - Confirmatory ELISA
 - Fusion
 - Clonal Expansion
 - Media Collection (Mab) and cell freeze-down
- Solution: Partial automation with separate workstations
 1. ELISA workstation
 2. Fusion workstation
 3. Expansion / Freezing

Fusion Process

Spleen Harvest and Splenocyte Preparation



Viable Cell Count (Splenocytes and SP-2)



Mix Splenocytes and SP-2



Wash cell suspension



Fusion with PEG at 37°C



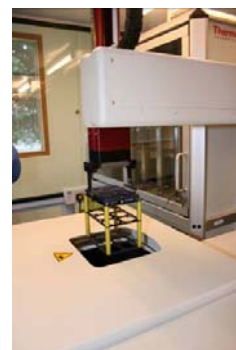
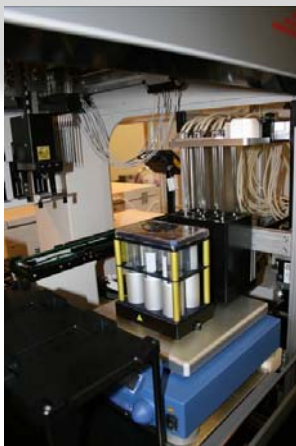
Plate fused cells at limiting dilution or in Semi-Solid Medium (HAT selection)

Mab Production: Fusion

- Biomek® NX^P Span-8 with Gripper
- Vi-CELL XR Cell Viability Analyzer to measure viable cell count
 - splenocytes and SP-2 cells are mixed in a predefined ratio
- IKA shaker (KS130)
 - thermal exchange unit for warming tubes
- Six Position Falcon Tube Rack
 - open tube but has a micro-titre plate lid
- Bulk Dispense / Wash System
- Automated Centrifuge
- Cytomat 6001 and Cytomat Linear 8 Hotel (for tube racks)
- Sterile enclosure (Bigneat)

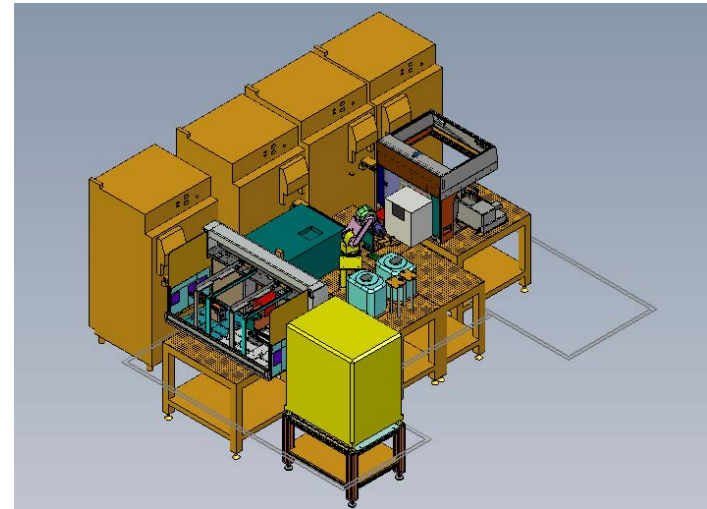


Mab Production: Fusion



Mab Production: Expansion / Freezing

- Expansion from 96 well through 24 / 48 well plates to 6 well plates / Autoflasks
- Hitpicking based upon ELISA data
- Media exchange
- Harvesting of cells from 6 well plates / Autoflasks
- Freezedown of cells and collection of Ab-rich media
- Biomek NX^P Span-8 with Gripper
 - expand cells through 96, 24 and 6 well plates
 - media sampling for confirmatory ELISA; hitpicking
- Motoman HP3JC (waist-based)
 - Sigma 6K15 Centrifuge
 - FluidX Xsd-96 screwcap decapper
 - Grant Asymptote EF600 controlled rate freezers (x2)
- Biomek FX^P Dual Hybrid
 - Span-8 bulk transfer system
 - Simultaneous transfers using four probes across two pieces of labware
- Cytomat 24 Incubators



Mab Production: ZONE 1 - Expansion

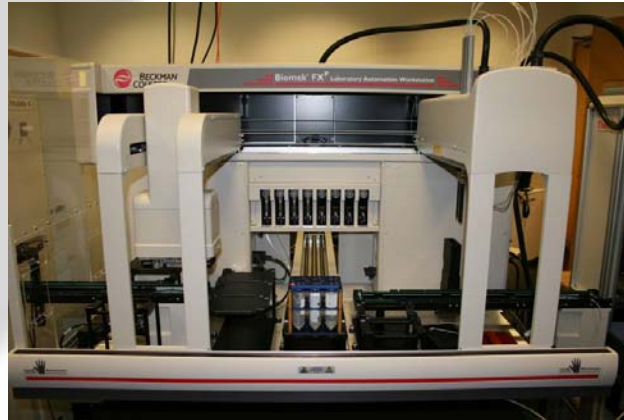
- Biomek NX^P Span-8 with Gripper
 - expand cells through 96, 24 and 6 well plates
 - media sampling for confirmatory ELISA; hitpicking
- Integrated Multidrop Combi
 - (custom Fillit software) for large volume media additions (24 and 6 well plates)
- Cytomat 24C incubators (x2)



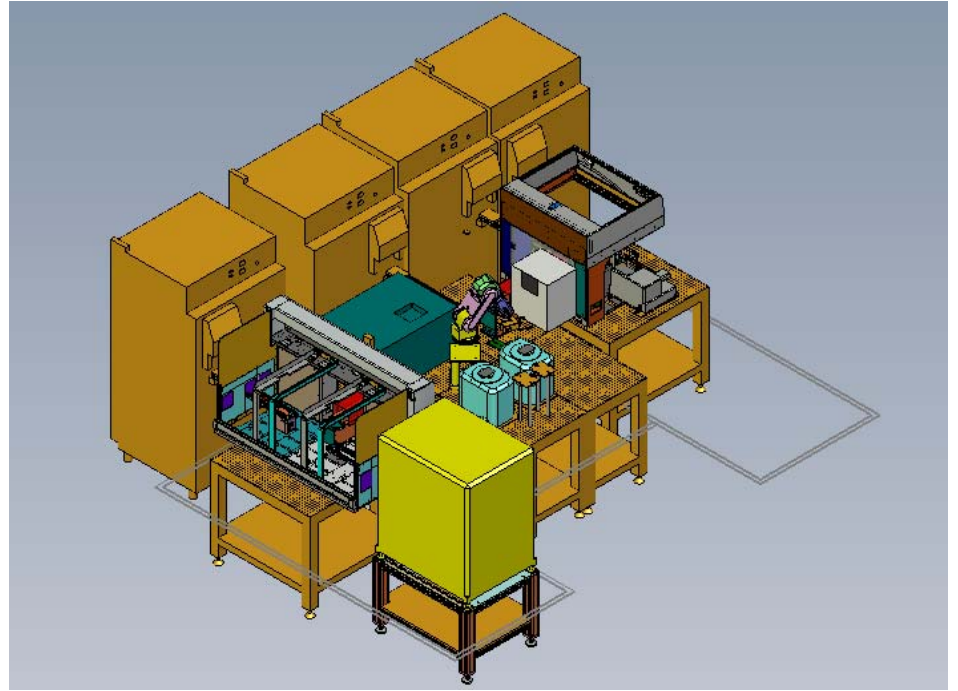
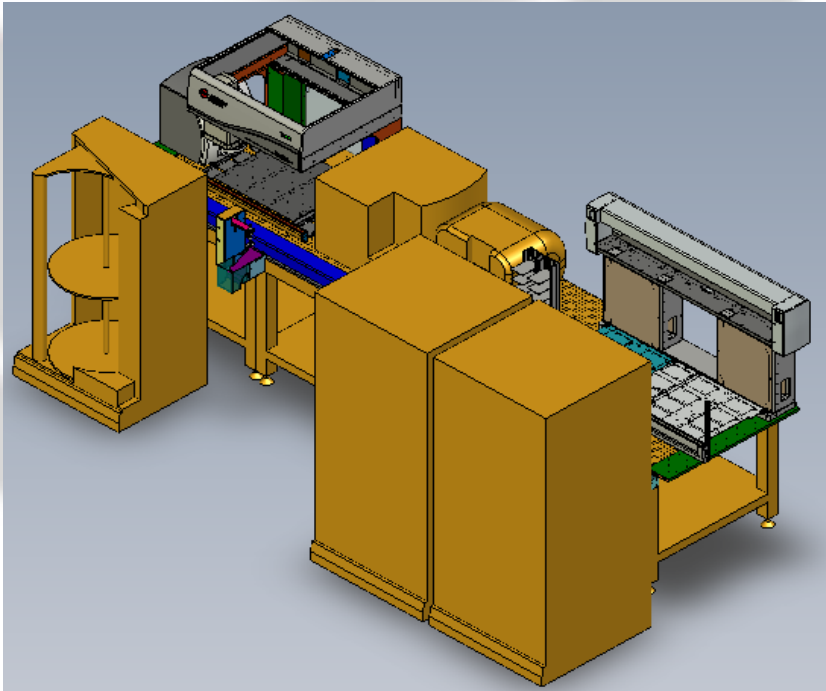
Mab Production:

ZONE 2 – Freezing & Media Harvest

- Biomek FX^P Dual Hybrid
 - Span-8 bulk transfer system
 - Transfers from 6 well plates to 50 ml Falcon tubes
 - Cell transfer to Nunc Cryovials
- Motoman HP3JC (waist-based)
 - Sigma 6K15 Centrifuge
 - FluidX Xsd-96 screwcap decapper
 - controlled rate freezers (x2)
- Cytomat 24 Incubators

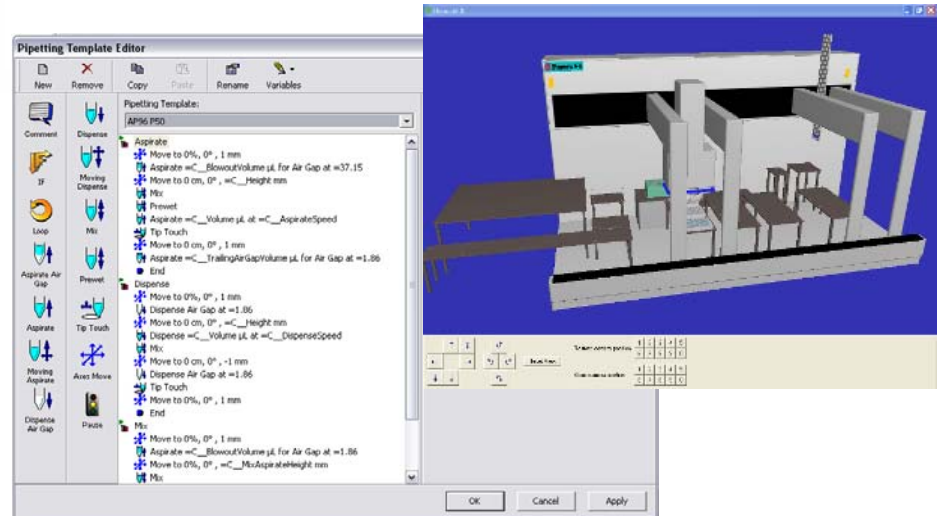


Software Control



Software Control

- Biomek Software
 - Common to Biomek FX^P and Biomek NX^P
 - Pipetting Template Editor controls pipetting operations (e.g. harvesting semi-adherent cells such as Hybridomas)
- SAMI[®] Workstation EX Scheduling Software
 - Optimising Scheduler interleaves complex processes
 - Draw the flow of plates through the process.
 - Configure liquid transfers (e.g. hitpicking), incubations (e.g. trypsinisation), device actions (e.g. multidrop media additions)
 - dynamic reschedule feature for data-driven liquid transfers (CloneSelect, Vi-CELL[®] XR data)

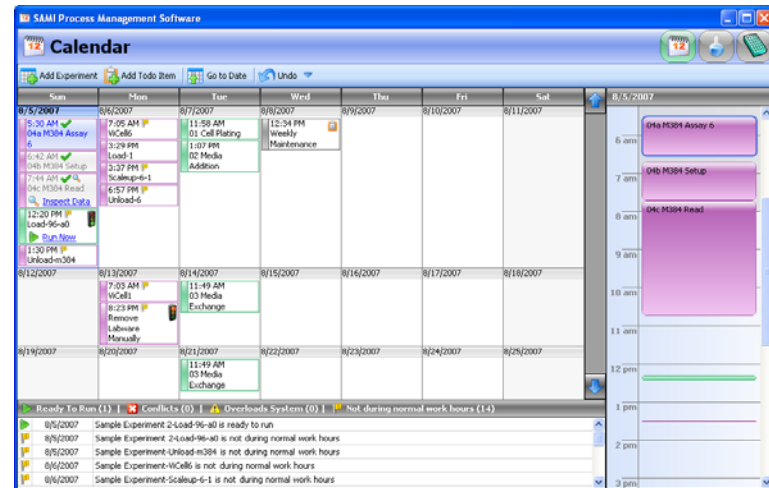
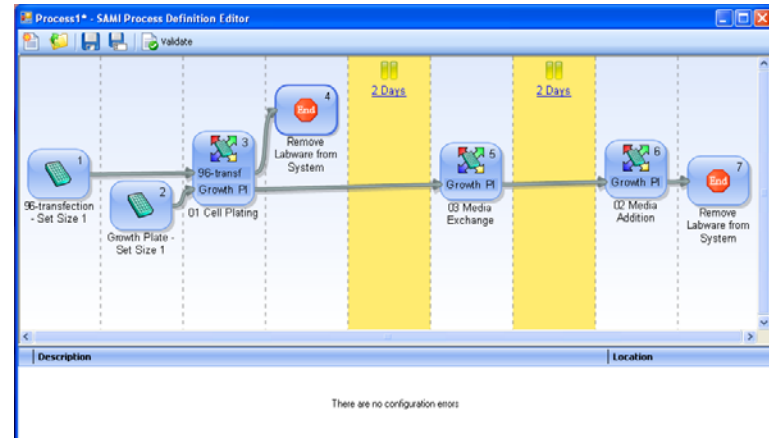


Process Management? - The Problem

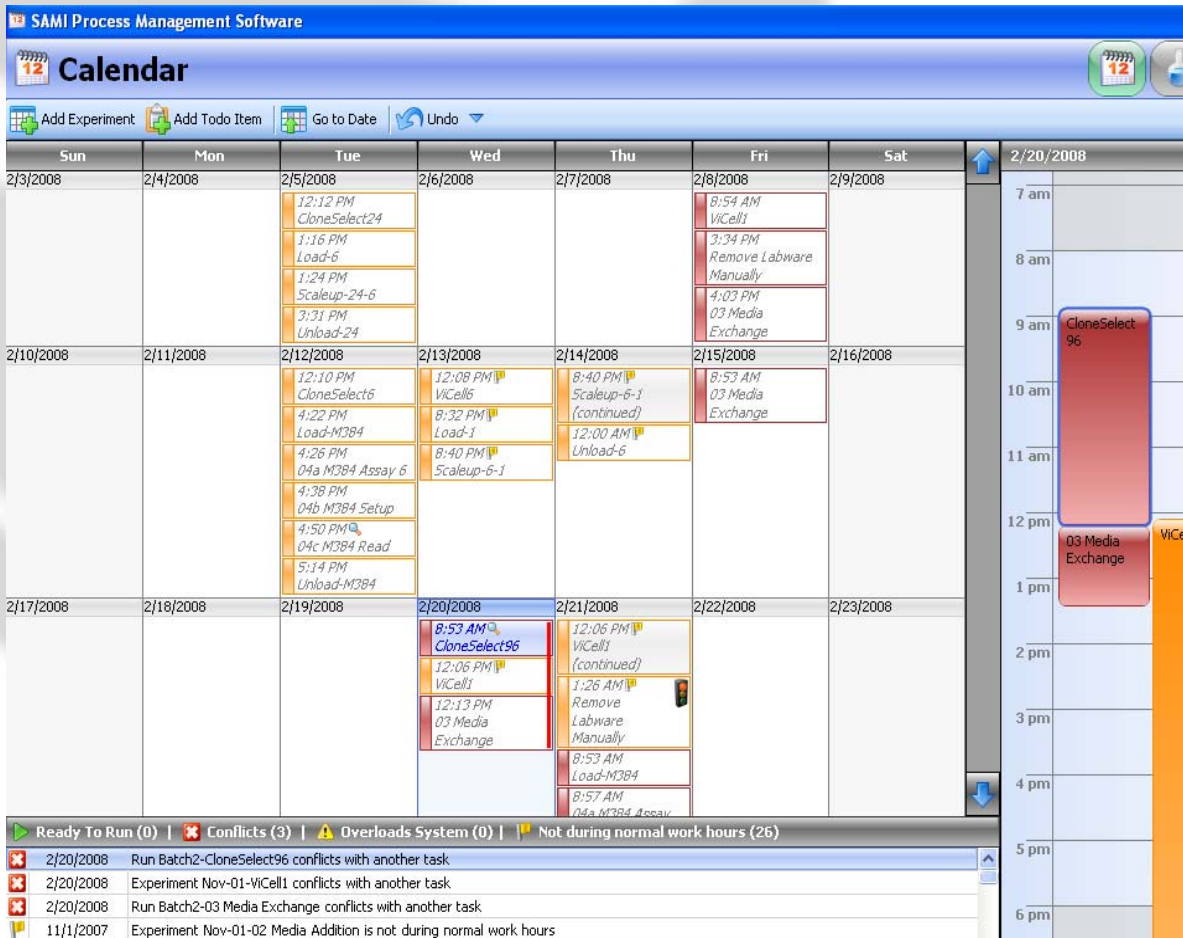
- How do I connect multiple SAMI EX methods?
- What happens to the data for these plates?
- How do I keep track of 4 batches X 6 methods per batch?
- Where are all my plates!?

SAMI[®] Process Management Software

- Process Editor allows to define processes that take weeks to months to complete
- Process Management allows multiple processes to move forward independently
- Calendar view for process tasks with conflict warning / resolving options
- Integrated Database allows data flow from process to process (Data Acquisition & Reporting Tool)
- SAMI PM will track plate location and data throughout the process



The Process Management Tool



- Generates conflict warnings:
 - Conflict with another task
 - Task not in normal working hours or weekends
 - System capacity (labware)
- Easy conflict resolution
 - Shift single tasks
 - Shift experiment
- Unattended start functionality for tasks occurring outside normal working hours

Summary

- Novel hardware solutions for automating complex cell culture automation
- First bespoke system for cell fusion from splenocyte preparation to Hybridoma
- First system for combined expansion and controlled rate freezing of cell lines
- SAMI Process Management is a unique software package allowing different cell culture processes and experiments to move forward independently
- DART enables data flow from process to process

Beckman Coulter: Your Partner in Cellular Analysis

