

Success Story

Leading Biopharma Uses Linkit[®] AX to Increase the Efficiency of Their Aliquoting Process

A Leading Biopharma was looking for a way to speed up their CAR-T media preparation. The fully manual and time intense media preparation process was not set up for GMP manufacturing. After 3 years of collaboration, Sartorius developed a semi-automated media blending and aliquoting solution – the Linkit[®] AX Aliquoting Solution – that has been implemented and is now part of a GMP CAR-T manufacturing process.

Customer Challenge

- Because the system was fully manual, it was a time intensive process for the customer
- Due to the amount of handling required, the risk of error and inaccuracy was high
- Unknown extractable profile
- There was no strategy to scale up the media preparation from lab scale to a GMP environment

Provided Solution

Standardized semi-automated solution that can be used to reduce intense time effort for media prep, minimize handling errors and provide high aliquoting accuracy – while allowing for sustainable growth for increasing demand.

Case Profile

Company Type:
Large Biopharma

Related Therapy:
CAR-T or T cells

Process Steps:
Cell Culture Media and Buffers



Before

- 2 days of media preparation for 50 L batch per media
- 2+ employees



After

- 4 h for media preparation and aliquoting for a 50 L batch per media
- <2 employees

Over 10× faster

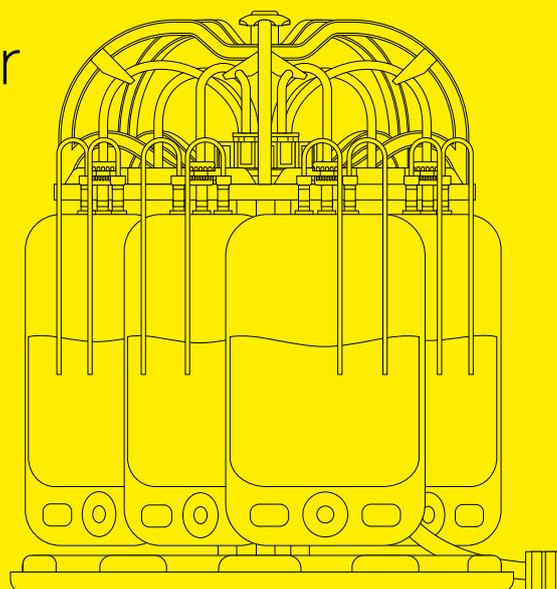
4 hours instead of 2 days.

Less errors

caused by operator handling.

Faster turnaround

of media preparation and aliquoting.



 **Contact a Sartorius Specialist or Visit**
www.sartorius.com/linkit