

Fraunhofer IZI Services for Antibodies

Fraunhofer IZI provides GMP Antibody Manufacturing Services with Sartorius Stedim Biotech single-use technologies



Success Story

Fraunhofer Institute for Cell Therapy and Immunology Department of Therapy Validation Leipzig, Germany

In recent years, the increasing number of therapeutic monoclonal antibody candidates under preclinical and clinical development require new flexible, efficient and economic opportunities for GMP production of therapeutic antibody candidates. Small-scale batch production of test samples for late preclinical GLP animal studies or for phase-1 and phase-2 clinical studies is often not appropriate for large-scale manufacturing facilities in the industry.

The Fraunhofer IZI, founded in 2005, investigates and develops solutions to such specific problems at the interfaces of medicine, life sciences and engineering.

State-of-the-art equipment

- Isolator | Incubators for inoculation from the master cell bank
- Sartorius Biostat® RM20 & RM50 for seeding
- Sartorius Biostat® STRplus 200 L for production
- Sartorius FlexAct® Controlling systems for media preparation, cell harvest and virus removal
- GE Healthcare ÄKTApilot chromatography system for protein purification
- Sartorius Sartoflow® Advanced for ultra and diafiltration
- Colanar semiautomatic filling machine for product formulation

Process development

- Transfer of the Research Cell Bank into a GMP Master Cell Bank with strategic partners for characterization
- Customized design for all parts of the manufacturing process
- Establishment of a labscale process and process optimization e.g. expression media screening, fed-batch development
- Identification of critical process parameters for cell growth and protein production
- Establishment and validation of product QC

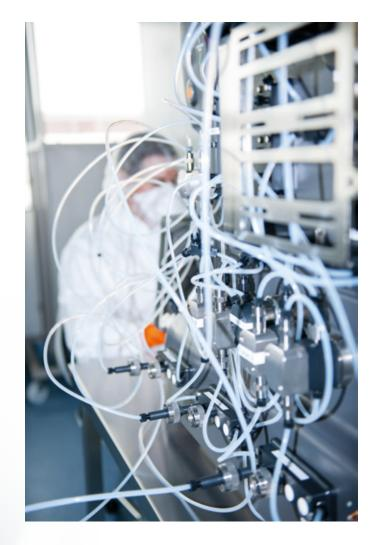


Upstream

- Transfer to production scale up to 200 L of total cultivation volume
- Based on single-use bag assemblies
- Equipment for aseptic liquid transfer e. g. Biowelder® TC, Biosealer® and Biosafe® Transfer Ports

Downstream

- Cell removal via depth filtration
- Capture, purification and polishing with three step affinity and ion exchange chromatography
- virus inactivation and removal
- Setting the final buffer for the product
- End up with a minimal volume of 200 mL



Aseptic filling

- Final product as bulk ware or in glass vials
- Filling in clean-room class A in class B environment
- Available glass vials in different sizes

Single-use Capabilities

- High flexibility
- Easy change-over to different products
- Fast implementation of technology changes
- Customized production
- Ideal for preclinical and early clinical trials
- Ready-to-use GMP compliant products by integrated sample filling



Services

- Transfer of your promising biopharmaceutical candidate from research to clinical development
- Manufacturing of Master and Working Cell Banks
- Process development from Upstream to Downstream
- Design of your user-specific process by single-use materials
- GMP-compliant production of e.g. human monoclonal antibodies in 200 L-scale









