



CLINICAL TRIAL AND COMMERCIAL BIOLOGICS MATERIAL

EUROGENTEC
BIOLOGICS

- GMP Recombinant proteins
 - GMP Plasmid DNA
- GMP Protein conjugation
 - US FDA inspected

Eurogentec is a GMP accredited manufacturer of parenteral biologics. We produce clinical trial material for all major markets according to FDA and EMA requirements. As experts in the manufacturing of biologics from bacterial and yeast sources, we offer significant know-how in high cell density fermentation, purification by refolding of inclusion bodies, isolation of periplasmic or extracellular secreted proteins, purification of intracellular soluble proteins as well as production of plasmid DNA.

Comprehensive GMP Experience

- ▣ GMP accredited since 1994
- ▣ FDA inspected 2011, 2013, 2014
- ▣ >100 custom GMP processes developed, > 500 GMP batches released
- ▣ Manufacturing to FDA 21 CFR Part 210 & 211, EU 2003/94/EC and Eudralex Vol 4

Experience in All Clinical Phases

- ▣ Manufacturing for Phase I, II, III and commercial
- ▣ Process development: USP, DSP, QC, preformulation
- ▣ QC qualification and validation plan
- ▣ Process characterization
- ▣ Process validation
- ▣ In-house QP release of DS and DP

Product Experience

- ▣ Recombinant proteins (eg enzymes, cytokines, antibody fragments, fusion proteins)
- ▣ Plasmid DNA, API and starting material
- ▣ PEGylated proteins
- ▣ Peptide-protein conjugates

Unique Manufacturing Platforms

- ▣ Strain specific high density fed-batch fermentation methods
- ▣ Continuous / non-batch based purification methods, highly scalable
- ▣ Off-patent *Pichia pastoris* systems
- ▣ Low O-glycosylation fermentation conditions for *Pichia pastoris*
- ▣ Plasmid manufacturing to 200 g scale

Comprehensive Service Offering

- ▣ Cell line development (*E.coli* and *P.pastoris*)
- ▣ GMP Cell banking
- ▣ USP, DSP and QC development
- ▣ Stress stability studies
- ▣ Preformulation development
- ▣ API Manufacturing
- ▣ Fill & Finish, Packaging, Labeling
- ▣ Tox batch manufacturing
- ▣ GMP Clinical trial manufacturing
- ▣ Process characterization & validation
- ▣ GMP Commercial manufacturing
- ▣ ICH Stability studies on drug substance and drug product





Host System Experience

Manufacturing with all the important microbial strains

- *E. coli*
- *P. pastoris*
- *H. polymorpha*
- *S. cerevisiae*
- Biosafety level 2 micro-organisms that are non-sporulating

Technical Expertise

- Fermentation development using a Design of Experiment approach with parallel 4x5L fermentors
- Purification development by parallel screening of resins for multiple process performance properties
- In-house development of QC tests, IPC & release tests incl cell based potency assays
- Formulation by Design of Experiment based Stress stability studies
- Scale-down model validation
- Statistical approach to process analysis and specification setting



Comprehensive Expression Experience

- Refolding of inclusion bodies
- Periplasmic expression
- Soluble cytoplasmic expression
- Extracellular secretion



Multi-Product Manufacturing Facility

- 3 GMP Fermentation suites (up to 500L)
- 2 GMP Purification suites
- 1 GMP Sterile Filtration suite
- FDA inspected 2011, 2013, 2014

FDA INSPECTED AND APPROVED FOR COMMERCIAL MANUFACTURING

→ FOR FURTHER INFORMATION, PLEASE CONTACT

PASCAL BOLON, Ph.D.
biologics@eurogentec.com
<http://biologics.eurogentec.com>
 TEL.: +32 4 366 6116
 FAX: +32 4 365 1604

