

Establishing human embryonic stem cells master cell banks under gmp conditions

Anders Aspegren, Katarina Andersson, Catharina Brandsten, Kerstin Dahlenborg, Maria Flood, Dan Lundin, and Kristina Runeberg



Introduction

The increased interest in using human pluripotent stem cells for advanced therapy medicinal products (ATMP) has revealed the urgent need for safe and GMP compliant established hESC Master Cell Banks. The culture conditions for human pluripotent stem cells have evolved since 1998, leaving the use of mouse feeders and animal components behind, and instead one employ defined, feeder free culture conditions. However, the absolute majority of the derived human ES cells are made with feeders and outside the GMP environment, with animal components. Further, the starting material, the blastocysts, are not sourced according to FDA's guidelines, lacking proper donor testing, and/or sourced in non-prion free regions of the world.

Tissue Establishment

- For Tissue Establishment we applied, according LFVS 2008:12 on quality and safety norms when handling human tissue and cells.
- We are a registered Tissue Establishment since Sept 2017.
- (US) IRB and local (Swedish) ethical permit
- Import license
- Document storage > 30 years
- Validated transport company
- Control of donors
- Controlled dry shippers
- Storage of donor material
- Tissue Establishment license

Sourcing of Blastocysts

- Starting material:
 - Human fertilized eggs has to be sourced from FDA approved regions ("prion free"), such as USA, Iceland, Israel, Australia or New Zealand)

- Ethical approval and donor consent
 - Donor consent including the right to commercialize products

- Recruitment of donor couples
 - Signed donor consent

- Selection of donor couples
 - Medical history interview
 - Donor testing for infectious diseases/viruses*
 - Physical examination
 - Fulfillment of all acceptance criteria

- Sourcing of material
 - Cryopreservation
 - Transfer to Takara Bio Europe AB

*Donor screening maximum 7 days before/after IVF donation of blastocysts/embryos by Clinical Laboratory Improvement Amendments (CLIA) certified laboratories according to FDA recommendations.

GMP Facility

- Cell therapy is classified as an Advanced Therapy Medicinal Product (ATMP).
- Manufacturing License, according to LVFS 2004:7 and inspection on Eudralex Volume 4, EU GMP guidelines.
- Manufacturing License for the manufacture and storage of Master Cell banks for use in Clinical trial material

- Eudralex Volume 4, EU GMP guidelines (Annex 1 and 15)
- Manufacturing License (Annex 16)
- Quality Risk Management according to International Conference on Harmonization, ICH Q9 (Part 3 EU GMP)

- Continuous Environmental Monitoring
- Aseptic Process Qualification
- Process Flow Validation
- Material flow
- Environmental control system (Air, Equipment)
- Continuous Gowning Qualification
- Life Cycle Management
 - validations, (IQ, OQ, PQ)
 - equipment,
 - maintenance

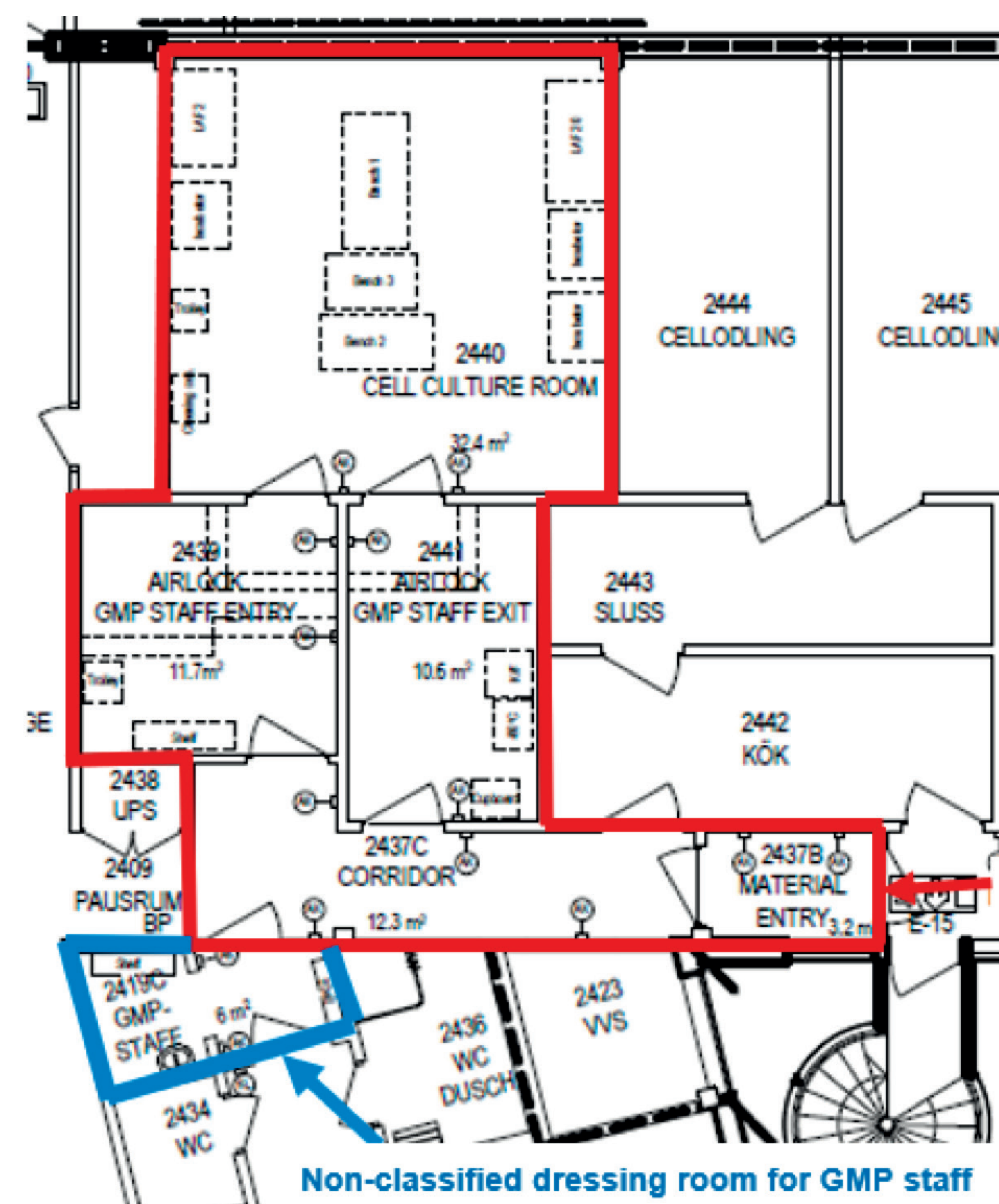
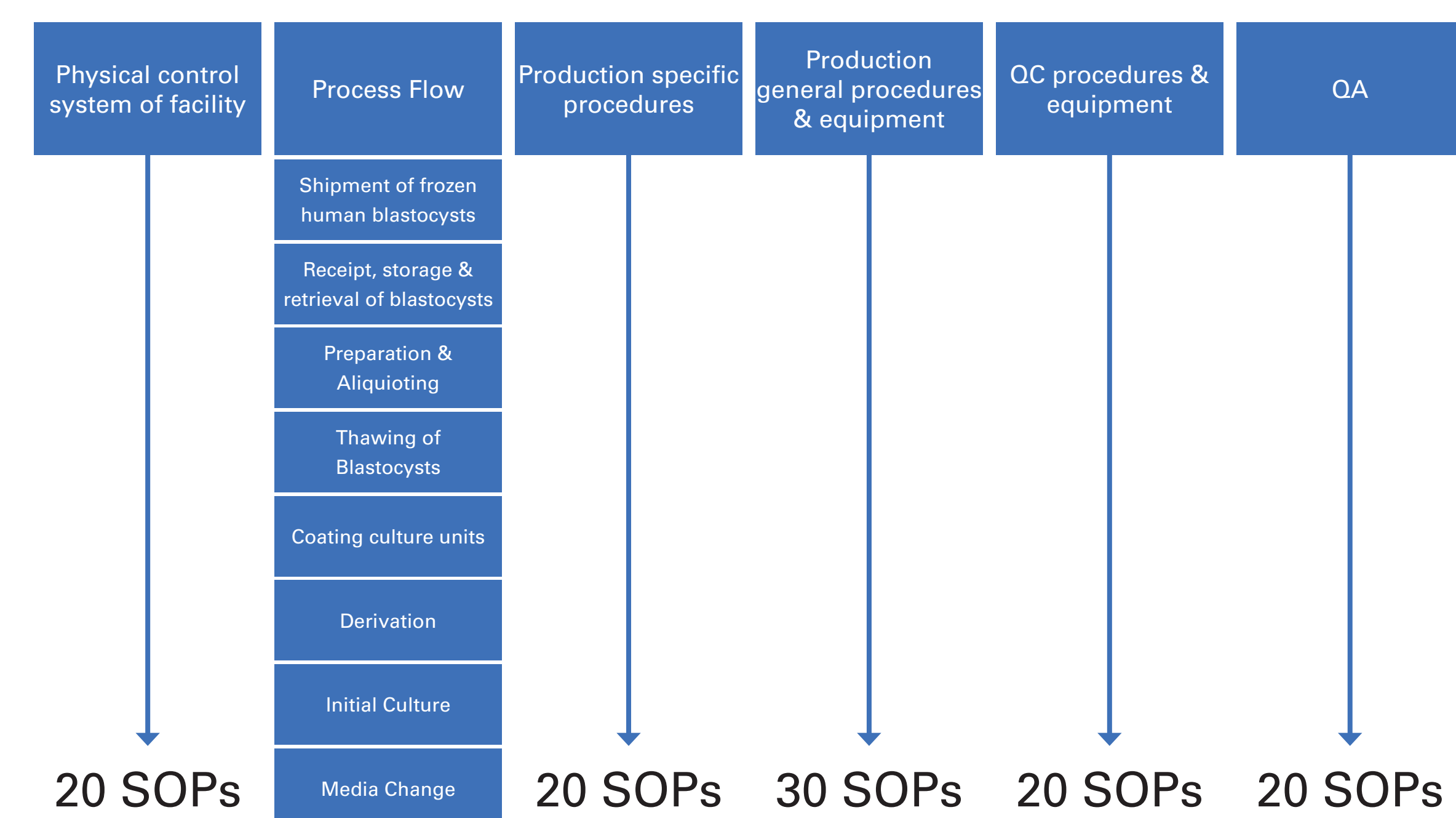
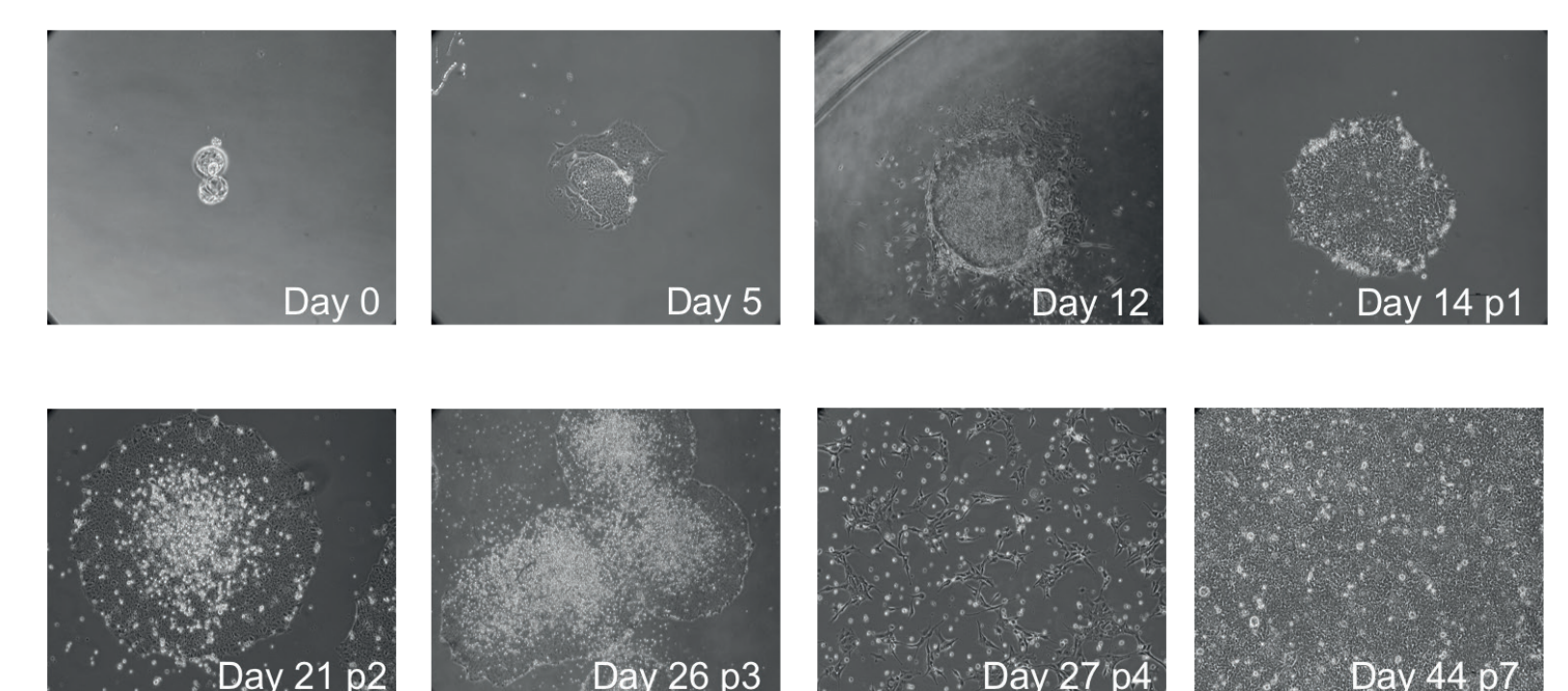


Figure 1. Detailed overview of GMP facility for ATMP. HEPA filtered air with regulated and controlled over-pressure (from Area B to D), together with control of strict gowning procedures, personell-and material flow and continuous Environmental Monitoring (static and dynamic) will assure the GMP set demands.

Processes surrounding the hESC establishment workflow



Previously Established hESC (non-GMP)



Quality Control and Release

Assessment of production data:

- Environmental control results from QC
- Lists and alarms from Production regarding Surveillance systems
- All Deviations from Manufacture and their Status
- Operational readiness document status
- Review of Batch records against the 'Batch tree' and GMP
- Review of QC assessment of analytical results against Product specification, CoA
- Release of SEED / MCB in relation to the reviewed documentation from Production and QC.
- Issue Certificates CoC, followed Manufacturing license, customer Quality agreement and GMP as expressed through TBEAB Quality system
- BSE/TSE according to EMA 410/01/revision 03 (current revision)
- Certify in a register telling where documents are stored and that released for further manufacturing of Clinical trial material