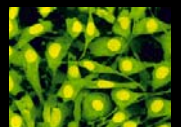


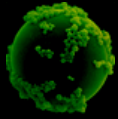


# Developing High Throughput Mammalian Expression Platforms

BPN Conference, Brisbane 2009

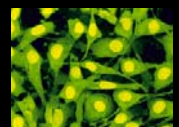
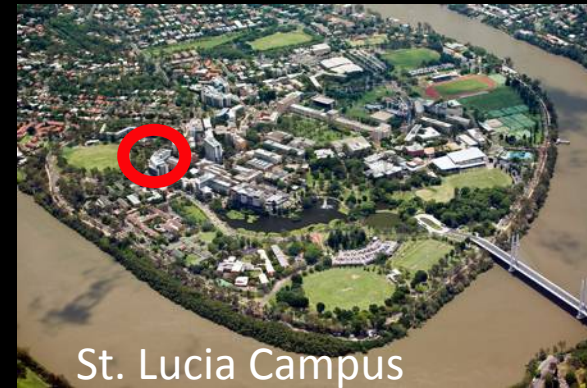
**Ben Hughes**  
([ben.hughes@uq.edu.au](mailto:ben.hughes@uq.edu.au))

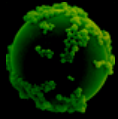




# AIBN, University of QLD

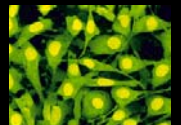
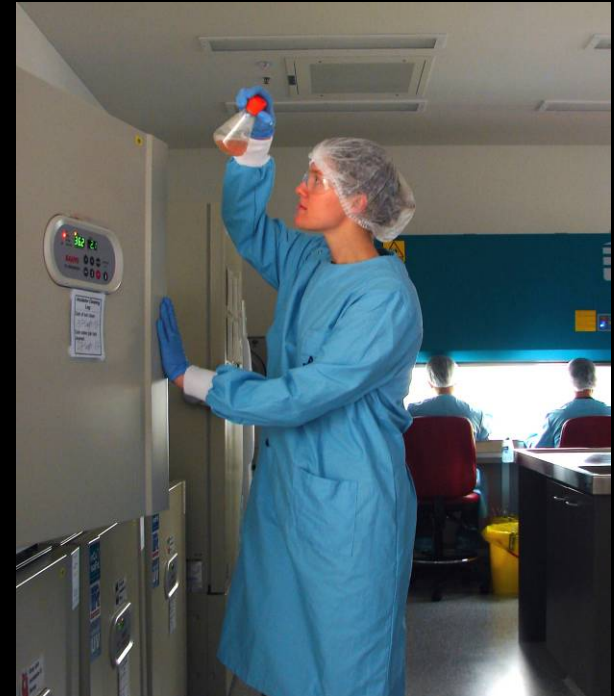
- **AIBN houses > 350 researchers**
  - Varied human health research areas
    - Biotherapeutic production
    - Vaccines and delivery systems
- **NCRIS Biologics Facility**
  - 200m<sup>2</sup> of dedicated clean rooms for mammalian cell culture and downstream processing
  - Bioreactors to 100-L scale
  - Protein analysis and characterisation
- **Acyte Biotech Pty. Ltd.**
  - Spin-off company (UNSW/UQ)
  - Novel IP for mammalian expression

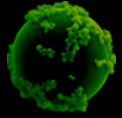




# Presentation Outline

- **Background and drivers**
  - Use of mammalian cells for biotherapeutic manufacture
    - mAbs and CHO cells
- **Developing high throughput platforms**
  - Transient protein production
    - *EpiCHO* system
  - Stable cell lines
    - Leveraging high throughput instruments
    - Fed-batch optimisation



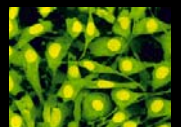


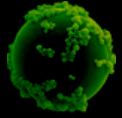
# Mammalian Biotherapeutic Manufacture

- **Widely used host system**
  - Account for ~ 50% of marketed recombinant proteins
    - > 70% produced in CHO<sup>1</sup>
  - Well characterised → platform processes to 20KL
- **Produce bioactive molecules**
  - Correct folding and desirable post-translational modifications
  - Proven track record / acceptance by regulatory bodies
- **Ideally suited for mAb production**
  - > 30% of biopharmaceuticals in current clinical trials<sup>2</sup>
    - Sustained double digit growth through 2007
    - Surpassed growth factors as the highest selling category of biologics

<sup>1</sup> Aggarwal S, 2007, Nat. Biotechnol

<sup>2</sup> Walsh G, 2006, Nat. Biotechnol.

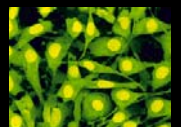


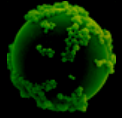


# Challenges with Mammalian Systems

---

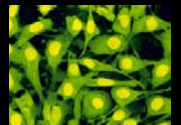
- **Time and resource-intensive**
  - Technically complex, lower cell density and yields
    - Higher COGs
  - Selection of high-producing, quality cell lines
    - Clone identification / verification ranges from 4-12 months
    - Cell line instability issues
- **Restricted access**
  - Major advances of the last decade are not 'freely' available
    - Heavy licensing costs for access to production cell lines / vectors
    - Proprietary media and fed-batch processes

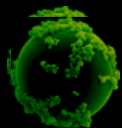




# Rapid Production of mAbs

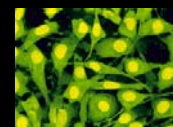
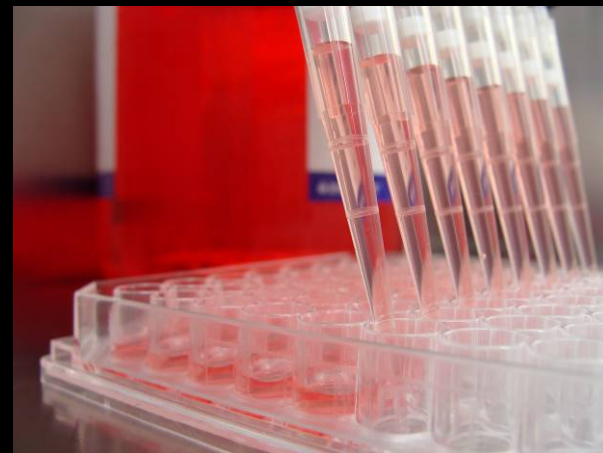
- **Recognising these challenges our group directs research towards:**
  - Improved cell lines (focus on CHO and mAbs)
  - Molecular engineering of expression platforms
  - High-throughput and intelligent single cell isolation
- **Consortium formed under Federal ICIP grant**
  - Low-cost mAb production (transient and stable mammalian cells)
    - Acyte Biotech Pty. Ltd
    - Agen Biomedical Ltd.
    - BioPharmaceuticals Australia
  - Demonstrate that this technology is available to Australian researchers and companies
    - There exist alternative options for mammalian expression
    - Adding value to your process

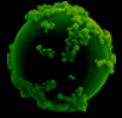




# Presentation Outline

- Background and drivers
  - Use of mammalian cells for biopharmaceutical manufacture
    - mAbs and CHO cells
- **Developing high throughput platforms**
  - Transient protein production
    - *EpiCHO* system
  - Stable cell lines
    - Leveraging of FACS and ClonePix instruments
    - Fed-batch optimisation

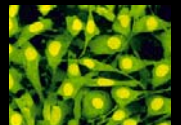




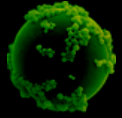
# Transient Expression Systems

---

- **Rapid generation of recombinant protein**
    - Allows for timely screening of pre-clinical drug candidates
      - Gram quantities in 1-3 weeks
    - Expression in a wide variety of mammalian systems
      - Transient systems in HEK continue to outperform CHO
  - **Challenges for transient systems:**
    - Rapid dilution of plasmid DNA as cell proliferates
    - Continuity of host cell line
      - Early stage material often manufactured in HEK, while final production host is CHO
- Ideally require a CHO-based transient system that maintains high expression levels over an extended period







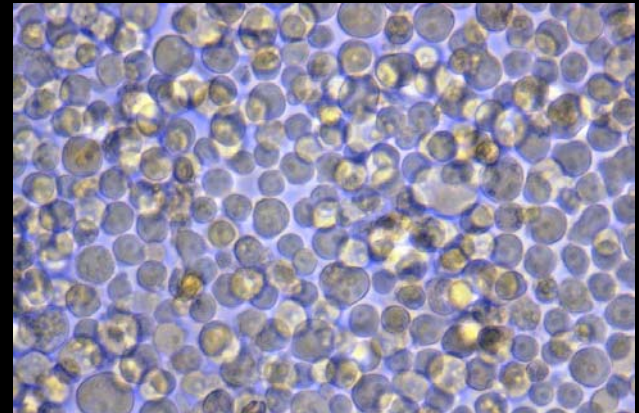
# The *Epi*CHO System

- *Epi*CHO is a transient expression system that provides for amplified and prolonged recombinant protein expression

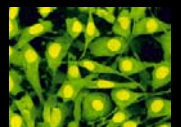
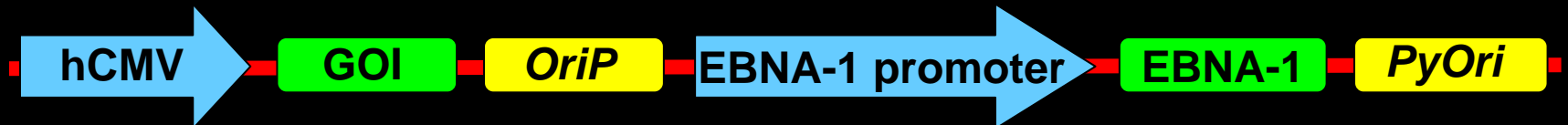
- Two key elements:

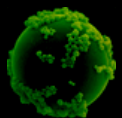
1. The host cell line **CHO-T**

- Suspension adapted to SFM
- Constitutively expressing PyLT

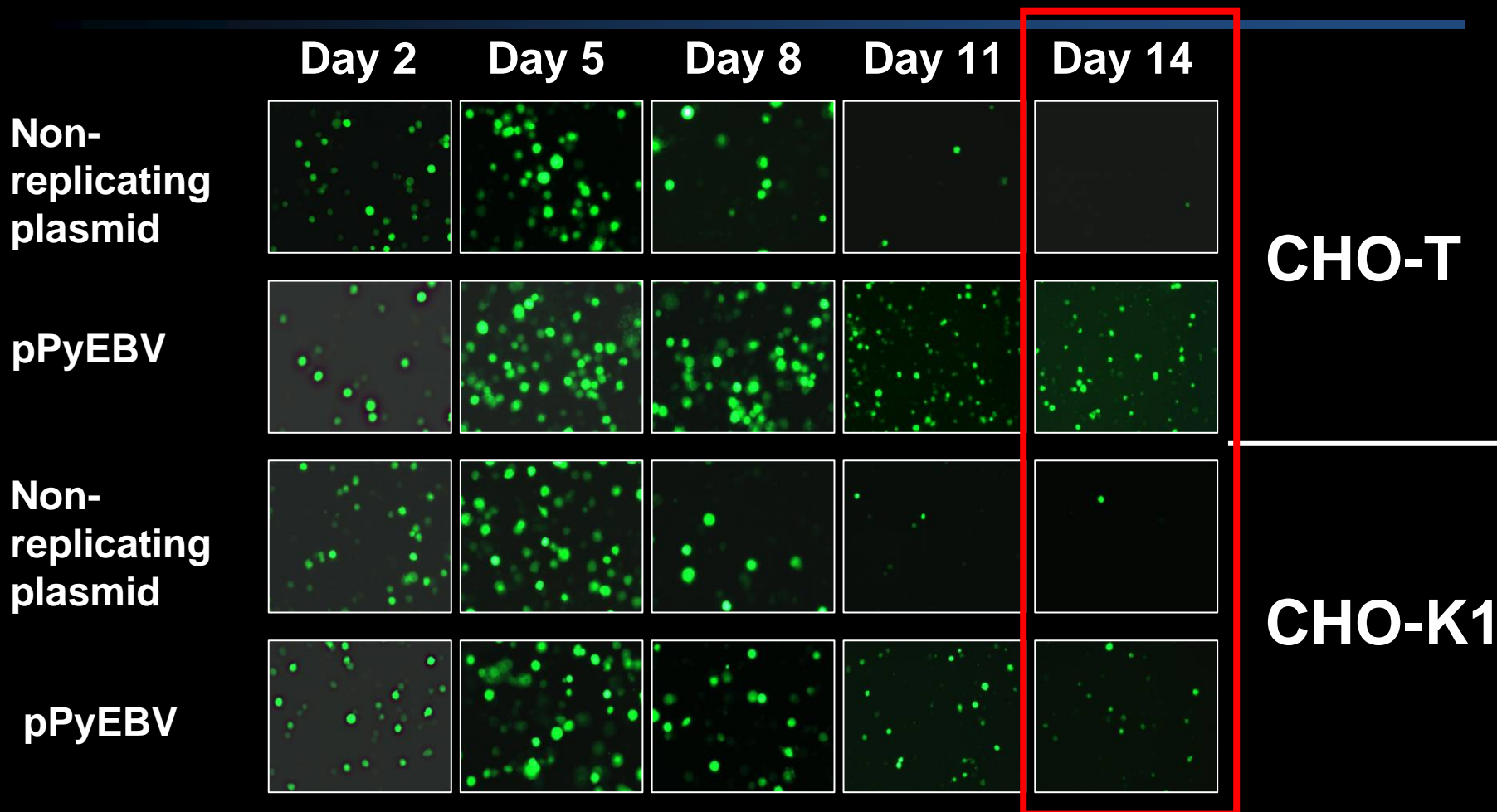


2. The expression vector **pPyEBV**

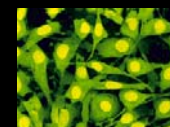


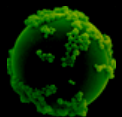


# Prolonged Expression with *EpiCHO*



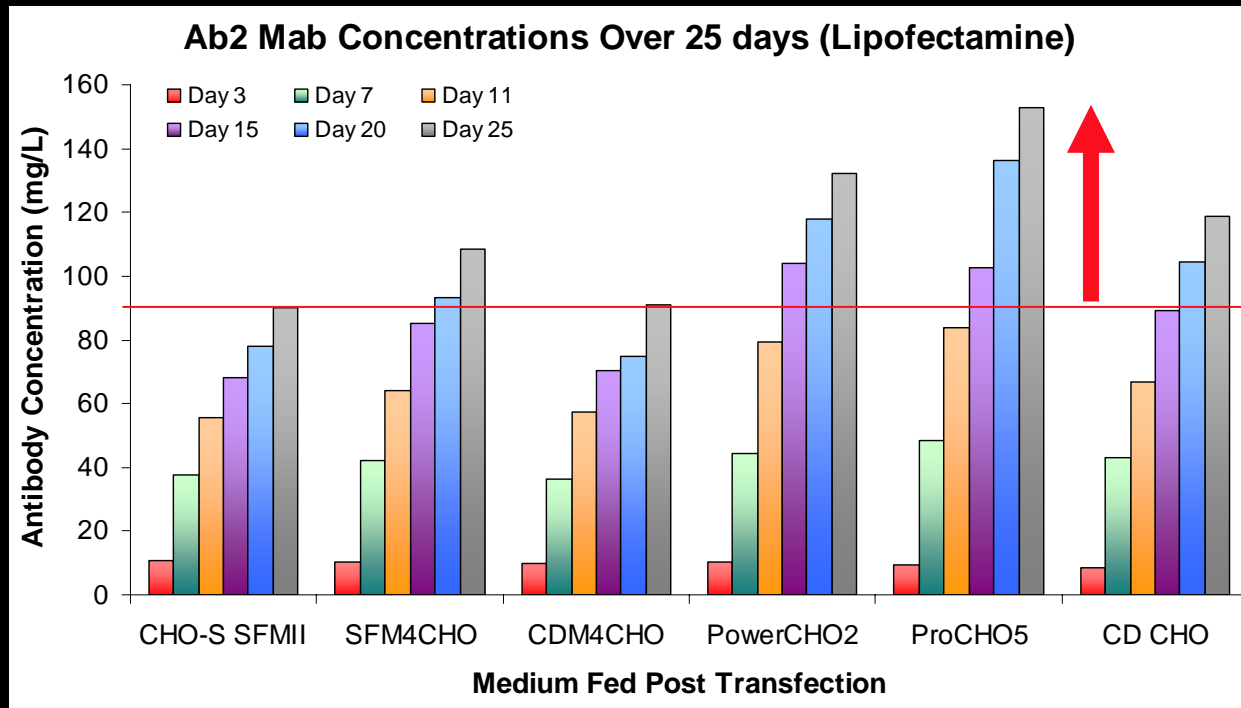
- Transient transfection with destabilised EGFP reporter



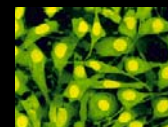


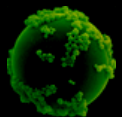
# Prolonged mAb expression with *Epi*CHO

- **CHO-T cells transfected with Agen Ab**
  - Evaluated productivity in different growth media
    - 50% media dilution 4 hours post transfection
    - Shift to hypothermic conditions (32°C)



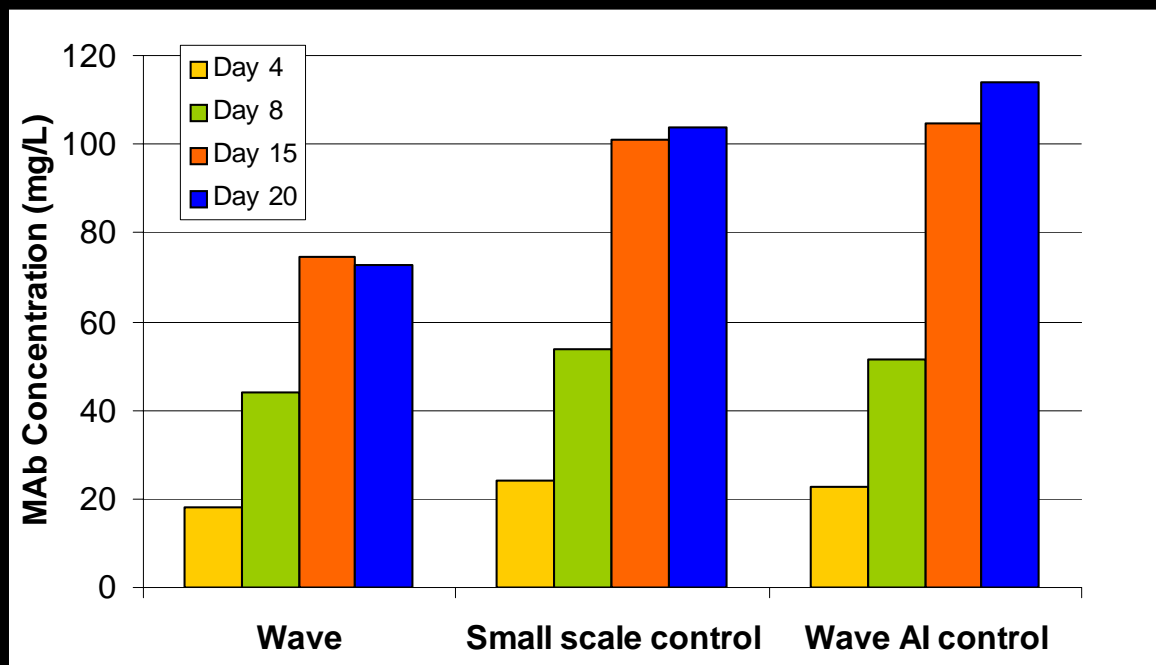
60% increase



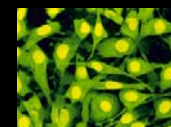


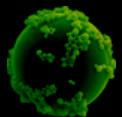
# Scalable Transient Production

- Traditionally challenging to maintain productivities and efficiency upon scale-up
  - Transfected Agen mAb at 7L volume in a disposable Wave reactor
    - Small scale controls (duplicate 125mL shake flasks)



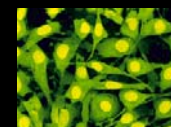
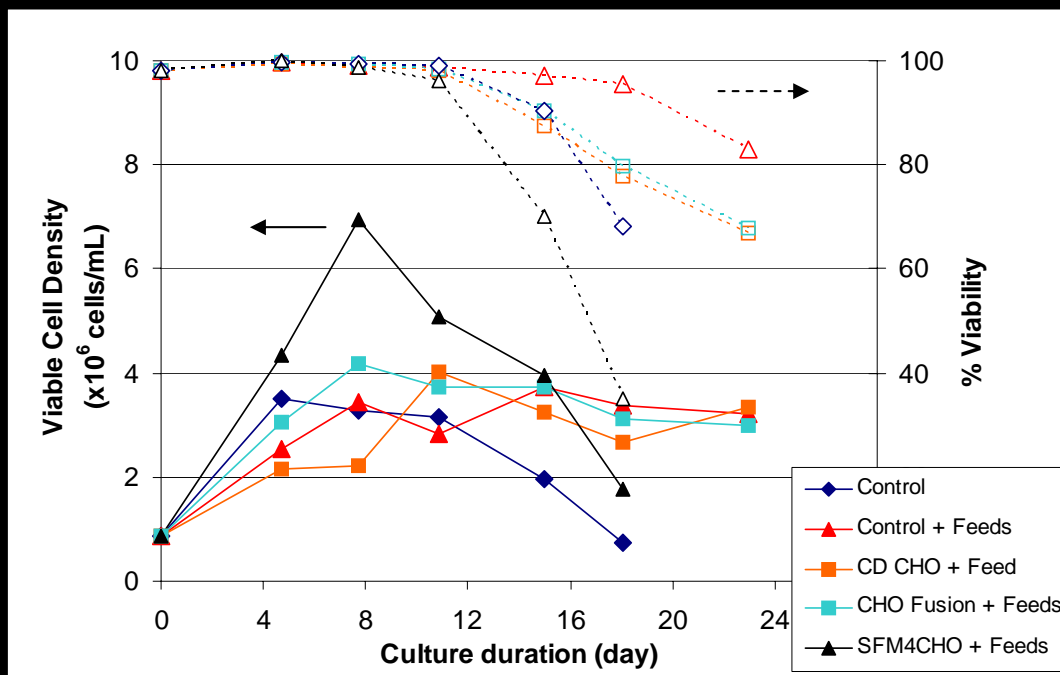
- Wave culture not linearly scaled
- Gram quantity from a 10L wave should be achievable

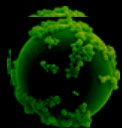




# EpiCHO Path Forward

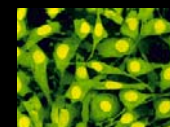
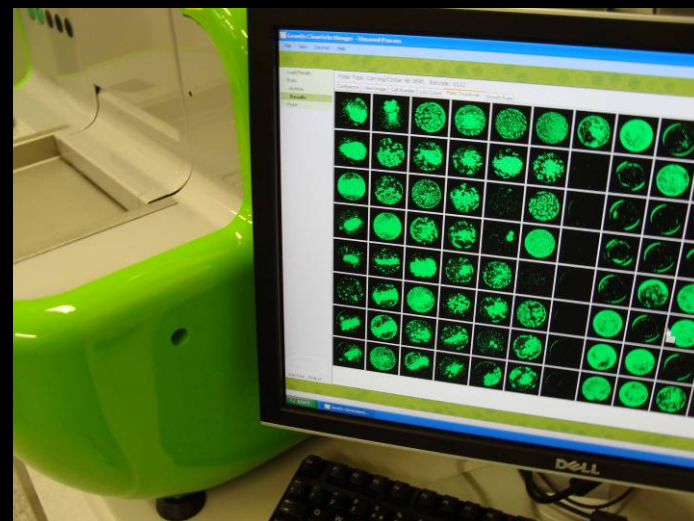
- **Continue large-scale transfections**
  - Optimise wave growth (base addition for pH control)
  - Evaluate alternative culture vessels (3-L disposable flasks)
- **Evaluation of new media (data for 32°C growth only)**
  - Fed-batch
    - Hydrolysates
  - Supplements
    - IGF-1
    - rTransferrin
  - Extended culture duration and viability
  - 50% increase in IVC

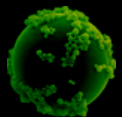




# Presentation Outline

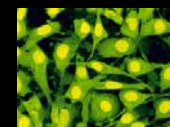
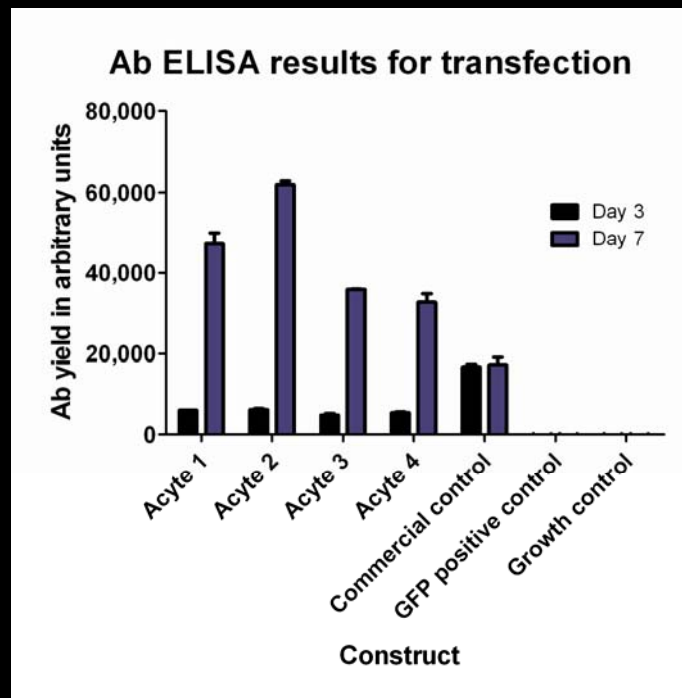
- Background and drivers
  - Use of mammalian cells for biopharmaceutical manufacture
    - mAbs and CHO cells
- Developing high throughput platforms
  - Transient protein production
    - *EpiCHO* system
  - Stable cell lines
    - Leveraging high throughput instruments
    - Fed-batch optimisation



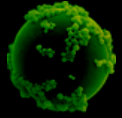


# Stable Expression Technology

- We have assembled a complete set of “in-house” expression vectors
  - Modular design
  - Proprietary high-level expression technology
  - Out-performs commercial mammalian vectors
  - Options for several expression units on a single plasmid
- Stable cell line
  - Based on a low-passage ATCC CHO-K1 variant
    - Suspension adapted
    - Maintained in fully defined conditions
  - Custom designed metal amplification system
  - Modified DHFR+ vectors also in use







# High Throughput Technologies for Cell Selection

- **Automated high-throughput instruments:**

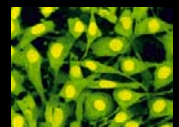
Genetix ClonePix FL



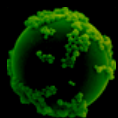
BD FACS AriaII



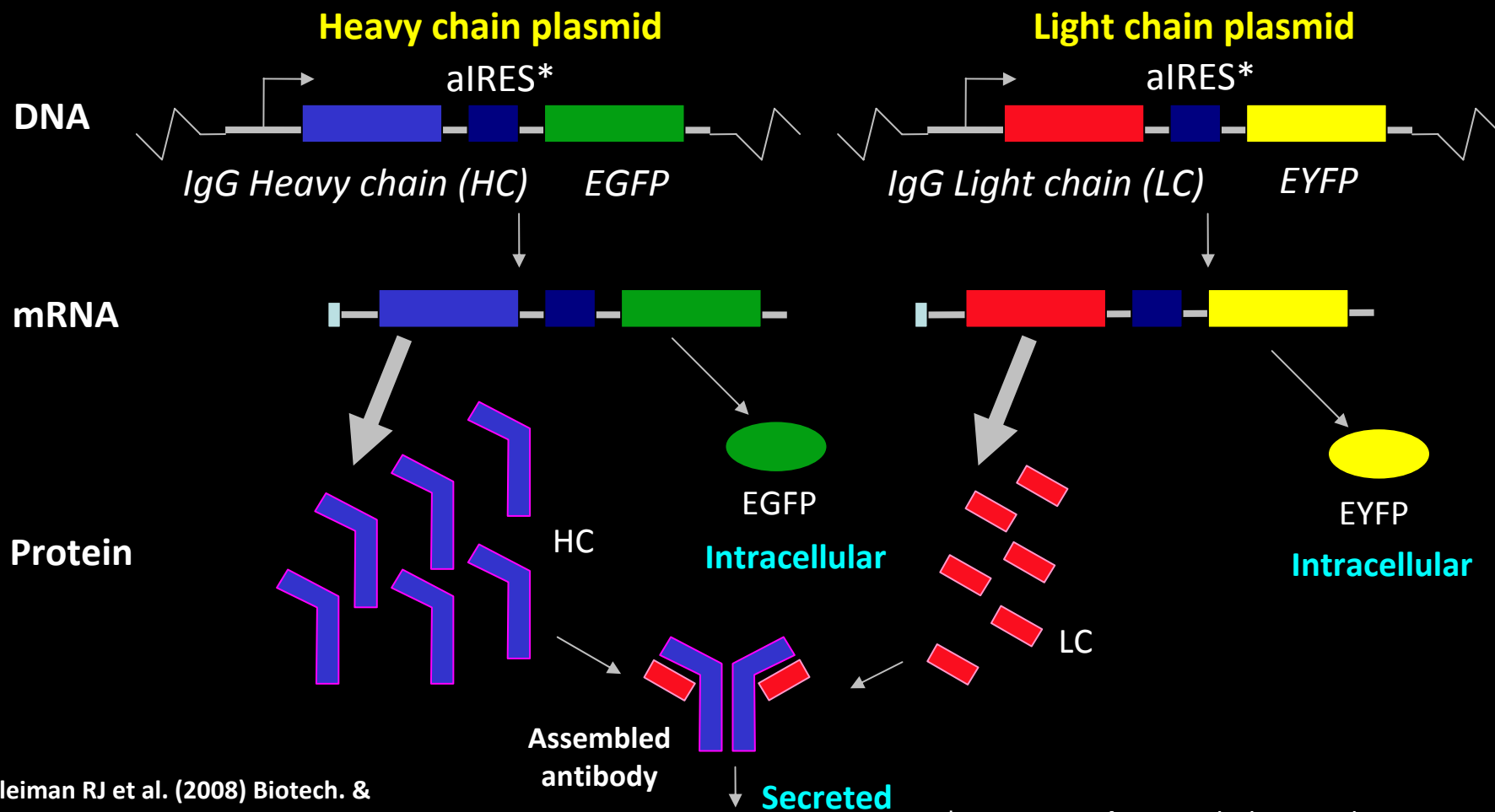
- **Facilitates rapid selection of high-expressing clones**
  - Utilises Acyte technology
    - Expression vectors with two-colour transcriptional reporters





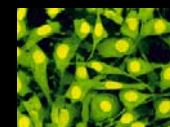


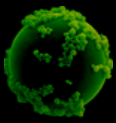
# Acyte Dual Bicistronic Reporter / Expression Vectors



Sleiman RJ et al. (2008) Biotech. & Bioeng.

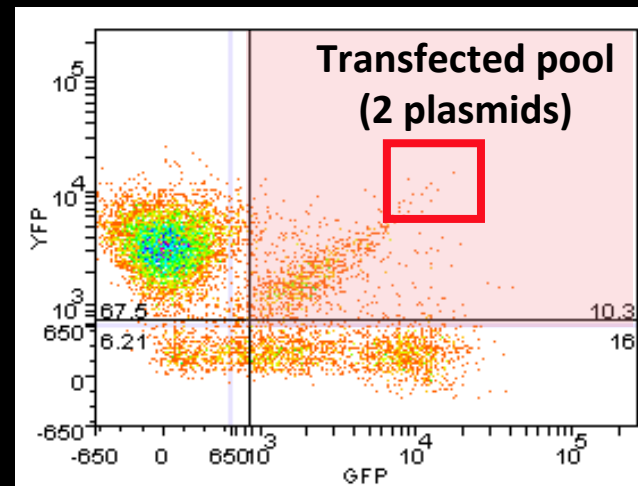
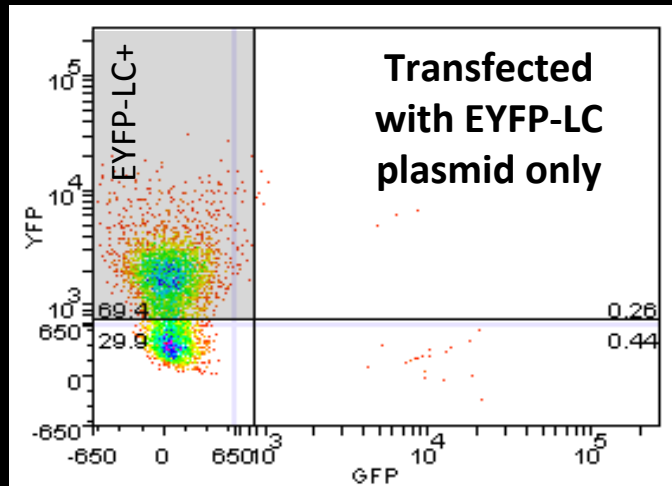
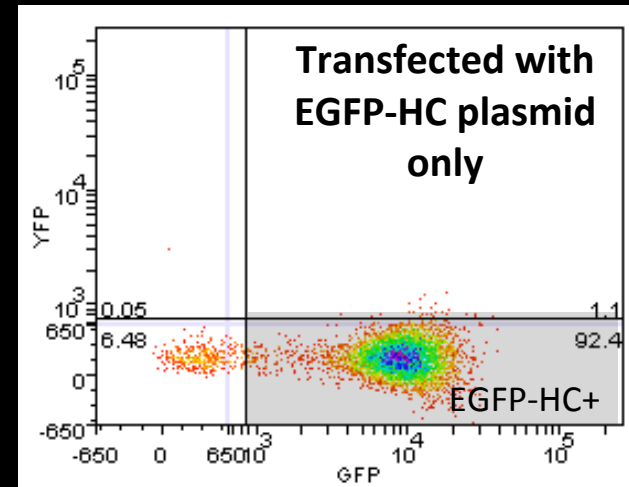
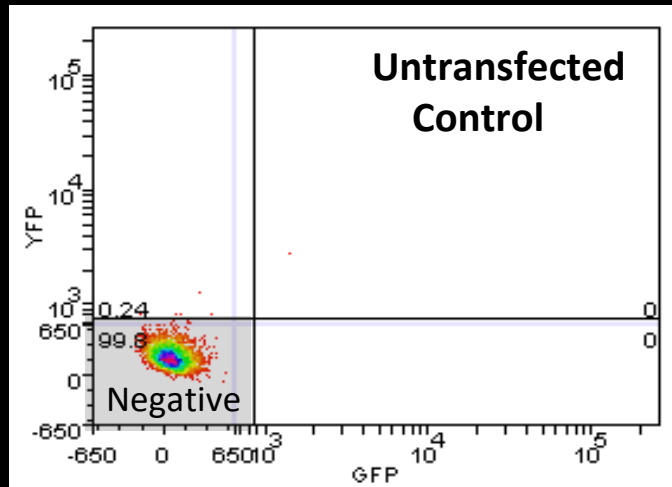
\* Attenuated internal ribosomal entry site





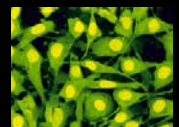
# FACS Single cell analysis of IgG LC/HC transcription in transfected pools

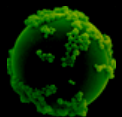
EYFP Fluorescence (IgG LC Transcription)



Antibody  
expressing cells  
(both chains)  
**High mAb  
expressers**

EGFP Fluorescence (IgG HC Transcription)





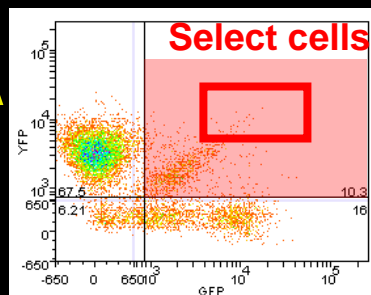
# Typical FACS Workflow

- BD FACSAriaII™: high speed data acquisition and cell sorting at rates up to **70,000 cells/sec** in up to 15 parameters, along with single cell deposition for clonal isolation

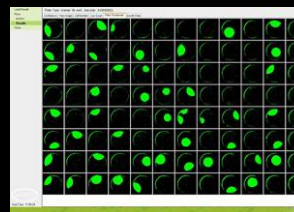


BD FACSAriaII

→  
EYFP (mAb LC)



→  
EGFP (mAb HC)



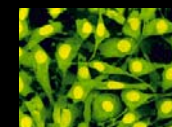
96-well plate culture. Verify clonality, screen for growth and productivity

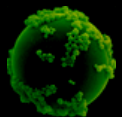


CloneSelect Imager



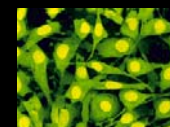
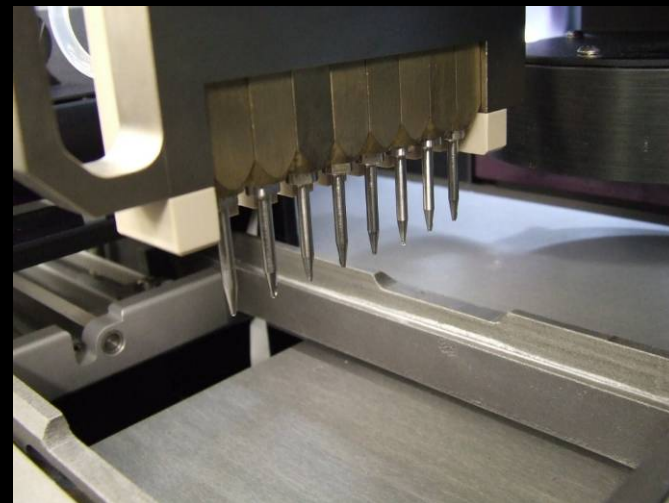
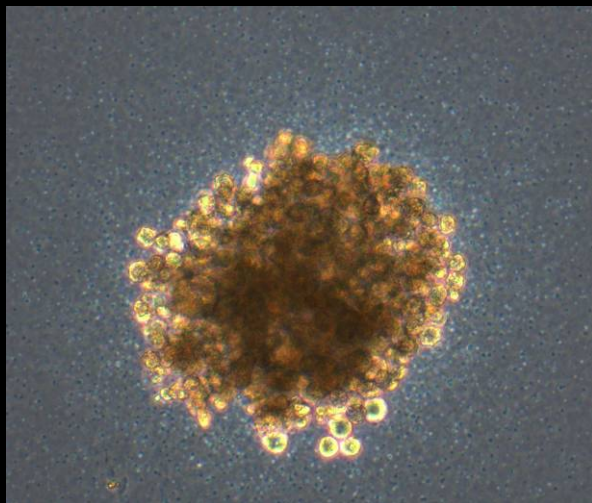
Expansion and Production

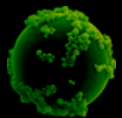




# Genetix ClonePix FL

- **Automated high resolution imager with sterile robotic picking head**
  - Selection of cells secreting recombinant proteins (e.g. mAbs).
  - Plate cells at low concentration in semi-solid media
    - Single cells grow into discrete colonies (7-14 days)
  - Secreted recombinant product detected by fluorescently tagged antibodies specific for the target protein.



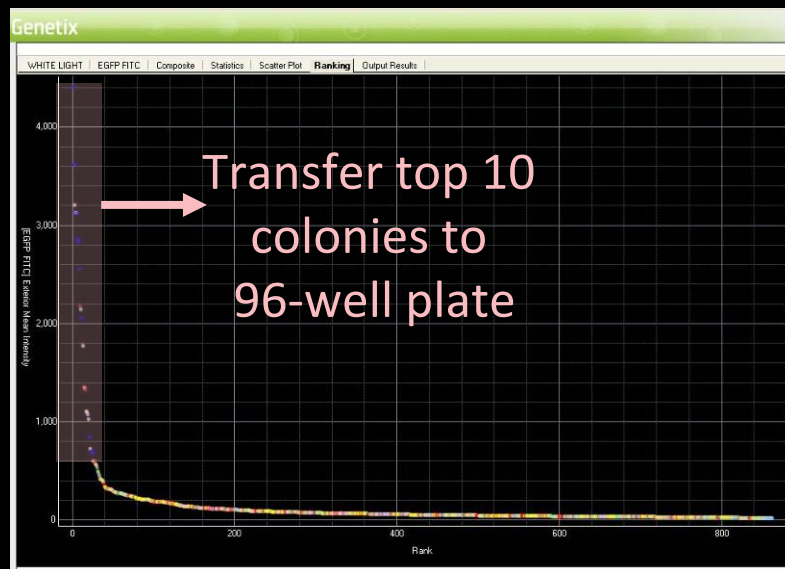


# Selecting mAb Producing Clones

- **Example composite image (white-light and florescent images)**
  - Clearly identifies secreting colonies of interest
  - Selection criteria can be customised to select high-producing, robustly growing colonies

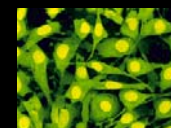


Fluorescence Intensity

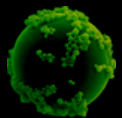


Colony #

Evaluate 1000's of cells / plate

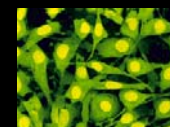
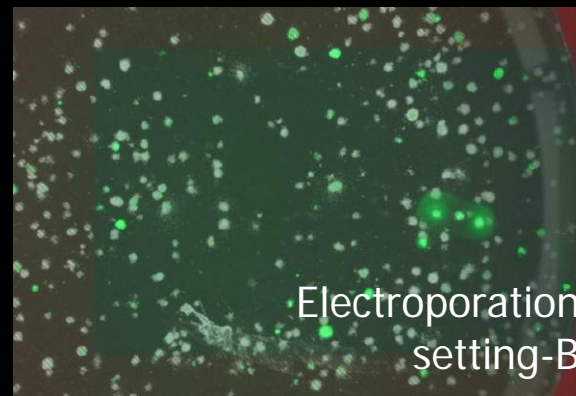
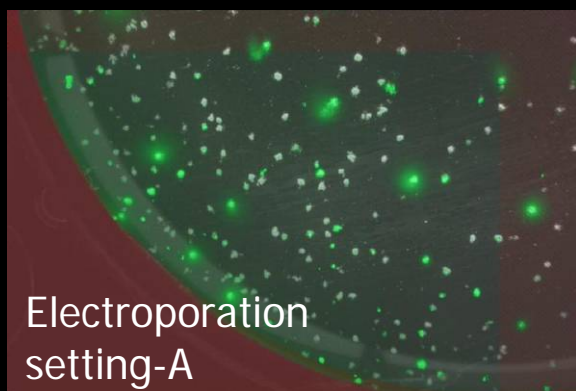
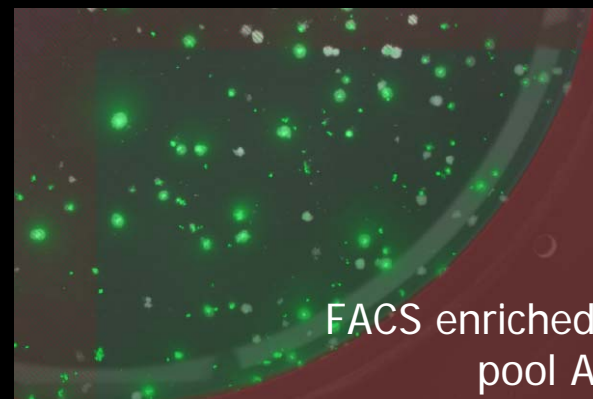
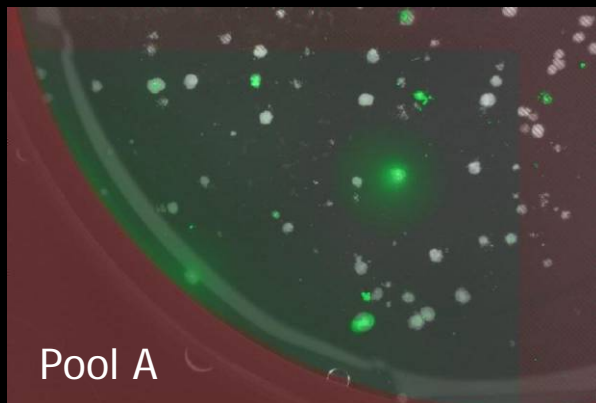


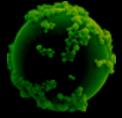




# ClonePix Advantages

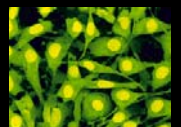
- Increased success rate for growth post picking
- Provides a visually intuitive feedback on cell line and transfection conditions

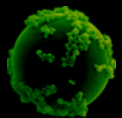




# Stable Expression of the Agen mAb

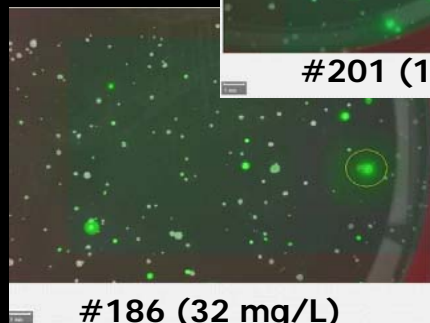
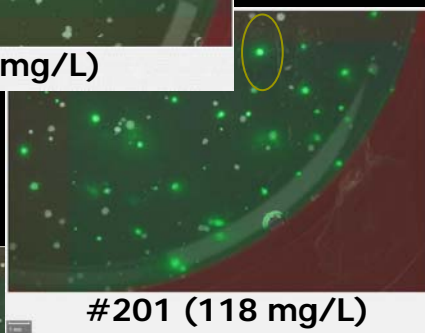
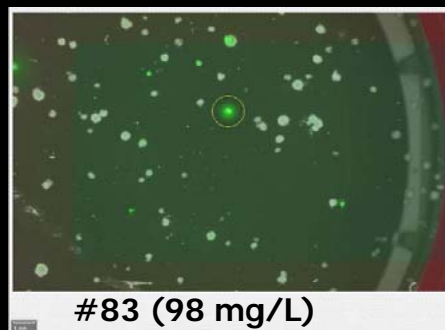
- **Stable pools generated with Agen antibody cDNA**
  - clones selected for scale-up in ~5 months
  - Project timeline:
    - cDNA resynthesis 3 weeks
    - Vector construction / verification 4 weeks
    - Electroporation and G418 selection 3 weeks
    - Clonal isolation:
      - FACS and ClonePix 2-4 weeks
    - Clonal screening / adaptation 2-4 weeks
    - Clone verification:
      - Fed-batch analysis / monitoring stability 6-8 weeks
  - Note: Pools were not amplified





# Analysis of Selected Clones

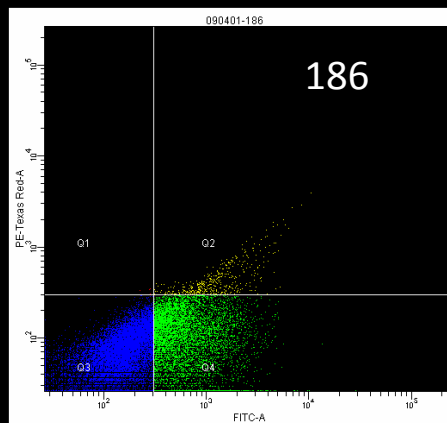
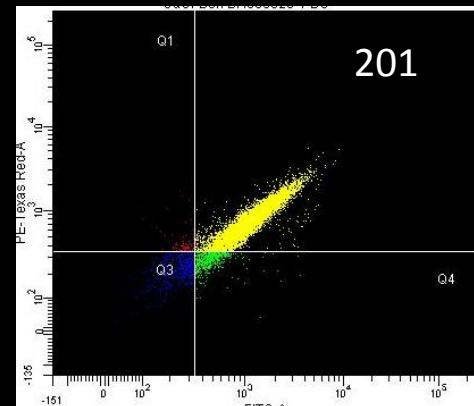
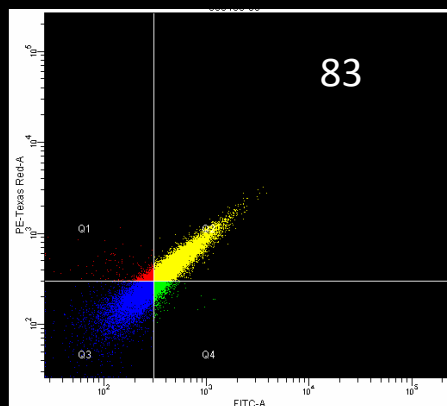
- ClonePix



(7-day batch accumulations)

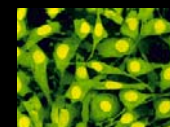
## FACS

mCherry Fluorescence (IgG LC Transcription)

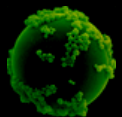


- Good correlation with ELISA productivity

EGFP Fluorescence (IgG HC Transcription)







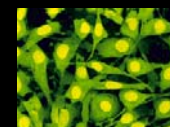
# Fed Batch Evaluation

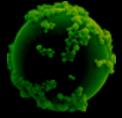
- Investigated commercially available media / supplements

- Spent media analysis
  - Glucose / glutamine
  - Amino acids
  - Trace minerals
- Hydrolysates
- Temperature shift
- CHO Efficient Feed

- Standard conditions:

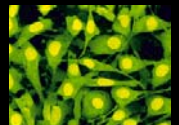
- ICD =  $0.2 \times 10^6$  cells/mL
  - 100% media exchange
- 125mL baffled shake flasks
  - 250mL square bottles
- 37°C, 7.5% CO<sub>2</sub>, 170-200 RPM

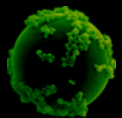




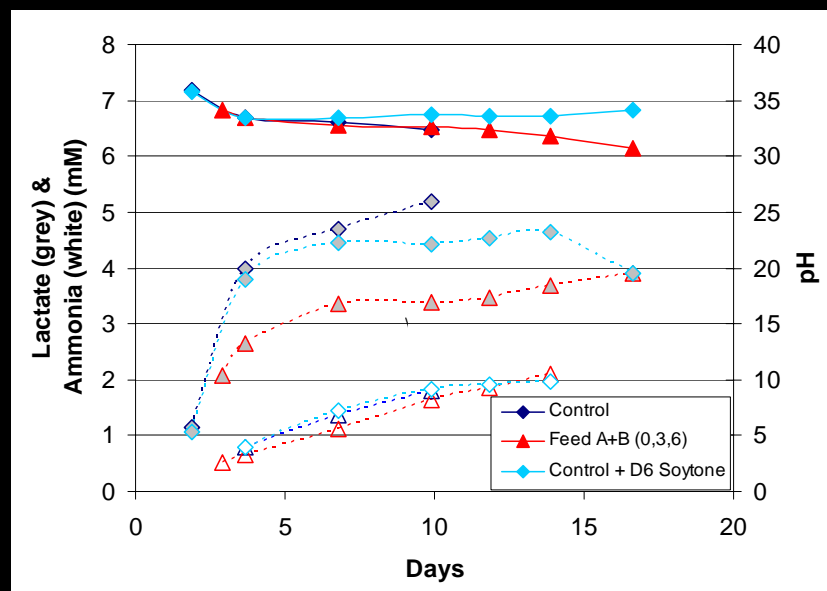
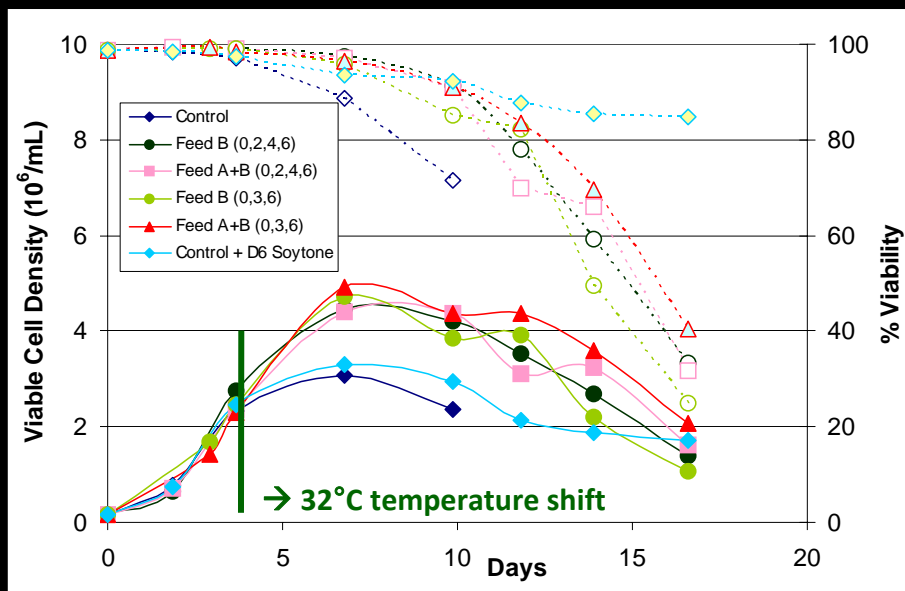
# Efficient Feed Supplements “A” and “B”

- **Chemically defined, protein free**
  - *Contains:* Carbon source, concentrated amino acids, vitamins, salts, trace minerals
  - *Does not contain:* Lipids, hydrolysates, or growth factors
- **Mini-DOE for multi-day feeding:**
  - 15% initial feed was beneficial
  - Adding B or A+B yielded highest titer
    - Noted altered cellular metabolism with reduced lactate / ammonia accumulation
    - Increased cell diameter, higher osmolality (~350 mOsm/L)
- **Follow-up experiment with most promising conditions:**
  - Additionally investigated addition of hydrolysates
  - Hypothermic temperature shift



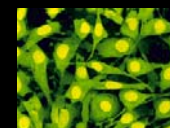


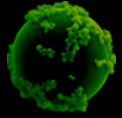
# Efficient Feed Optimisation



- **Feed A+B (day 0,3,6) yields best cell growth and mAb titer**
  - Positive effect of Day 6 Soytone addition (cell growth and pH)

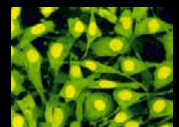
→ Identified optimal feeding conditions

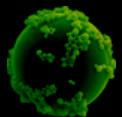




# Scale-up of Fed-batch process

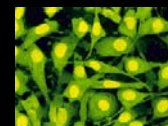
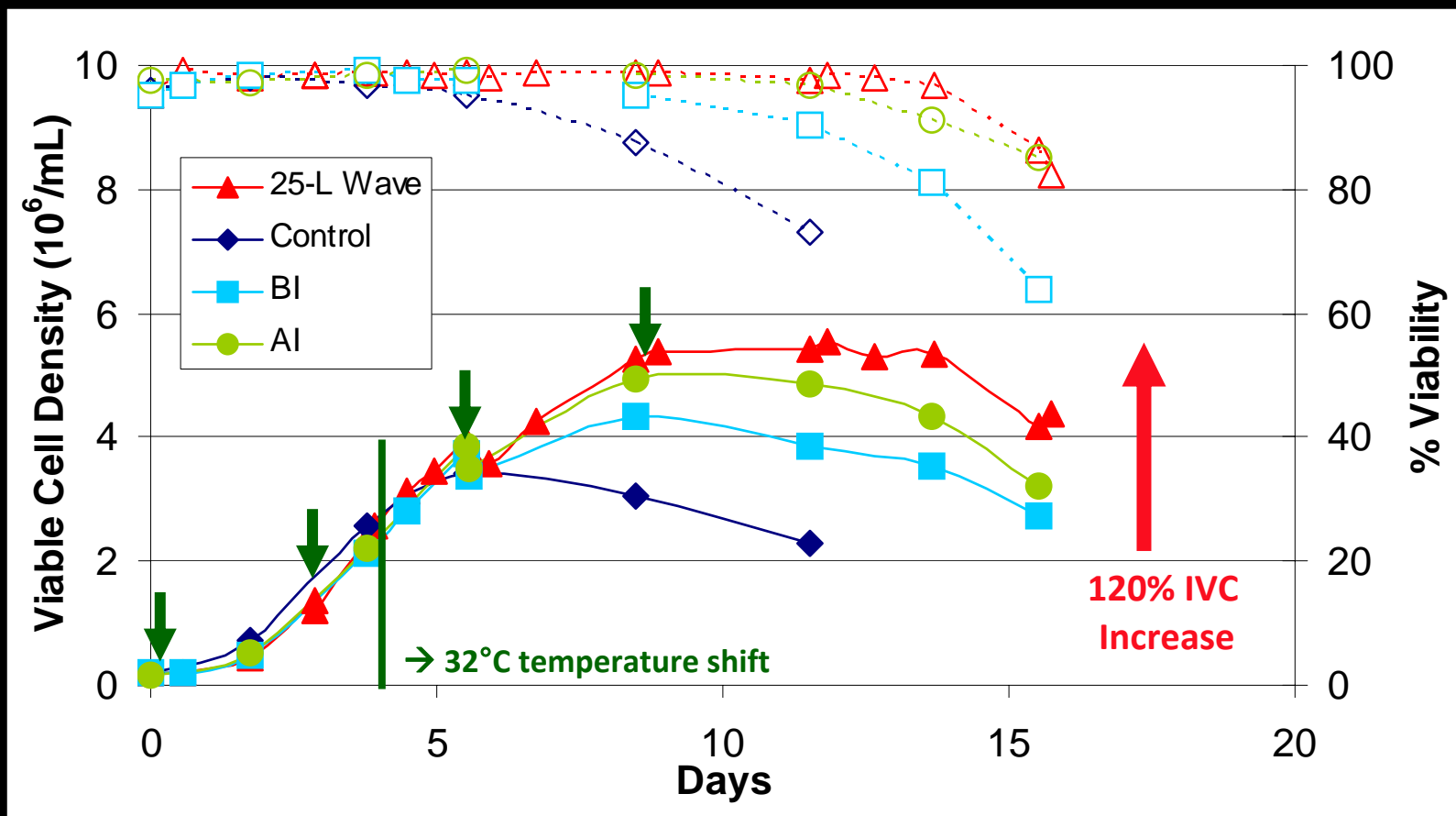
- Performed 25-L fed-batch process (50-L Cultibag)
  - pH manually adjusted (%CO<sub>2</sub> and 1M Na<sub>2</sub>CO<sub>3</sub> addition)
  - Culture Feeds:
    - Efficient Feed A+B
      - 15% Day 0, 3, 10% Day 6
    - L-glutamine
      - 8mM Day 0, 5mM Day 6
    - Soytone Hydrolysate
      - 5g/L Day 9

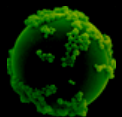




# 25-L mAb Production Run

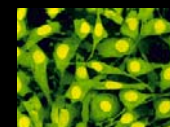
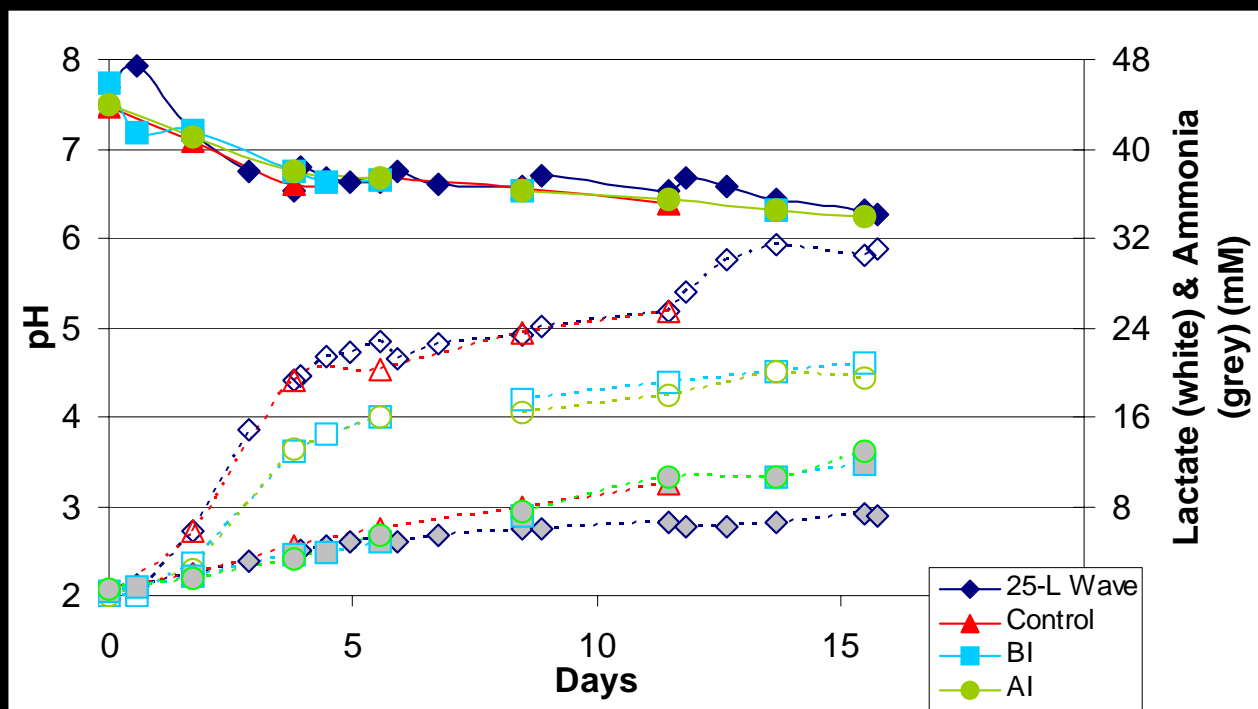
- Cell growth and viability:

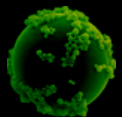




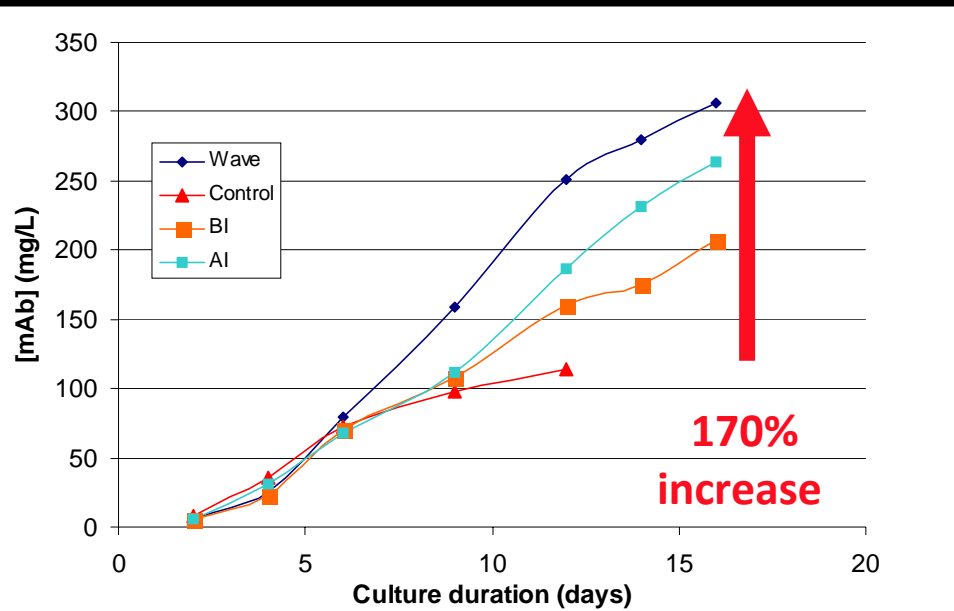
# Metabolites and pH profiles:

- Wave pH maintained at  $\sim 6.7$ 
  - No base addition to small scale cultures
- Reduced lactate in Efficient Feed cultures
  - Wave suffered elevated pH (7.9) for initial 12 hours





# 25-L mAb Productivity Results

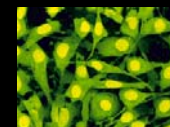


- Significant mAb increase over batch control
  - Product quality retained as compared to Agen reference mAb
- Acceptable productivity for non-amplified clone ( $q_p \approx 7$  pg/cell.day)

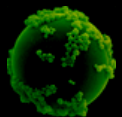
- Continued fed-batch evaluations

- Re-adaptation of above clone to alternative media and feeds:

Media	Feed	IVC increase	Titer increase
CD CHO (control)	15% A+B	-	-
SFM4CHO	15% A+B	2%	-54%
CHO Fusion	None	19%	76%
CHO Fusion	Fusion Hydrolysate	24%	18%

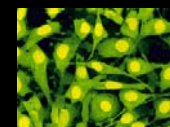




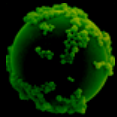


# Summary

- **Successfully demonstrated mAb production both transient and stable expressions systems**
  - This technology is readily available to the Australian biotech industry
- ***EpiCHO* is a transient system that promotes elevated and prolonged expression**
  - Platform for rapid scalable production
- **High-throughput equipment are powerful tools for efficient and intelligent clonal isolation**
  - Provide continuous feedback for streamlining clonal selection
- **Significant process improvements can be realised with fed-batch optimisation**
  - “Off the shelf” solutions (+ a liberal amount of sweat)



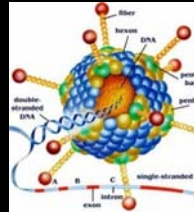




# Acknowledgements

- **AIBN – Mammalian Cell Engineering Group**

- Peter Gray
- Trent Munro
- Warren Pilbrough
- Joe Codamo
- Stefanie Dietmair
- Jeff Hou



Images courtesy of <http://en.ec.cn>,  
[www.theanimalfiles.com](http://www.theanimalfiles.com)

- **AIBN – NCRIS Biologics Facility**

- David Chin
- Karen Hughes
- Steve Mahler
- Michael Song
- Ashleigh Linville
- Jongwei Wooh
- Kristin Raphaelli
- Kym Hoger
- Cherrine Chan



- **Agen**

- Michael Gerometta
- Katey Whitham

- **BioPharmaceuticals Australia**

- David Hughes
- Brett Whitecross

- **QLD and Federal Governments**

- **Suppliers**

- InVitro
- SAFC
- Invitrogen
- BD
- Millipore
- Novozymes
- GE
- Sartorius-Stedim
- Genetix

