

Development of a customized quantification of engineered proteins with the Octet

Hervé Dubois, Sanofi R&D

Contents

- INTRODUCTION
 - Vitry Site, Bio-Manufacturing Analysis
- OBJECTIVES
 - Context
 - Assay requirements
- ASSAY DEVELOPMENT
 - Assay strategy, preliminary test with BLtz
 - Ligand loading
 - Regeneration buffer selection
 - Assay optimisation
- CONCLUSIONS
- ACKNOWLEDGEMENT

Personal presentation

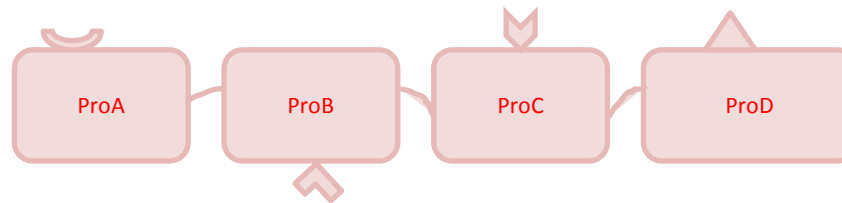
-
- Rhone-Poulenc Rorer R&D Vitry Site in 1992
 - Technician in Solid Phase Peptide Synthesis and Combinatory Chemistry
 - Gencell/Centelion, Aventis Group
 - Technician in Quality Control Department for release lot of Recombinant Adenovirus and Plasmidic DNA
 - Sanofi Industrial Affairs, Process Development Biochemistry
 - Analyst of recombinant protein (IgG)
 - Sanofi R&D, Bio-Therapeutic
 - Analyst of recombinant protein



- **Bio-Process Analytics:**
 - **Analytical support to Bioprocess Development: Cell Engineering, Up-Stream Processing at different production scales, Down Stream Processing**
 - **Analysis: Product titration at different stages of process, quality of product, contaminants titration, media composition during process...**

- Method to be developed for clone selection: High Throughput Method

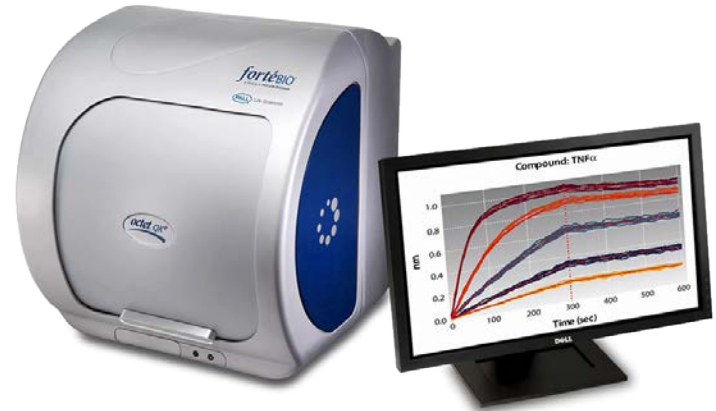
- Molecule:



- Tetrameric protein where each monomere is specific for a target
- Monomeres are separated by a peptidic spacer
- No specificity :
 - No tag (Like poly-Histidine or streptavidin...)
 - No fragment easy to detect (Fc fragment for example)

Requirement

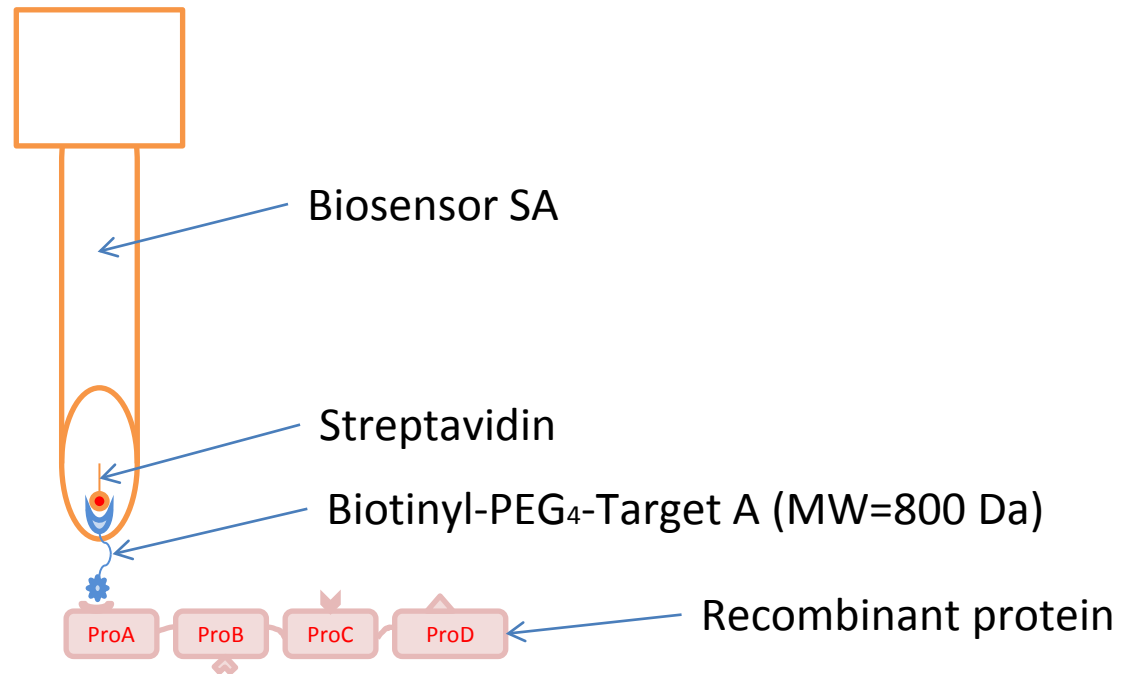
- For first screening, protein assay needs to be:
 - **Rapid to develop** because of short delay
 - **Fast:** 1000 samples should be generated
 - **Fast:** Because of cell passage, results are needed in 2 days
 - **Easy:** Without sample preparation if possible
 - **Sensitive:** 1st screening is tested at Day 4 of culture; range between 0 to 100 $\mu\text{g/ml}$.
 - **Specific:** samples are assayed in supernatant. Some cell proteins are present
 - **Low cost:** Because of the large number of samples
 - **Simple** data reprocessing



Preliminary Tests



Preliminary tests and titrations were realized with BLitz system by Frankfurt Analytical Team of Bio-Innovation, Sanofi R&D



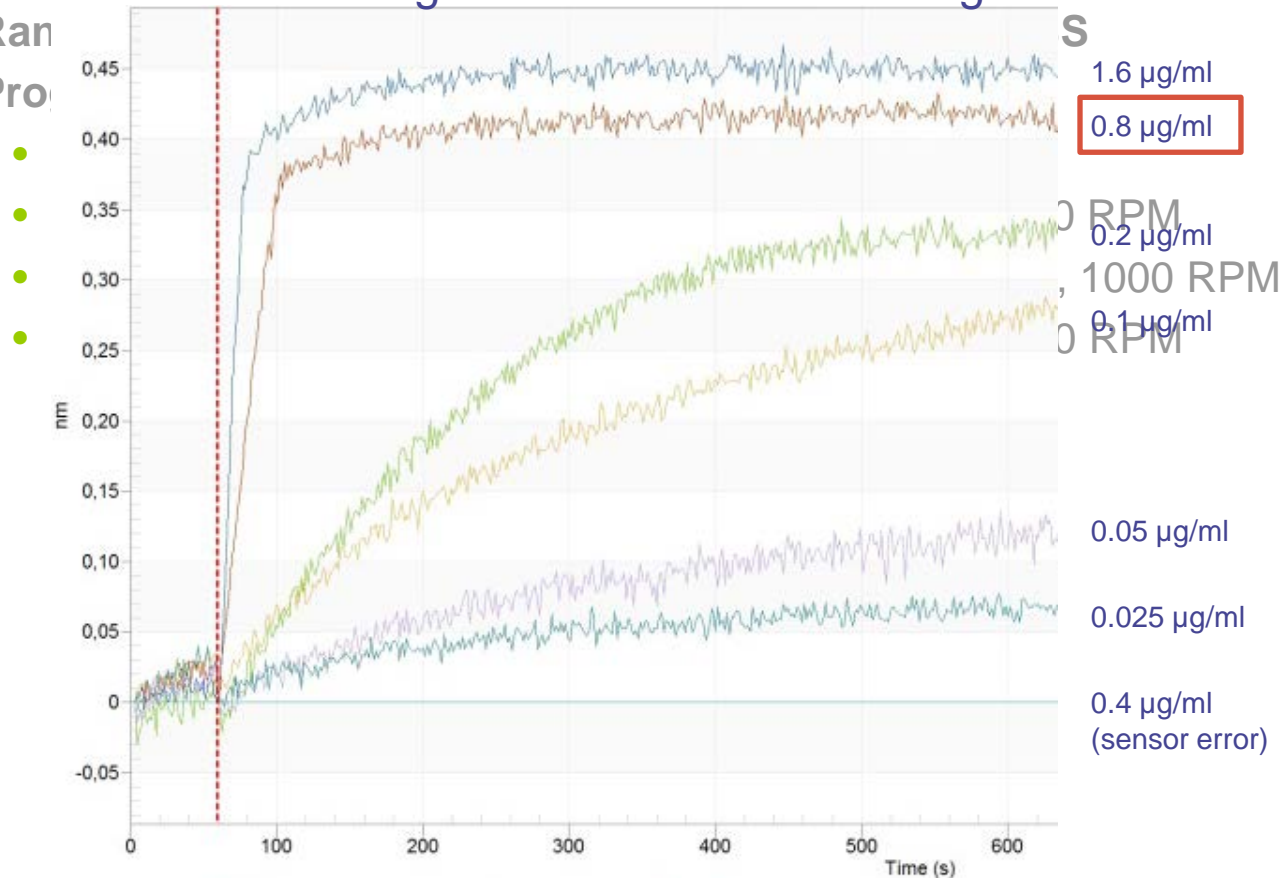
Assay developed by colleagues is functional but Biosensors aren't reusable

Ligand Loading

- Principle:

- Protein concentration was constant: 100 $\mu\text{g/ml}$ in Cell Culture media (CCM)

- Ran
- Pro



● Principle

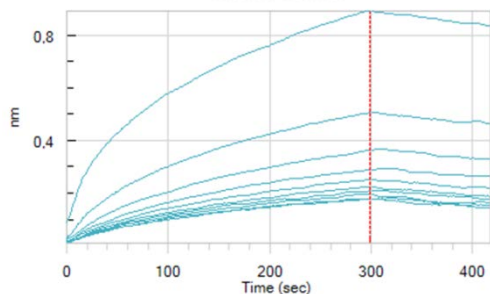
● Test of a large number of buffers:

- Glycine 10 mM pH1.7,
- Glycine 10 mM pH1.7, NaCl 5M
- Glycine 10 mM pH1.7, Tween 20 0.5%
- Glycine 10 mM pH1.7, Triton X100 0.5%
- NaOH 0.01M
- NaOH 0.05M
- NaOH 0.1M
- NaCl 5M
- HCl 0.1M
- H3PO4 0.5M
- DPBS
- DPBS, Triton X100 0.1%
- DPBS, Triton X100 0.5%
- DPBS, EDTA50mM
- DPBS NaCl 5M
- Acetic Acid 0.1%
- Urea 8 M, pH=7
- Urea 4M, pH=7
- Urea 8M in Glycine buffer (pH=3)
- Urea 6M in Glycine buffer (pH=2)
- Urea 8M in NaOH 0.05M (pH=14)
- Urea 6M in NaOH 0.05M (pH=14)
- Urea 8M in Acetic Acid 0.1% (pH=5)
- Urea 6M in Acetic Acid 0.1% (pH=4.5)
- Guanidin 8 M, pH=7
- Guanidin 6M, pH=7
- Guanidin 8M in NaOH 0.05M (pH=14)
- Guanidin 6M in NaOH 0.05M (pH=14)
- Guanidin 8M in NaOH 0.1M (pH=14)
- Guanidin 6M in NaOH 0.1M (pH=14)
- Guanidin 8M in Acetic Acid 0.1% (pH=4)
- Guanidin 6M in Acetic Acid 0.1% (pH=4)

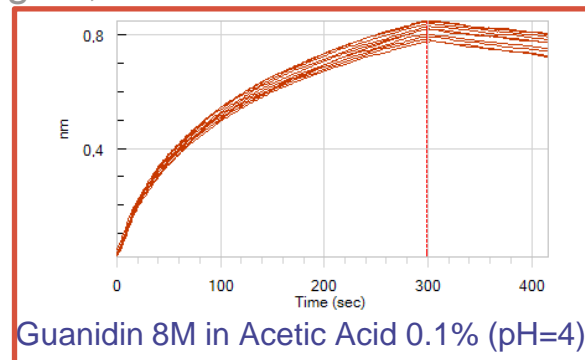
Regeneration Buffer Selection

● Principle:

- Constant protein concentration : 100 µg/ml in CCM
- Constant Biot-Target A: 0.8 µg/ml in PBS
- Program in Basic Kinetic Mode:
 - Baseline in PBS, 60 seconds, 1000 RPM
 - Loading in Biot-Target A in PBS, 120 seconds, 1000 RPM
 - Custom Baseline: Cell Culture media, 60 seconds, 1000 RPM
 - Regeneration: tested buffer, 5 seconds, 400 RPM
 - Neutratisation: Cell Culture Media, 5 seconds, 400 RPM
 - Association: Tetrameric protein solution 100µg/ml, 300 seconds 1000 RPM



Glycine buffer 10mM, pH=1.7



Guanidin 8M in Acetic Acid 0.1% (pH=4)

- Best result is obtained with Guanidin 8M with 0.1% of Acetic Acid

Assay Optimization

- Experiment:

- Read 3 times this plate
- Same program

	Standard	Known concentration samples (µg/ml)							Regeneration			
	1	2	3	4	5	6	7	8	9	10	11	12
A	100.0	5	10	20	30	40	60				Regeneration: Guanidin 8M 0.1% Acetic Acid	Neutralisation: Cell Culture Media
B	50.0	5	10	20	30	40	60					
C	25.0	5	10	20	30	40	60					
H	12.5	5	10	20	30	40	60					
E	6.3	5	10	20	30	40	60					
F	3.1	5	10	20	30	40	60					
G	1.5	5	10	20	30	40	60					
H	0.0	5	10	20	30	40	60					

- Calcul model: R-Equilibrium

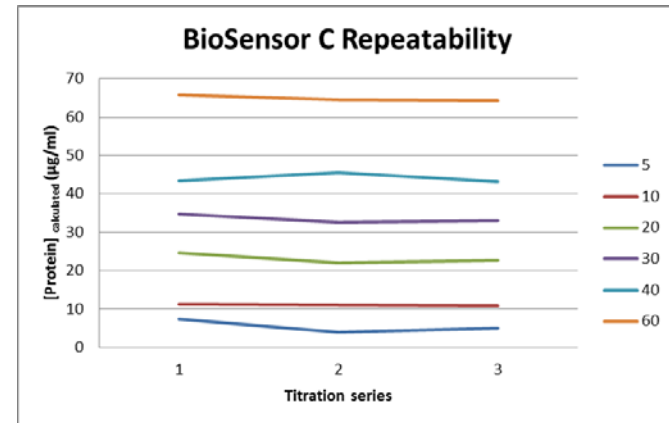
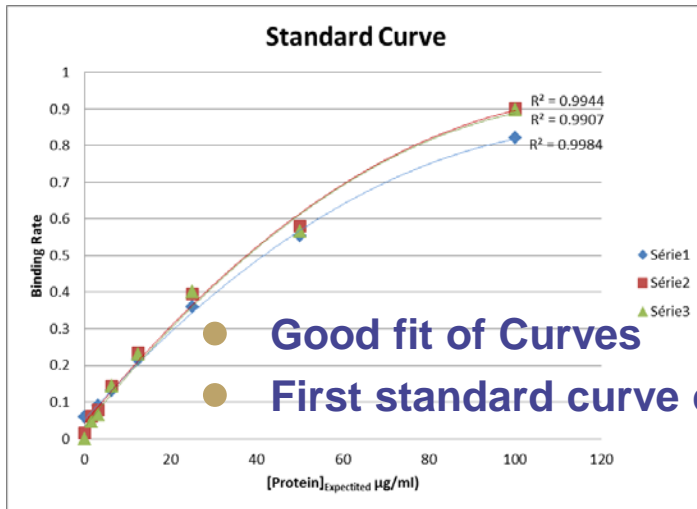
- Initial slope model is not suitable, association too slow

- Integration Time: 150 seconds is a good compromise

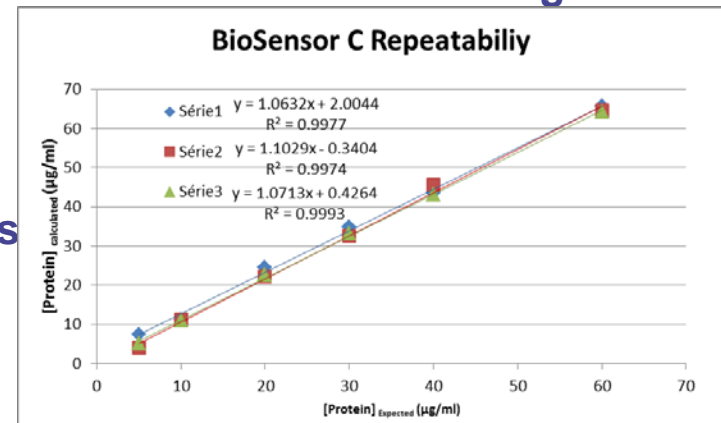
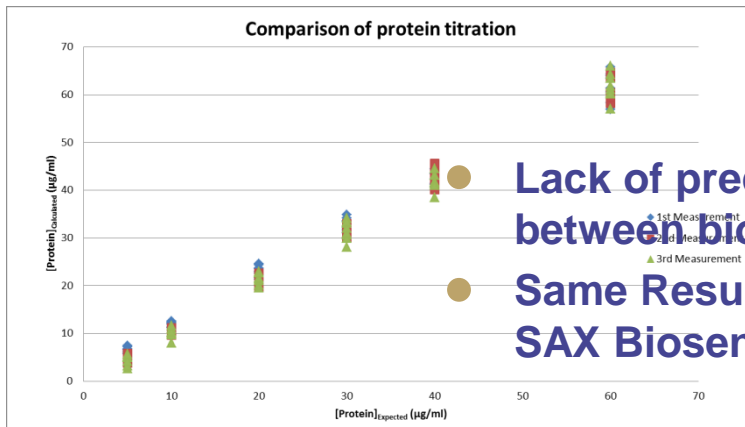
- 100 Sec: Too much variation (VC > 10%) for concentrations <20µg/ml
- 300 Sec: Increase analysis time

- Range between 1.5 to 100 µg/ml:

Result observations



- **Good repeatability of Biosensors**
- **No-decrease between regeneration**



Conclusions

Final assay conditions

Biosensor
SA

Temperature: 30°C

Step 1: Hydration, 5 min in DPBS

Step 2: Loading Target

Biotinyl-PEG4-Target A: 0.8µg/ml in DPBS, 180 sec, 1000RPM

Neutralisation: CCM, 120 sec, 1000 RPM

Step 3: Regeneration/neutralization

Guanidin 8M + Acetic Acid 0.1%, 5 sec, 400 RPM

CCM, 5 sec, 400 RPM

} 3 times

Step 4: Association

Titration: 0-100µg/ml of protein in CCM, Integration time 150 sec, 1000 RPM

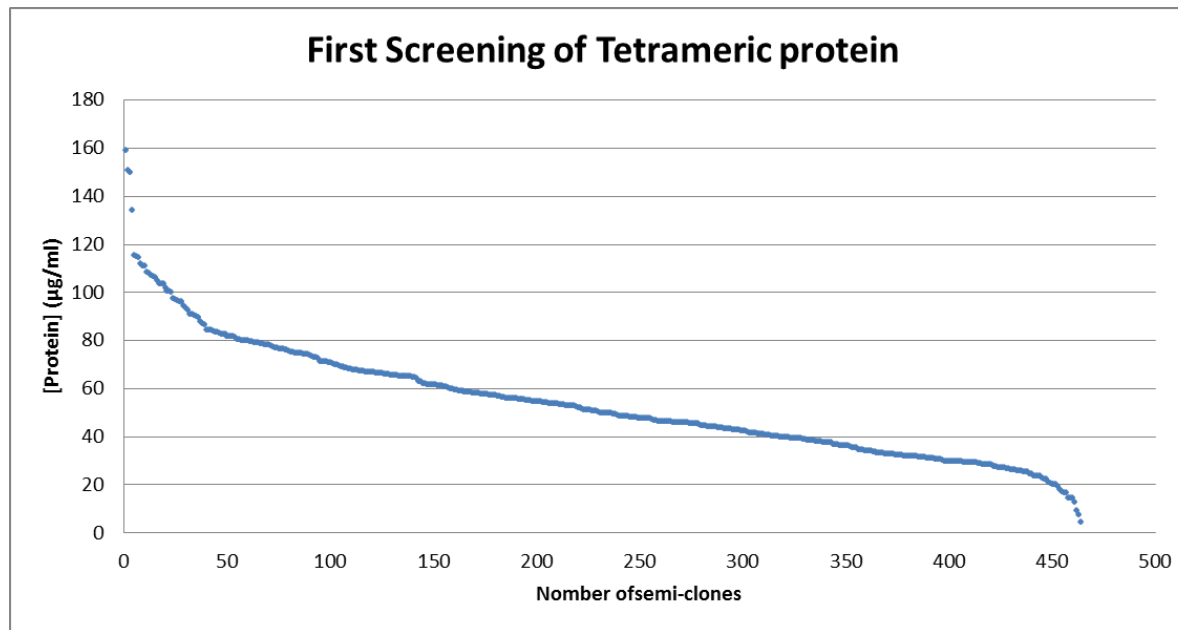
} 10 times

- Limitations of assay:
 - Lake of precision
 - Repeatability
- Advantages of assay
 - Easy to use
 - Fast



Conclusions

- Optimization and validation of method have been beneficial
- Method was efficient to carry out the first screening of the project:
 - Productivity of 450 clones measured at D4 of culture in a day
 - 45 best producers selected for the next step
 - Final selection of cell line is in progress



ACKNOWLEDGEMENT

- Cell Line Engineering (Vitry):
 - Corinne Gillard
 - Séverine Durant
 - Jennifer Di Carlo
 - Laetitia Maçon
- Up Stream Processing (Vitry)
 - Valérie Besset
- Chemist (LGCR, Vitry)
 - Nathalie Karst
- Frankfurt R&D Team
 - Christian Lange
 - Carsten Corvey
 - Ingo Focken
- Pall-FortéBio
 - Arnaud Vonarburg

