

# Host Cell Protein Analysis on the ForteBio Octet Platform

ForteBio Workshop  
Bio-Processing Summit Conference  
August 23

Martha Jackson  
Bioassay and Impurities Testing  
Pfizer, Andover



# Topics

- **Introduction to Host Cell Proteins (HCP)**
- **HCP Analysis**
- **HCP Method**
- **ForteBio HCP Application for the Octet**
- **Results from two Pfizer sites**
- **Conclusions**

# What are Host Cell Proteins?

- **Process related impurities**
  - They are housekeeping proteins that co-express with the protein of choice.
- **Heterogeneous population of proteins that represents the four quadrants of a two dimensional gel**
  - CHO has 10,000's of individual proteins
  - E.coli has 1,000's of individual proteins
- **Host Cell Proteins are not expressed at constant concentrations nor does each species give a similar response in an ELISA**

# Host Cell Proteins Analysis

- **HCP analysis is mentioned in ICH guideline 6B**
- **HCP can be a marker during purification development**
- **HCP can be a marker to monitor batch consistency, column performance and process validation**
- **HCP can be a specification for product quality**

# Method Development

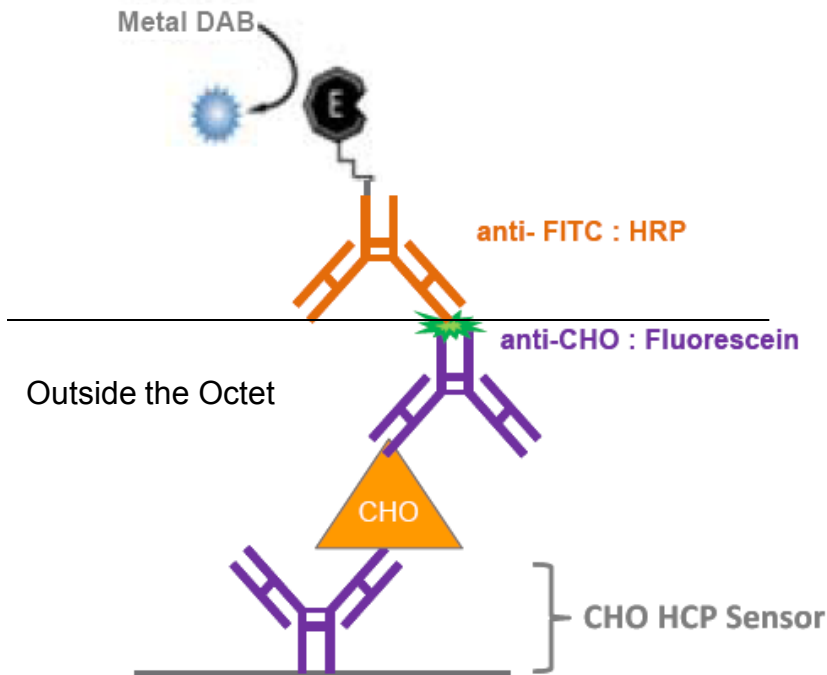
- **Immunological method**
  - Immunogen and standards
    - A set of HCPs collected typically from a null transfected cell line
      - Broad spectrum
      - Enrichment of a subset of HCPs
  - Anti-HCPs
    - Polyclonal
      - IgG cut
      - Affinity Purified against the immunogen or standard
    - Capture reagent
    - Detector reagent
  - Method Output
    - Provide a numeric value for HCP

# Method Comparison

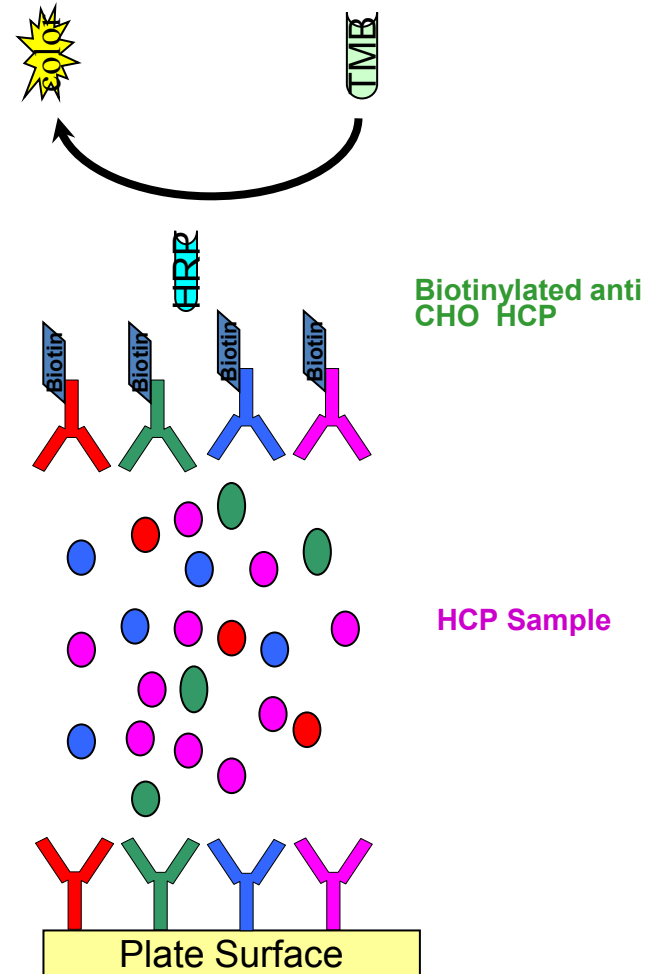
## ForteBio Method

Assay Protocol:

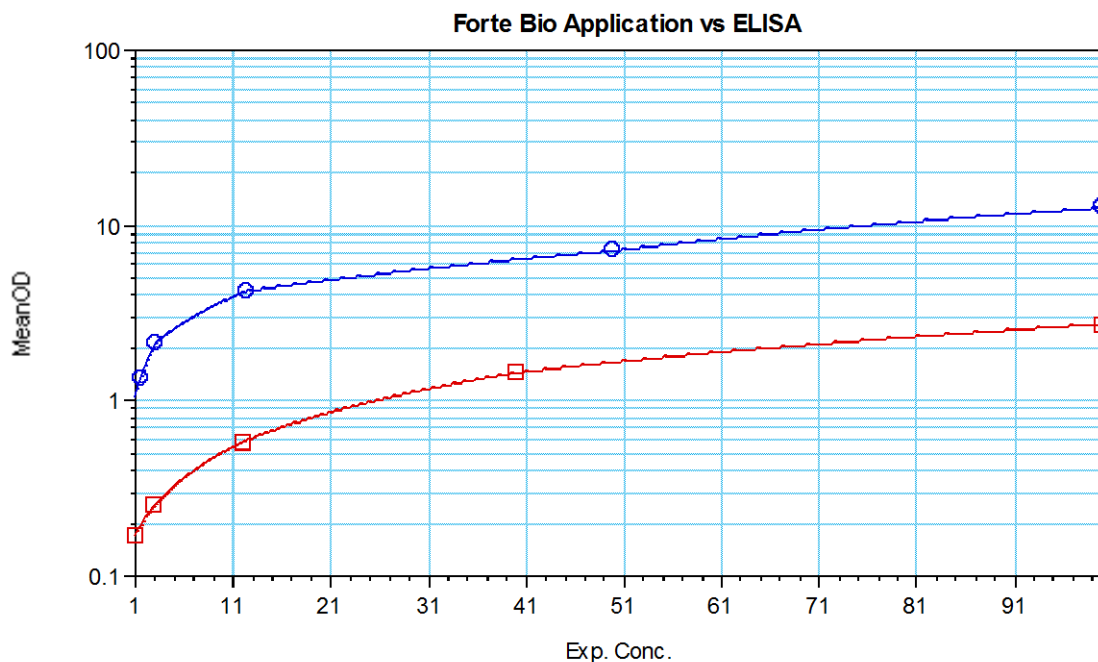
Assay Configuration:



## HCP ELISA



# CHO HCP Standard Curves



Point-To-Point Fit

- Forte Bio (Forte Bio: Conc. vs MeanOD)
- CygnusELISA (Cygnus ELISA Standards: Concentration vs MeanOD)



# Beta Testing

- **Two-Site Blinded Study**
- **Performance Parameters**
  - Specificity
  - Dilutional Linearity
  - Robustness
  - Accuracy and Recovery
  - Sample Analysis



# Dilutional Linearity

- ForteBio HCP standard was spiked into sample diluent at 500 or 200 ng/mL. The samples were tested in a three fold dilution series for a total of eight dilutions

Dilution	Dilution Factor	Site 1 HCP (ng/mL)	Site 1 Adjusted Result (ng/mL)	Site 2 HCP (ng/mL)	Site 2 Adjusted Result (ng/mL)
1	1	173	173	168	168
2	3	137	411	80.2	241
3	9	111	999	22.9	206
4	27	29.3	791	8.16	220
5	81	11.5	931	2.93	238
6	243	4.5	1093	0.772	188
7	729	2.02	1473	0.341	248
8	2187	0.49	1081	BLOQ	BLOQ
Mean			1057		223
SD			257		23.3
CV			24		10



Dilutional Linearity is observed starting at 100ng/mL  
 HCP results are 211% and 112% from expected concentration

## Control Performance

- Intra assay Precision: Cell Culture Sample Tested in a Dilution Series

	Site 1 Ctrl 01 (ng/mL)	Site 1 Ctrl 02 (ng/mL)	Site 1 Ctrl 03 (ng/mL)	Site 1 Ctrl 04 (ng/mL)	Site 2 ELISA (ng/mL)	Forte Bio (ng/mL)	Site 2 ELISA (ng/mL)	Forte Bio (ng/mL)
	252	83.4	35.1	14.7	21275	7199	258	111
	211	96.5	49.9	22.7	15118	7516	208	98.0
		54.1	30.9	9.22	10879	5757	185	84.3
		63.4	40.1	12.2	10191	6382	173	68.2
Mean	232	74.4	39	14.7	12063	6713	206	90.3
SD	28.6	19.2	8.2	5.8	2668	797	37.5	18.2
CV	12.3	25.8	21.0	39.3	22.1	11.9	18.2	20.2
Min	115	65.8	36.3	15.6				
Max	193	104	57	35.8				



Assay Precision is similar to microtiter plate ELISA's

# Intermediate Precision

- Two process samples were analyzed with the ForteBio assay on three different days to determine intermediate precision for the assay.

Sample	Column 1 pool (ng/mL)	Final filter pool (ng/mL)
<b>Exp 1</b>	<b>6713</b>	<b>90.3</b>
<b>Exp 2</b>	<b>9270</b>	<b>96.2</b>
<b>Exp 3</b>	<b>15000</b>	<b>138</b>
Average	10328	108
Std Dev	4243	26.1
%CV	41	24

Intermediate precision for an ELISA is  $\leq 10\%$

# Assay Performance: Recovery and Precision

- HCP spiked into ForteBio sample diluent

	0.1ng HCP/mL Site 2	1 ng HCP/mL Site 2	1 ng HCP/ mL Site 1	5 ng HCP/ mL Site 1	10 ng HCP/mL Site 2	10 ng HCP/ mL Site 1
	0.321	0.982	0.894	4.11	11.5	7.66
	BLOQ	1.05	0.850	5.15	13.1	12
	BLOQ	0.911	0.883	4.13	11.0	13.5
	0.330	0.985	0.841	4.95	10.8	12.8
	BLOQ	1.01	0.761	5.48	12.0	11.3
	BLOQ	0.896	0.749	5.07	10.2	10.7
	BLOQ	0.851	0.894	6.29	10.4	15.1
	0.376	0.835	0.806	3.83	9.72	8.83
Mean		0.94	0.83	4.88	11.1	11.5
SD		0.8	0.06	0.82	1.1	2.43
CV		8.3	6.9	16.8	9.8	21.2
% Recovery		94.0	83.5	97.5	111	115



% Recovery and precision at low concentrations of HCP is acceptable

# Assay Performance: Accuracy

- HCP spiked into 1mg/mL Ab matrix

	1 ng HCP/ mg Ab 1ppm	2 ng HCP/ mg Ab 2ppm	5 ng HCP/ mg Ab 5ppm	10 ng HCP/ mg Ab 10ppm
	0.757	1.84	4.37	11.6
	0.711	1.57	4.17	10.8
	0.809	1.63	6.63	11.2
	0.739	1.64	3.91	10.2
<b>Mean</b>	<b>0.754</b>	<b>1.67</b>	<b>4.77</b>	<b>11.0</b>
<b>SD</b>	<b>0.04</b>	<b>0.12</b>	<b>1.25</b>	<b>0.6</b>
<b>CV</b>	<b>5.5</b>	<b>7.0</b>	<b>26.3</b>	<b>5.5</b>
<b>% Accuracy</b>	<b>75.4</b>	<b>83.5</b>	<b>95.4</b>	<b>109.5</b>



% Accuracy was demonstrated at 1ng HCP/mg Ab matrix (1ppm)

# BioTherapeutic Research and Development

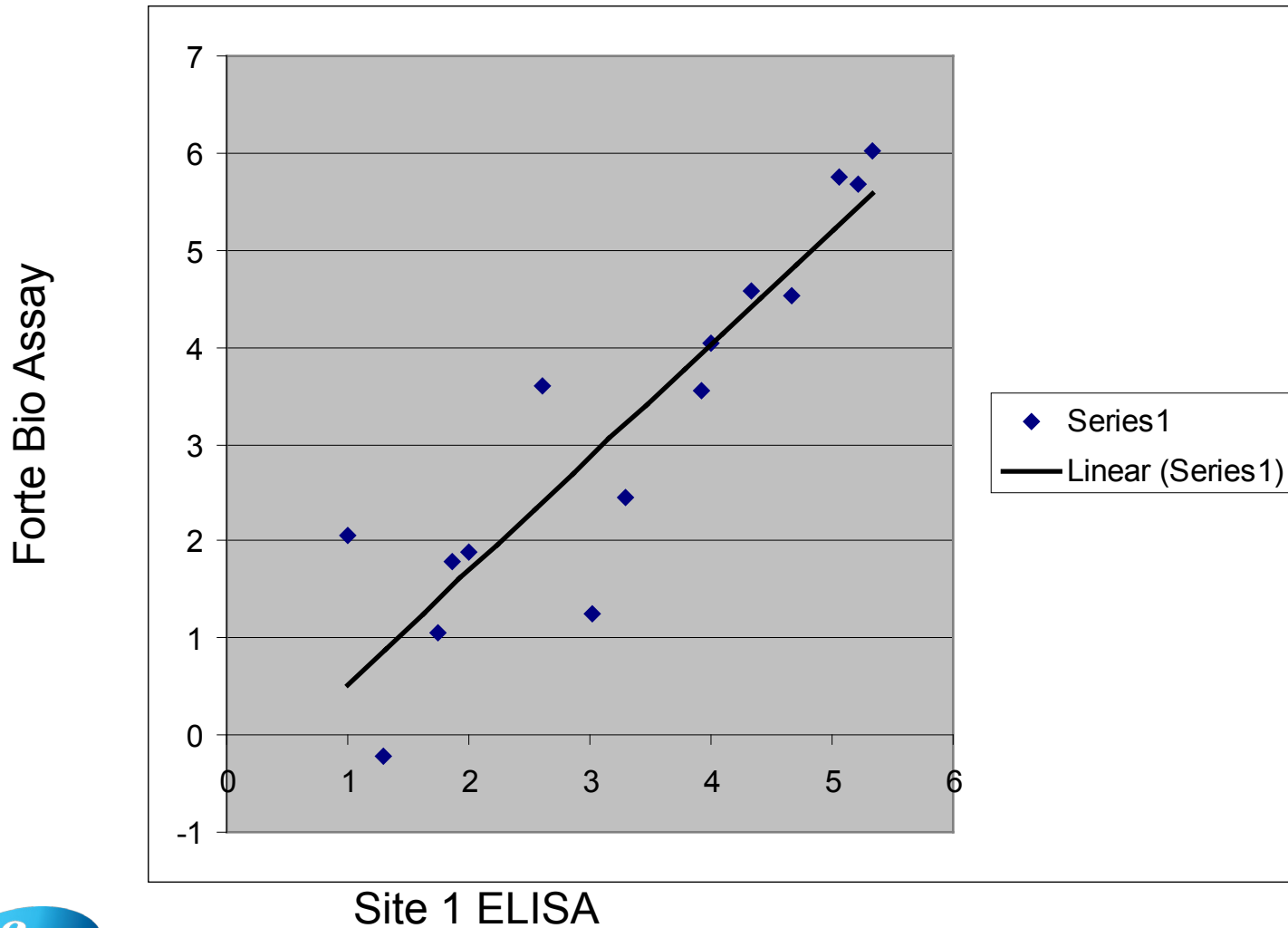
## Sample Performance Site 1

- Three antibodies and two proteins were monitored in the ForteBio HCP Assay

Sample Description	Site 1ELISA Results (ng/mL)	ForteBio Results (ng/mL)
Ab 1 Harvest	114860	562000
Ab 1 Column 1 Load	215209	1048000
Ab 1 Column 2 Load	1046	18.3
Ab 1 Column 2 Peak	<20	0.612
Ab1 Final Pool	<10	114
Protein 1 Column 2 Load	46921	34280
Protein 1 Column 3 Load	21074	37000
Protein 1 Column 3 Peak	8205	3496
Protein 1 Column 3 Wash	100	76
Ab 2 Column 2 Load	72	60
Ab 2 Column 2 Peak	56	11
Ab 2 Column 2 Final Pool	10046	10900
Ab 3 Final Pool	414	4000
Protein 2 Column 1 Load	166411	472500
Protein 2 Column 2 Load	2013	280
	<b>R<sup>2</sup></b>	<b>0.96</b>
	<b>Slope</b>	<b>4.27</b>
	<b>Intercept</b>	<b>-21992</b>



## HCP results of the 15 samples plotted on a log, log scale



## Sample Performance

- ForteBio HCP Analysis Across Purification Steps

<b>Protein</b>				
<b>Step</b>	<b>Batch 1 ng/mL</b>	<b>Fold Reduction</b>	<b>Batch 2 ng/mL</b>	<b>Fold Reduction</b>
<b>Cell Culture</b>	<b>525000</b>		<b>1739500</b>	
<b>Column 1 Load</b>	<b>320400</b>		<b>1583100</b>	
<b>Column 1 Peak</b>	<b>8.43</b>	<b>38007</b>	<b>49.5</b>	<b>31981</b>
<b>Column 2 Peak</b>	<b>11.4</b>	<b>0.74</b>	<b>302.4</b>	<b>0.16</b>
<b>Antibody</b>				
<b>Column 1 Load</b>	<b>305500</b>			
<b>Column 1 Peak</b>	<b>129500</b>	<b>2.4</b>		
<b>Column 2 Peak</b>	<b>1918</b>	<b>67.5</b>		
<b>Final Pool</b>	<b>2639</b>	<b>0.7</b>		



## Comparison of ForteBio and ELISA for HCP Clearance Site 2

	CHO HCP ForteBio Assay		PS CHO HCP ELISA	
	HCP (ng/mg)	Fold Removal	HCP (ng/mg)	Fold Removal
<b>mAb 1</b>				
Harvest	68163		95476	
Column 1 Pool	744	92	599	159
Column 2 Pool	96.2	7.7	61.8	10
Column 3 Pool	21.6	4.5	47.9	1.3
Final Filter Pool	19.7	1.1	29.3	1.6
<b>mAb 2</b>				
Column 1 Pool	300		169	
Column 2 Load	40.2	7.4	18.4	9.2
Column 3 Load	32.9	1.2	14.6	1.3
Final Filter Pool	1.0	33	1.1	14
<b>mAb 3</b>				
Column 1 Load	258,611		229,326	
Column 2 Load	475	544	287	800
Column 3 Load	109	4.4	33.5	8.6
Column 3 Eluate	10.1	11	10.2	3.3

# Conclusions

- **The ForteBio HCP assay range is 1 to 100 ng HCP/mL**
- **Dilutional Linearity is observed from 0.4 to 100 ng HCP/mL**
- **Recovery was demonstrated at 1ng HCP/mL**
- **Several of the assay parameters are equivalent to established HCP ELISA**
- **The specificity of the reagents demonstrated adequate HCP removal across purification processes**
- **This application would be applicable for measuring HCP in down stream samples**
- **The ForteBio HCP application offers a sample turn around time of 90 minutes for 7 samples tested as duplicates or 14 samples tested in single wells**

# Thanks

- **ForteBio**
  - John Proctor
- **Pfizer**
  - Phoebe Baldus
  - Keith Davis
  - Laura Bass
  - Ned Mozier
  - William Flaherty
  - Martha Jackson
  - Mark Hardy

