

# Application of the Octet Red96 system to characterize protein interactions

Alan Wong  
(James Rini Lab)

ForteBio Workshop and User Meeting,  
Toronto  
Dec 11, 2014

# Protein interaction analysis

Proteins are fundamental molecules in biology

Interactions are essential for protein functions

Protein-protein  
(Ab-Ag), (ligand-receptor)

Protein-carbohydrate  
(lectin-sialic acid)

Protein-lipid  
(protein-phospholipid)

## **Early stage characterization**

Rapid, easy to use

Relatively crude preparation

Qualitative

Rough ranking

## **Later stage characterization**

Rapid, easy to use

Purified samples

Quantitative

Publishable results

# Octet Red96 biolayer interferometry



# Rapid qualitative antigen-antibody interaction characterization

1. Hybridoma produced IgG
2. Stably expressed IgG of phage display Ab

IgGs are secreted in tissue culture media

For early stages:

Use crude tissue culture media containing IgG

Quantify IgG production?

Does the IgG bind the antigen?

# Rapid qualitative antigen-antibody interaction characterization

IgG quantification (reproducible test)

Ab: Human phage display IgG1

IgG produced from serum free media (no purification)

Probe: protein A (ProA) biosensor

Well: IgG containing serum free media

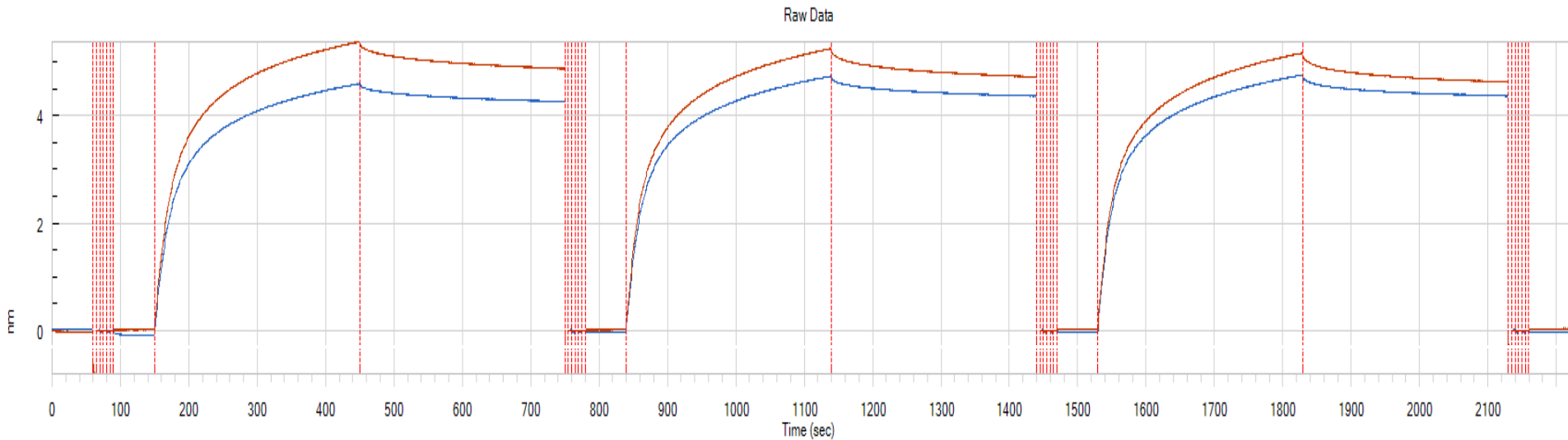
# Repeated measurement of IgG using proA probe after subsequent acid washes (100mM citric acid pH3)

	1	2	3	4	Probe
A	Hyb Med	IgGZ	100mM citric acid pH3	100mM HEPES pH7.5	proA
B	..	..	..	..	..

Baseline

Regenerate

Neutralization



Both lanes generated curves of similar level

Proceed with further characterization

# Rapid qualitative antigen-antibody interaction characterization

## Hybridoma produced IgG-Ag interaction

Ab: mouse IgG1, mouse IgG2

IgG produced from serum free hybridoma media (no purification)

Ag: untagged purified soluble glycoproteins

gel filtration purified (>95%) antigens expressed in HEK293 cells

biotin label Ag on primary amines

Probe: streptavidin (SA) biosensor (externally loaded biotin-Ag)

Well: crude hybridoma media

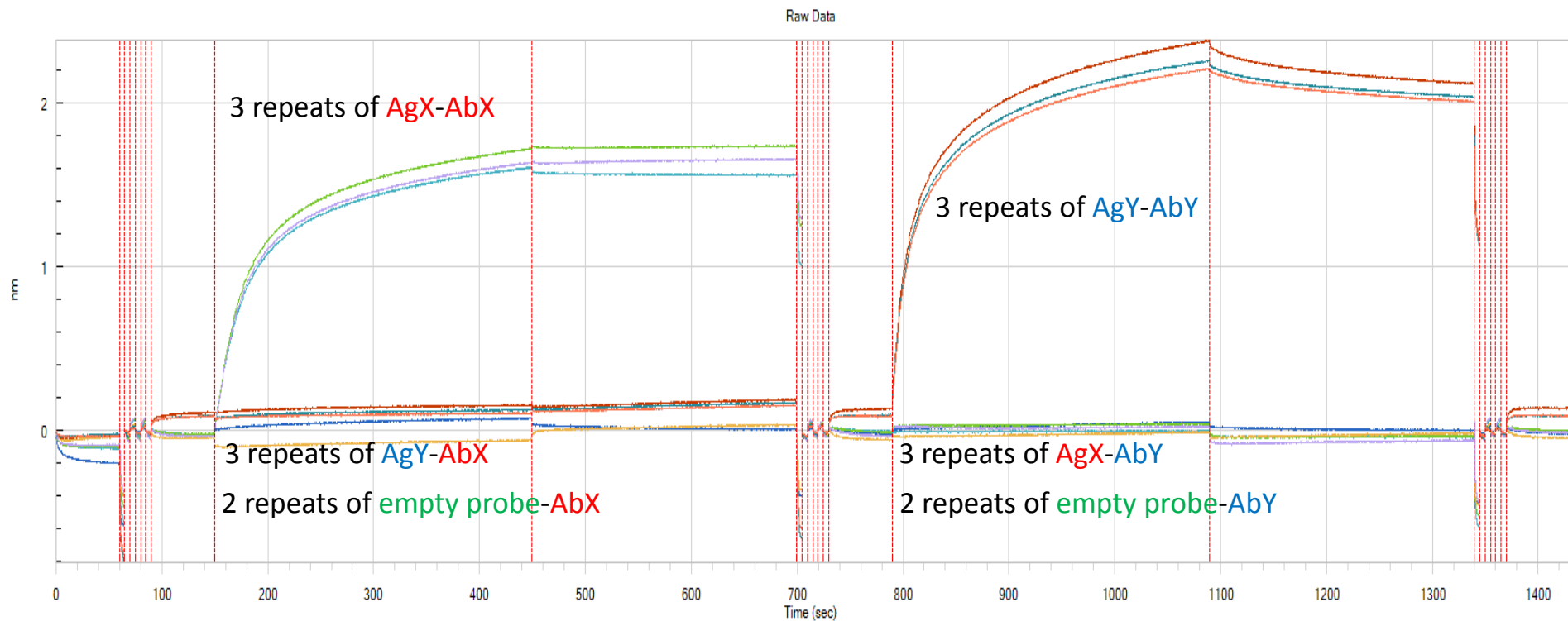


	1	2	3	4	5	Probe
A	Hyb Med	IgGX	IgGY	50mM Acetic acid pH 4.5	100mM HEPES pH7.5	SA: Empty
B	..	..	..	..	..	SA: AgX
C	..	..	..	..	..	SA: AgY
D	..	..	..	..	..	SA: AgY
E	..	..	..	..	..	SA: Empty
F	..	..	..	..	..	SA: AgY
G	..	..	..	..	..	SA: AgX
H	..	..	..	..	..	SA: AgX

Baseline

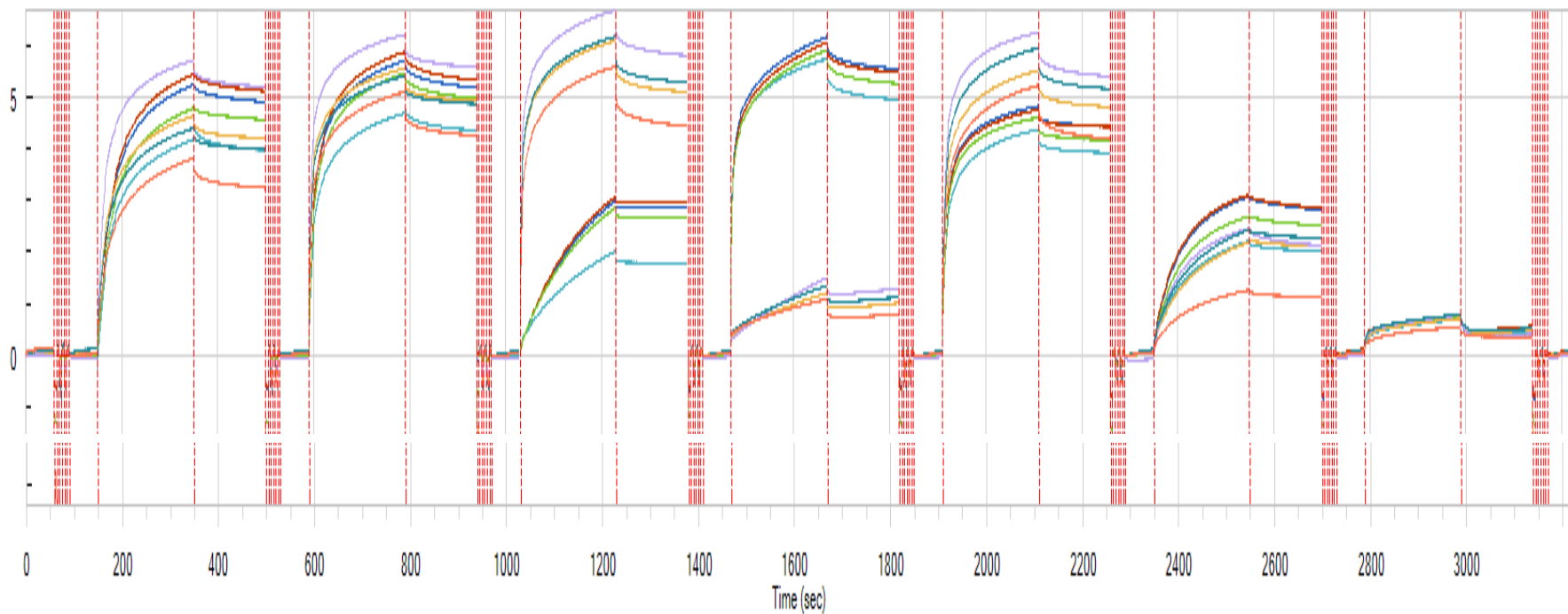
Regeneration

Neutralization



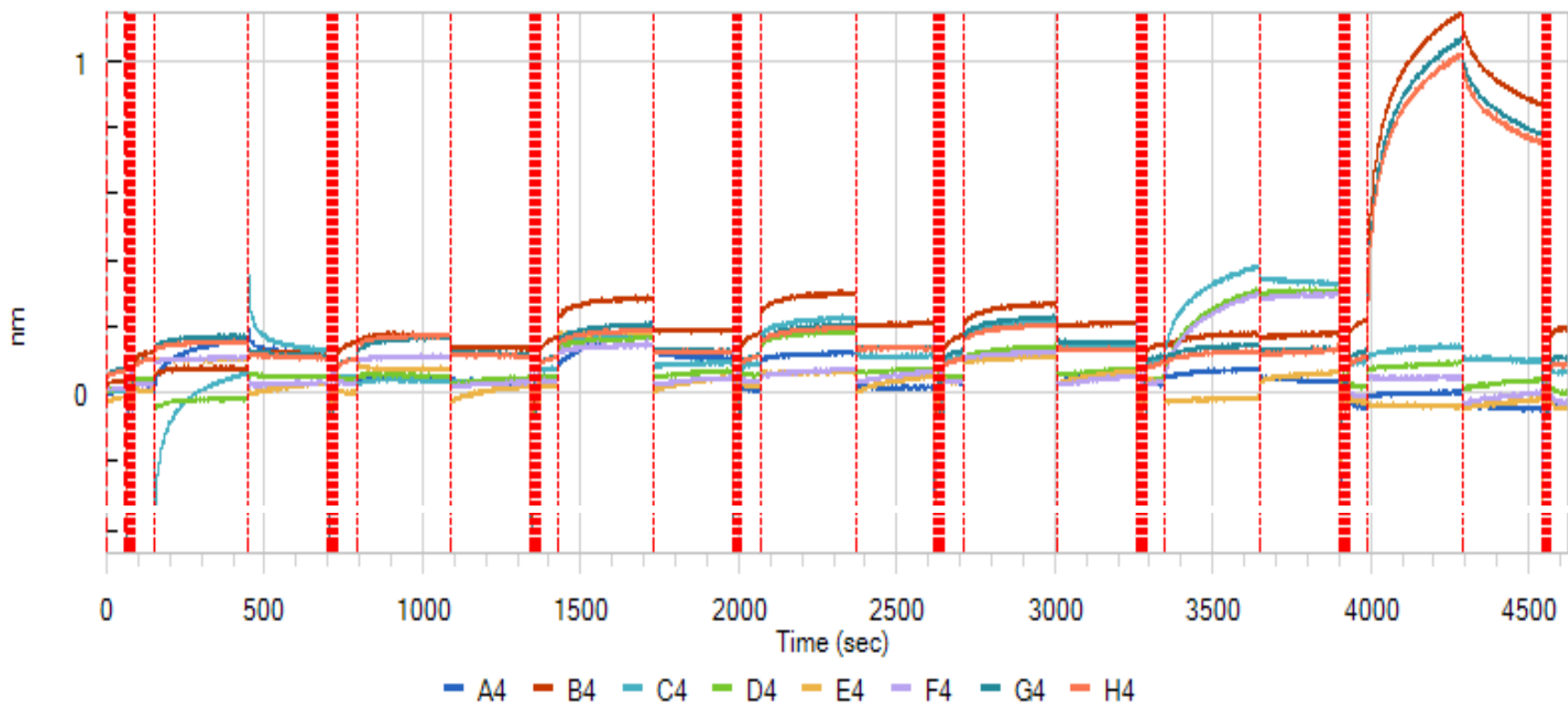
	1	2	3	4	5	6	7	8	9	10	Probe
A	100mM Citric Acid pH3	100mM HEPES pH7.5	SFM	IgGA	IgGD	IgGF	IgGH	IgGJ	IgGL	IgGM	proA
B	..	..	..	..	..	..	..	..	..	..	..
C	..	..	..	..	..	..	..	..	..	..	..
D	..	..	..	..	..	..	..	..	..	..	..
E	..	..	..	IgGB	IgGE	IgGG	IgGI	IgGK	..	..	..
F	..	..	..	..	..	..	..	..	..	..	..
G	..	..	..	..	..	..	..	..	..	..	..
H	..	..	..	..	..	..	..	..	..	..	..

Raw Data



	1	2	3	4	5	6	7	8	9	10	Probe
A	100mM Acetic Acid pH4.5	100mM HEPES pH7.5	SFM	IgGA	IgGD	IgGF	IgGH	IgGJ	IgGL	IgGM	SA: Empty
B	..	..	..	..	..	..	..	..	..	..	SA: AgX
C	..	..	..	..	..	..	..	..	..	..	SA: AgY
D	..	..	..	..	..	..	..	..	..	..	SA: AgY
E	..	..	..	IgGB	IgGE	IgGG	IgGI	IgGK	..	..	SA: Empty
F	..	..	..	..	..	..	..	..	..	..	SA: AgY
G	..	..	..	..	..	..	..	..	..	..	SA: AgX
H	..	..	..	..	..	..	..	..	..	..	SA: AgX

Raw Data (Sensor Location)



# Quantitative protein-protein interaction characterization

## Protein-protein interaction

Can we obtain binding constants of protein-protein interaction using purified proteins?

# Quantitative protein-protein interaction characterization

## Mammalian cell line produced protein-protein interaction

Protein 1: untagged purified soluble glycoprotein  
gel filtration purified (>95%) proteins expressed in HEK293 cells

Protein 2: untagged purified soluble glycoproteins  
gel filtration purified (>95%) proteins expressed in HEK293 cells  
biotin label Ag on primary amines

Probe: streptavidin (SA) biosensor (externally loaded biotin-protein 2)

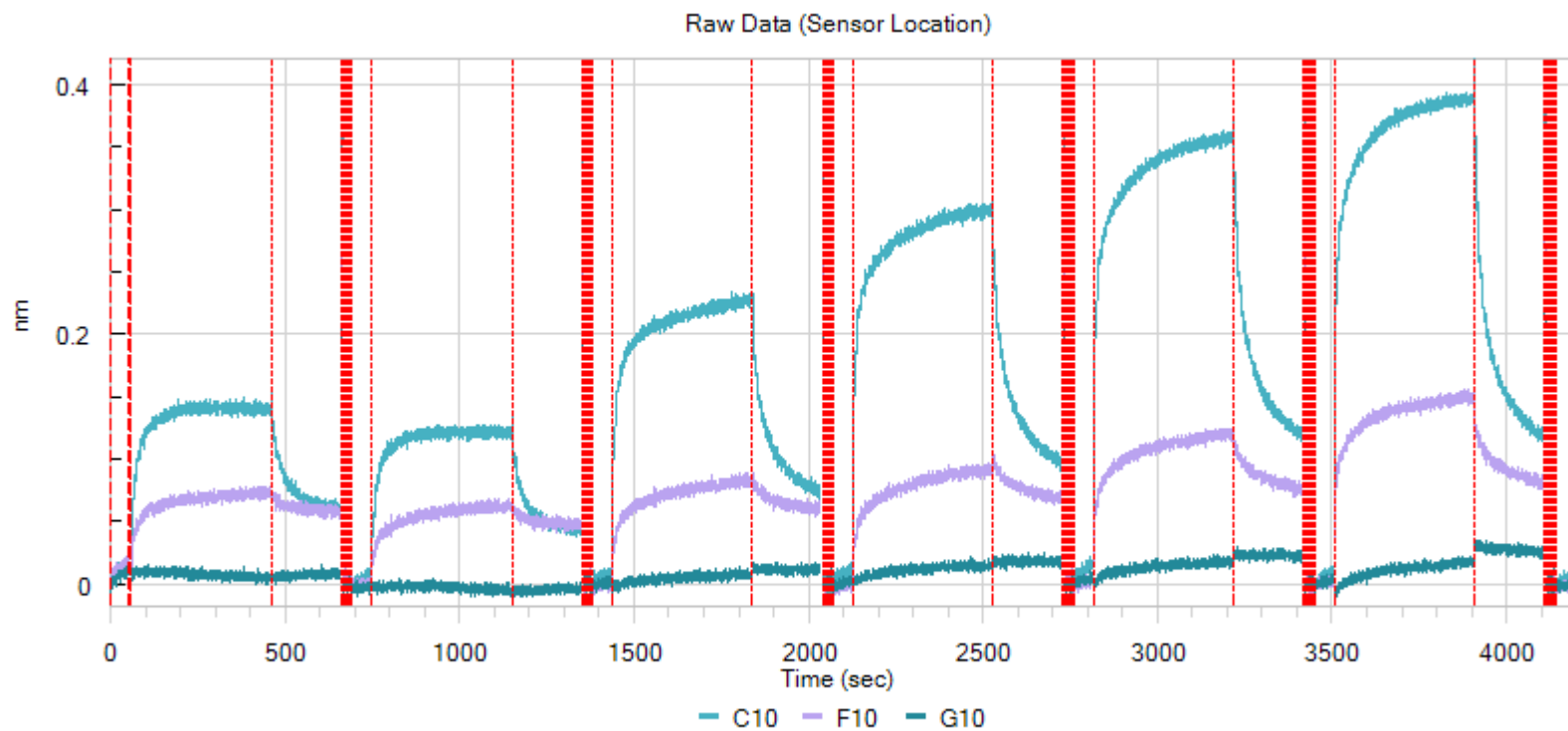
Negative control: Biotin-non-interacting protein 3

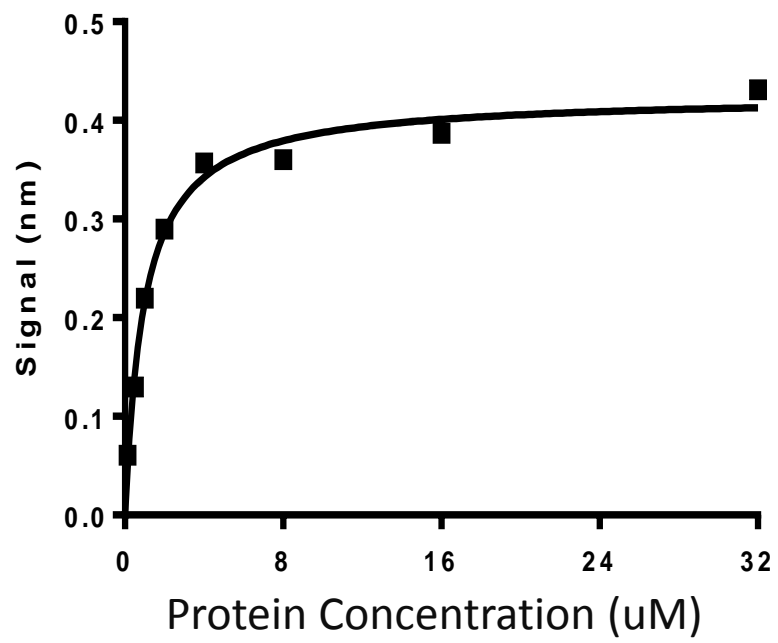
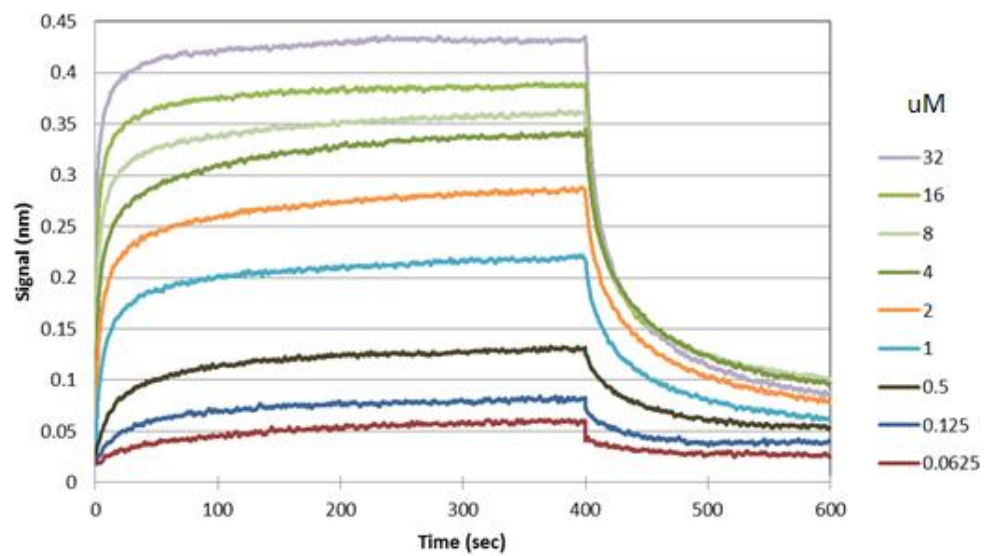
Well: titration of purified protein 1

	1	2	3	4	5	6	7	8	9	10	11	12	Probe
C	50mM Citric acid pH 4	100mM HEPES pH7.5	100mM HEPES 100mM NaCl pH7.5	0.5uM protein 1	100mM HEPES 100mM NaCl pH7.5	1uM protein 1	100mM HEPES 100mM NaCl pH7.5	2uM protein 1	100mM HEPES 100mM NaCl pH7.5	4uM protein 1	100mM HEPES 100mM NaCl pH7.5	8uM protein 1	SA: Protein 2
F	..	..	..	..	..	..	..	..	..	..	..	..	SA: Non- interacting protein 3
G	..	..	..	..	..	..	..	..	..	..	..	..	SA: Empty

Regeneration

Neutralization





$K_d = 0.98 \text{ uM}$

# Quantitative lectin-carbohydrate interaction characterization

## Lectin-carbohydrate interaction

Can we obtain binding constants of describing protein-carbohydrate interaction using purified lectins?



# Quantitative lectin-carbohydrate interaction characterization

## Lectin-sialic acid interaction

Lectin: sialic acid binding soluble glycoprotein

Gel filtration purified (>95%) lectin expressed in HEK293 GnT1<sup>-</sup> cells  
(no sialic acid on the lectin)

Carbohydrate: soluble sialic acid containing glycoproteins

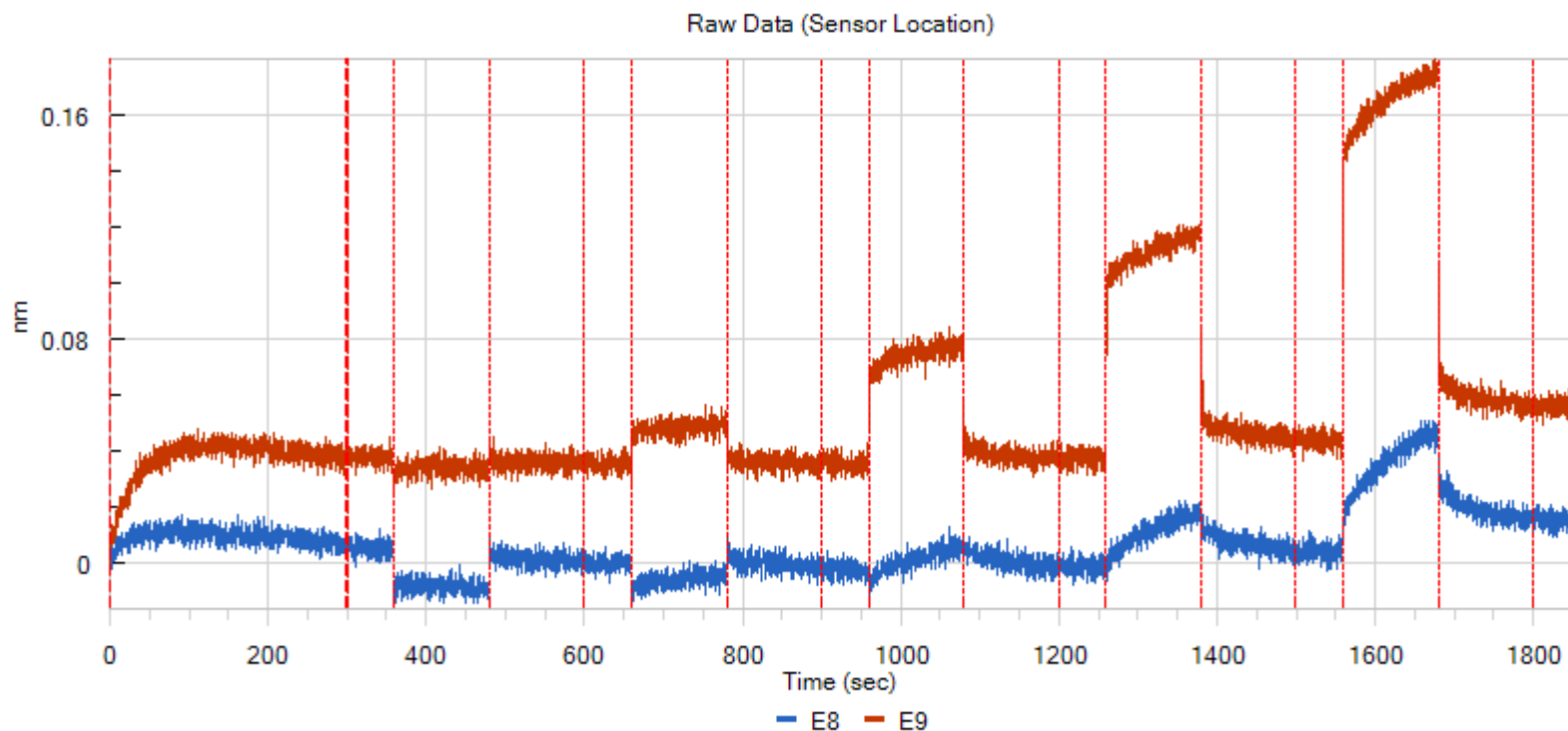
Biotin label sialic acid containing glycoproteins on primary amines

Probe: streptavidin (SA) biosensor (externally loaded biotin- sialic acid containing glycoprotein)

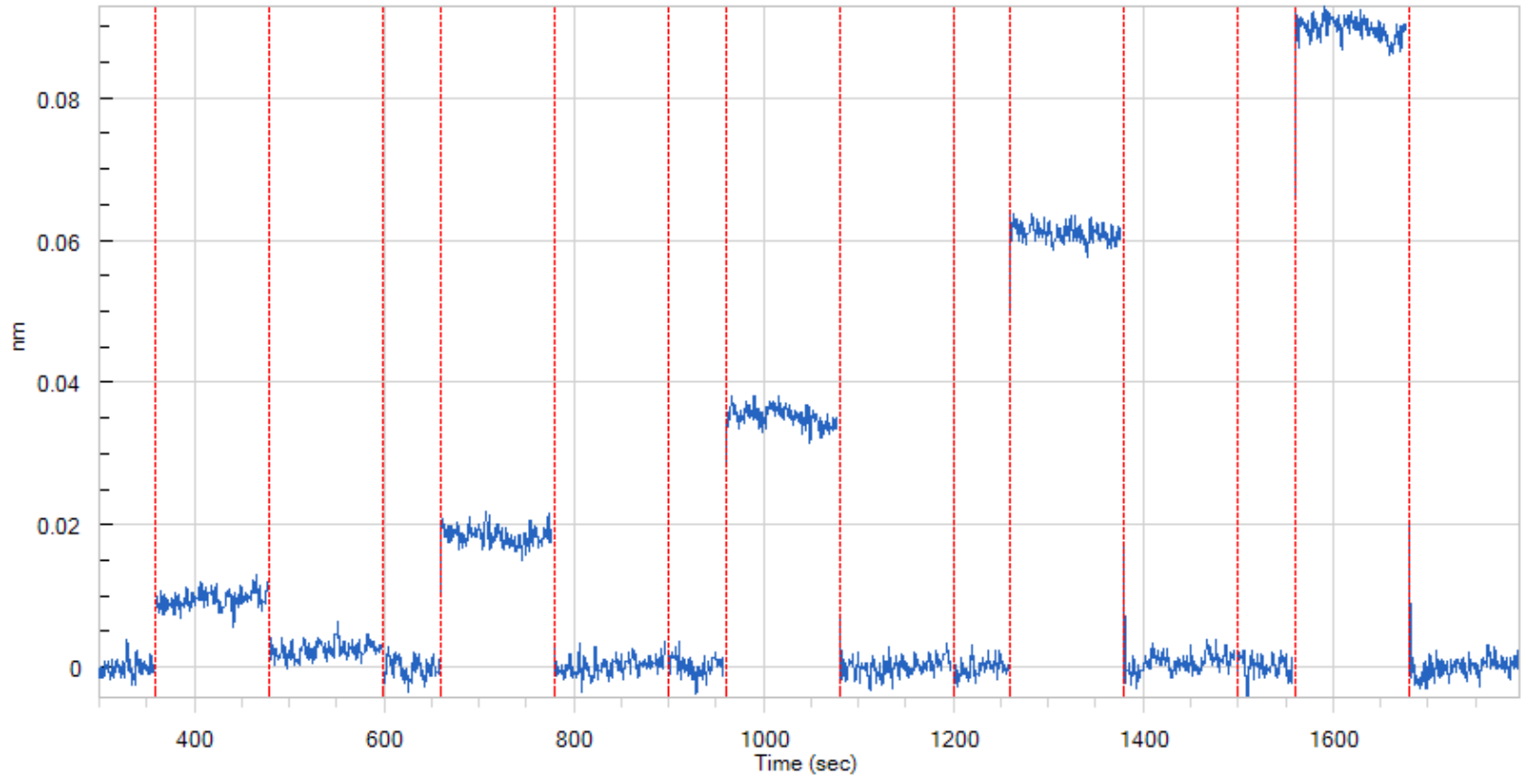
Negative control: probe incubated with neuraminidase (external)

Well: titration of lectin

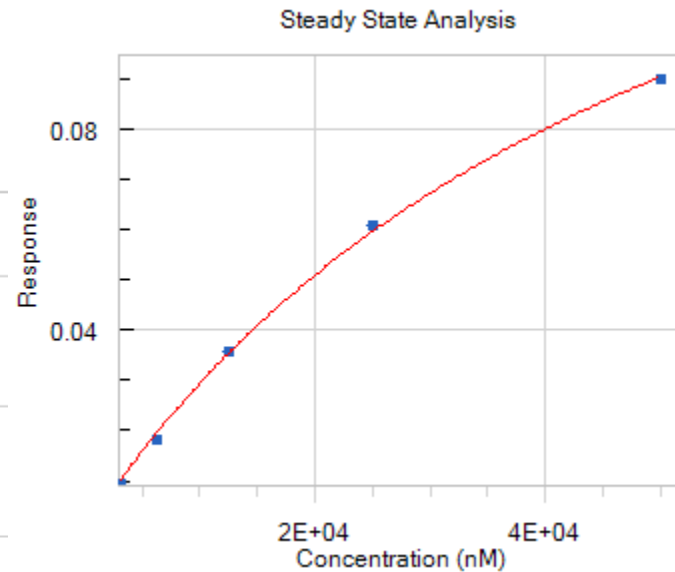
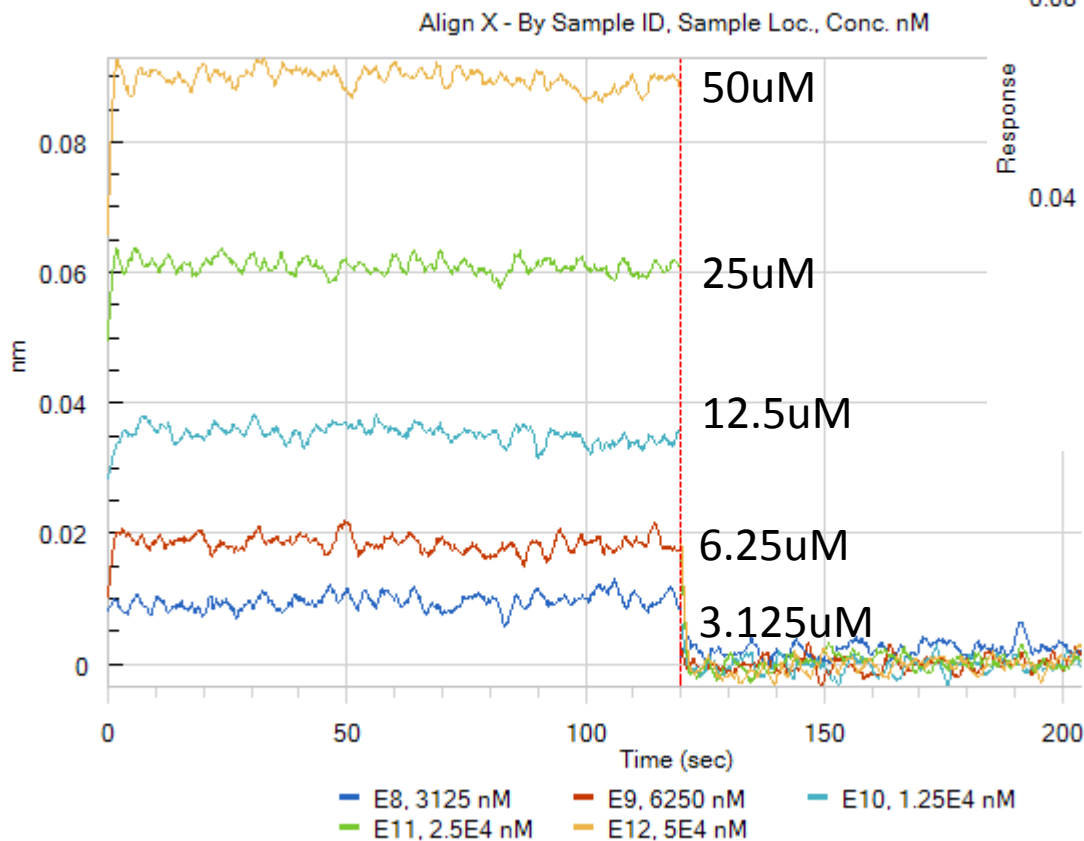
	1	2	3	4	5	6	7	8	9	10	11	12	Probe
E9	100mM HEPES 100mM NaCl pH7.5	3.125uM Lectin	100mM HEPES 100mM NaCl pH7.5	6.25uM Lectin	100mM HEPES 100mM NaCl pH7.5	3.125uM Lectin	100mM HEPES 100mM NaCl pH7.5	12.5uM Lectin	100mM HEPES 100mM NaCl pH7.5	25uM Lectin	100mM HEPES 100mM NaCl pH7.5	50uM Lectin	SA: Glycoprotein
E8	..	..	..	..	..	..	..	..	..	..	..	..	SA: Neuraminidase treated glycoprotein



Align Y



Take response value at 20 to 100 sec



Data Used	Response	
Y Weighting		
Chi <sup>2</sup> /DoF	1.47857E-06	
R <sup>2</sup>	0.998973223	
RMax	0.187001208	0.007974
KD	5.30E-05 ±3.7E-06M	

KD from BLI = 53uM

KD from SPR = 45 uM

# Qualitative protein-lipid interaction characterization

## Protein-lipid interaction

Can we obtain detect binding of protein with phospholipid in a liposome?

# Qualitative protein-lipid interaction characterization

## Protein-liposome interaction

Protein: Phospholipid binding protein, multiple lipid binding sites  
Gel filtration purified (>95%) protein expressed in E.coli

Liposome: biotin labeled liposome

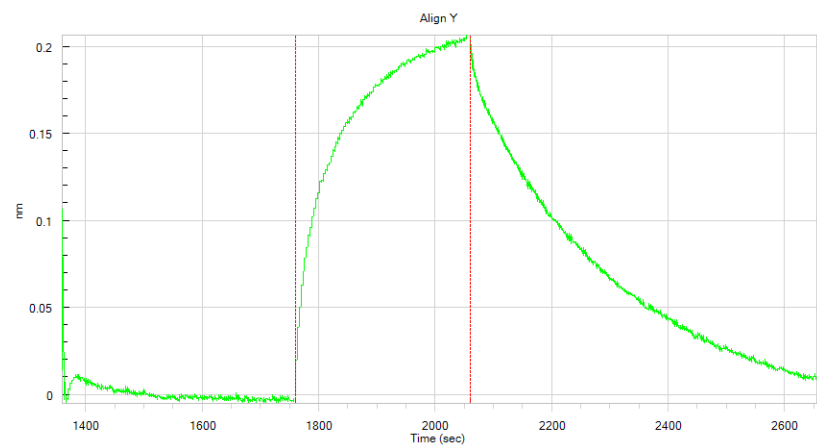
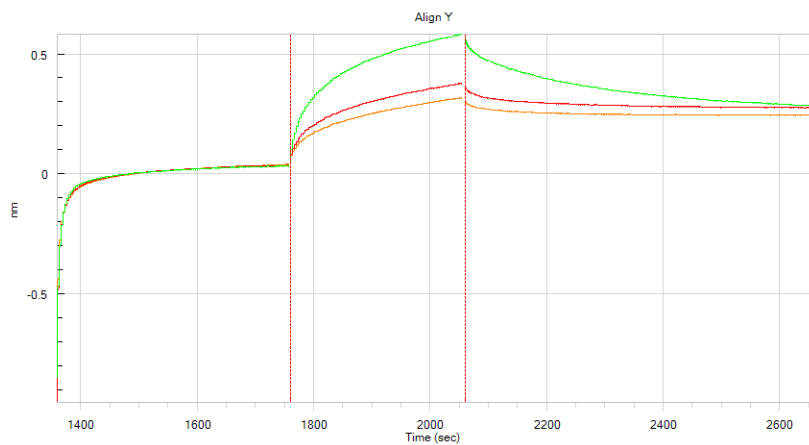
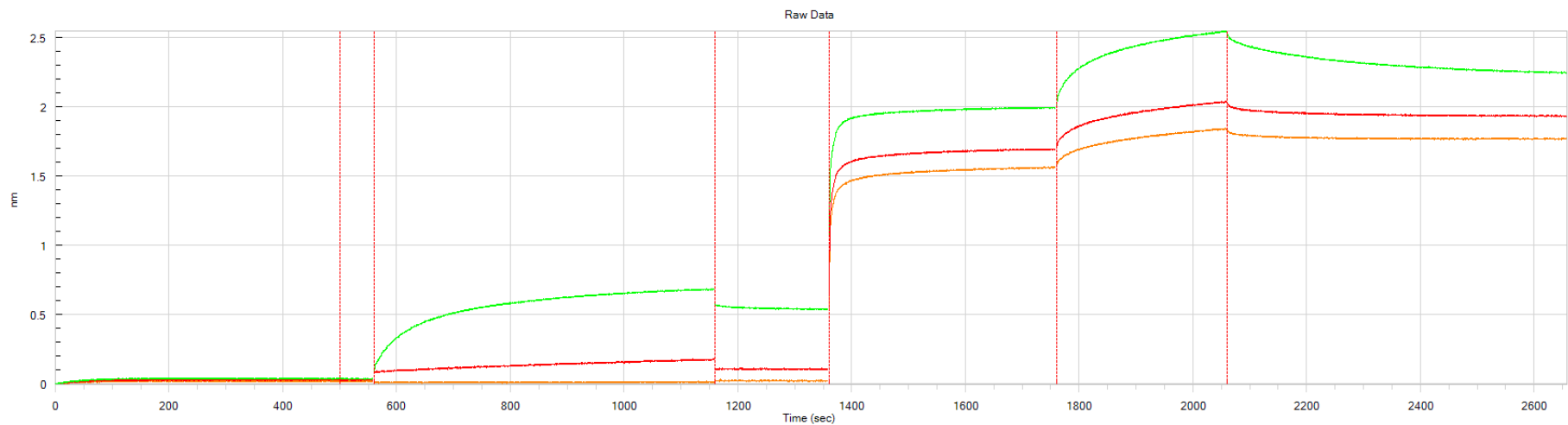
Probe: streptavidin (SA) biosensor

Load: liposome with biotin labeled PE

Negative control: liposome with normal PE

Well: titration of phospholipid binding protein

	1	2	3	4	Probe
A	100mM HEPES 100mM NaCl pH7.5	67%PC / 30%PS / 3% Biotin-PE	Skim Milk	2uM Protein in skim milk	SA
B	..	67%PC / 30%PS / 3%PE	..	..	..
C	..	100mM HEPES 100mM NaCl pH7.5	..	..	..



# Conclusion

Octet RED 96 system can be used for:

Detection of Ab production and rapid identification of Ag binding at early stage  
Alternatives: Anti-Mouse IgG quantification and anti-mouse Fc capture probes

Determine protein-protein and protein-carbohydrate binding properties

Possible to characterize protein-lipid interaction  
Alternatives: Aminopropylsilane probes



# Acknowledgments

## Supervisor

James Rini

## Rini lab

Kristina Han  
Xuyao Li  
Zhijie Li  
Malathy Satkunarajah  
Dongxia Zhou

## BLI

Trevor Moraes

## Liposome

Sharon Eisenberg  
Sergio Grinstein

