



Assessment of Monoclonal Antibody/Fc Receptor Interactions Using Octet:

Establishment of a Toolbox Panel for Characterization of Therapeutic mAbs and mAb Biosimilars

Michael Sadick, PhD
Senior Manager, Large Molecule Analytical Chemistry
Catalent Pharma Solutions, Kansas City, MO



DEVELOPMENT



DELIVERY



SUPPLY

more products. better treatments. reliably supplied.™

Introduction

- **Assessment of biopharmaceutical therapeutic molecules, or their associated derivative biosimilar molecules, requires an orthogonal analytical approach.**
 - No one particular analytical technique can fully define the physical and biological characteristics.
 - Combination of physicochemical and biological analyses
 - Physicochemical analyses
 - Mass spectrometry,
 - HPLC (SEC, IEX, HIC, etc),
 - SDS-electrophoresis (gel or capillary)
 - Isoelectric focusing (gel or capillary)
 - Glycoanalyses
 - Biological analyses
 - Cell-based bioassay
 - ELISA
 - Binding assays
 - Surface plasmon resonance
 - Bio-layer interferometry).
 - Complete and integrated understanding of the biotherapeutic molecule's form and function.

Antibody Binding Properties and Functions

- **IgG monoclonal antibody (mAb) biotherapeutics,**
 - Assays needed both for MOA characterization/definition as well as release (CMC)
 - Multiple functions for one molecule
 - Interactions between “fab” complementarity determining region (CDR) and the target
 - Specific biological function (neutralization, inhibition, etc)
 - Interactions between the Fc region with all of the potential human IgG-Fc-binding cell-surface receptors.
 - Low affinity CD16a and CD16b receptors (FcγRIIIA and FcγRIIIB, respectively),
 - Responsible for the majority of effector function — antibody-dependent cellular cytotoxicity, or ADCC — associated with mAb therapeutics (chiefly IgG1)
 - Low affinity receptors, CD32a and CD32b/c (FcγRIIA and FcγRIIB/C, respectively)
 - Associated with antibody-dependent cellular phagocytosis, or ADCP
 - High affinity receptor CD64 (FcγRI)
 - Responsible for mediating multiple immune responses .
 - ‘Neonatal’ FcRN,
 - Great impact on the in vivo half-life of a therapeutic mAb .

Analyses of mAb Structure and Function

- **There are many alterations to the mAb that can impact its activities**
 - Amino acid substitution(s)
 - Oxidation (met)
 - Glycosylation
 - Etc
- **Biological impact of the changes may be assessed with functional assays**
 - Loss of inhibition (bioassay or ELISA)
 - Loss of effector function (bioassay)
- **Exact changes may be identified via physicochemical means**
- **However, measurements of binding, either CDR/Target and/or Fc/FcR provide crucial understanding of the underlying changes or consistency of mAb biotherapeutics**

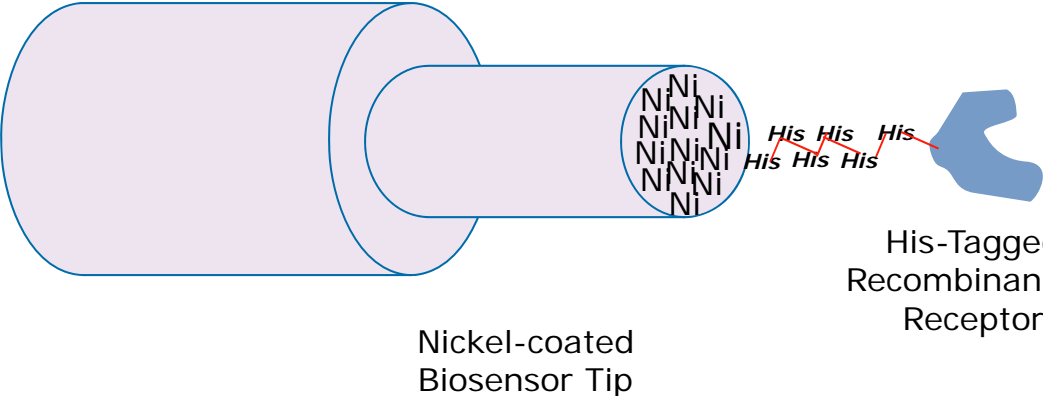
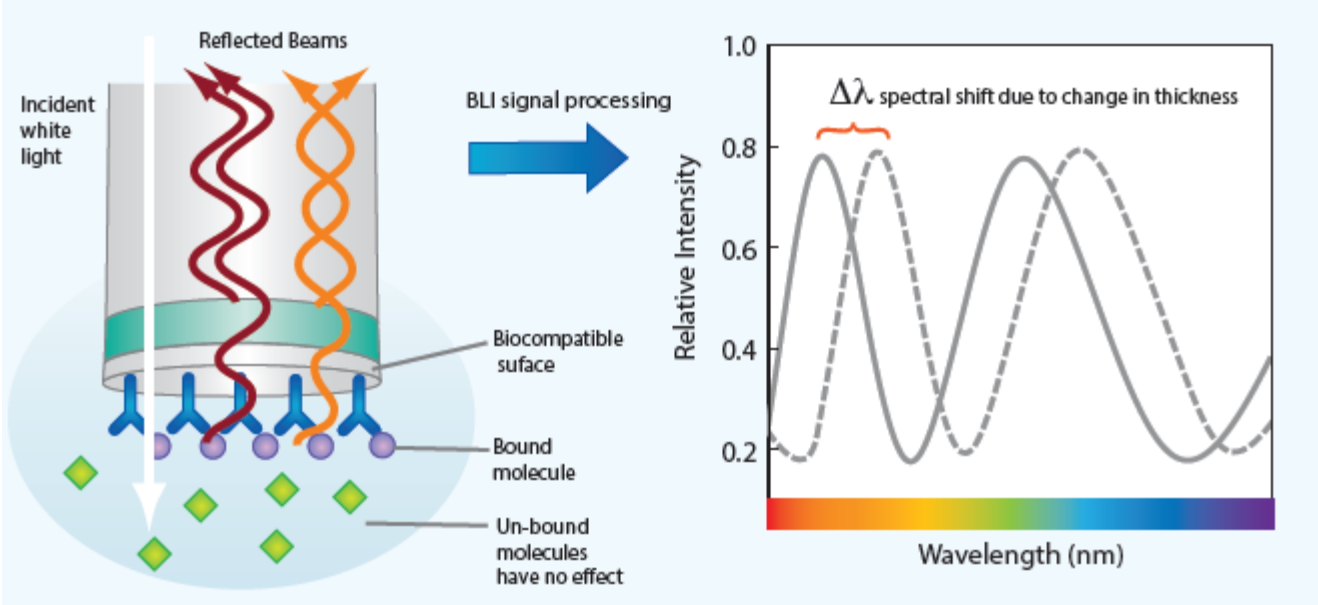
Driver Behind Development of an FcR Binding Panel

- **Important to test therapeutic mAb for binding with the complete panel of human IgG-Fc-binding receptors for characterization of original therapeutics and/or testing of biosimilar mAbs .**
 - If the mAb is an effector mAb, then it is necessary to show consistency in binding to the high and low affinity receptors CD64, CD32 and CD16.
 - If the mAb is a neutralizing moiety only (no effector function), one might then expect minimal binding, if any, to the Fc γ receptors.
 - However, all antibodies will bind in some fashion to the Fc γ receptors.
 - More importantly, all antibodies will bind to FcRN, which despite the name 'neonatal' is present in adult bone marrow-derived cells, as well as other tissue.
 - When an antibody binds to the FcRN, the Ab/FcRN complex is endocytosed and ends up in an endosomal vesicle with a low pH.
 - In the low pH environment of the endosome, the FcRN is in its high affinity state and as such retains a very tight hold on the antibody, essentially sequestering it from the blood stream.
 - When the endosome recycles back to the cell surface, the Ab/FcRN complex is exposed to neutral pH serum/plasma, and the FcRN shifts to its low affinity state, allowing the Ab to detach .
 - It is in this way that the FcRN is thought to play an important role in the PK of mAb therapeutics .

Toolbox Panel Approach

- **We (Catalent Pharma Solutions, Large Molecule Analytical Chemistry, Kansas City, Mo), set out to develop essentially a toolbox approach for assessing mAb/Fc receptor binding kinetics.**
 - The intent was to eventually use the analytical panel both for characterization of biosimilar molecules and comparison to originator molecules, as well as for CMC-related release and stability testing.
 - Use of SPR was one possible approach
 - We chose to proceed with the more robust, user friendly and more economical method of BLI, with an OctetRed96™.
 - A kinetics binding panel was developed on the Octet RED96 platform with recombinant
 - CD16a (FcγRIIIA), CD16b (FcγRIIIB),
 - CD32a (FcγRIIA), CD32b/c (FcγRIIB/C),
 - CD64 (FcγRI)
 - FcRN.

Basic Strategy For FcR Panel Development



- Full Panel:
- His-tagged (Commercially Available)*
 - huFcγRI
 - huFcγRIIA
 - huFcγRIIB/C
 - huFcγRIIIA
 - huFcγRIIIB
 - raFcRN

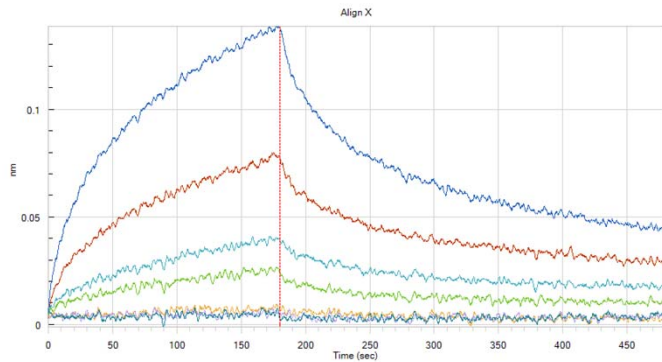
Case #1: FcR K_D Values for an Effector IgG₁

First Attempt at Establishing Fc Receptor Panel

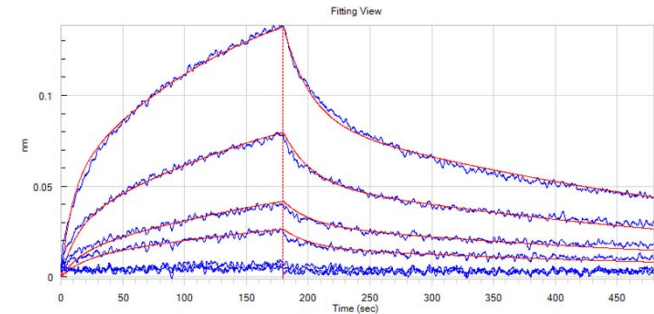
- **mAb with known effector activity**
- **Goal: to characterize/measure the K_D for Fc binding to each of the Fc Receptors**
- **Optimized load conditions for each receptor type**
- **Tested multiple association/dissociation conditions for each mAb/FcR interaction**
- **Project was cancelled before anything beyond optimizing reactions for the initial mAb/FcR K_D determinations**

Case #1: FcR K_D Values for an Effector IgG₁

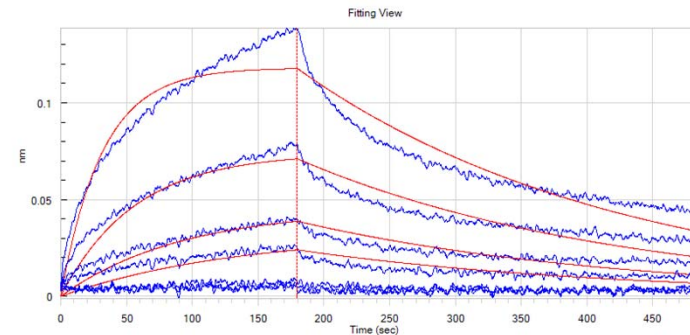
First Attempt at Establishing Fc Receptor Panel



2:1 Global
Fitting Model



1:1 Global Fitting
Model
(not as pretty,
but statistically
more accurate)



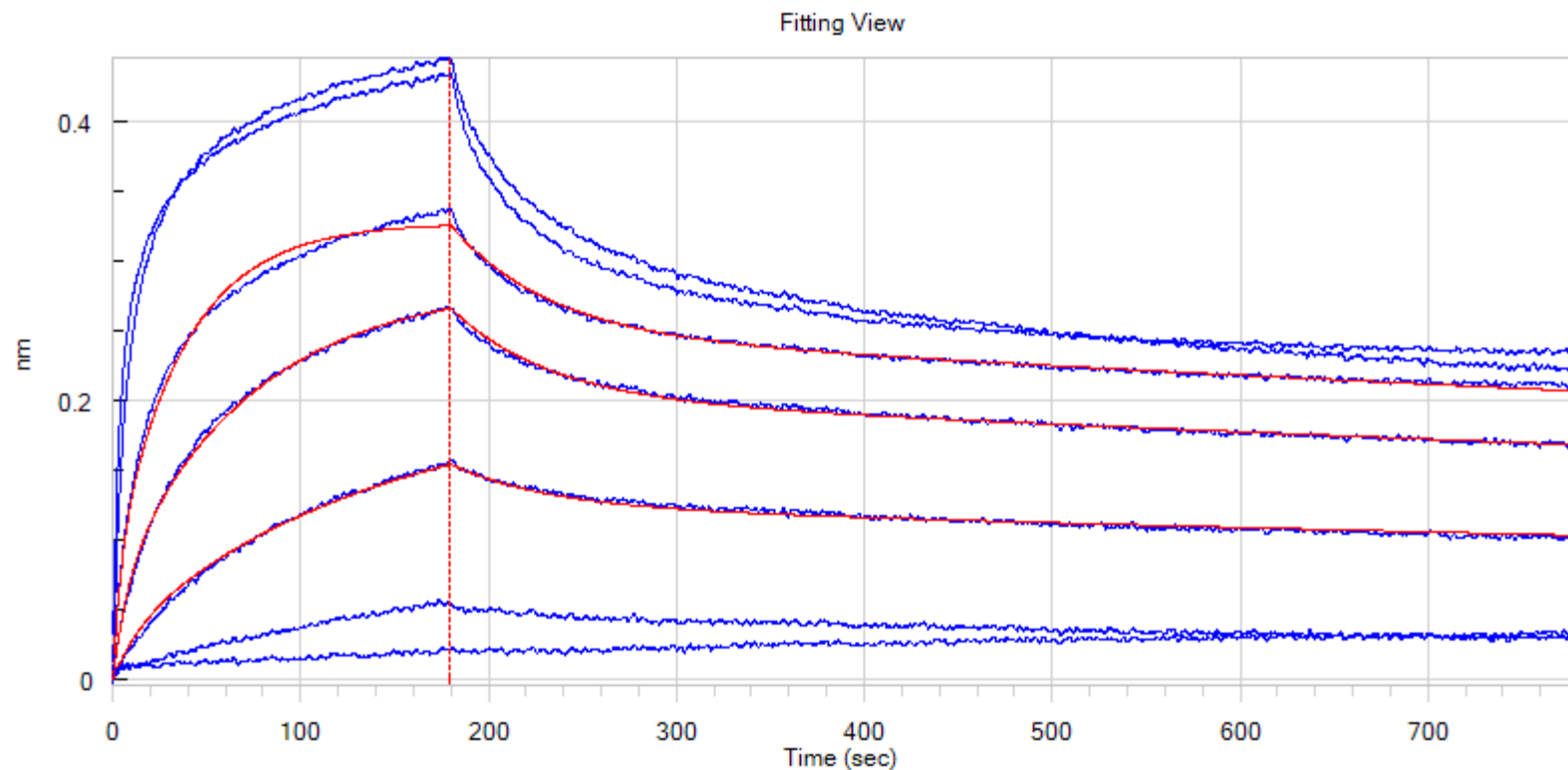
Case #1: FcR K_D Values for an Effector IgG₁

First Attempt at Establishing Fc Receptor Panel

FcR Type	CD nomenclature	Octet-determined mean K_D values (from 5-6 different conditions) (M)	Averaged K_D Error	Averaged K_D Error (% of K_D)	Octet-determined K_D value variability (n = 5 - 6) (%CV)
FcγRI	CD64	2.2×10^{-9}	8.3×10^{-11}	4%	18%
FcγRIIA	CD32a	1.4×10^{-7}	3.6×10^{-9}	3%	11%
FcγRIIB/C	CD32b/c	2.4×10^{-7}	1.4×10^{-8}	6%	37%
FcγRIIIA	CD16a	1.8×10^{-7}	4.7×10^{-9}	3%	30%
FcγRIIIB	CD16b	1.0×10^{-7}	1.4×10^{-9}	1%	23%
FcRN (@pH7.4) FcRN (@pH 6.0)	NA	2.1×10^{-7} 1.6×10^{-9}	5.1×10^{-9} 4.3×10^{-11}	2% 3%	53% 34%

Case #1: FcR K_D Values for an Effector IgG₁ First Attempt at Establishing Fc Receptor Panel

Analysis of FcRN Binding, High Affinity (Acidic pH) Condition



Keep this in mind for later

Case #2: Recent Project using Fc Receptor Panel

- **IgG₁ mAb with no known effector activity**
 - MOA: Neutralization of a soluble factor
- **Goal: to characterize/measure the K_D for Fc binding to each of the Fc Receptors**
- **Establish assay for eventual use as part of release and stability testing**
- **Started with receptor load conditions for each receptor type based upon optimization from prior project**
 - Required minimal reoptimization
 - Primarily around selection of buffer with minimal non-specific signal
- **Tested multiple association/dissociation conditions for each mAb/FcR interaction**
 - This did require reoptimization (load time and dissociation time)
- Final execution done with 1 reference standard and 2 lots of drug substance
 - Each sample run as independent duplicates

Case #2: FcR K_D Values for a Neutralizing IgG1

FcγRI (CD64)

Reference Standard						Drug Substance #1					Drug Substance #2						
	KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²
Rep1	2.102E-09	1.466E-10	7.0%	0.451	0.999	Rep1	9.676E-09	2.076E-10	2.1%	1.283	0.991	Rep1	9.267E-09	1.925E-10	2.1%	1.136	0.994
Rep2	3.991E-09	2.082E-10	5.2%	0.667	0.997	Rep2	8.176E-09	1.685E-10	2.1%	13.477	0.991	Rep2	8.081E-09	1.693E-10	2.1%	1.488	0.992
Ave	3.047E-09	1.801E-10				Ave	8.926E-09	1.891E-10				Ave	8.674E-09	1.813E-10			
RD	1.336E-09					RD	1.061E-09					RD	8.386E-10				
%RD	44%					%RD	12%		% of RS K _n	293%		%RD	10%		% of RS K _n	285%	

FcγRIIA (CD32a)

Reference Standard						Drug Substance #1					Drug Substance #2						
	KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²
Rep1	3.572E-08	1.078E-09	3.0%	0.025	0.999	Rep1	4.768E-08	1.320E-09	2.8%	0.028	0.999	Rep1	4.951E-08	1.643E-09	3.3%	0.036	0.998
Rep2	NA	NA	NA	NA	NA	Rep2	4.704E-08	1.365E-09	2.9%	0.025	0.999	Rep2	4.036E-08	1.810E-09	4.5%	0.047	0.996
Ave	NA	NA				Ave	4.736E-08	1.343E-09				Ave	4.494E-08	1.729E-09			
RD	NA					RD	4.525E-10					RD	6.470E-09				
%RD	NA					%RD	1%		% of RS K _n	133%		%RD	14%		% of RS K _n	126%	

FcγRIIB/C (CD32b/c)

Reference Standard						Drug Substance #1					Drug Substance #2						
	KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²		KD	Error	% Error	χ ²	R ²
Rep1	6.319E-08	2.651E-09	4.2%	0.074	0.994	Rep1	8.908E-08	3.937E-09	4.4%	0.068	0.993	Rep1	8.565E-08	4.610E-09	5.4%	0.071	0.986
Rep2	5.586E-08	2.743E-09	4.9%	0.056	0.993	Rep2	8.676E-08	4.631E-09	5.3%	0.075	0.990	Rep2	7.210E-08	4.053E-09	5.6%	0.096	0.981
Ave	5.953E-08	2.697E-09				Ave	8.792E-08	4.298E-09				Ave	7.888E-08	4.340E-09			
RD	5.183E-09					RD	1.640E-09					RD	9.581E-09				
%RD	9%					%RD	2%		% of RS K _n	148%		%RD	12%		% of RS K _n	133%	

Case #2: FcR K_D Values for a Neutralizing IgG1

Fc γ RIIIA (CD16a)

Reference Standard						Drug Substance #1						Drug Substance #2					
	KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²
Rep1	1.746E-07	1.542E-08	8.8%	0.105	0.979	Rep1	6.056E-07	3.300E-08	5.4%	0.122	0.979	Rep1	3.842E-07	2.316E-08	6.0%	0.098	0.979
Rep2	2.145E-07	1.316E-08	6.1%	0.049	0.957	Rep2	4.191E-07	2.252E-08	5.4%	0.122	0.973	Rep2	2.314E-07	1.743E-08	7.5%	0.250	0.989
Ave	1.946E-07	1.433E-08				Ave	5.124E-07	2.825E-08				Ave	3.078E-07	2.050E-08			
RD	2.821E-08					RD	1.319E-07					RD	1.080E-07				
%RD	15%					%RD	26%		% of RS K_D	263%		%RD	35%		% of RS K_D	158%	

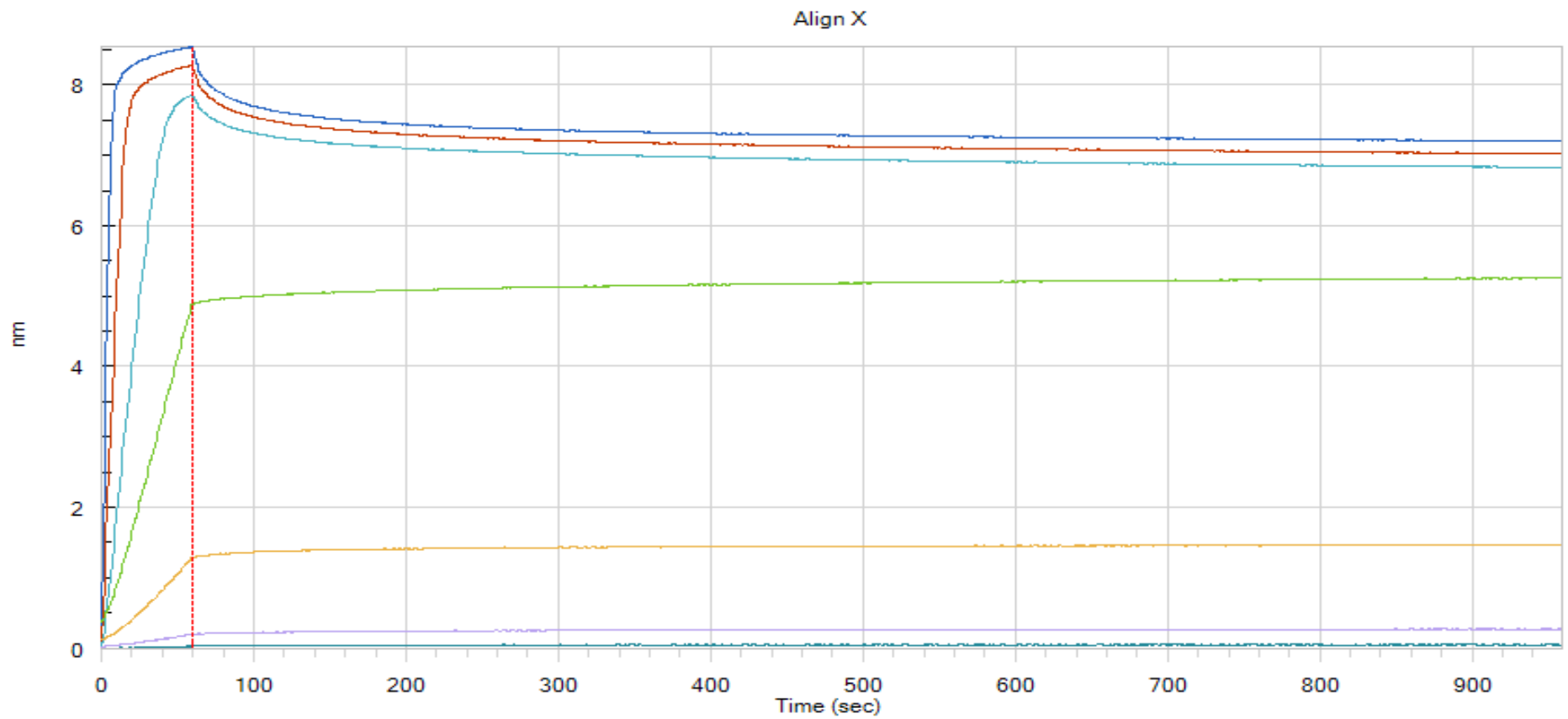
Fc γ RIIIB (CD16b)

Reference Standard						Drug Substance #1						Drug Substance #2					
	KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²
Rep1	1.231E-07	1.210E-08	9.8%	0.050	0.882	Rep1	3.159E-07	1.644E-08	5.2%	0.142	0.986	Rep1	3.936E-07	2.059E-08	5.2%	0.215	0.979
Rep2	9.526E-08	8.266E-09	8.7%	0.028	0.969	Rep2	2.686E-07	1.820E-08	6.8%	0.082	0.977	Rep2	1.964E-07	1.208E-08	6.2%	0.258	0.984
Ave	1.092E-07	1.036E-08				Ave	2.923E-07	1.734E-08				Ave	2.950E-07	1.688E-08			
RD	1.969E-08					RD	3.345E-08					RD	1.394E-07				
%RD	18%					%RD	11%		% of RS K_D	268%		%RD	47%		% of RS K_D	270%	

Fc γ RN, pH 7.5 (Low Affinity Condition)

Reference Standard						Drug Substance #1						Drug Substance #2					
	KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²		KD	Error	% Error	X ²	R ²
Rep1	1.490E-07	1.192E-08	8.0%	0.018	0.984	Rep1	3.963E-07	1.668E-08	4.2%	0.067	0.993	Rep1	3.113E-07	1.446E-08	4.6%	0.137	0.993
Rep2	1.000E-07	8.228E-09	8.2%	0.010	0.980	Rep2	4.665E-07	2.294E-08	4.9%	0.075	0.990	Rep2	2.991E-07	1.397E-08	4.7%	0.135	0.992
Ave	1.245E-07	1.024E-08				Ave	4.314E-07	2.006E-08				Ave	3.052E-07	1.422E-08			
RD	3.465E-08					RD	4.964E-08					RD	8.627E-09				
%RD	28%					%RD	12%		% of RS K_D	347%		%RD	3%		% of RS K_D	245%	

FcRN Binding at Acidic (High Affinity) pH



Summary of mAb/Fc Receptor Binding Panel

- **Results derived for the K_D 's of IgG for the different Fc receptors fall in line with expectations.**
- **Fc γ RII and Fc γ RIII show low to medium binding kinetics in the ' μ M' range (0.01 – 1.0).**
- **Fc γ RI is high affinity with KD values in the nM range.**
- **FcRN has μ M range binding at neutral pH (low affinity condition), but, when measurable, nM range binding at acidic pH (high affinity condition)**
 - The FcRN binding kinetics at the two different pH conditions is consistent with the current understanding of the function of the FcRN receptor, and how it impacts antibody half-life in vivo.
 - Extremely high affinity for some FcRN binding (in acidic conditions) are more problematic to derive an actual K_D value.
- **Replicate analyses of the same sample demonstrate excellent reproducibility/precision.**

Conclusions

- **Successfully demonstrated development and establishment of a platform for characterization of IgG/Fc receptor binding at Catalent**
- **Utilizing full panel of appropriate Fc receptors and easy-to-use Octet instrumentation.**
- **As the immobilized moieties for this panel are the Fc receptors, the analyte of interest (the mAb) is in no way derivatized**
 - Allows objective analysis of the receptor panel with any mAb.
- **With minimal additional optimization, the Fc receptor panel established here can be quickly adapted for any mAb.**
- **In the process of finalizing validation of the Octet RED96 system, and will be further developing and qualifying/validating the Fc receptor panel testing for use not only in characterization, but in cGMP release and stability assessments when appropriate .**

A few other insights

Use of Octet in Determining Biosimilarity

CDR/Target Binding

Batch Number	K_D (M)	K_D Average (M)
1	4.8×10^{-8}	4.7×10^{-8}
2	3.5×10^{-8}	
3	2.9×10^{-8}	
4	2.8×10^{-8}	
5	7.5×10^{-8}	
6	6.7×10^{-8}	

K_D determination of originator mAb CDR/Ligand interaction

Biosimilar Candidate Number	K_D (M)
1	5.9×10^{-8}
2	4.6×10^{-8}
3	4.8×10^{-8}

K_D determination of mAb CDR/Ligand interaction for potential biosimilar candidates

Use of Octet in Determining Biosimilarity

Fc/FcR γ IIIa Binding

Batch Number	K _D (M)	K _D Average (M)
1	5.1x10 ⁻⁶	4.4x10 ⁻⁶
2	4.7x10 ⁻⁶	
3	6.2x10 ⁻⁶	
4	5.7x10 ⁻⁶	
5	2.3x10 ⁻⁶	
6	2.3x10 ⁻⁶	

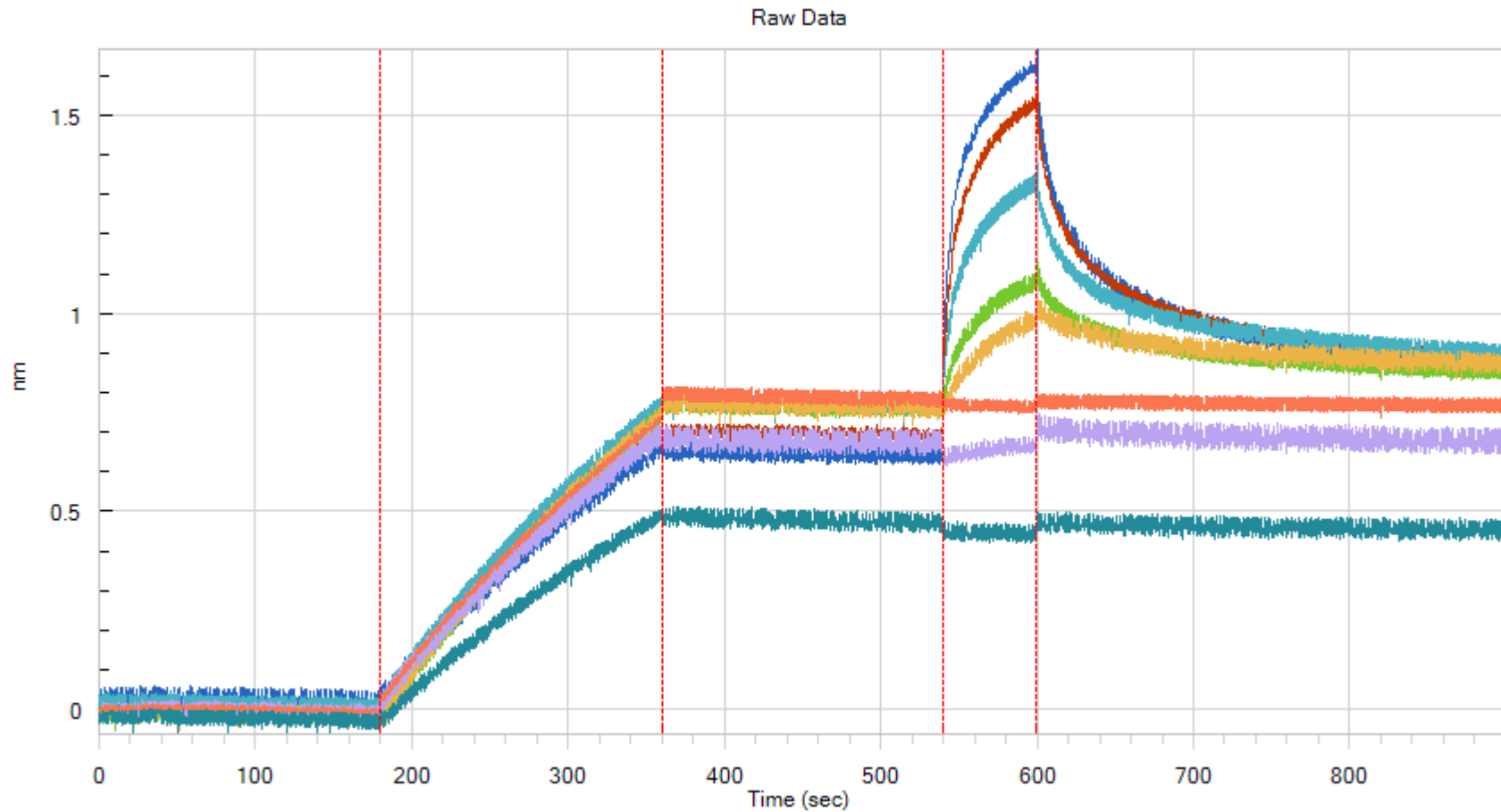
K_D determination of originator mAb Fc/FcR γ IIIa interaction

Biosimilar Candidate Number	K _D (M)
1	1.9x10⁻⁵
2	2.3x10 ⁻⁶
3	4.6x10 ⁻⁶

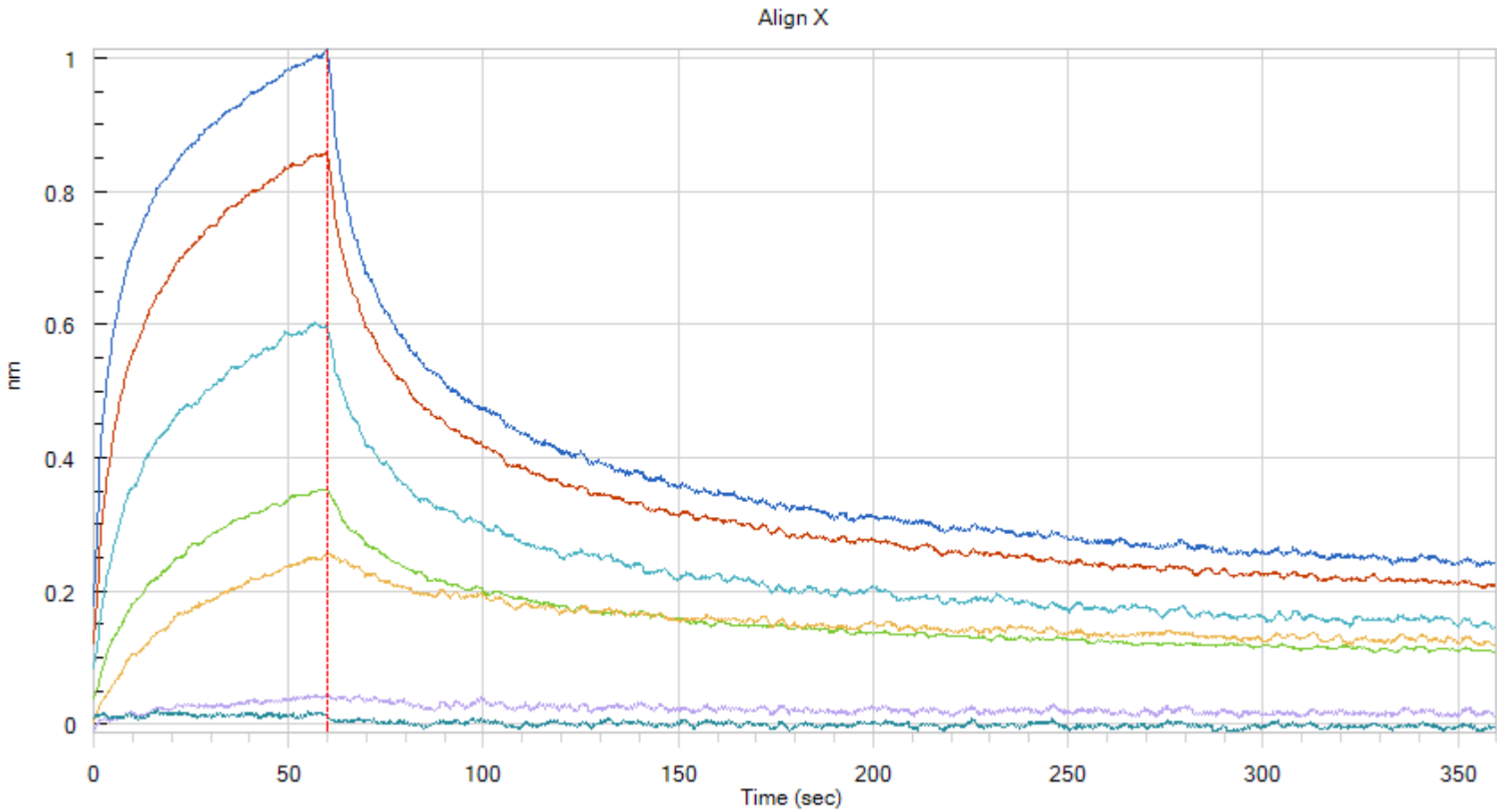
K_D determination of mAb Fc/FcR γ IIIa interaction for potential biosimilar candidates

A Funny Thing Happened on the Way to the Binding....

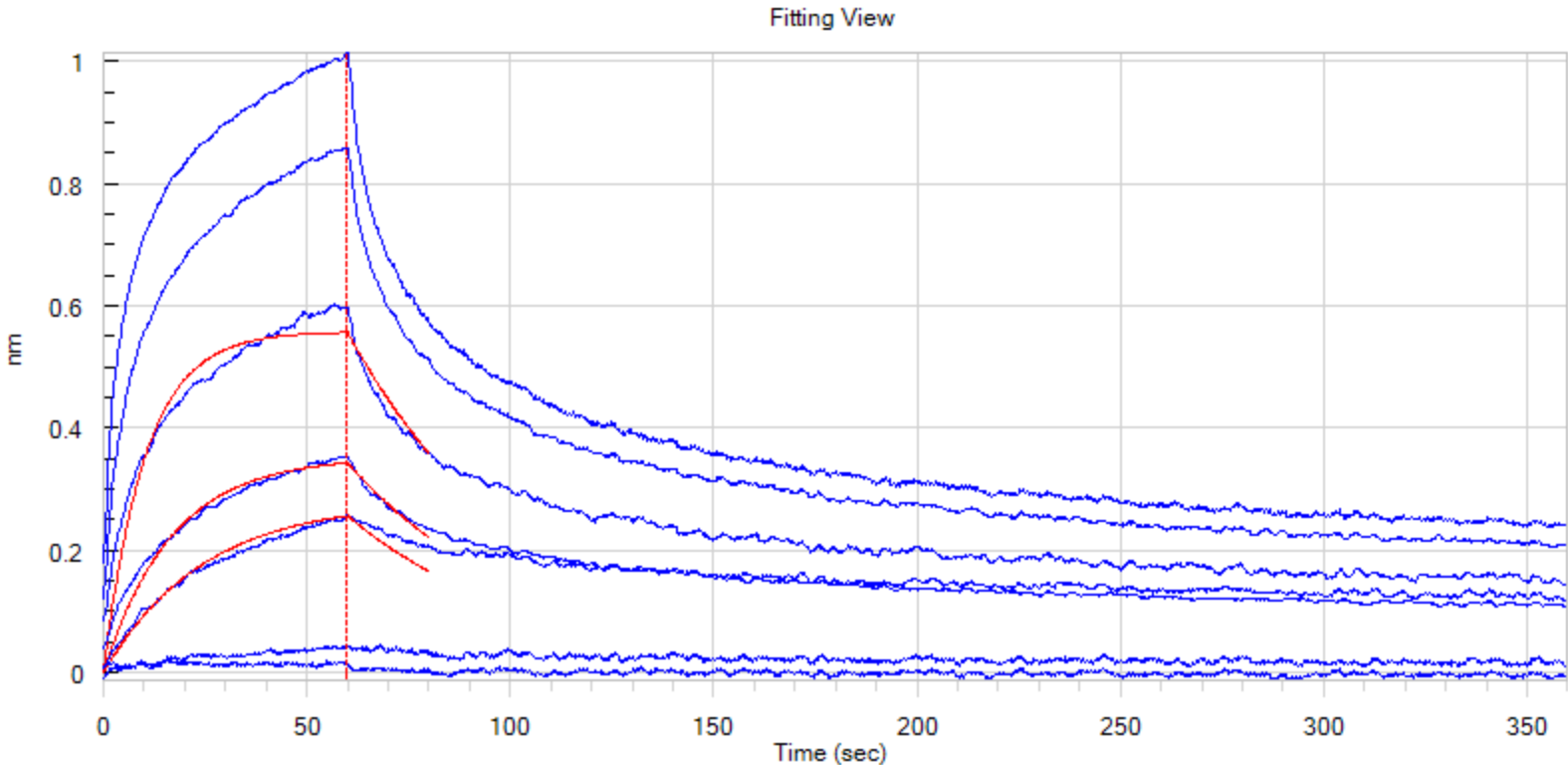
CD32b/c (Fc γ RIIB/C) Analysis



CD32b/c (Fc γ RIIB/C) Analysis



CD32b/c (Fc γ RIIB/C) Analysis

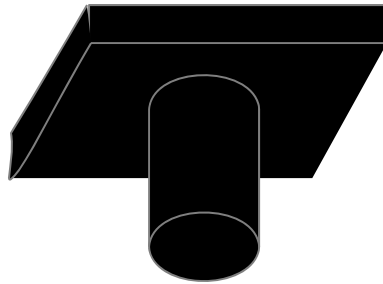


CD32b/c (FcγRIIB/C) Analysis

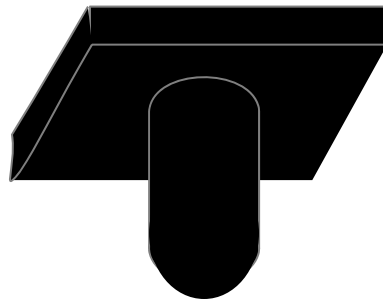
Loading Sample ID	Conc. (nM)	Response	KD (M)	KD Error	Full X ²	Full R ²	kon(1/Ms)	kon Error	kdis(1/s)	kdis Error	Rmax	Rmax Error
FcgRIIB/C	1000	0.9796										
FcgRIIB/C	500	0.8321										
FcgRIIB/C	250	0.5802	7.210E-08	4.053E-09	0.096064	0.981154	3.04E+05	1.25E+04	2.20E-02	8.43E-04	0.7187	0.015
FcgRIIB/C	125	0.3363	7.210E-08	4.053E-09	0.096064	0.981154	3.04E+05	1.25E+04	2.20E-02	8.43E-04	0.5557	0.0176
FcgRIIB/C	62.5	0.2356	7.210E-08	4.053E-09	0.096064	0.981154	3.04E+05	1.25E+04	2.20E-02	8.43E-04	0.6042	0.0251
FcgRIIB/C	31.3	0.0369										
FcgRIIB/C	15.6	0.0125										

Answer to the Mystery.....

Look carefully at the plate you use!!!



Use this



Not this

Acknowledgements

Dan Papa, PhD
Tiffany Walker

*Special Thanks to
David Apiyo, PhD*



discover more.

CATALENT PHARMA SOLUTIONS
14 SCHOOLHOUSE ROAD
SOMERSET, NJ 08873

+ 1 866 720 3148

www.catalent.com



DEVELOPMENT



DELIVERY



SUPPLY

more products. better treatments. reliably supplied.™