

Design and Validation of a Non Cell-based Receptor Binding Assay for the Detection of Neutralizing Antibodies to a Biological Therapeutic

Shalini Gupta

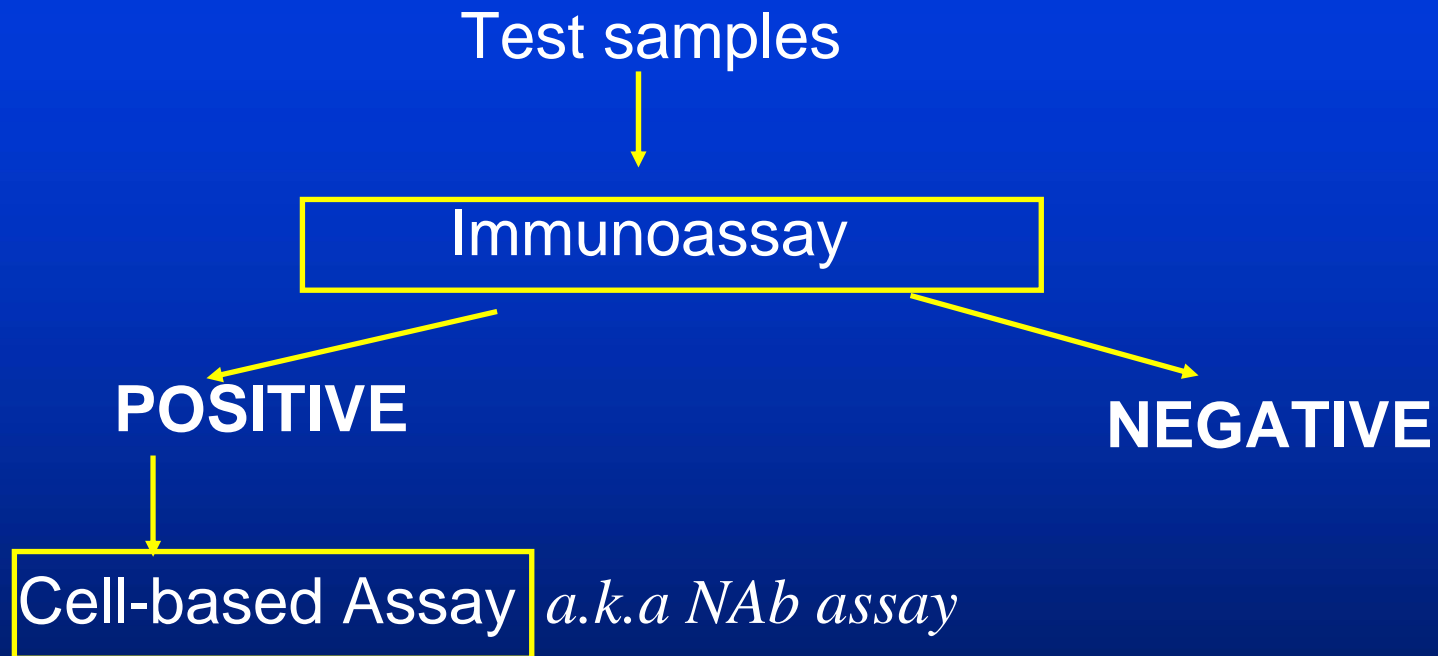
Associate Director

Amgen, Thousand Oaks, CA

shaliniq@amgen.com

May 24th, 2005

Immunogenicity Testing



Provide a functional biological system to assess if the Abs detected by the immunoassay have neutralizing capability

Cell-based NAb Assay Designs

- Utilize intracellular signaling events triggered within a cell line that responds to :
 - * the drug product (eg. cytokines)
 - * the ligand inhibited by the drug product (eg. MAbs)
- May also utilize study of extracellular binding events at the cell surface (eg. MAbs)

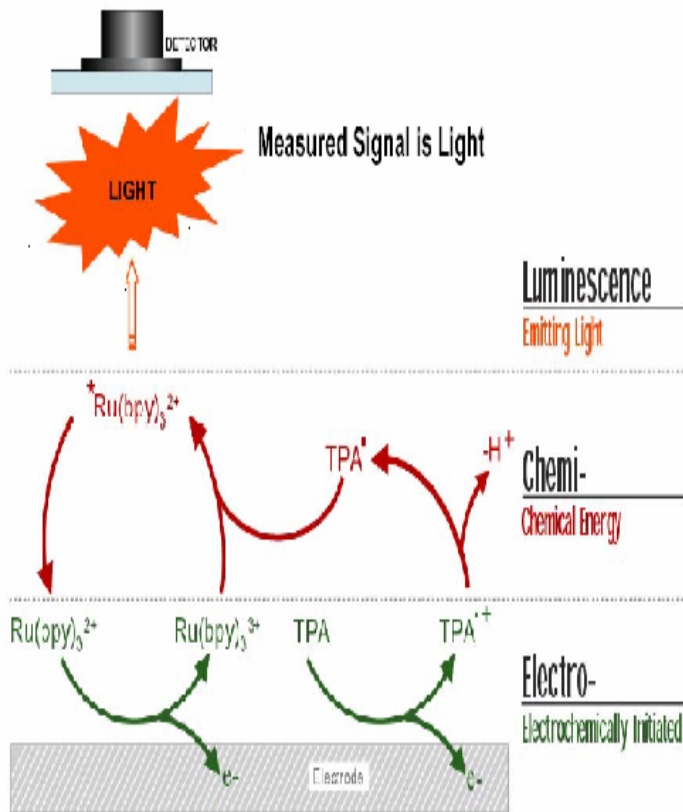
Requirements for Cell-based NAb Assays

- Stable cell line responsive to drug
- Measurable Readout (cpm, OD, conc of secreted protein)
- Robust signal to noise ratio in serum matrix
- Availability of a positive control antibody to the drug
- Medium to High Throughput
- Adaptable to Automation

Justification For a Non-cell Based NAb Assay

- Drug product : blocked ligand activity
- Limited choice of cell lines expressing the receptor for ligand that was target for drug product
- Only cell line available yielded at best 2-fold signal when treated with 100-200 ng/mL ligand
- Cell line had finite life span
- Difficult cells to culture
- Low throughput (24-well plates)
- Irreproducible bioassay results

Meso-Scale Discovery (MSD) Technology



Meso-Scale Discovery (MSD) technology employs microtitre plates fitted with a series of electrodes associated with the bottom of each well.

Using an MSD Sector PR™ plate reader, an electrical current is placed across the plate-associated electrodes in the presence of a Tripropylamine (TPA) containing buffer.

The result is a series of electrically induced oxidation-reduction reactions involving Ruthenium (from the captured complex) and TPA leading to a luminescent signal. The consequent electrochemiluminescent (ECL) signal is measured by photodiodes and is quantified as a relative unit (RU).

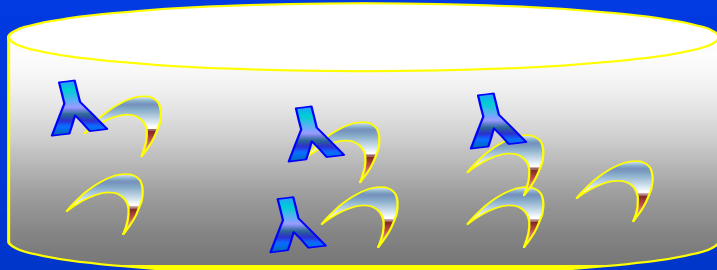
Assay Reagents

- **Purified Soluble receptor (ruthenylated)**
- **Purified Ligand (biotinylated)**
- **Drug product (TP)**
- **Streptavidin-coated MSD Plates**
- **Cynomolgus monkey serum**
- **Affinity purified rabbit anti-TP antibody**

Assay Methodology

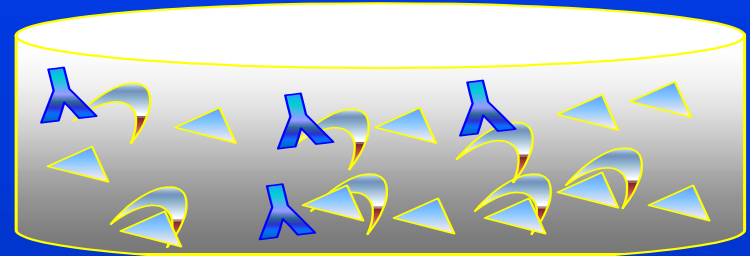
Assay Format

1.



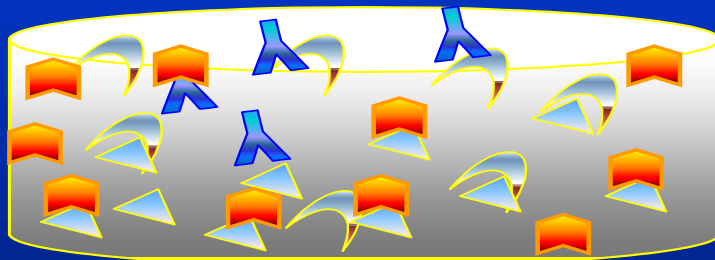
Incubate serum sample with TP - 15-30 Min

2.



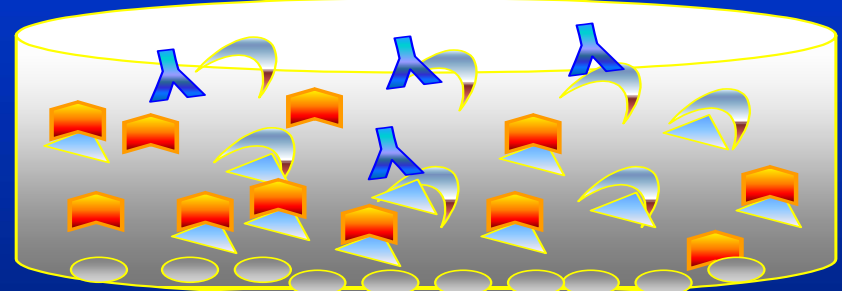
Add L-Biot - 15-30 Min

3.



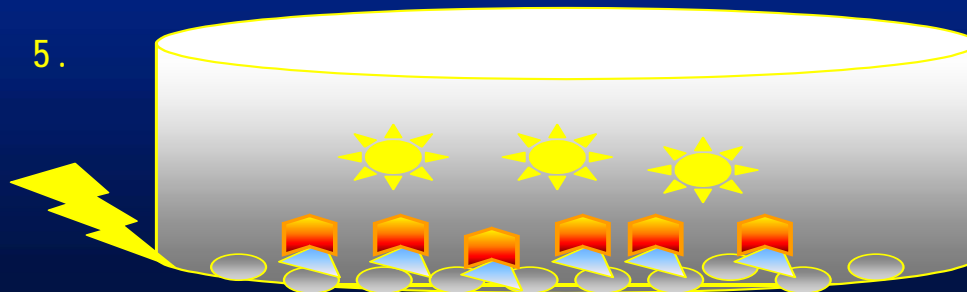
Add R-Ru - 15-30 Min

4.

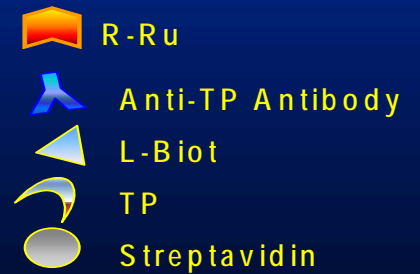


Add Mixture to blocked SA-Coated Multiarray Plate - 30-60 Min

5.



Wash 2X and add TPA
Containing read buffer
Read on MSD reader.



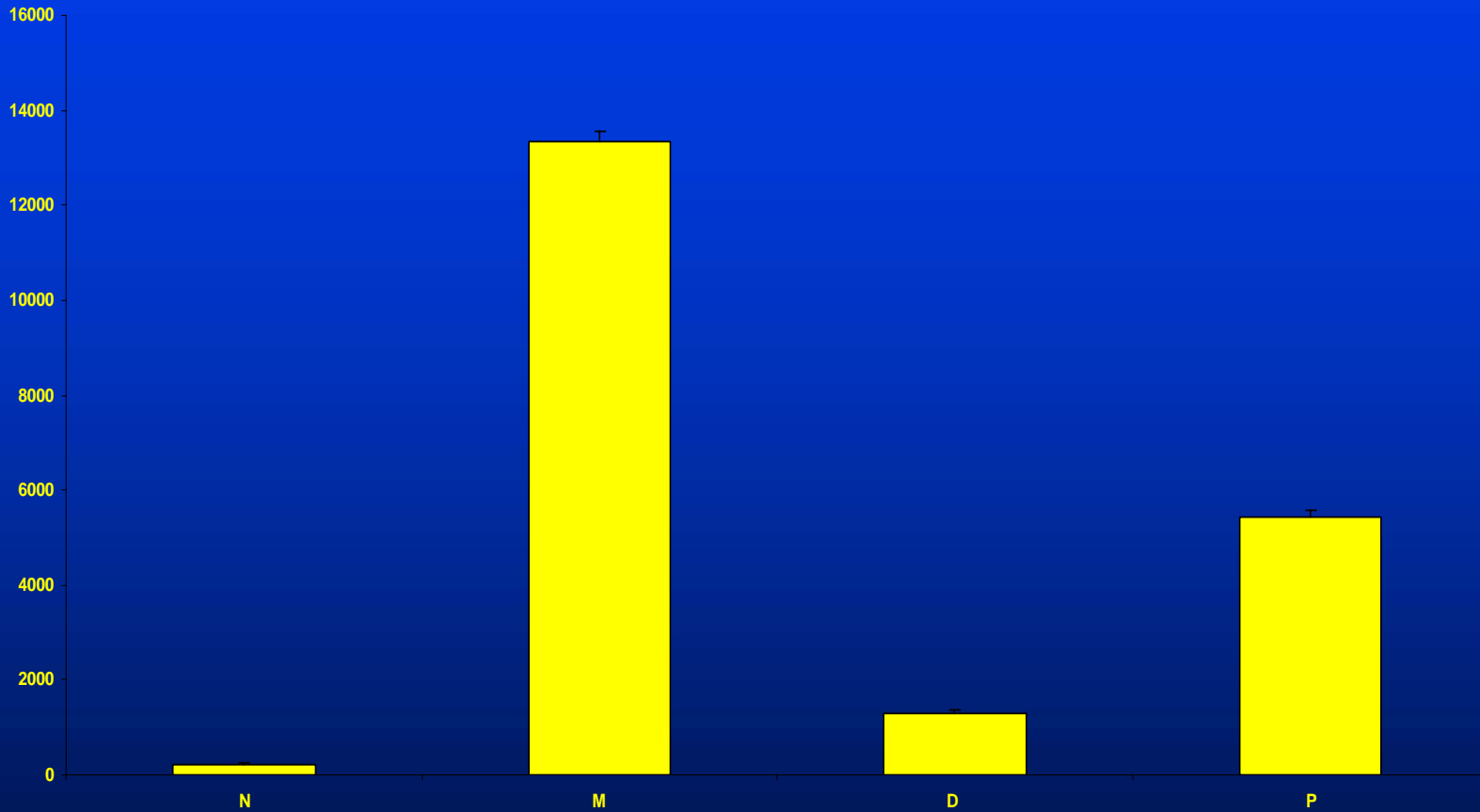
Assay Development

- **Optimized ratio between Ru-R and Biot-L for optimal signal**
- **Used 50% cyno serum for initial development expts**
- **Due to donor (n=10) variability serum was reduced to 15%**
- **Optimized order of mixing of reagents (capture vs. homogenous)**
- **Improved precision and inter-day repeatability in homogenous format with assay controls**

Optimized Assay Conditions

- 15% cynomolgus monkey serum
- 250 ng/mL Biot-Ligand
- 900 ng/mL Ru-R
- 80 ng/mL TP

Assay Controls



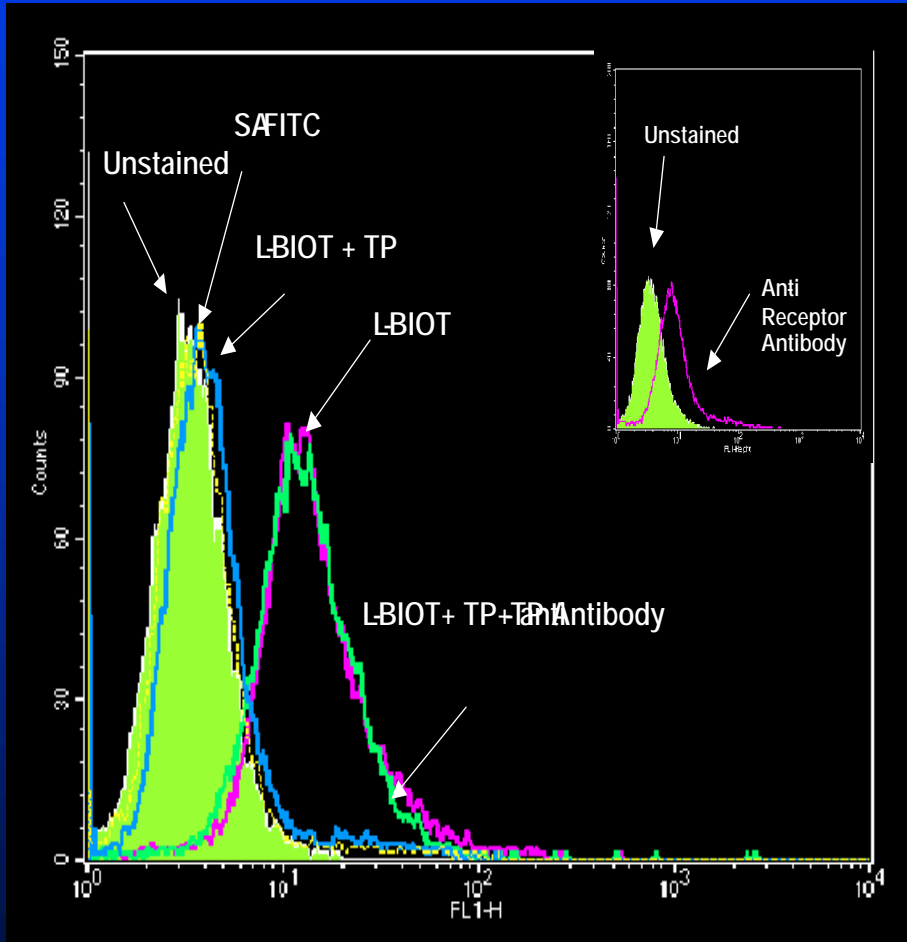
N = background (15% PCS with R-Ru)

M = maximum binding (15% PCS with R-Ru and L-biot)

D = drug control (15% PCS with TP, r-RU and L-biot)

P = positive control (D with anti-TP antibody)

FACS Analysis of a Cell Line Naturally Expressing the Target Receptor



- A cell line naturally expressing the receptor was used to demonstrate:

- L-Biot binding, inhibition of binding by TP
- Restoration of binding by the anti-TP NAb.

- Binding of L-Biot to the receptor was detected using a SA-FITC conjugate.

- Receptor expression was confirmed using a biotinylated goat polyclonal anti-receptor Ab and detected with SA-FITC.

- Binding of L-Biot to the target receptor was inhibited by the TP and restored by the anti-TP NAb.

NAb Testing Strategy

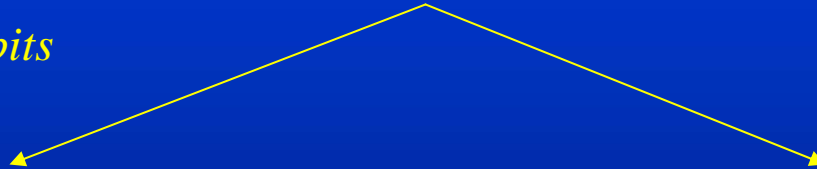
IM-Positive Samples



NAb Receptor Binding Assay

*Sample inhibits
TP activity*

*Sample does not inhibit
TP activity*



Specificity Assay

(in absence of TP)

*(sample shows non-specific
activity)*

NEGATIVE



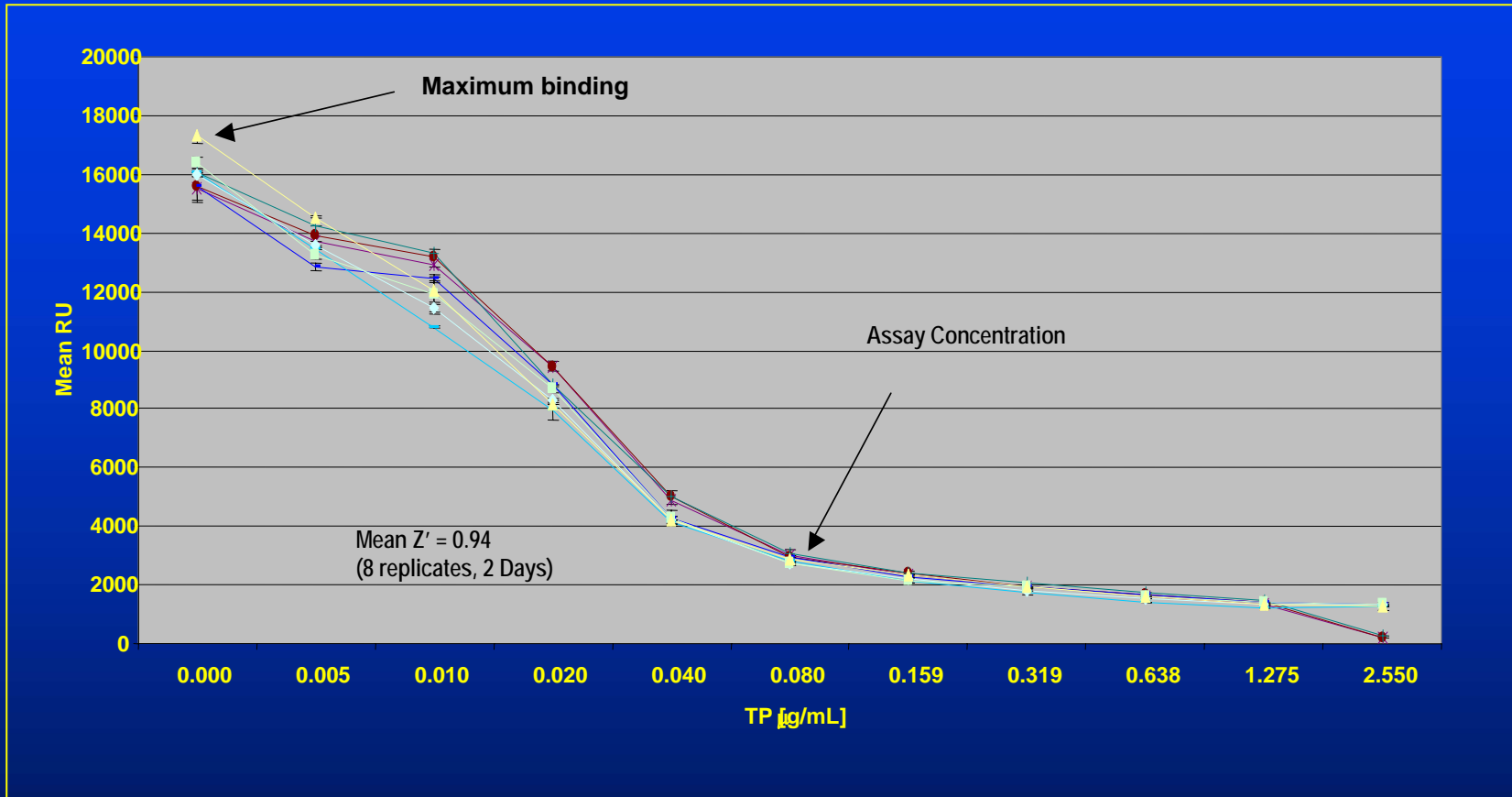
*Sample does not show non-
Specific activity)*

POSITIVE

Validation Experiments

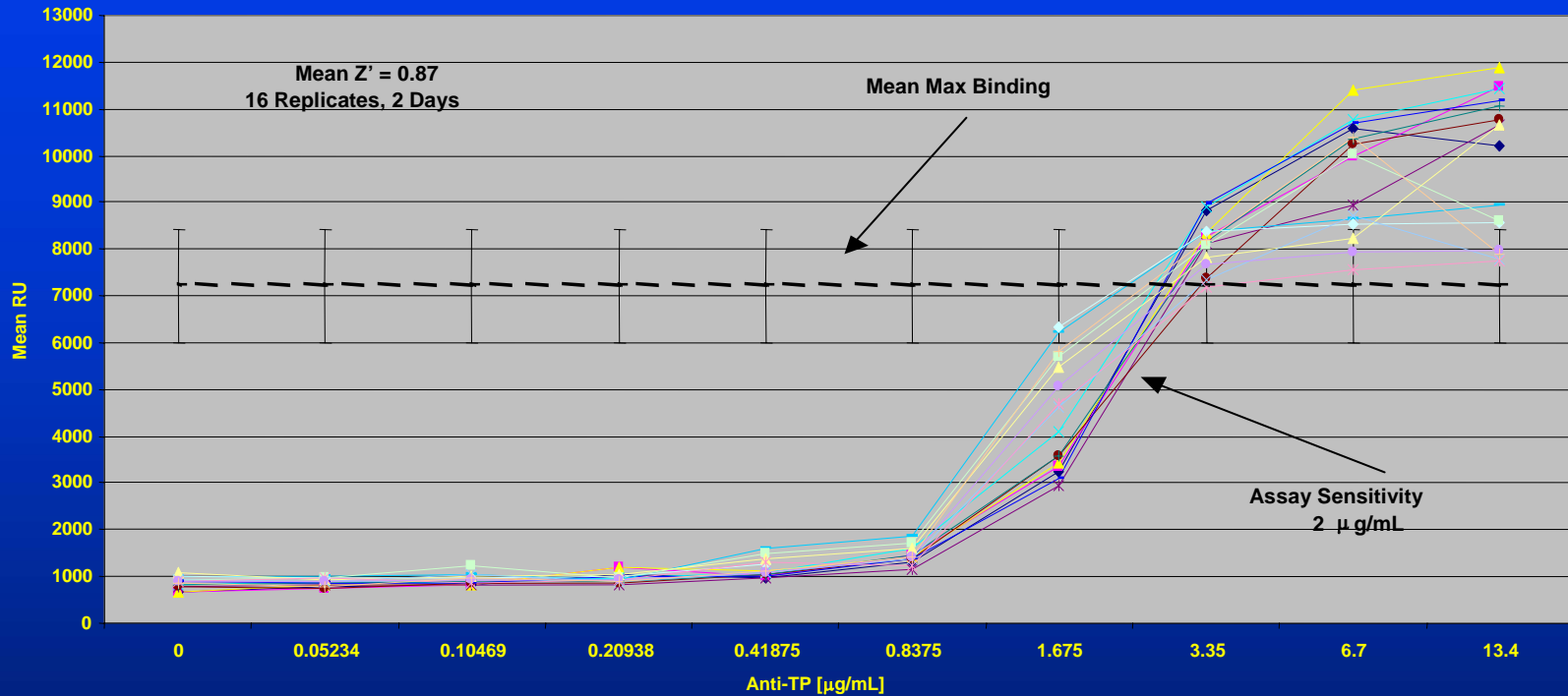
- **TP dose response curves**
- **Positive Control Ab curves**
- **NAb assay cutoff**
- **Specificity Assay Cutoff**
- **Interference by drug product**
- **Freeze-thaw stability**

Inhibition of L-Biot Binding to R-Ru By TP in 15% Pooled Cynomolgus Monkey Serum



- TP was incubated at increasing concentrations with 250 ng/mL L-Biot followed by incubation with 900 ng/mL R-Ru in 15% PCS.
- TP showed a dose dependent inhibition of L-Biot/R-Ru binding.
- A concentration of 80 ng/mL TP was chosen to be used for the Screening Assay.

Neutralization of TP and Restoration of Ligand Binding by Anti-TP Neutralizing Antibody in Pooled Cynomolgus Monkey Serum

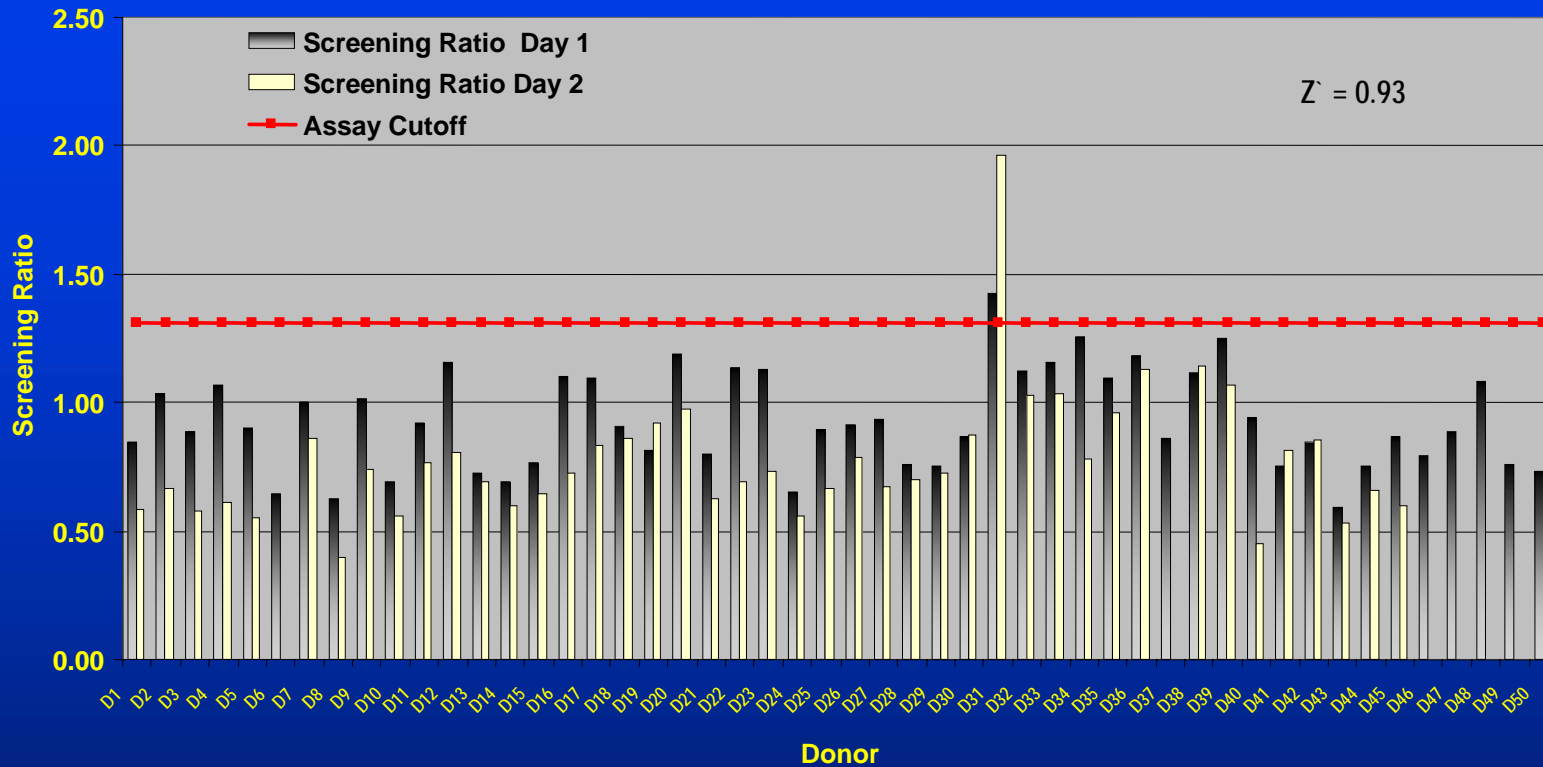


- Rabbit Polyclonal anti-TP Antibody (NAb) was added to neat PCS.
- The Ab containing PCS was diluted to 15% and preincubated with 80 ng/mL TP followed by sequential incubations with 250 ng/mL L-Biot and 900 ng/mL R-Ru.
- Anti-TP antibody showed a dose dependent restoration of L-Biot/R-Ru binding to maximal binding signal (above curves represent 2 days testing)
- The lowest concentration of Anti-TP Nab to restore binding by twofold was 1 µg/mL

Precision and Analytical Recovery of the PAb

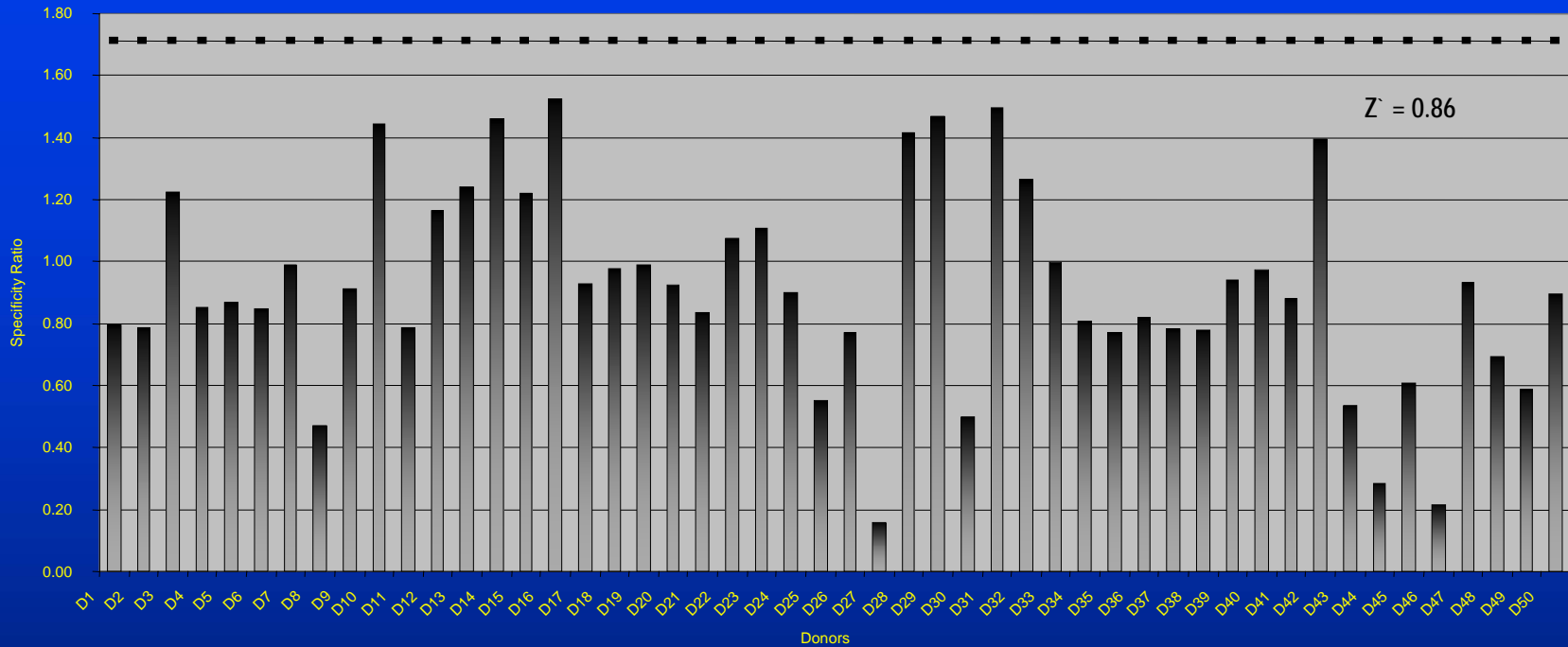
Ab (ug/ mL)	0.055	0.11	0.22	0.44	0.88	1.75	3.5	7	14
AR	356%	228%	117%	138%	98%	99%	98%	87%	68%
%CV									
Inter-day	19%	29%	17%	8%	11%	2%	4%	25%	77%
Inter-assay	35%	40%	44%	16%	8%	3%	7%	0.00	49%
Intra-assay	14%	20%	14%	13%	10%	2%	9%	26%	104%

Screening Assay Cutoff



- Serum from 50 individual cynomolgus monkeys was tested at a final dilution of 15% with 80 ng/mL TP, 250 ng/mL L-Biot and 900 ng/mL R-Ru.
- Samples results were compared with control results in PCS to by dividing Sample Results/Control Results (Screening Ratio).
- Samples were tested on two separate days and the Ratio 1 values were combined to establish an assay cutoff using the upper bound of the 95% prediction limit.
- The assay cutoff (Screening Ratio) was determined to be 1.26.

Specificity Assay Cutoff



- Serum from 50 individual cynomolgus monkeys was tested at a final dilution of 15%, 250 ng/mL L-Biot and 900 ng/mL R-Ru in the absence of the TP to assess for nonspecific enhancement of ligand binding.

- Samples results were compared with control results in PCS to by dividing Sample Results/Control Results (Ratio 3).

- The Ratio 3 values were used to establish an assay cutoff using the upper bound of the 95% prediction limit.

- The assay cutoff (Specificity Ratio) was determined to be 1.71.

AP Criteria for Positive/Negative

- **Assay Acceptance Criteria**
 - $M/N \geq 10$
 - $M/D \geq 2.0$
 - $P/D \geq 1.26$

Pre-dose: Ratio 1 ≥ 1.26

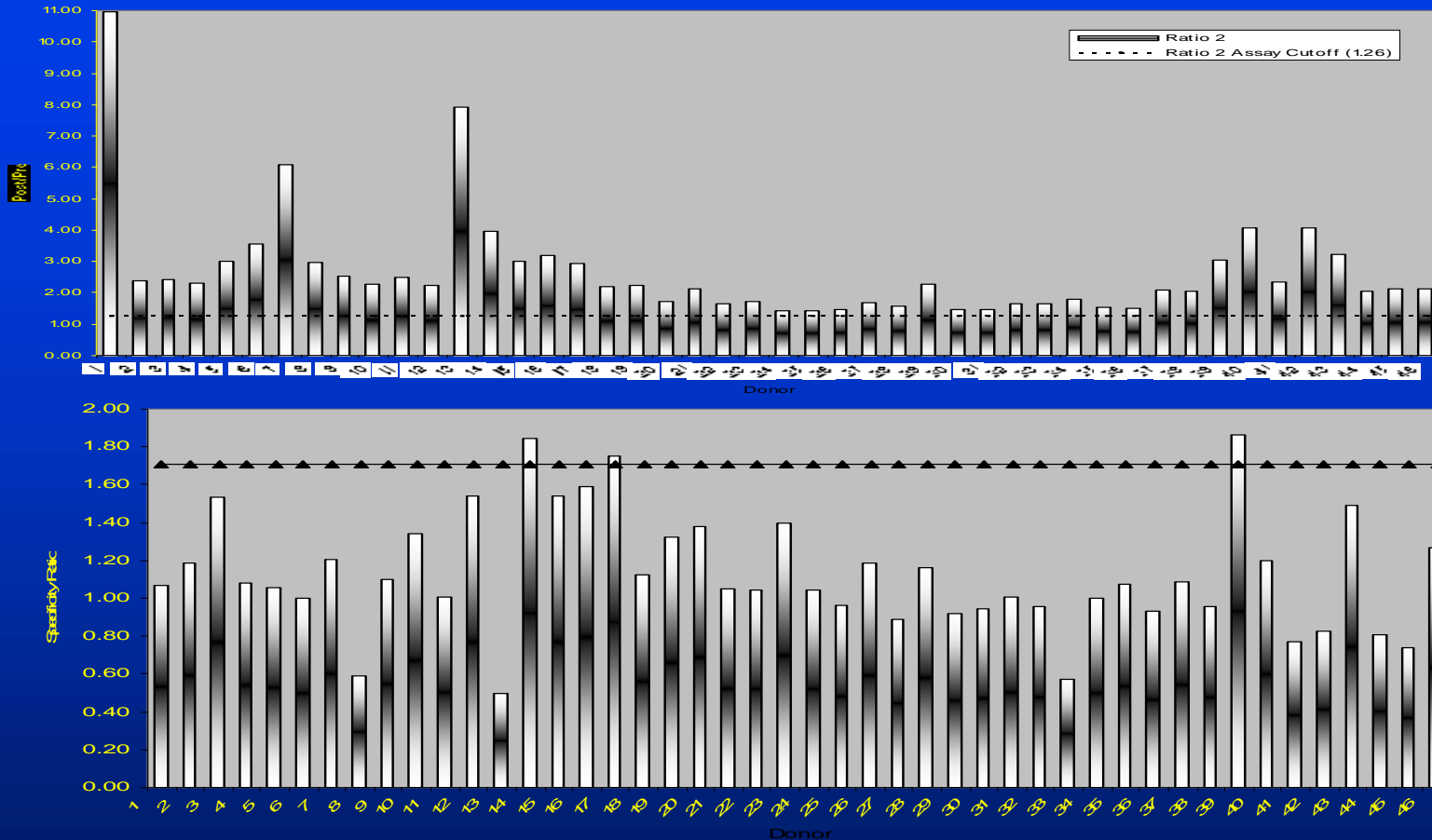
Ratio 3 < 1.71

Post-dose: Ratio 2 ≥ 1.26

Ratio 3 < 1.71

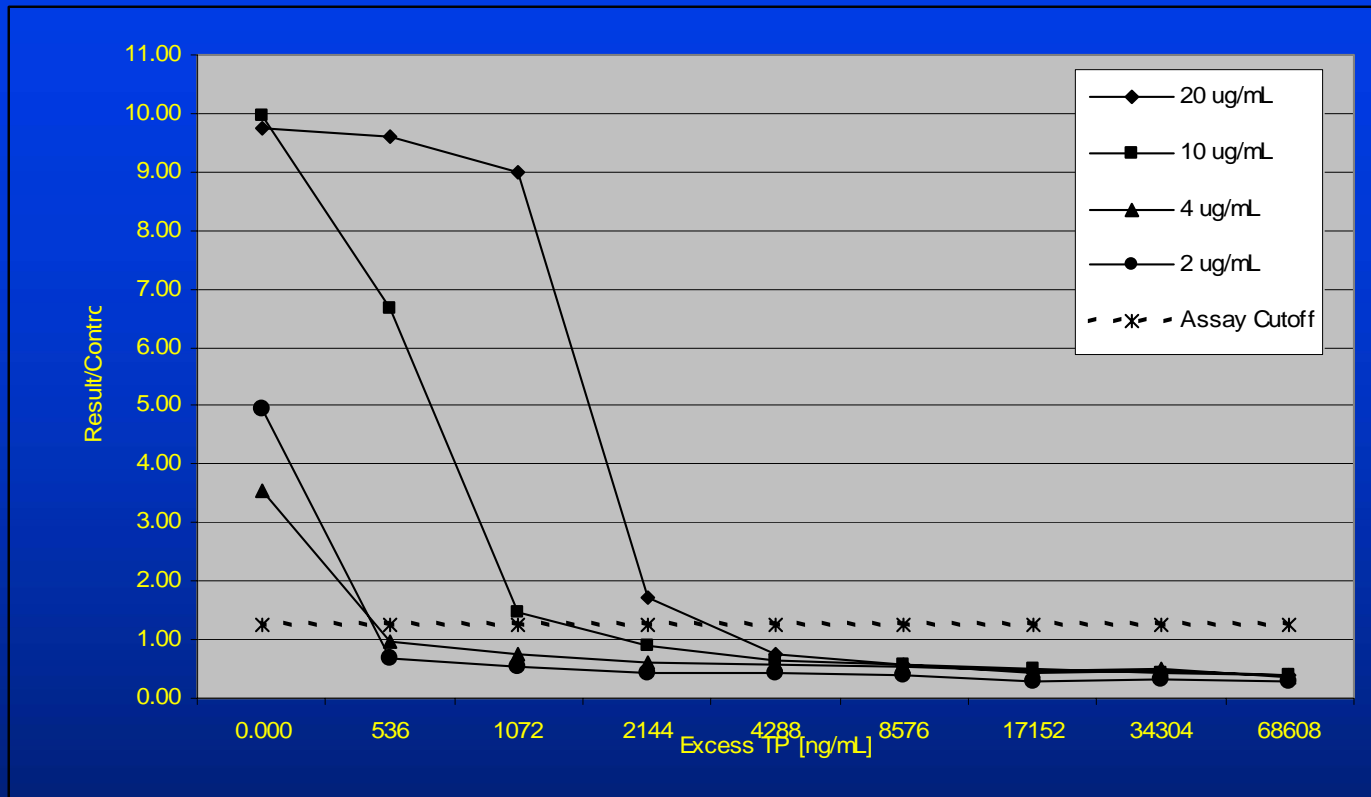
Post/Pre used because some individual animals yielded extremely low Ratio 1 values

Antibody Detection in Individual Monkeys



- Serum from 50 individual cynomolgus monkeys was tested treated and untreated with 2 $\mu\text{g}/\text{mL}$ anti-TP NAb at a final dilution of 15% in the Screening or Specificity Assay.
- A Ratio of Treated/Untreated samples (Post/Pre) was calculated and the Screening Ratio cutoff value was applied (1.26).
- Samples results were compared with Post/Pre and Specificity Ratio cutoff values.
- All samples treated with anti-TP NAb were above the Post/Pre Ratio cutoff.
- All samples treated with anti-TP NAb were below the Specificity Ratio cutoff except for three. These represent false negative samples.
- All untreated samples were below the Screening and Specificity Ratio cutoff (data not shown).

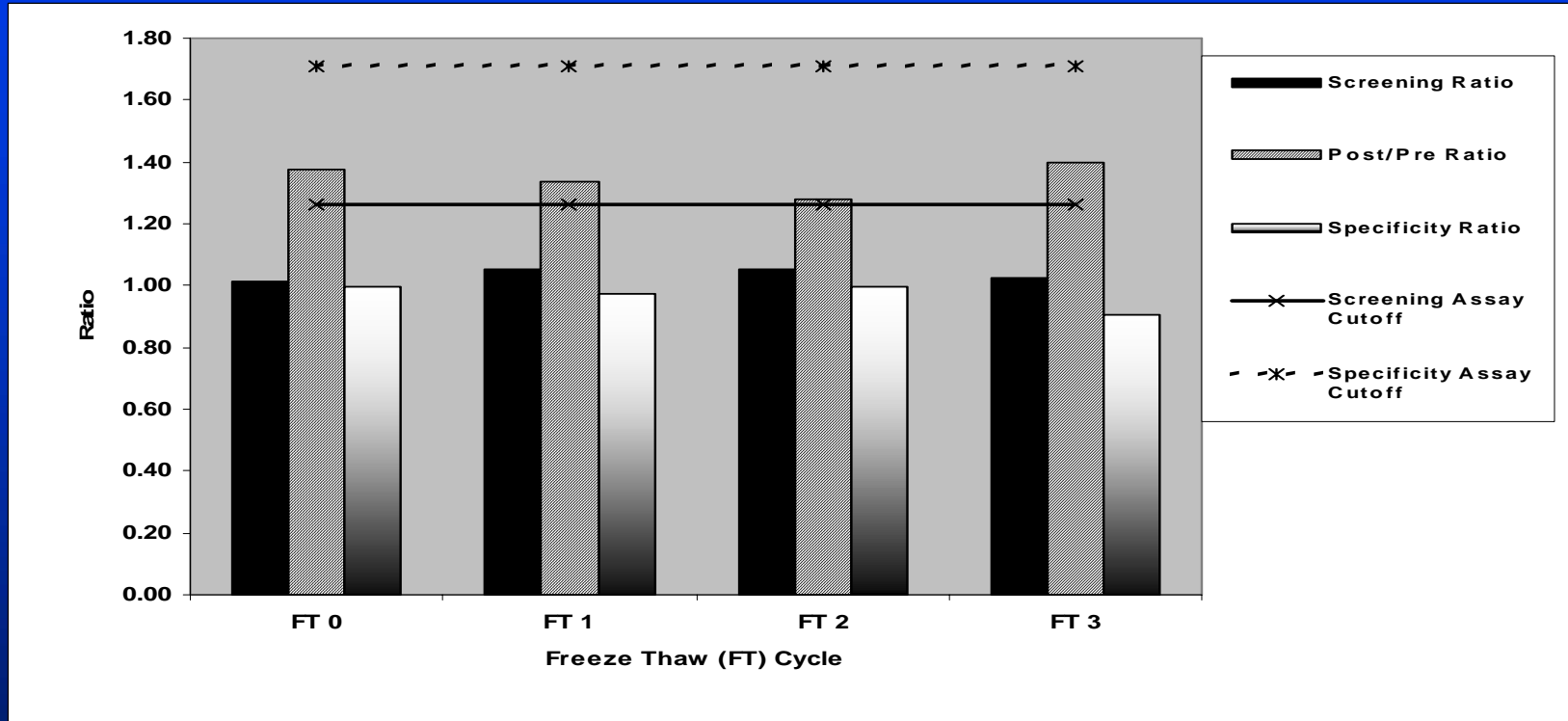
Excess Drug Interference



•Rabbit polyclonal anti-TP antibody was prepared in neat PCS at 20, 10, 4 and 2 $\mu\text{g}/\text{mL}$. At each concentration of antibody, the TP was titred from 68.6, 34.3, 17.15, 8.576, 4.288, 2.144, 1.072, 0.536 and 0 $\mu\text{g}/\text{mL}$ $\mu\text{g}/\text{mL}$ in neat PCS .

•The ability of the assay to detect 2 and 4 $\mu\text{g}/\text{mL}$ anti-TP antibody was inhibited by 80 ng/mL excess TP (536 ng/mL in neat serum). The assay is able to detect 10 $\mu\text{g}/\text{mL}$ anti-TP antibody in the presence of up to 160 ng/mL (1.07 $\mu\text{g}/\text{mL}$ in neat serum) excess TP and 20 $\mu\text{g}/\text{mL}$ in the presence of up to 320 ng/mL (2.14 $\mu\text{g}/\text{mL}$ in neat serum) excess TP.

Effects of Freezing and Thawing on Antibody Detection



- Positive control anti-TP antibody was added to PCS at the LOD of the assay (2 $\mu\text{g}/\text{mL}$).
- All samples containing antibody were positive up to 3 freeze thaw cycles while all untreated samples remained negative

Study Results

Animal	Dose Group	Immunoassay Result	Neutralizing Antibody Result
1	Low	Negative	Not Analyzed
2	Low	Negative	Not Analyzed
3	Low	Negative	Not Analyzed
4	Mid	Negative	Not Analyzed
5	Mid	Negative	Not Analyzed
6	Mid	Negative	Negative
7	High	Positive	Positive
8	High	Negative	Not Analyzed
9	High	Negative	Not Analyzed

•Serum Samples from Cynomolgus Monkeys dose with TP that were found to be Reactive in a Screening Immunoassay were tested in the Receptor-Binding Assay

•Of the samples tested one sample was found to be positive for neutralizing antibodies which correlated with the Immunoassay result.

Conclusions

A robust and sensitive receptor binding assay was developed for the detection of neutralizing antibodies to a TP in cynomolgus monkey serum

The assay was able to detect 2 $\mu\text{g}/\text{mL}$ rabbit polyclonal anti-TP antibody in monkey serum in the presence of 80 ng/mL TP.

The successful implementation of the assay detected antibodies in 1 animal from a preclinical pharmacokinetic study (11% incidence) .

Concluding Comments

- **Dependent upon quality of reagents**
- **Biological activity of reagents should be evaluated**
- **Sequence of assay steps extremely important**
- **Serum still an important factor**
- **Quick readout**
- **Assay development requires careful consideration**
- **Ratio of ligand/therapeutic important**
- **Not a slam dunk!**

Acknowledgements

- **Jason Pennucci**
- **Mike Moxness**
- **Gene Koren**
- **Jad Zoghbi**
- **Darin Asbury (Protein Sciences)**
- **John Thomas (Flow)**
- **Ling Wang (Research)**