

# APPROACHES AND LESSONS LEARNED IN ESTABLISHING HT-ADME ASSAYS

---

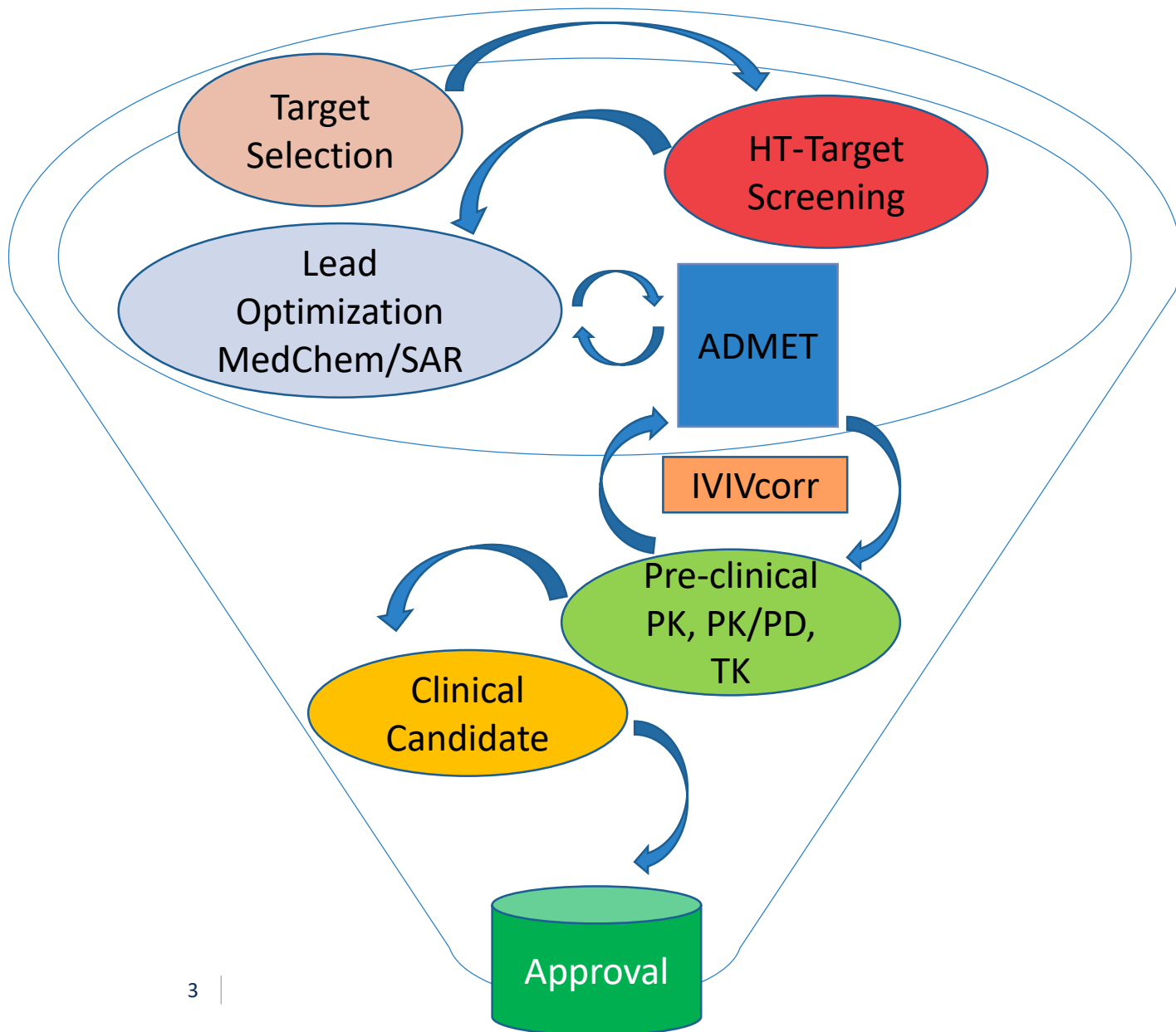
Solvo Meet-the-Experts Transporter Conference  
Adrian Sheldon, Charles River Labs, Worcester MA  
05-Sept-2019

EVERY STEP OF THE WAY

# PRESENTATION OUTLINE

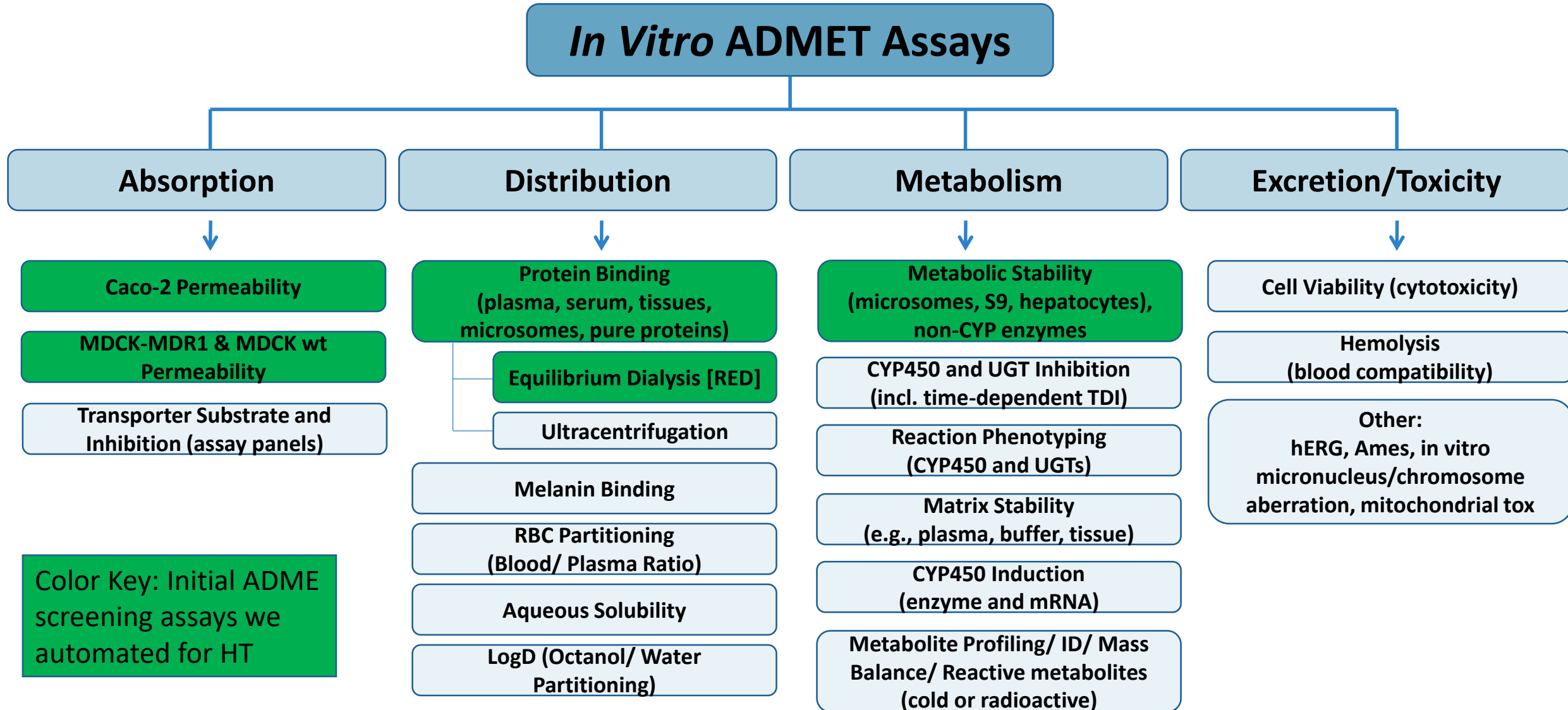
- WHY **ADME** AND WHY **HT**?
- APPROACH FOR SETTING UP HT-ADME PLATFORM
- QUALIFICATION ASSAY DATA / STATISTICAL ANALYSES TO DEMONSTRATE PERFORMANCE
- LESSONS LEARNED / SUMMARY

# IMPORTANCE OF ADME FOR DRUG DISCOVERY & DEVELOPMENT



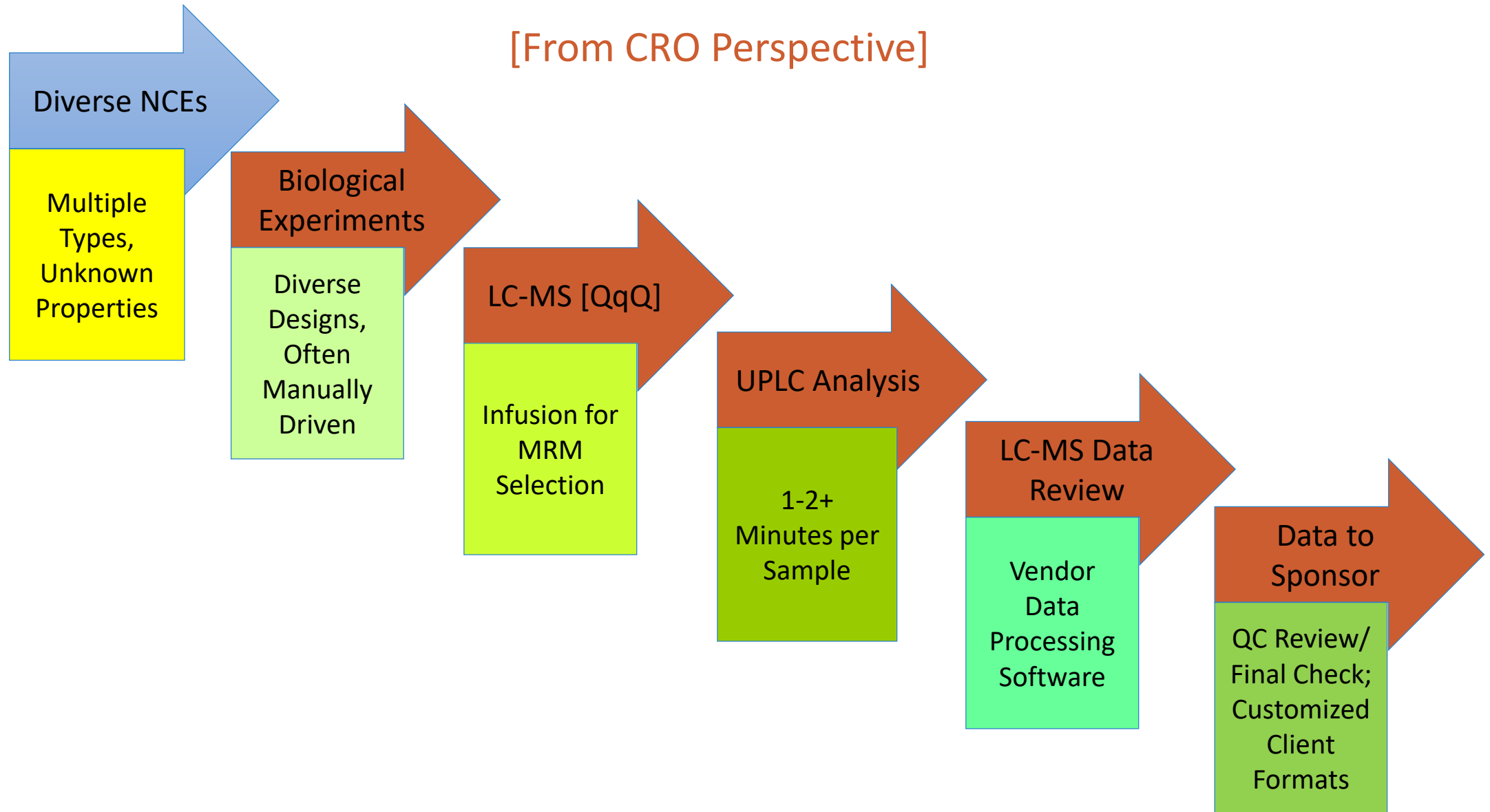
- *In vitro* ADME data are important for triaging compounds in early discovery to enable *in vivo* studies: PK, PK/PD, efficacy and TK studies, before IND selection
- Critical to optimize DMPK properties to enable successful outcomes in preclinical and clinical studies
- Saves time and money to establish an IV-IV-correlation during the discovery phase

# VARIETY OF IN VITRO ADME ASSAYS



# BOTTLE-NECKS WITH HIGH-VOLUME ADME NEEDS

[From CRO Perspective]



# HT-ADME PLAN TO ADDRESS BOTTLE-NECKS

- Life science industry needs capacity and rapid TAT to drive drug discovery programs
- Target objectives (based on clients' requests):
  - Focus on first tier assays:
    - Liver microsomal metabolic stability
    - Plasma protein binding (RED; Rapid Equilibrium Dialysis)
    - Permeability (Caco-2, MDCK-MDR1)
  - Fast data TAT:  $\leq 5$  Days from compound receipt to delivery of results
  - Capacity: ~500 compounds/week for primary screening assays
  - Reformatting: ability to cherry-pick compounds and to handle a variety of inputs (a challenge in CRO environment)
- Implement assay automation and HT-LC-MS/MS workflow (within ~8 month period)
- LC-MS/MS optimization and data
  - Facile MRM development and optimization (DQ)
  - Efficient data processing and review (leverage Gubbs GMSU)
- ADME LIMS
  - Efficient data storage, flexible calculation and reporting

# RESOURCE INVESTMENTS TO ENABLE HT-ADME

- Investments initiated in 2017 to enable HT-ADME capability to support both LT- and HT-ADME demands (capacity and flexible experimental designs)
- Liquid handling automation:
  - Evaluated 3 vendors in detail (hardware and software)
  - Purchased 2 Hamilton Vantage (2m; 8- and 96-channel; temp-control, shaking)
- HT-LC-MS/MS
  - Evaluated LC systems and visited labs with ADDA's; considered QE accurate mass
  - Purchased 2 Apricot ADDA with Agilent 1260s pumps and Sciex 5500 MS/MS
- 2 systems provide redundancy along with capacity (to minimize downtime)
- A small dedicated team with internal and external expertise to enable HT set-up and qualification/validation
- LIMS: Evaluated 3 vendors in detail, selected Edge BioRails/Morphit (UK)

# HT AUTOMATION (TO GENERATE AND ANALYZE SAMPLES)



## Hamilton Vantage robot [2m deck; N=2]

- Compound cherry-picking
- Heating, cooling, shaking, timed incubations
- Barcoding

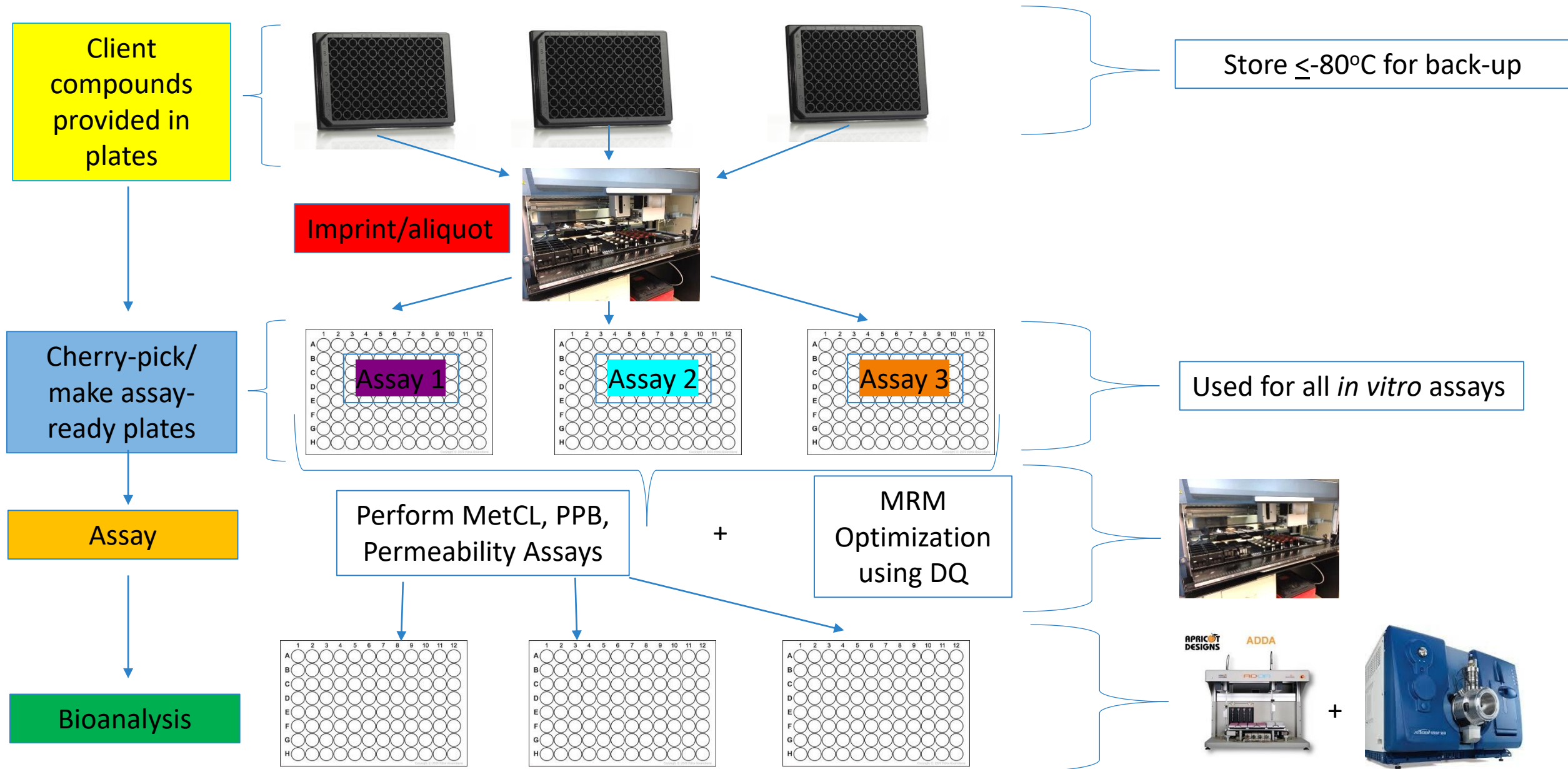
## ADDA LC-MS/MS [N=2]

- 96/384-well plates
- Automated MRM method development
- T/E and gradient modes
- Flexible LC column and mobile phase selections

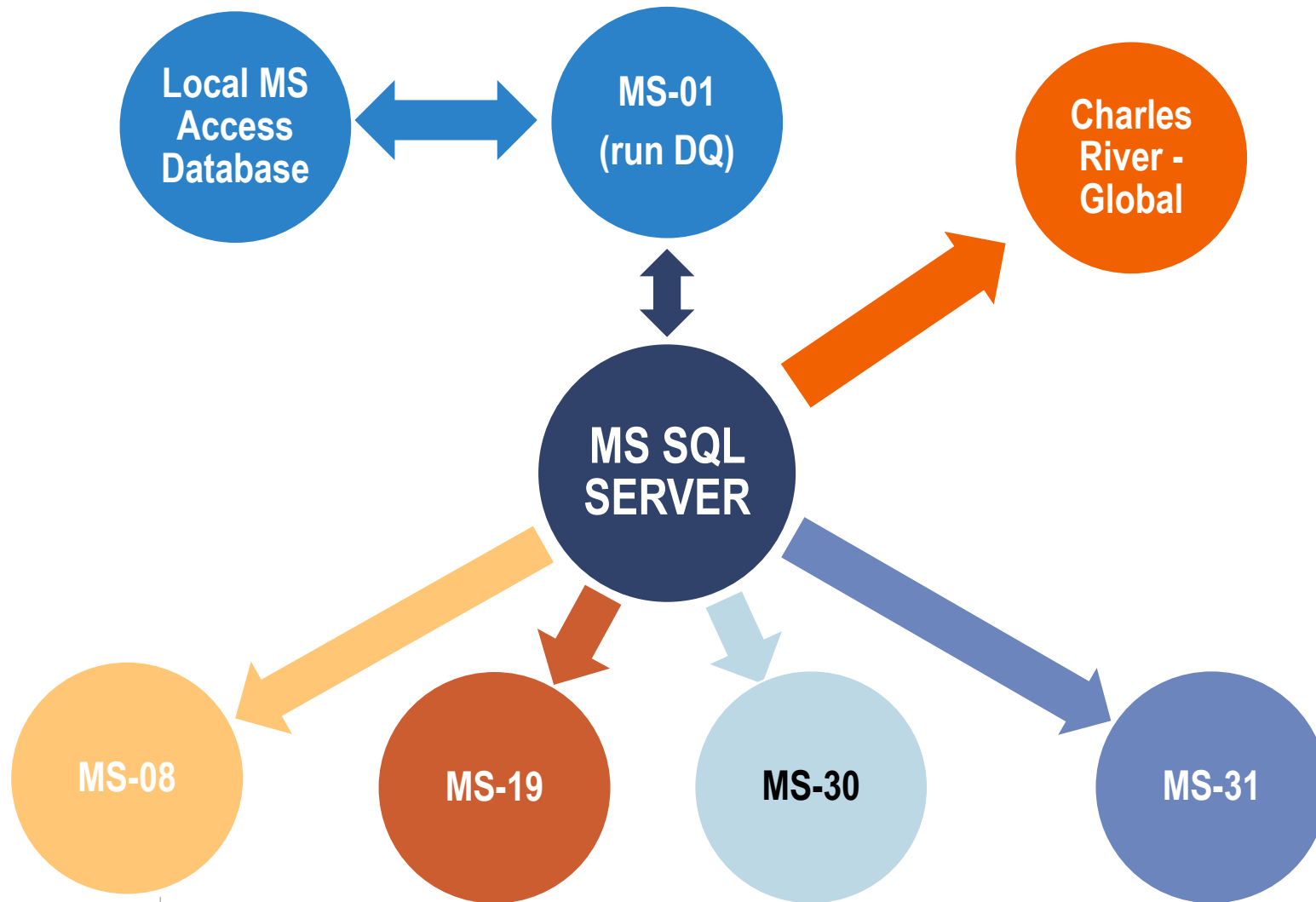




# HT-ADME WORKFLOW: SOURCE → ASSAY → BIOANALYSIS



# DISCOVERYQUANT<sup>®</sup>: MRM OPTIMIZATION



- Single database which serves as central repository of all compounds assayed within the *in vitro* ADME group
- Optimize once on a single MS (4000)
  - Run compound plates overnight
  - ~1 min/compound (fast tuning)
  - ~2 min/compound (fine tuning)
- Utilize the central database without having to repeat manual tuning and eliminate redundancies
- Global MS SQL server
  - Share MRM conditions
  - Review spectra

# ADDA HT-LC/MS/MS SYSTEMS

➤ Dual-arm, high-speed multiplexing autosampler

- ~10 seconds/injection

➤ Two modes of operation:

- Trap and elute
  - TIS for “clean” samples
  - APCI for plasma samples
- Gradient

➤ Software: Sound Analytics

- DiscoveryQuant (DQ) for automated MRM optimization
- LeadScape for acquisition

APRICOT  
DESIGNS

ADDA



+

UFLC Pumps



ABI 5500 MS

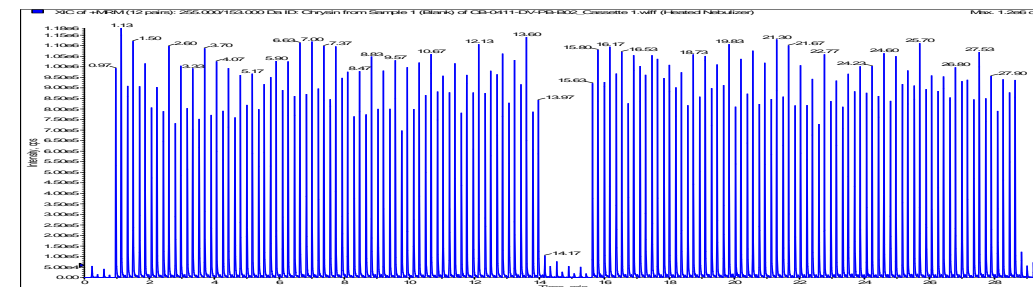


# ADDA HT-LC-MS/MS PLATFORM



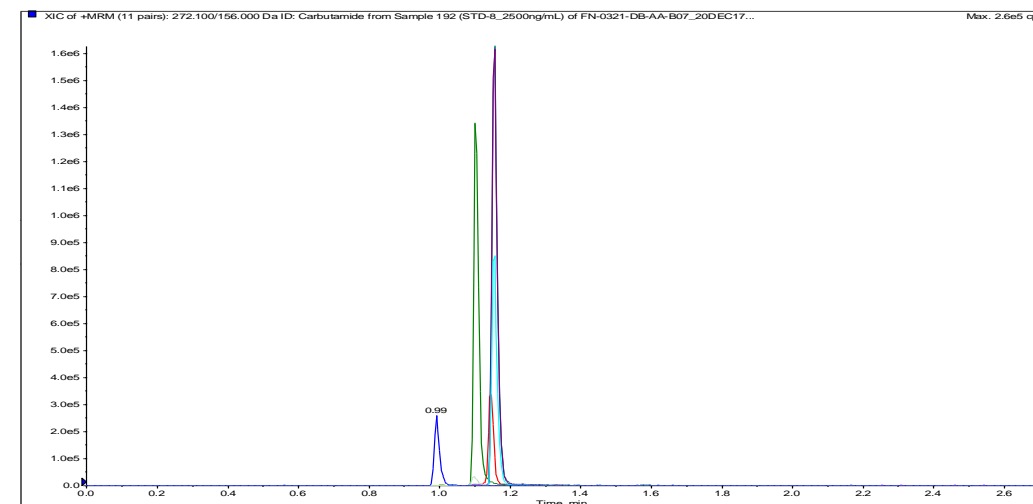
T/E

- Performs T/E and gradient modes on the fly
  - 384-well, 96-well, etc.
  - High-quality peak shape
- Diverse column chemistries [T/E]
  - RP C<sub>18</sub>: 10 mm to 30 mm, 3 μM
  - HILIC for polar molecules
- ADDA proven to be very robust and reliable
- More flexible than RapidFire
- Need data processing software to be able to integrate multiple peaks and analytes in a single file



HT-Permeability Assay  
[160 Samples Analyzed in 30 minutes]

Gradient



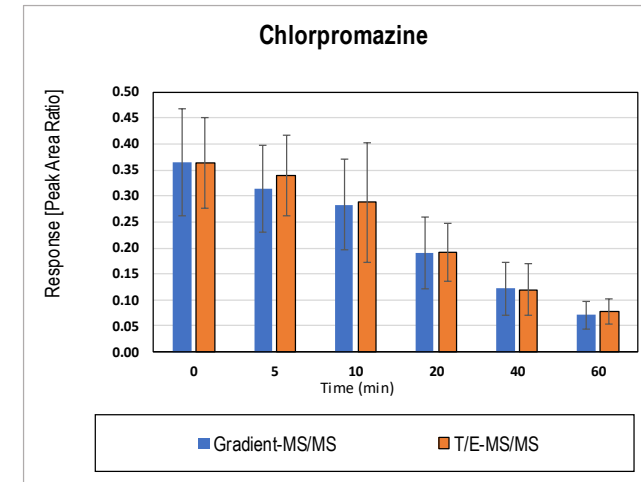
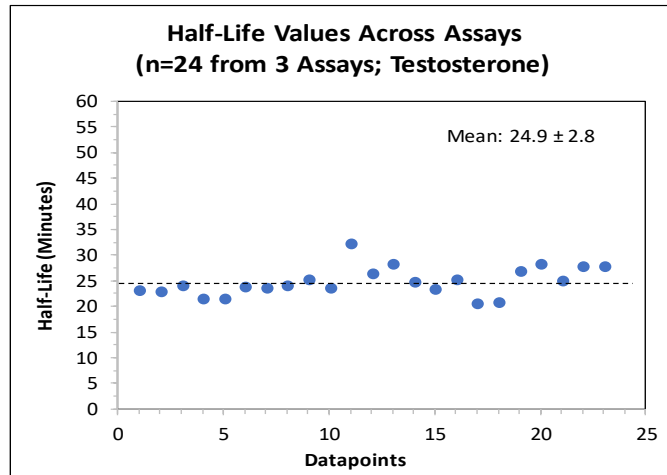
# STEPS FOR PROCESS QUALIFICATION

- Define final assay procedure
- Optimize pipetting (liquid classes, liquid into dry plate or solution, lab temp/humidity)
- Optimize workflow efficiency (deck layout, minimize back&forth movements)
- Run through procedure in simulation mode, then blank reagents, then real assay
- Demonstrate results agree with expected values (literature and internal historical)
- Assess data consistency across runs, time
- Demonstrate agreement between manual vs. robotic, and gradient vs. T/E
- Assess performance using known drug compounds, plus test “real world” discovery compounds

# QUALIFICATION: METABOLIC STABILITY ASSAY

# HT-ADME ASSAY VALIDATION RESULTS FOR METABOLIC STABILITY ASSAY

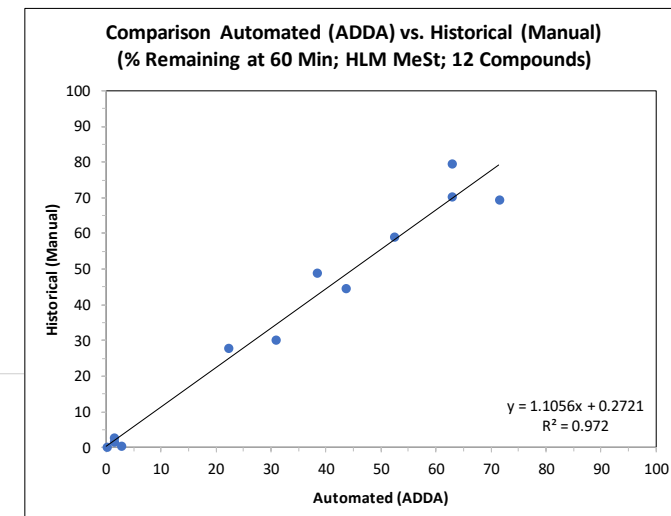
➤ Demonstrated consistent robot and HT-LC/MS performance:



T/E vs. Gradient:  
1  $\mu$ M compound  
0.5 mg/mL HLM  
37°C  
Time course

➤ Demonstrated equivalent results between automated assay and manual (historical) MeSt assay:

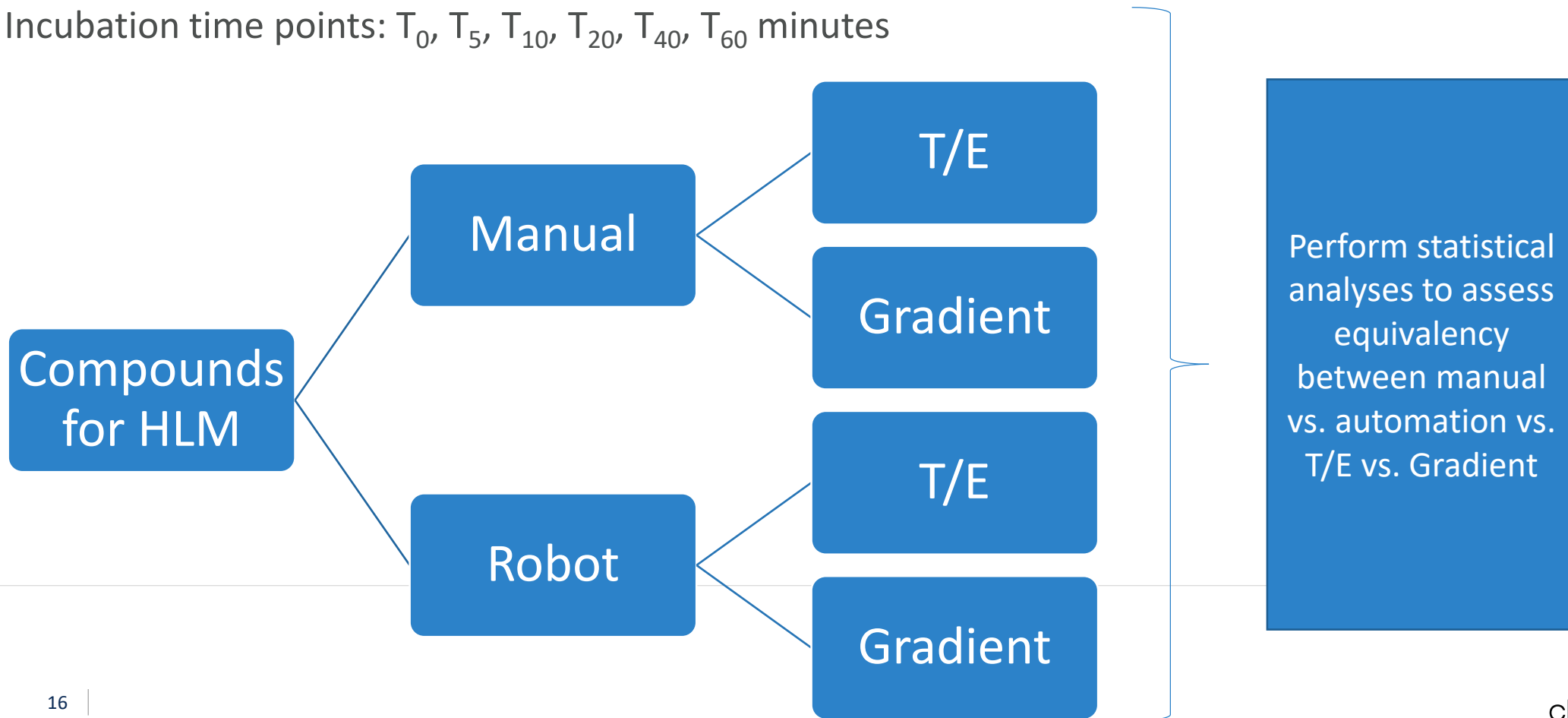
Compound ID	Mean % Remaining at T <sub>60min</sub>		Mean T 1/2 (min)		Mean CL <sub>int</sub> (mL/min/kg)	
	Automated	Manual	Automated	Manual	Automated	Manual
7-EC	2.8 ± 2.5	1.8 ± 0.2	9.7 ± 1.4	9.1 ± 0.3	131 ± 15	137 ± 4
Testosterone	22.3 ± 6.3	24.0 ± 4.6	24.9 ± 2.8	29.0 ± 2.7	50.7 ± 5.5	43.3 ± 3.9
Imipramine	52.4 ± 4.2	57.8 ± 6.6	66.7 ± 10.9	79.0 ± 17.0	19.2 ± 3.0	16.4 ± 3.2
Terfenadine	1.5 ± 0.8	1.5 ± 1.1	11.3 ± 1.7	10.4 ± 1.8	113 ± 17	123 ± 20
Quinidine	62.9 ± 5.9	70.3 ± 5.7	103 ± 31	108 ± 25	13.3 ± 4.0	12.1 ± 3.1



[Equivalency also demonstrated with 40 sponsor discovery compounds]

# QUALIFICATION USING UNKNOWN NCE'S

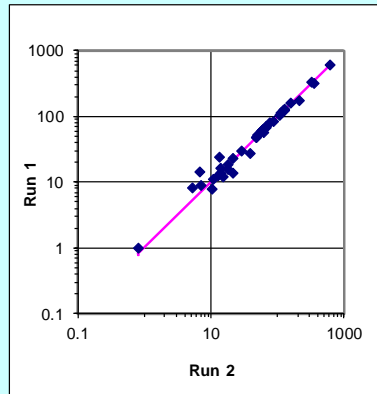
- 40 discovery compounds submitted by Sponsor for HLM
- Concentration: 1  $\mu$ M
- Incubation time points:  $T_0$ ,  $T_5$ ,  $T_{10}$ ,  $T_{20}$ ,  $T_{40}$ ,  $T_{60}$  minutes





# HLM: MANUAL VS AUTOMATION FOR NCEs [N=40] T/E & GRADIENT [CL<sub>INT</sub>]

T/E vs. GRADIENT with 45-degree Line

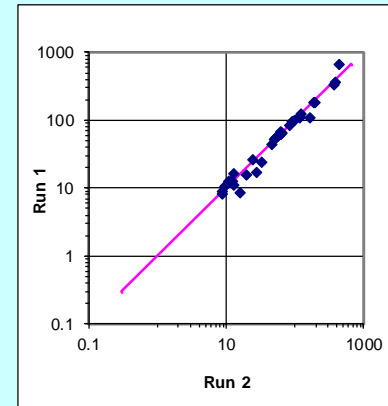


Correlation Coefficient of **LOG-LOG** Plot  
= 0.988

15 points below the 45-degree line  
23 points above the 45-degree line

Manual  
Experiment

T/E vs. GRADIENT with 45-degree Line

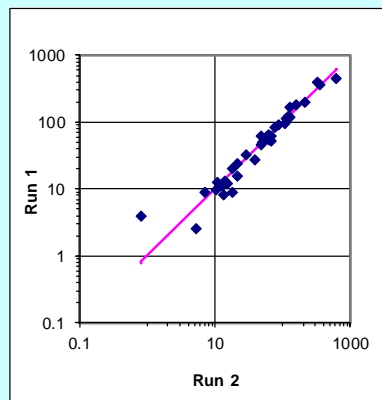


Correlation Coefficient of **LOG-LOG** Plot  
= 0.987

16 points below the 45-degree line  
19 points above the 45-degree line

Automation  
Experiment

Robot vs. Manual with 45-degree Line

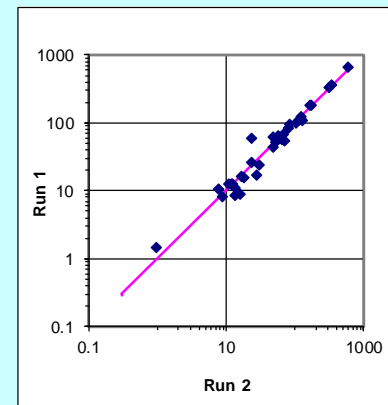


Correlation Coefficient of **LOG-LOG** Plot  
= 0.962

19 points below the 45-degree line  
19 points above the 45-degree line

Gradient Analysis

Robot vs. Manual with 45-degree Line



Correlation Coefficient of **LOG-LOG** Plot  
= 0.980

17 points below the 45-degree line  
20 points above the 45-degree line

T/E Analysis

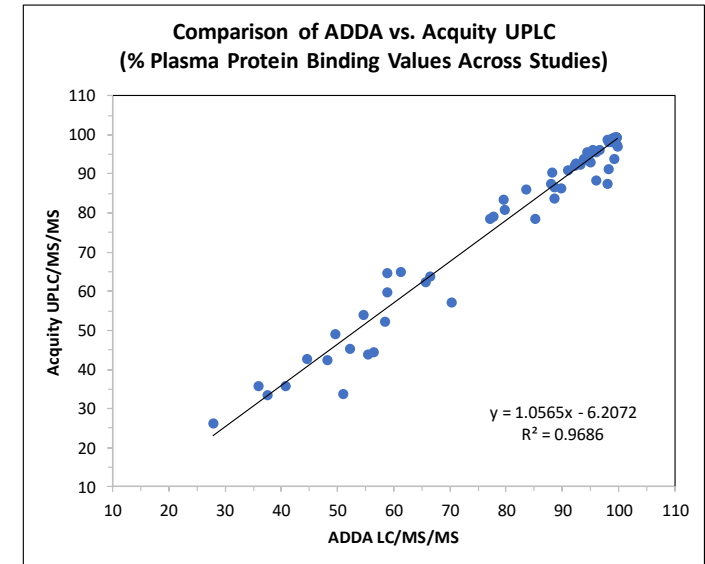
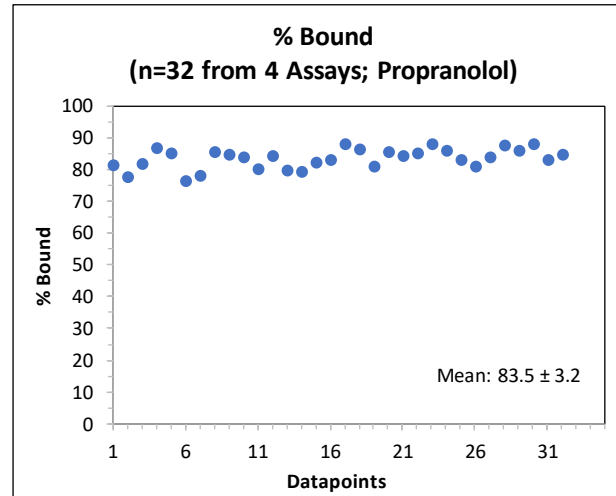
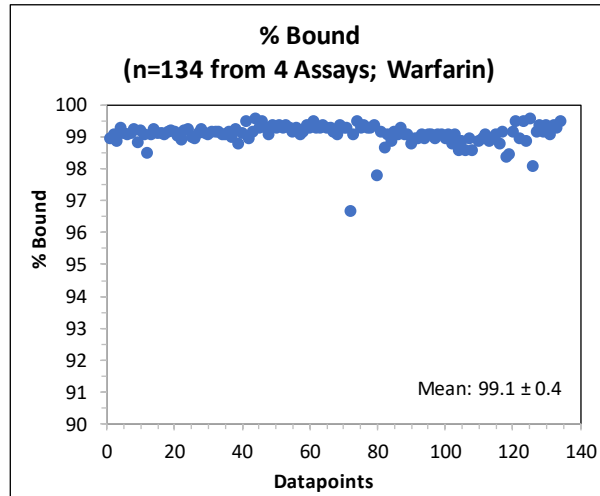
# QUALIFICATION: PROTEIN BINDING ASSAY (RED)

# HT-RED PLASMA PROTEIN BINDING

- Typically test 1-2  $\mu\text{M}$  compound (from 1000X DMSO stock)
- Replicates: N=1 to 3
- Incubation time: 6 hrs with gentle mixing in 5%  $\text{CO}_2$  at 37°C (to provide longer time to reach equilibrium, and to maintain pH)
- Matrices: Mouse, rat, dog, monkey, human plasma
  - Human lots are pre-screened (warfarin % binding, common control compounds, since we have noticed more unacceptable lots in recent years)
  - May be valuable to pre-screen with AGP binders too (plasticizers etc. can affect)
- Often include a concurrent matrix stability and recovery control ( $T_0$  vs  $T_6$ )

# HT-ADME ASSAY VALIDATION RESULTS: RED PPB ASSAY

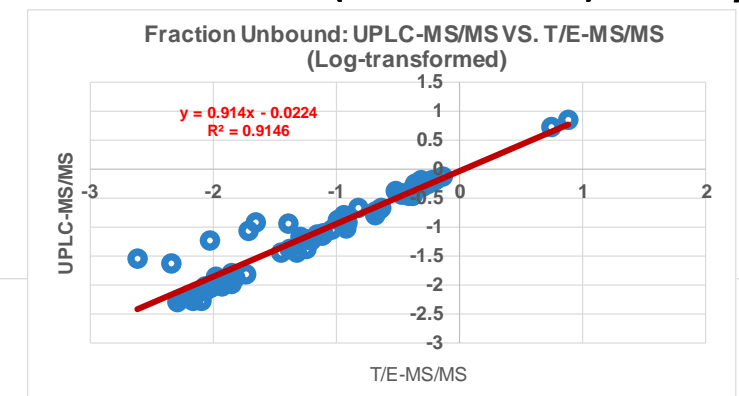
➤ Demonstrated consistent robot performance:



➤ Demonstrated equivalent results between automated assay and manual (historical) assay:

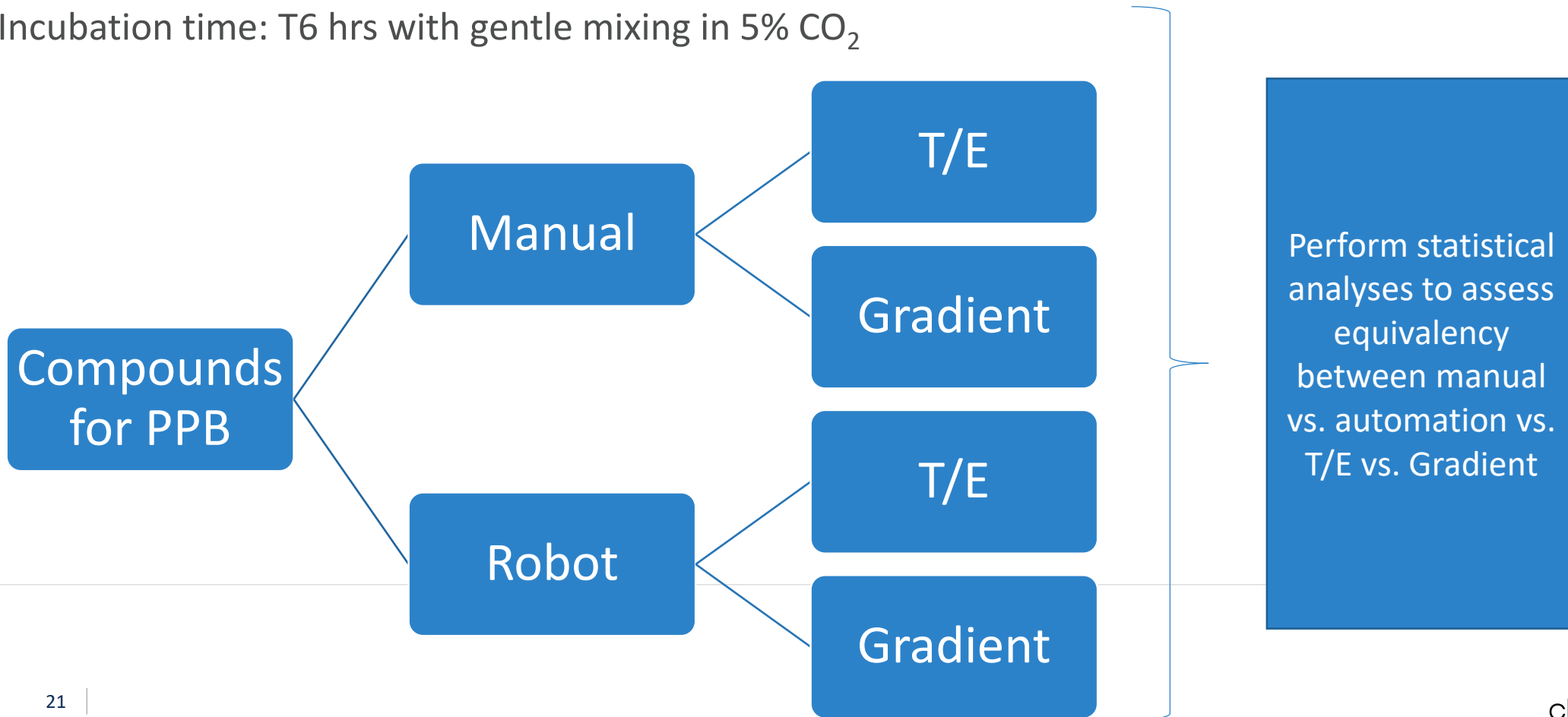
Compound ID	Mean % Bound	
	Automated	Manual
Chlorpromazine	$97.3 \pm 1.4$	$95.8 \pm 5.1$
Digoxin	$47.7 \pm 18.3$	$45.6 \pm 12.2$
Propranolol	$83.5 \pm 3.2$	$83.8 \pm 0.6$
Verapamil	$87.4 \pm 0.8$	$87.7 \pm 3.6$
Warfarin	$99.1 \pm 0.4$	$99.1 \pm 0.3$

Equivalency also demonstrated with 68 sponsor discovery compounds:



# PPB [RED] VALIDATION USING UNKNOWN NCES

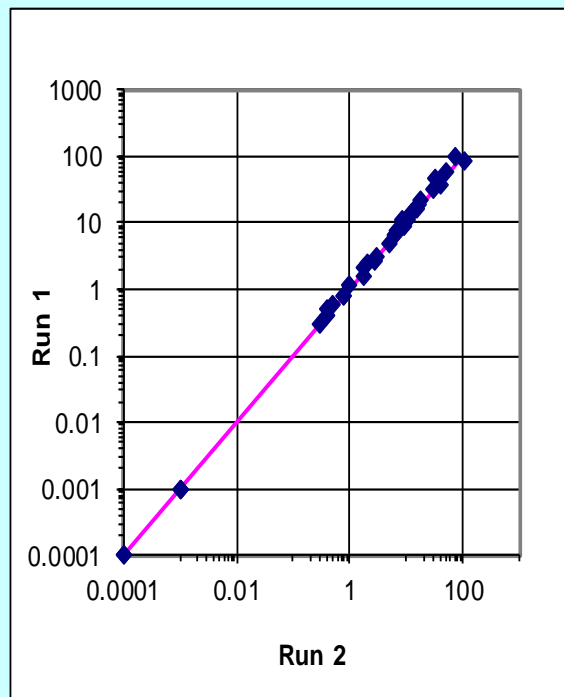
- 40 discovery compounds submitted by Sponsor for PPB
- Concentration: 1  $\mu\text{M}$
- Incubation time: T6 hrs with gentle mixing in 5%  $\text{CO}_2$



# PPB [RED] USING AUTOMATION: T/E AND GRADIENT ANALYSIS

## Fraction Unbound

T/E vs. Gradient with 45-degree Line

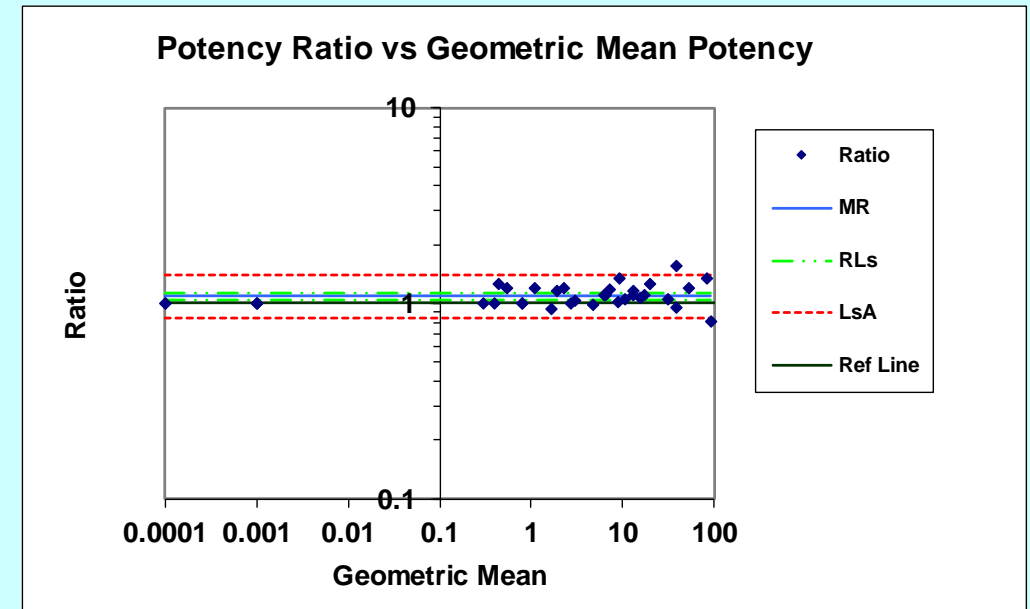


Correlation Coefficient of **LOG-LOG** Plot  
= 1.000

4 points below the 45-degree line  
21 points above the 45-degree line

Automation

PPB [RED] AUTOMATION Assay T/E & Gradient

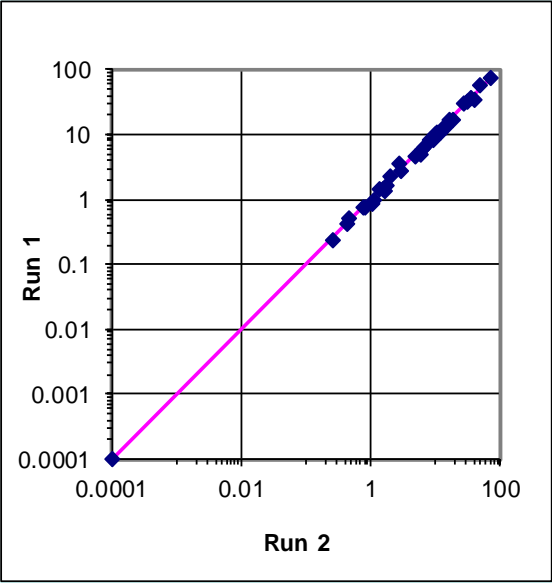


Statistical Analysis of "Bland-Altman" Plot			
Equivalence Test Results		Reproducibility Test Results	
N	35	MSR (Within-Run)	1.28
MR	1.09	LsA	0.85
RLs	1.04		1.39
	1.13	SD	0.0380
Sig Diff Between Runs Test, p =	0.0004		

# PPB [RED] USING MANUAL: T/E AND GRADIENT ANALYSIS

## Fraction Unbound

T/E vs. Gradient with 45-degree Line

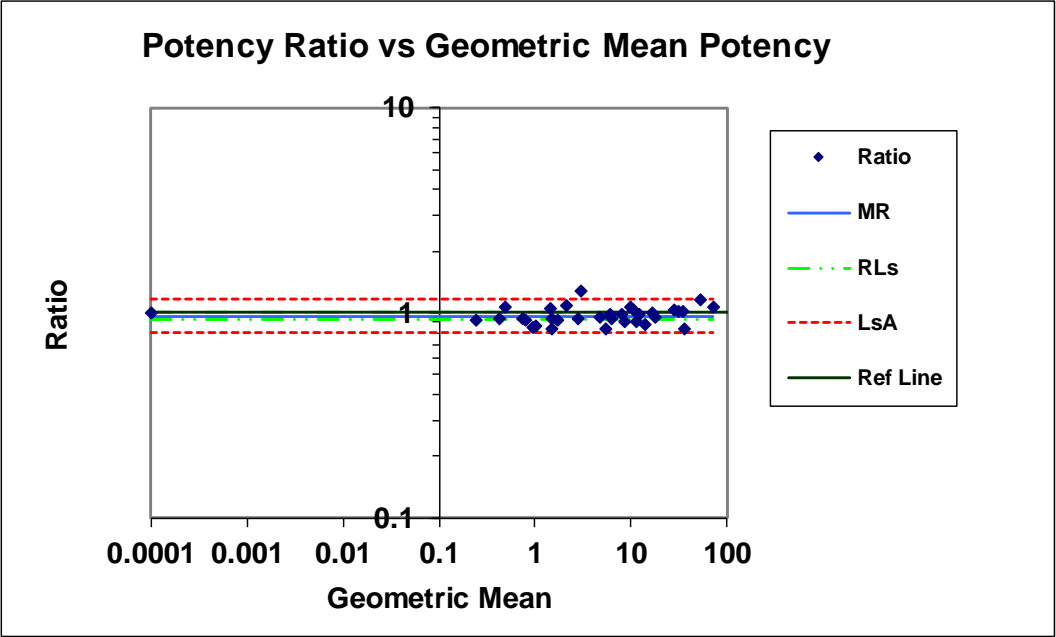


Correlation Coefficient of **LOG-LOG** Plot = 0.999

21 points below the 45-degree line  
12 points above the 45-degree line

Manual

PPB [RED] MANUAL Assay T/E & Gradient



T/E and Gradient Analysis Demonstrated Equivalent Performance By Deploying ApCI for T/E analysis

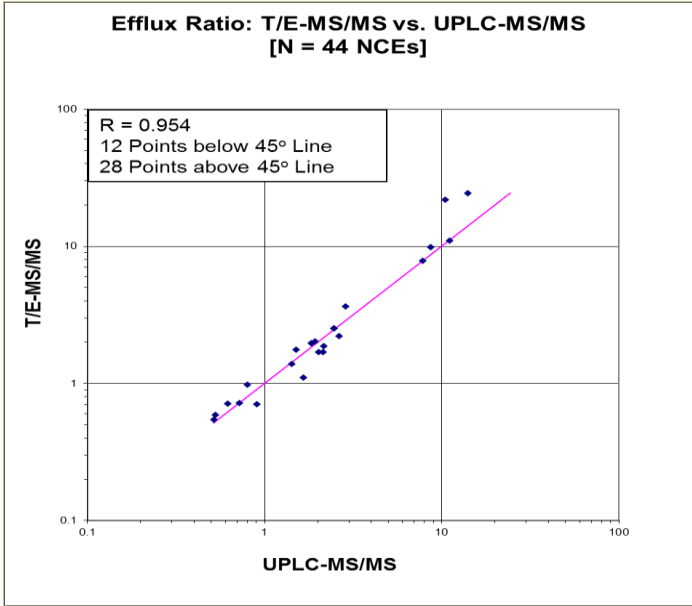
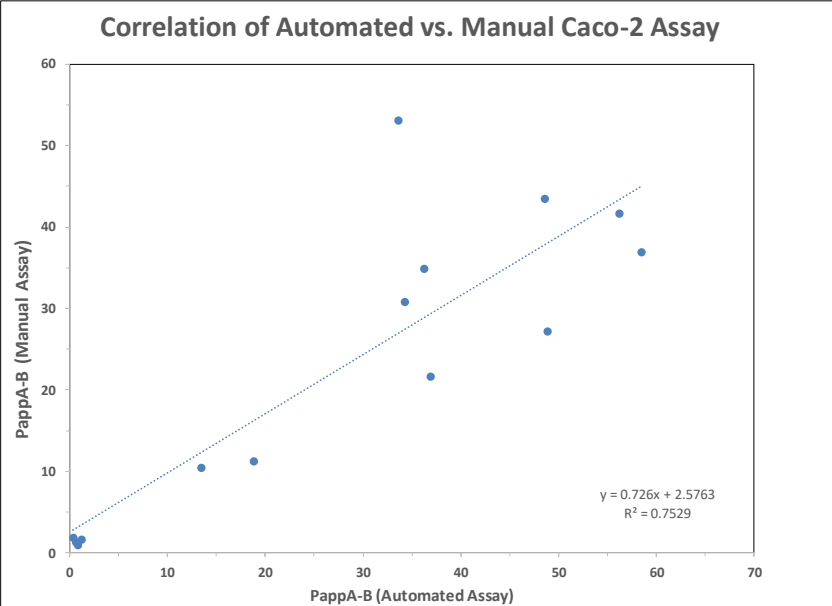
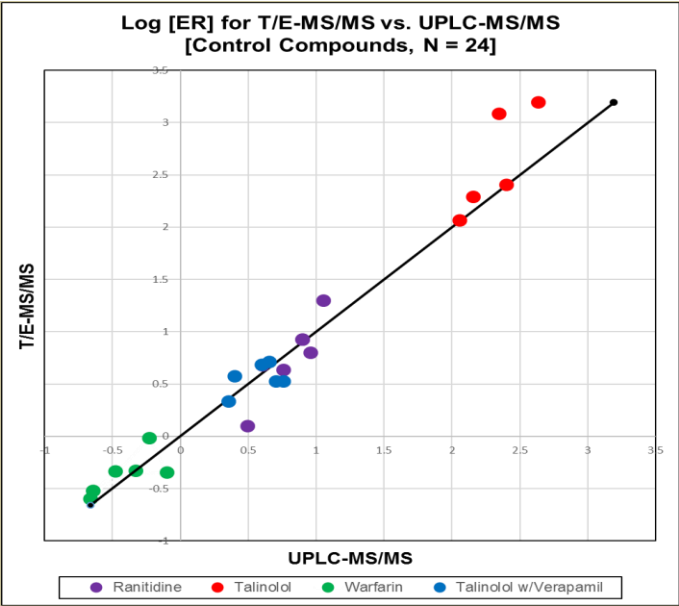
Statistical Analysis of "Bland-Altman" Plot			
Equivalence Test Results		Reproducibility Test Results	
N	34	MSR (Within-Run)	1.21
MR	0.97	LsA	0.80
RLs	0.94		1.17
	1.00	SD	0.0294
Sig Diff Between Runs Test, p =	0.0487		

# QUALIFICATION: PERMEABILITY ASSAY



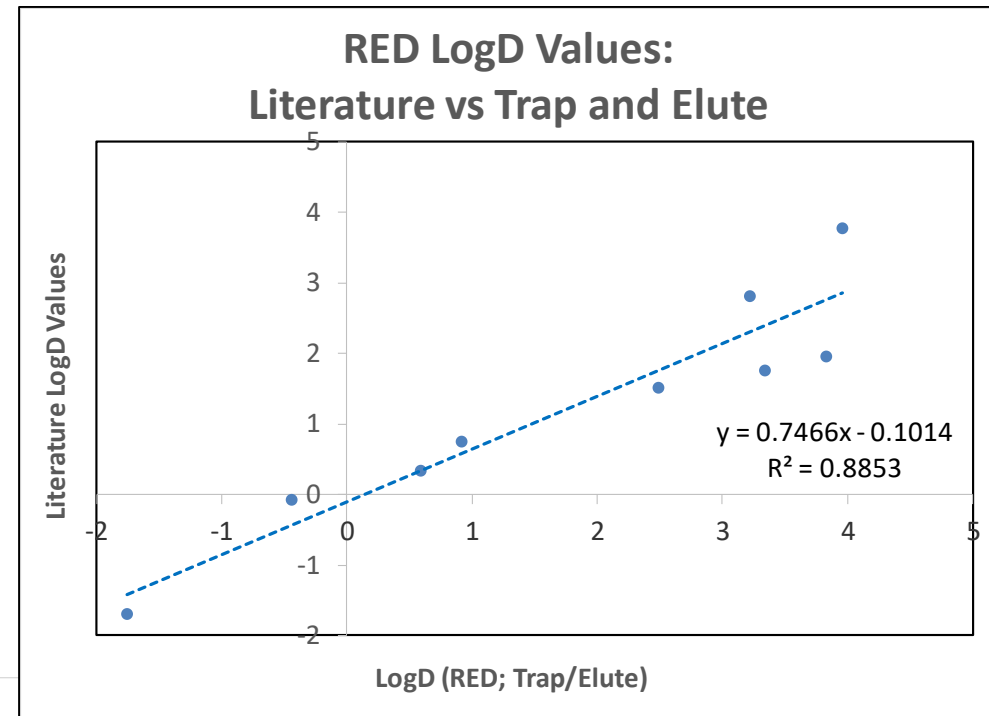
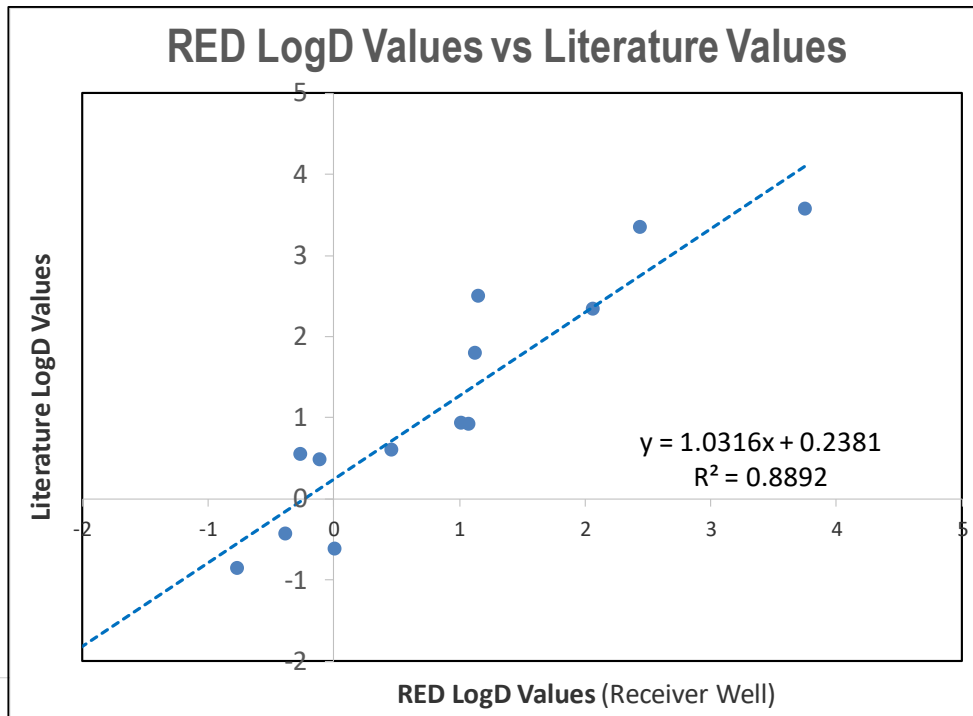
# HT-ADME ASSAY VALIDATION RESULTS FOR PERMEABILITY ASSAY

- Typically 10 μM, n=3, 2 hours incubation
- Demonstrated consistent performance (gradient vs T/E):

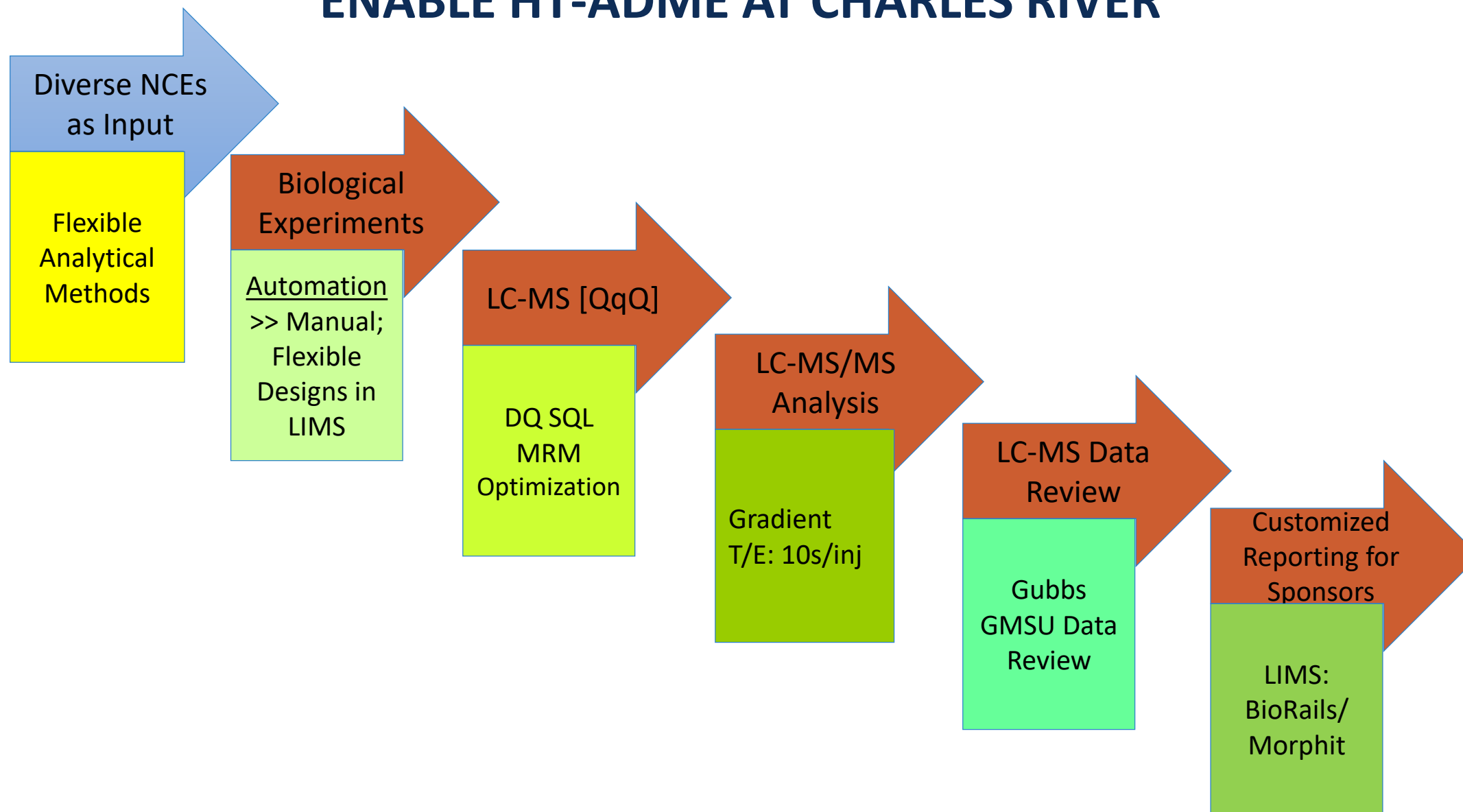


# OTHER HT-ADME ASSAY PROJECTS

- Have added additional assay capabilities (CYP inhibition, aqueous solubility, etc.)
- Working on logD (octanol/water) assay using RED units (spike PBS at 100  $\mu$ M, dialyze vs. octanol, sample PBS after 18 hours); format easier to automate than shake-flask.



# SUMMARY: BOTTLE NECKS ADDRESSED TO ENABLE HT-ADME AT CHARLES RIVER



# LESSONS LEARNED

## Liquid Handling Automation

- 1) Hamilton hardware is solid, dependable (following some tweaks after installation)
- 2) Should have looked into variety and details of available hardware options in more detail (some accessories were more useful than others)
- 3) Hamilton Venus (VoV) software is very powerful but was more complex and longer learning curve than expected (initial vendor assistance was very important; need internal expert; found different programmers have different styles)
- 4) Pipetting accuracy is very dependent on parameters (e.g., liquid class, speed, height above well or cells, etc.); use real reagents vs. water for testing; in hindsight would have been helpful to define optimal pipetting parameters and plasticware earlier in process

## HT-LC-MS/MS; Data Management

- 1) ADDA hardware observed to be solid, reliable
- 2) Slight stream differences observed, but is mitigated by running a complete set of samples for each compound within a single stream only
- 3) Fast data processing is key to reduce bottle-neck
- 4) Evaluated MultiQuant, LeadScape and GMSU, etc. (often using GMSU)
- 5) Evaluated diverse ADME LIMS options; critical to be track compounds in/data out; decide how much flexibility is required (decided to implement BioRails/Morphit; very happy with choice)

# ACKNOWLEDGMENTS

## Charles River:

- HT-*In Vitro* ADME team
  - David Plourde
  - Patty Walton
  - Allison Lewia
  - Jakal Amin
  - Guofeng Ye
  - Sarah Meloche
- *In Vitro* ADME team
- Charles River-Data Science
  - Viswanath Devanarayan
- Many other colleagues (IT)

## Software/ Equipment Vendors:

- Apricot
- SoundAnalytics
- Gubbs, Inc. (Larry Elvebak)
- Edge/BioRails (Andrew Lemon)

Any further questions:

[adrian.sheldon@crl.com](mailto:adrian.sheldon@crl.com)

508-471-5527