



The Next Frontier in ADME Science: Predicting and Verifying Tissue Drug Exposure

Jashvant (Jash) Unadkat

Milo Gibaldi Endowed Professor

Dept. of Pharmaceutics

School of Pharmacy

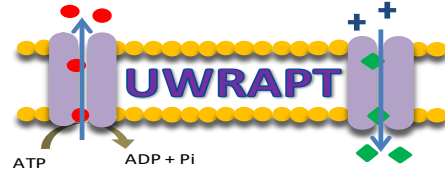
University of Washington

Seattle, WA

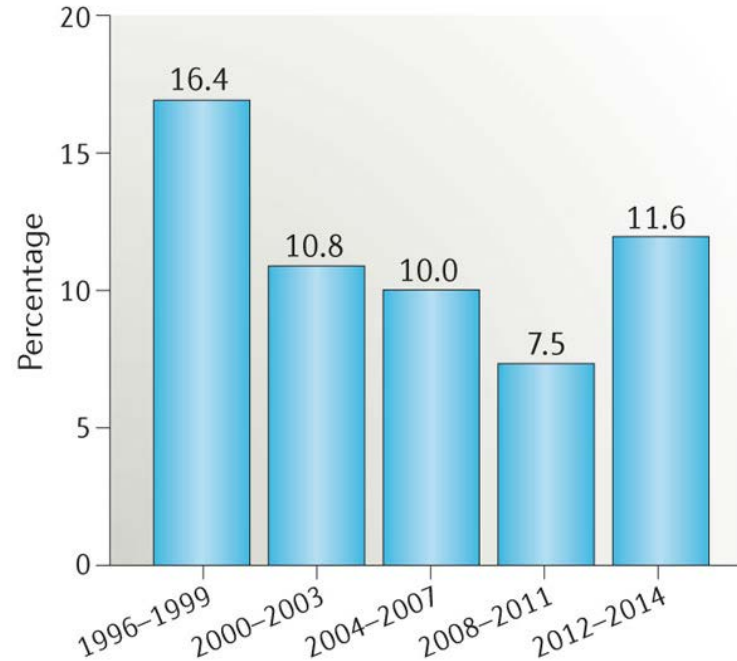
jash@uw.edu



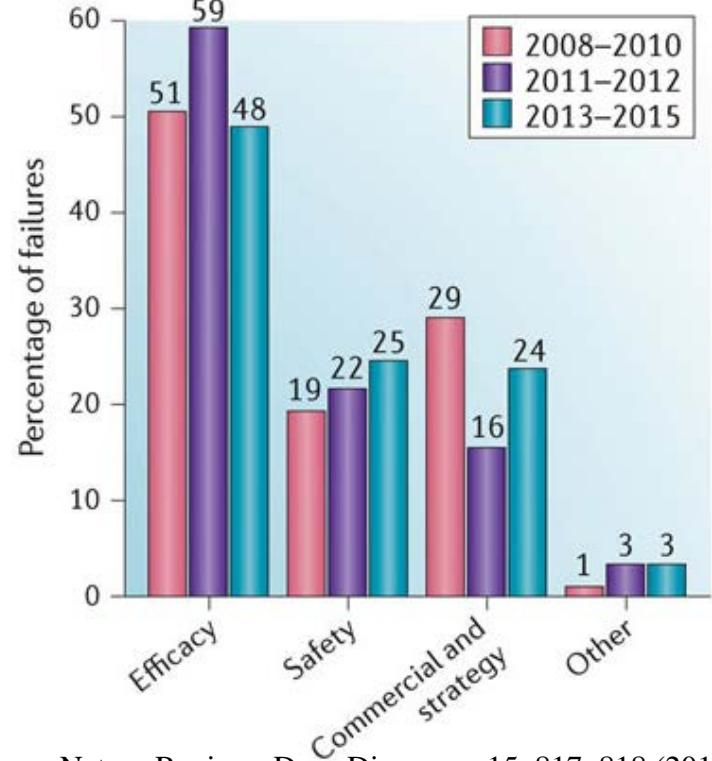
GAPS in ADME: Drug Development Failure Rate and Reasons



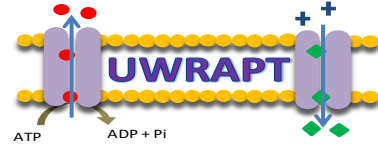
b Cumulative success rate Phase I to launch
Percentage likelihood of moving from Phase I to launch



b Reason for failure in phase II

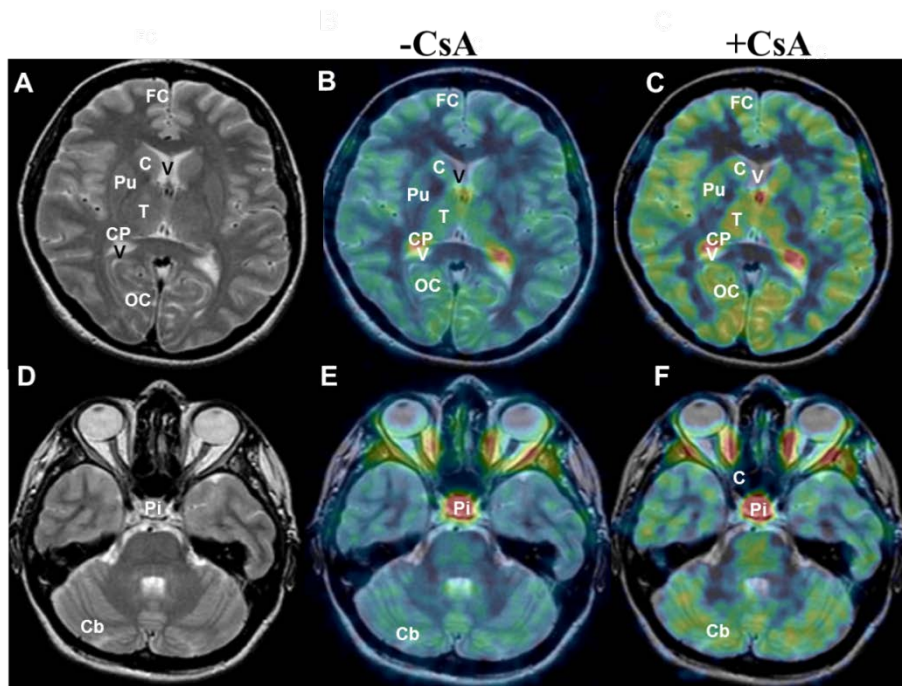


One Possible Reason for Lack of Drug Efficacy & Safety



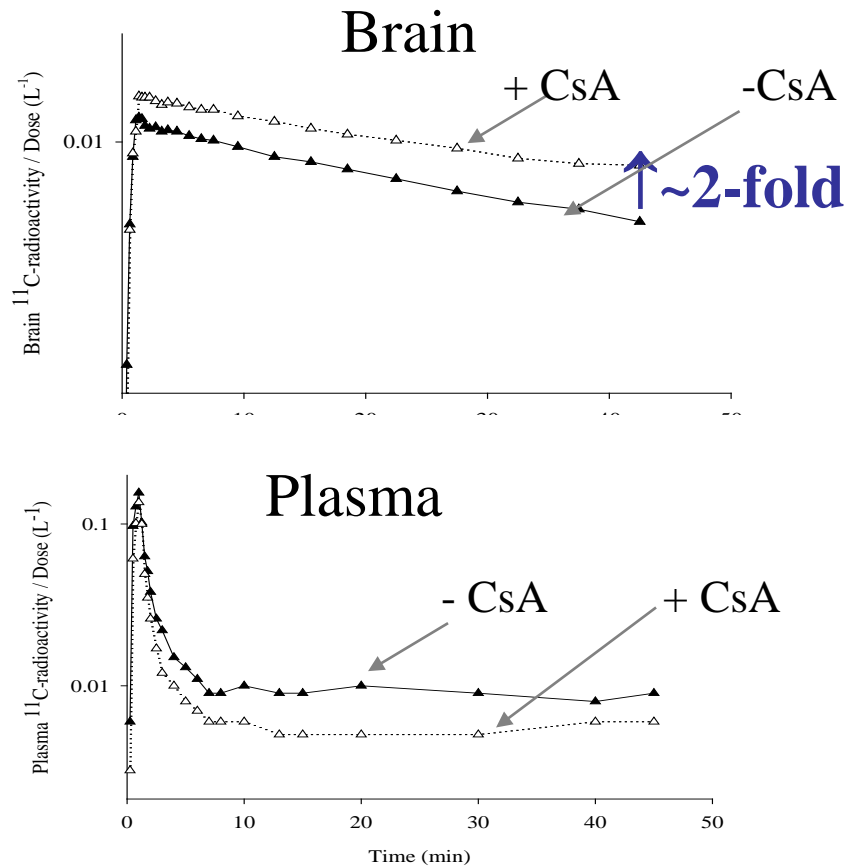
- Unable to routinely measure (or accurately predict) tissue drug conc.-time profile for drugs that are **transported** across tissue:blood barrier
- Transporters at the tissue:blood barrier (e.g. blood:brain barrier, BBB, liver:blood barrier)
 - Unbound tissue conc \neq unbound plasma conc. i.e. **asymmetry** in tissue:blood conc.
 - Impact differs between non-eliminating (e.g. brain) and eliminating organ (e.g. liver)

Asymmetry in Drug Conc. at the Human Brain:Blood Barrier: P-gp Efflux of ^{11}C -Verapamil

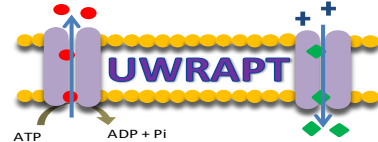


^{11}C -verapamil $\text{AUC}_{\text{brain}}:\text{AUC}_{\text{blood}}$ (20 min) - 0.42 ± 0.04

Eyal et al., *Clin Pharmacol Ther.* 2010

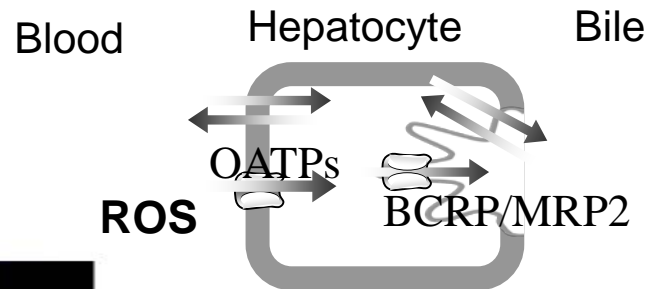
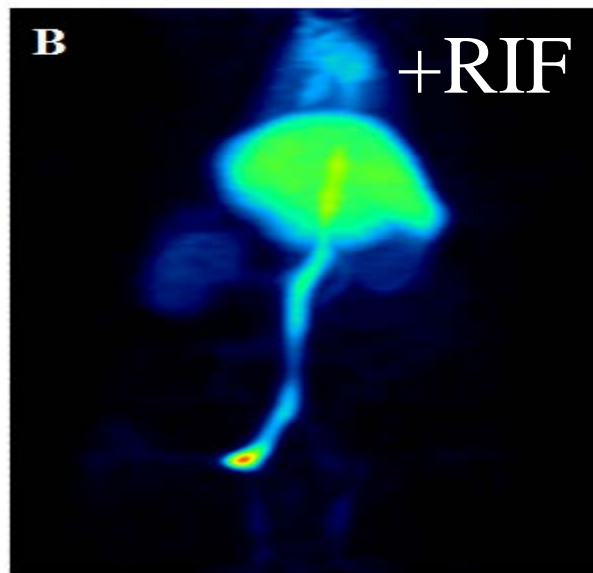
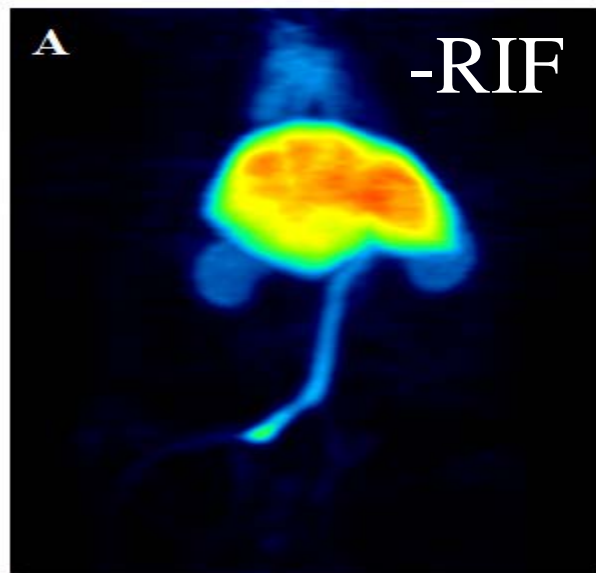


Asymmetry in Hepatic:Blood Conc. of ^{11}C -Rosuvastatin in the Rat

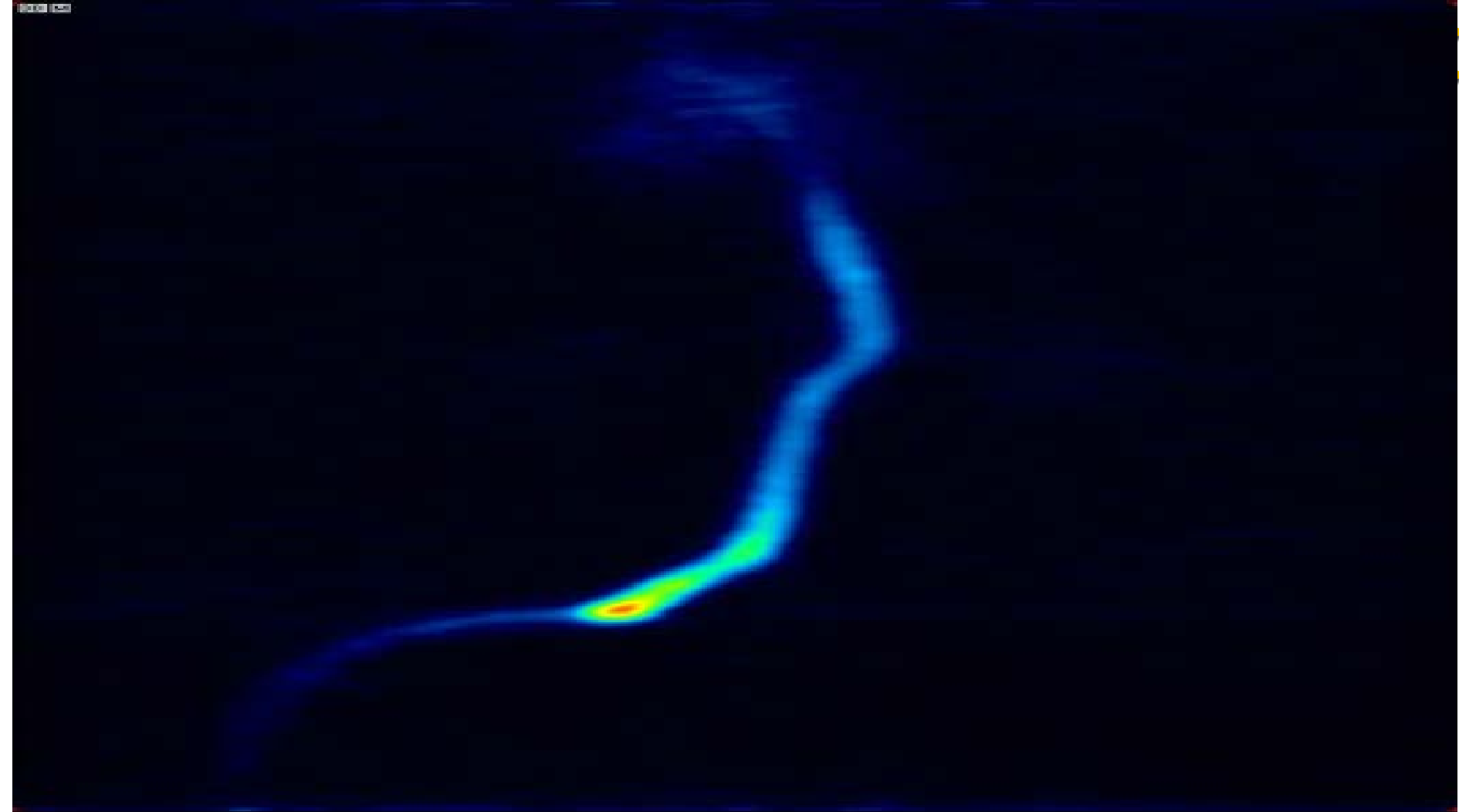


Coronal 2 min SUV images of ^{11}C -Rosuvastatin

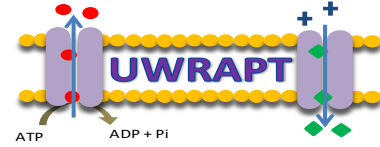
0  10^4



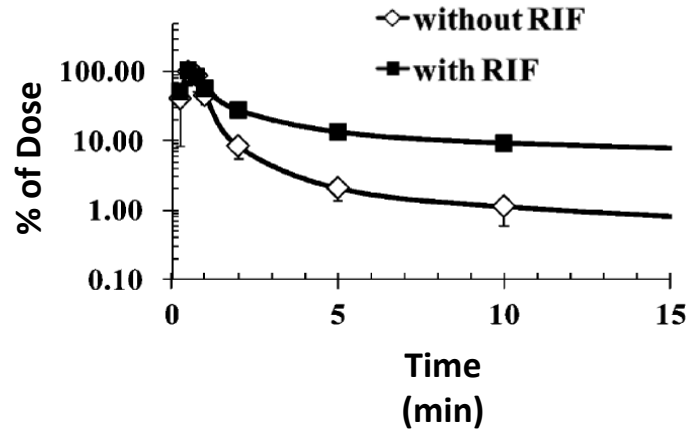
He et al., Mol Pharm., '14



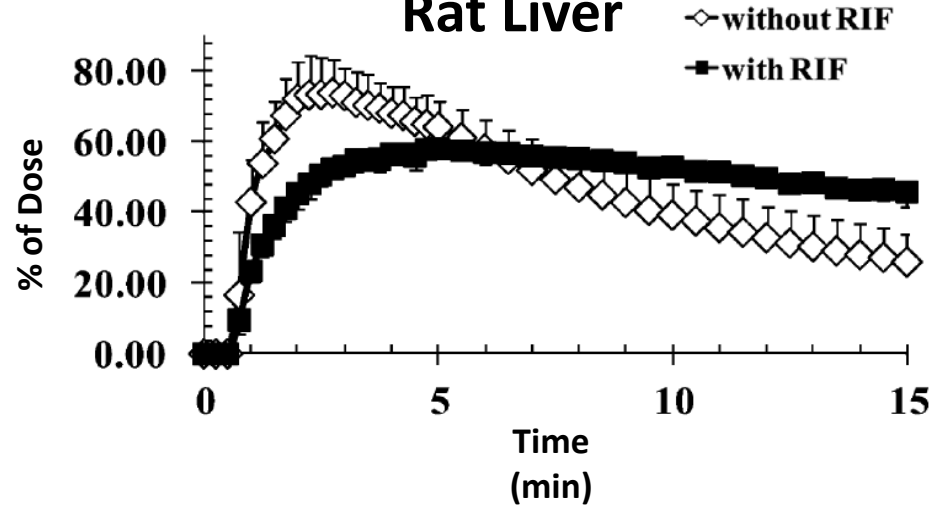
Asymmetrical change in rosuvastatin conc. in liver and blood in the presence of rifampin



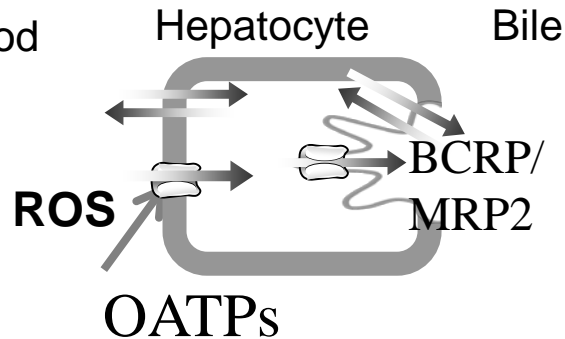
Rat Blood



Rat Liver

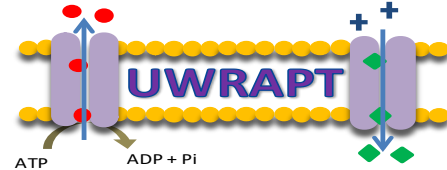


2.3-fold increase in plasma AUC
but no significant increase in liver
AUC





Hepatic Uptake and Biliary Excretion of $[^{11}\text{C}]$ Rosuvastatin \pm CsA



1 min

5 min

10 min

30 min

0

RSV

Liver

Liver

Liver

Liver

GallBladder

RSV +
CSA

Liver

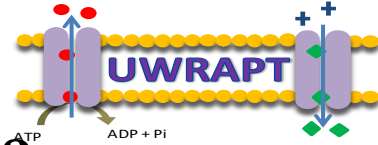
Liver

Liver

Liver

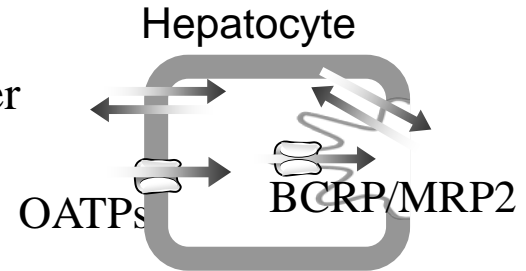
GallBladder

How Can we Predict/Measure Tissue Drug Conc. and Tissue:Blood Asymmetry in Humans?

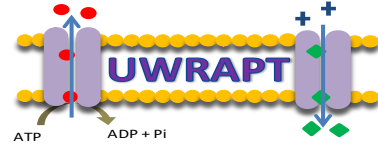


One could actually measure tissue drug conc:

- PET imaging (MRI and other imaging modalities do not have the required sensitivity):
 - Requires sophisticated equipment and radiochemistry
 - Costly (about \$20-40K/experiment/subject)
- Therefore we need alternative methods that will allow us to predict tissue conc. of drugs in humans
 - Predict all clearances associated with the tissue:blood barrier
 - Possible using REF but not RAF approach



REF Method to Predict Tissue Drug Conc.



Hypothesis: Predict tissue drug conc. by scaling in vitro CL in transporter expressing cells to in vivo using relative expression factor (REF)

In vitro CL

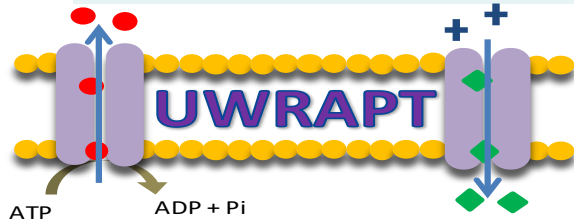
CL via transport of interest
in cell line expressing the
transporter

Relative Expression Factor (REF)

1. Transporter expression/g of
tissue
2. Tissue weight

In vivo CL

Contribution of individual
transporter in tissue
uptake/efflux



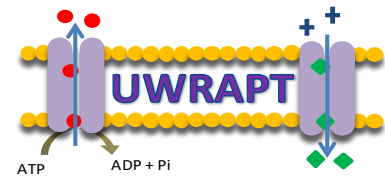
$$REF = \frac{[T]_{\text{ex vivo in organ}}}{[T]_{\text{in vitro}}}$$

- <https://sop.washington.edu/departments-of-pharmaceutics/research/research-affiliate-program-on-transporters-uwrapt/>

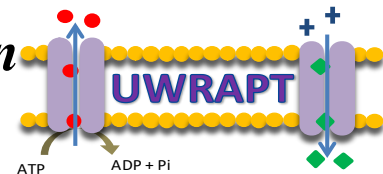
Can the use of these Quantitative Proteomics Data Successfully Predict Transporter-Mediated CL and Tissue Conc. of Drugs?

Criteria of success: within 2-fold of observed value

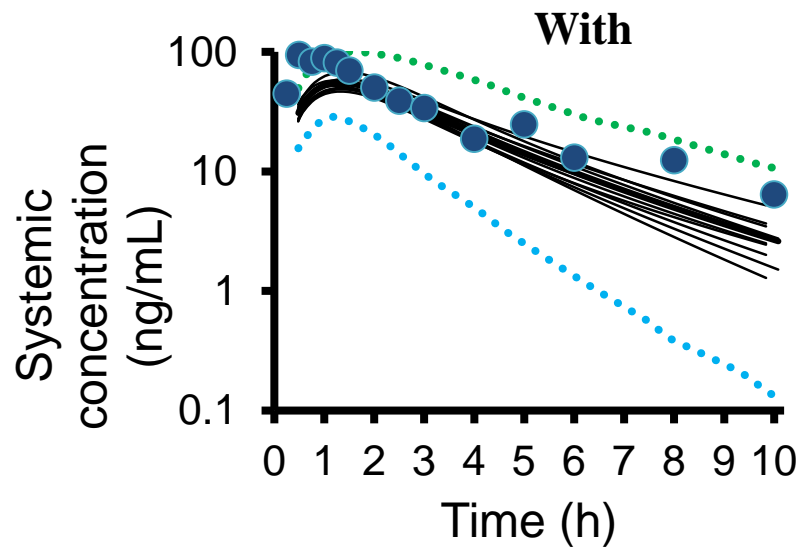
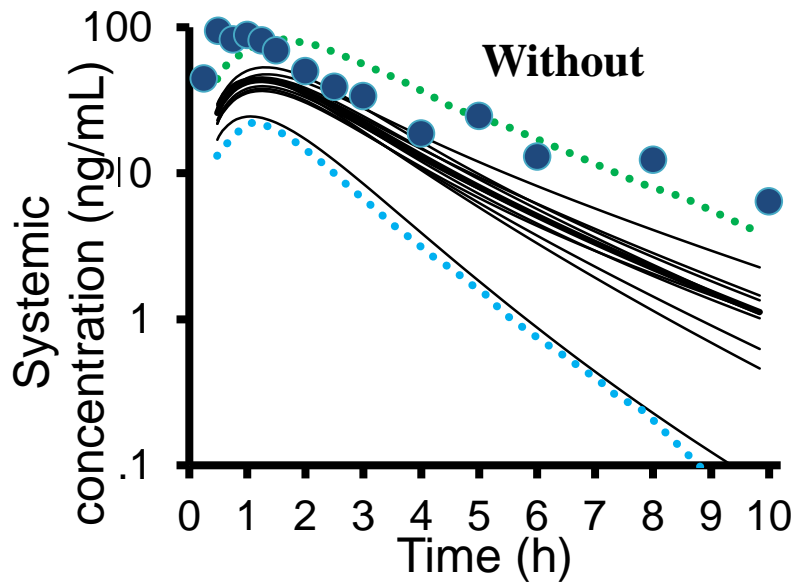
- Changes due to hepatic impairment (e.g. cirrhosis)
- Renal secretory CL of drugs
- Hepatic CL and tissue conc. of drugs – rat and human



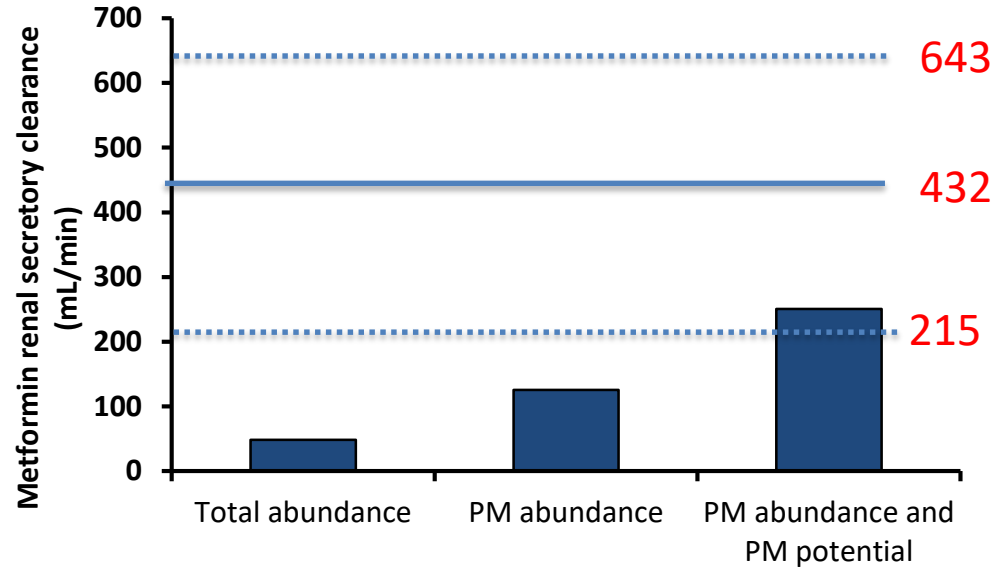
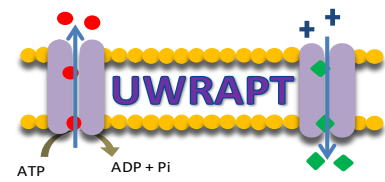
Improved PBPK Prediction of Repaglinide Pharmacokinetics in Liver Cirrhosis Patients When the Effect of Cirrhosis on OATP1B1 Abundance is Incorporated



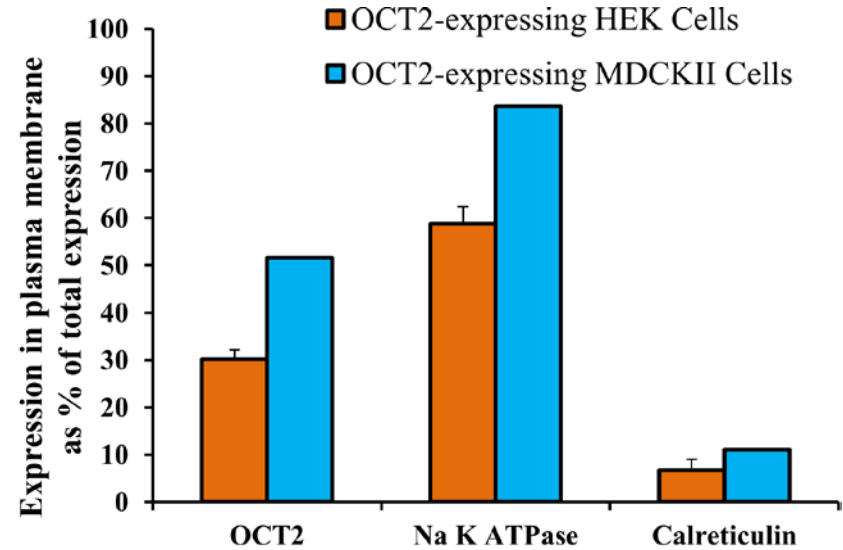
Virtual population (10 trials and n=12 for each trial)



Metformin renal clearance is reasonably well-predicted using OCT2 expressing cells

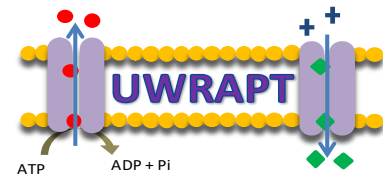


Observed metformin renal secretory clearance in humans: 432 (range 215-643) mL/min

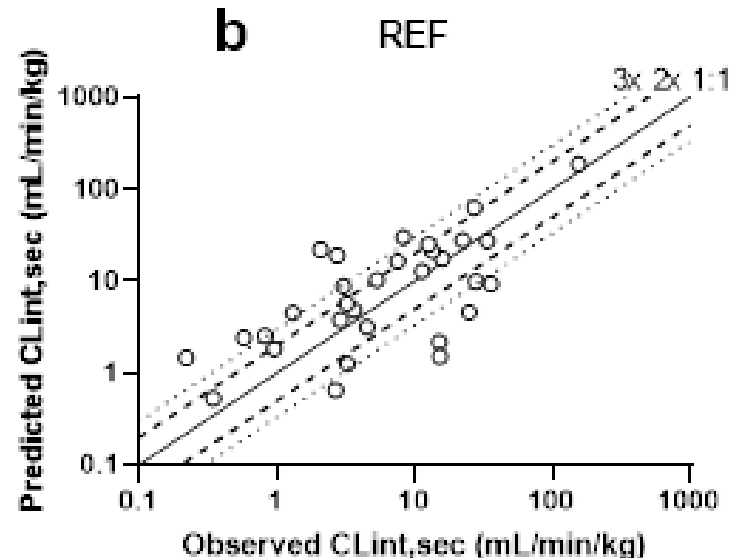
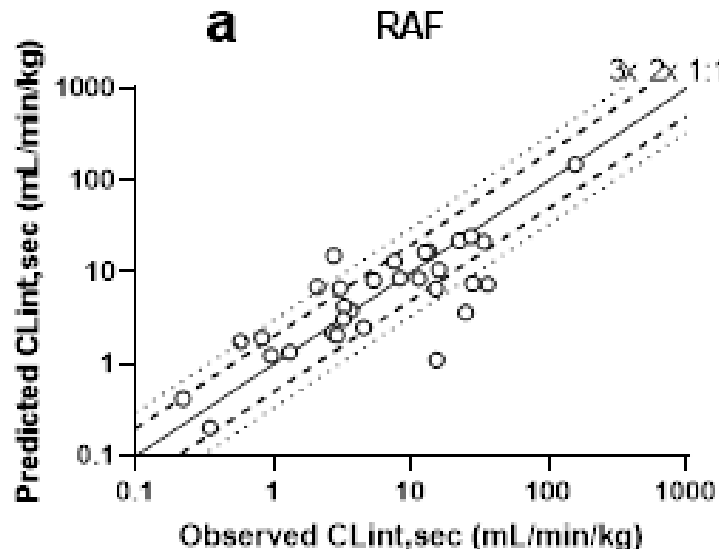


Kumar et al., DMD 2018

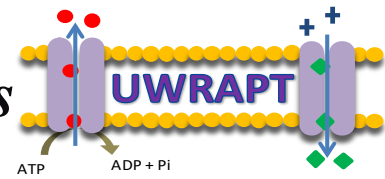
Renal Secretory CL by OATs: RAF vs. REF



- 31 drugs secreted by renal OATs
- RAF used tenofovir (OAT1), acyclovir and ganciclovir (OAT2), and benzylpenicillin/oseltamivir acid (OAT3) as probe substrates
- REF used quantitative proteomics data on renal OATs in human kidney cortex and transfected cells which were also used to determine *in vitro* uptake CL of the drugs.



Hypothesis: Predict Transporter-Mediated In-Vivo Hepatobiliary CL and Hepatic Concentrations of Drugs in Rats/Humans using REF

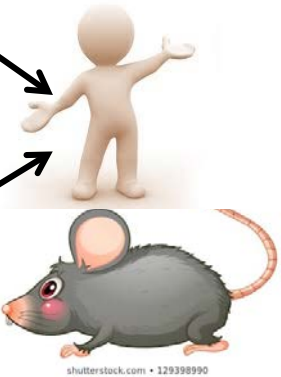


Transporter-expressing cells

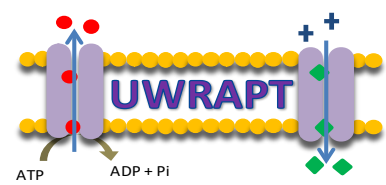
1. transporter-mediated drug CL
2. transporter abundance using quantitative proteomics
3. Obtain REF

Predict *in-vivo* hepatobiliary CL using REF and traditional approach

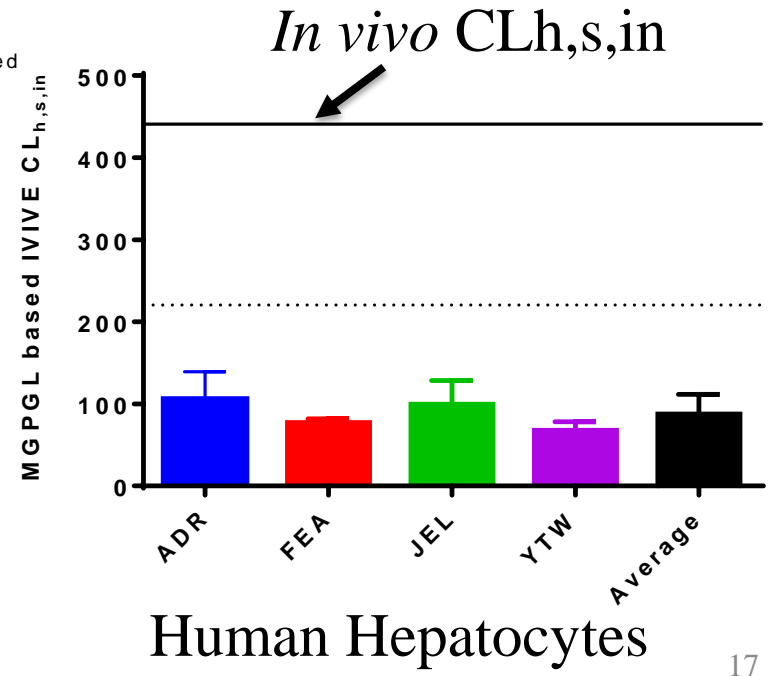
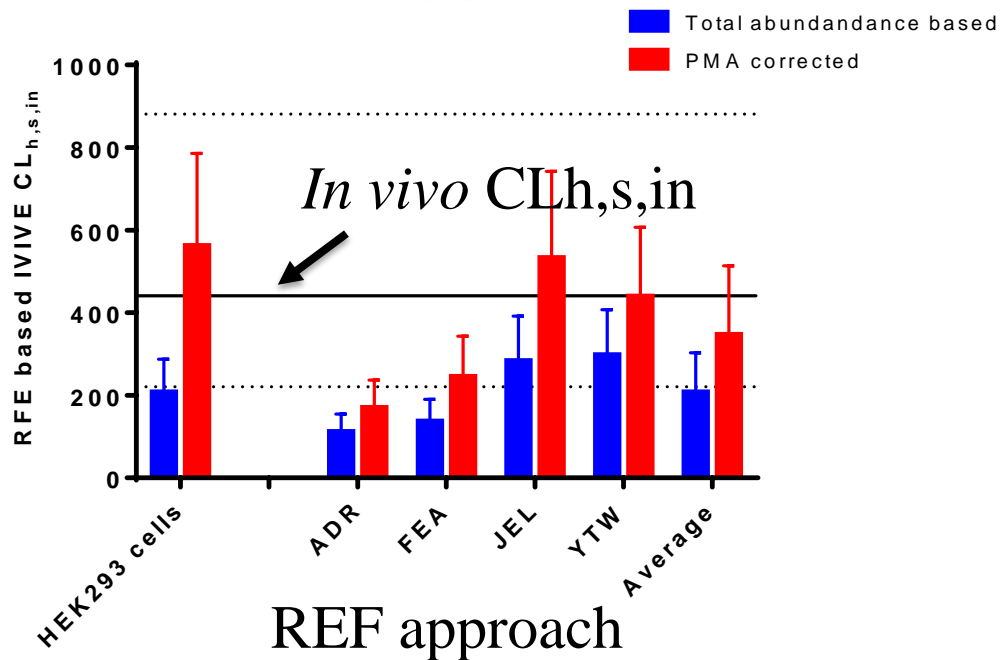
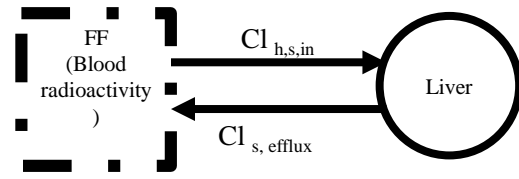
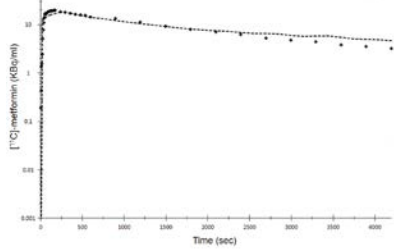
Verify predictions using PET imaging



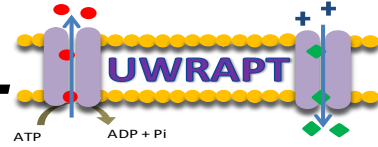
REF (using HEK293 cells or PHH) better predicts Hepatic Uptake CL of metformin (as measured by PET imaging) vs. PHH



Gormsen, et al., *J Nucl Med* 57(12): 1920-1926.

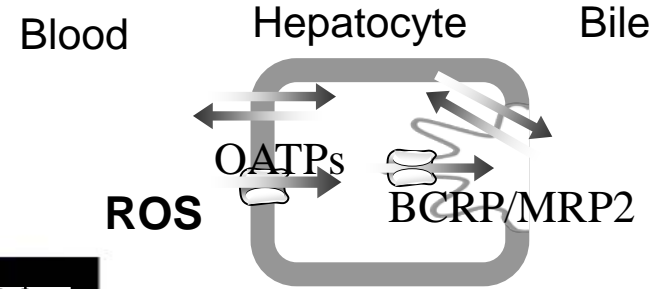
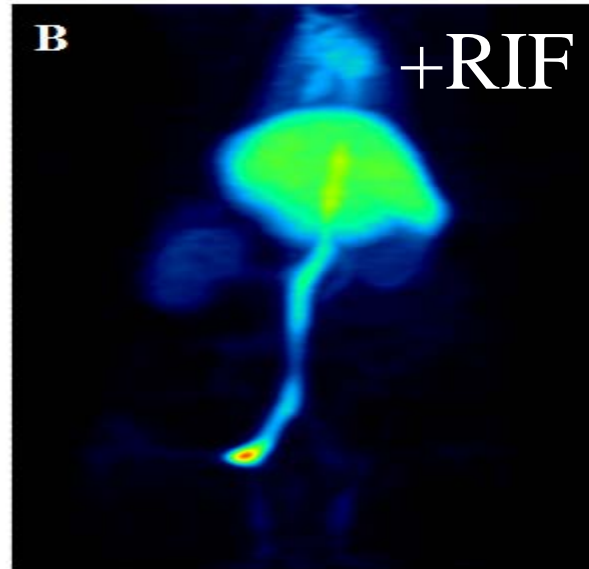
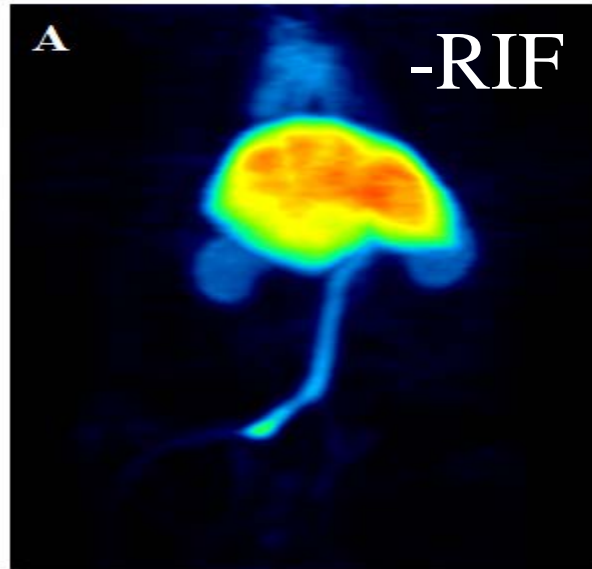


Hepatic Uptake and Biliary Excretion of ^{11}C -Rosuvastatin in the Rat



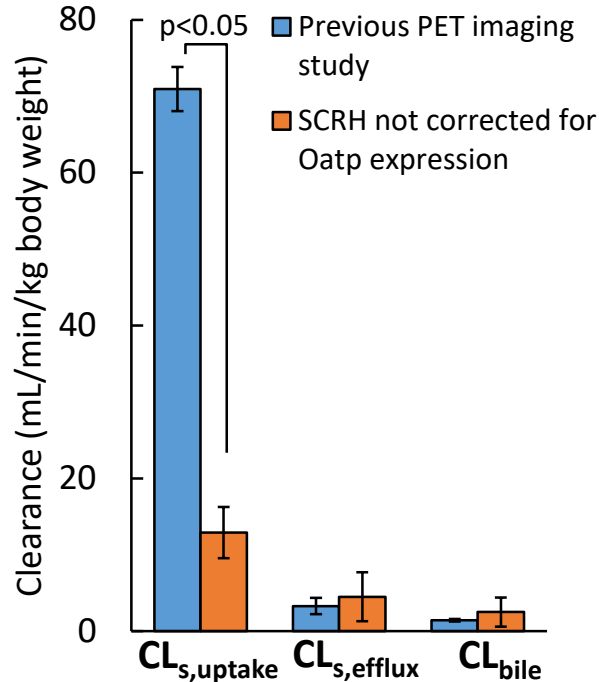
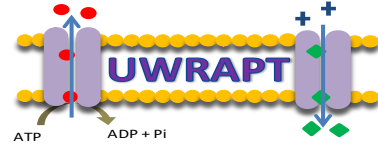
Coronal 2 min SUV images of ^{11}C -Rosuvastatin

0  10^4

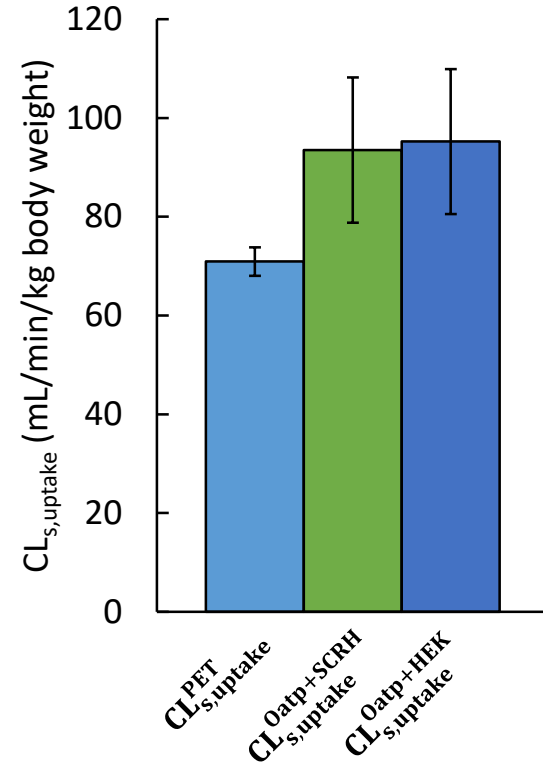
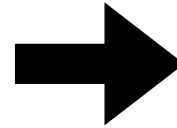
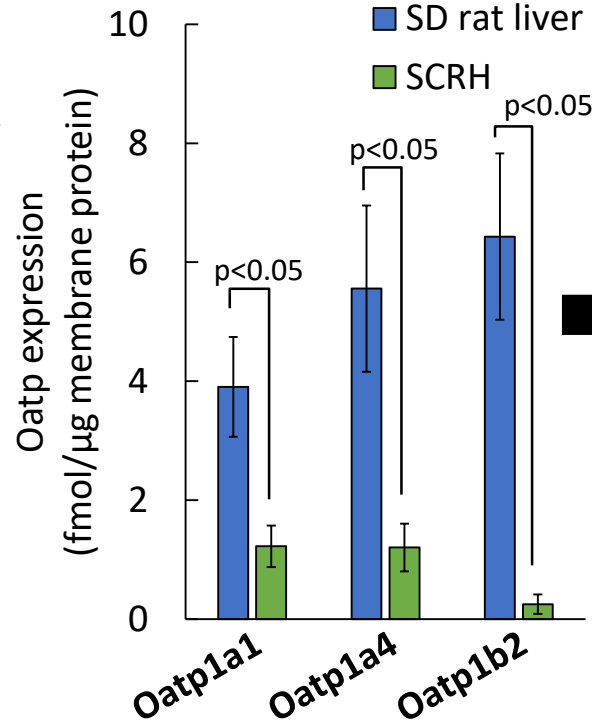


He et al., Mol Pharm., '14

Successful prediction of the hepatobiliary clearance of rosuvastatin using cell lines, sandwich-cultured rat hepatocytes and quantitative proteomics

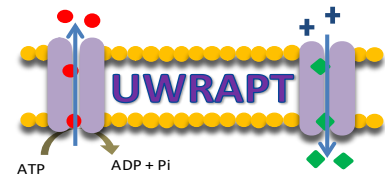
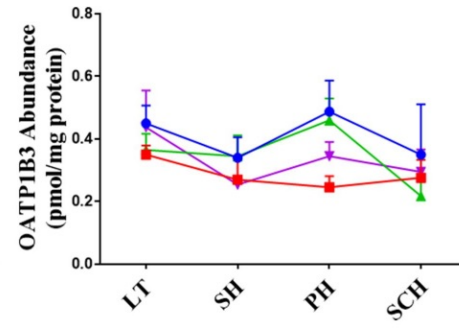
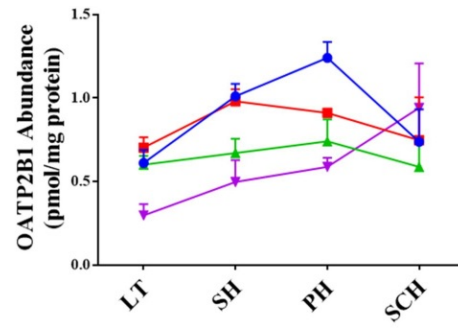
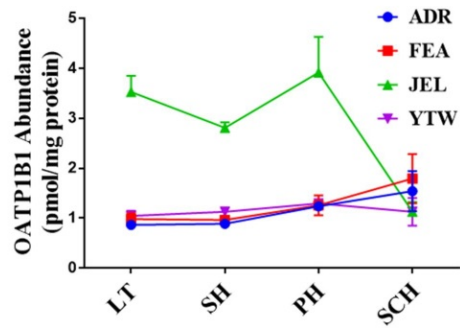


CL_{s,uptake}; sinusoidal uptake
 CL_{s,efflux}; sinusoidal efflux
 CL_{bile}; canalicular efflux

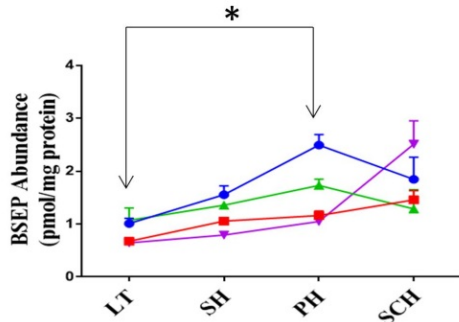
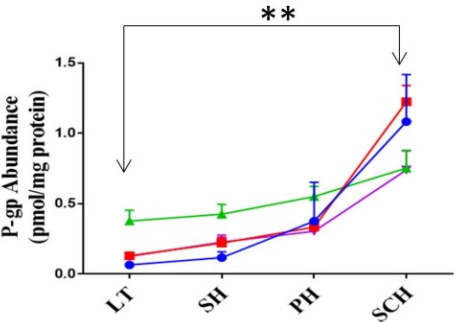
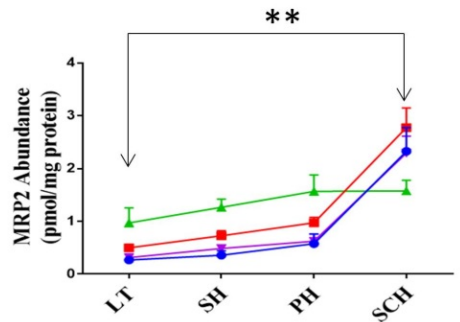
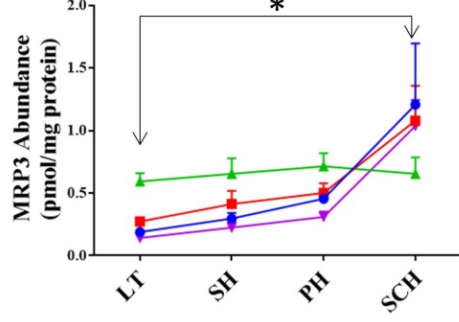
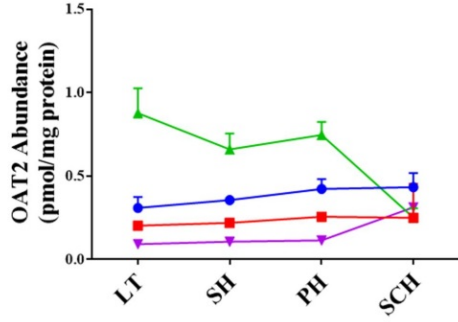
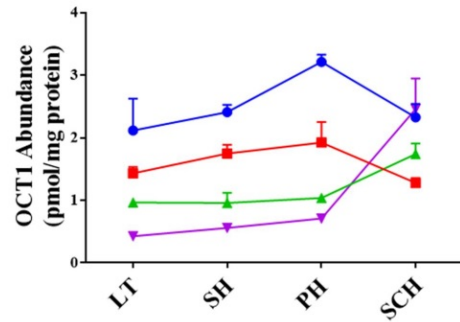




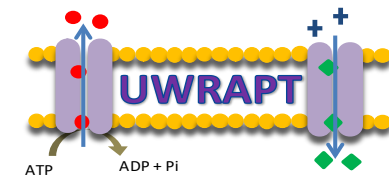
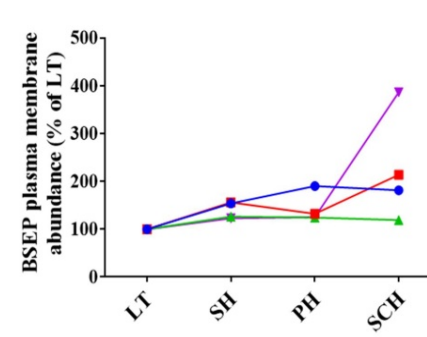
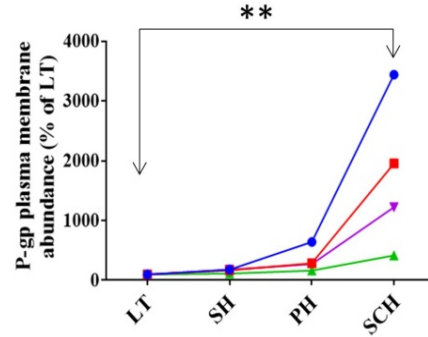
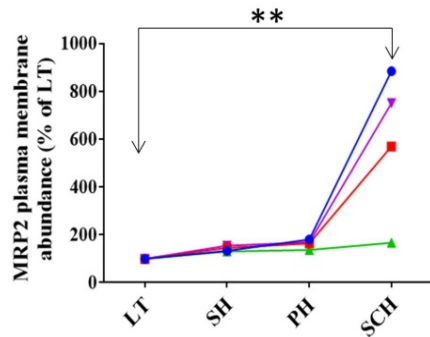
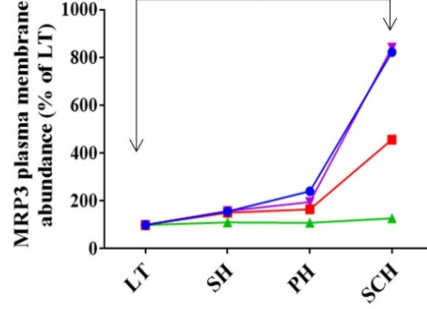
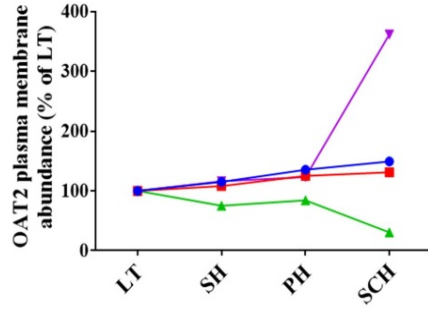
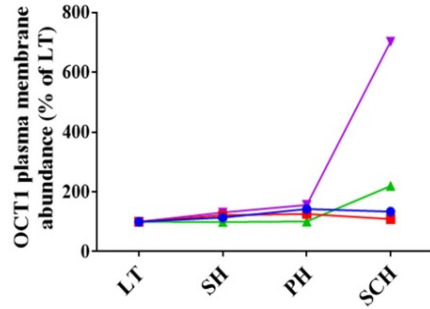
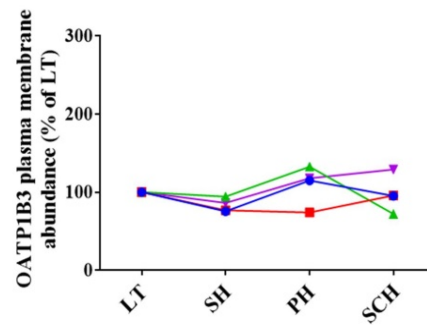
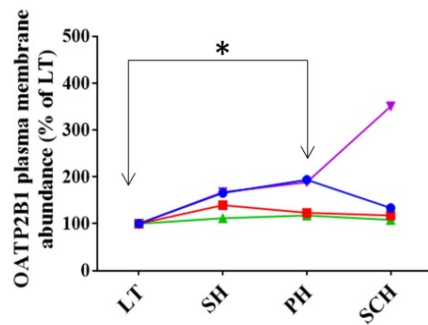
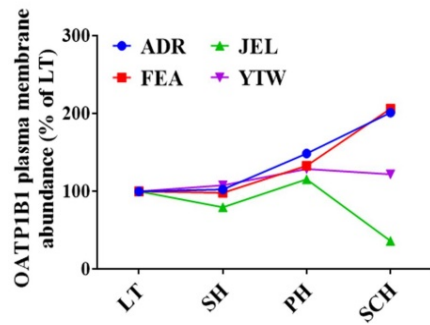
***Can Rosuvastatin Hepatobiliary CL and Hepatic
Conc. be Predicted in Humans?***



Total transporter abundance in suspended (SH), plated (PH), sandwich-cultured (SCH) hepatocytes and liver tissue



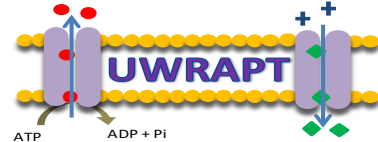
Kumar et al.,
DMD 2019



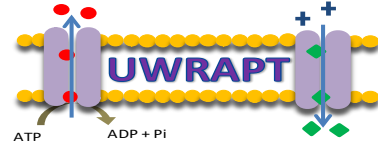
Plasma membrane transporter abundance in suspended (SH), plated (PH), sandwich-cultured (SCH) hepatocytes cf liver tissue

Kumar et al.,
DMD 2019

Summary



- Predicting transporter-mediated CL & tissue concentrations and therefore efficacy and toxicity of a drug is the next frontier in ADME research
- The hepatic ECL model clarifies when transporters will or will not affect the systemic and tissue PK of a drug
- Tissue conc. measurement is possible using PET. However, this method cannot be routinely applied
- IVIVE using transfected cells and quantitative transporter proteomics is a promising technique to predict tissue drug conc
- These predictions should be validated using PET imaging probes that interrogate multiple drug transporters



Major Contributors



Gabriela Patilea-Vrana



YuYang

Jiake He



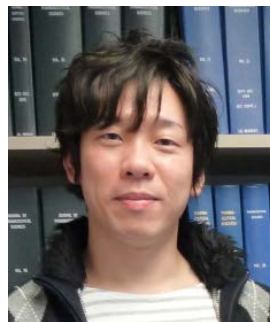
Li Wang



Sarah Billington



Vineet Kumar



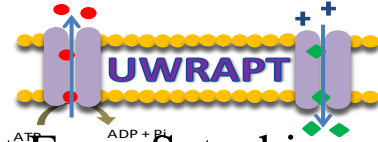
Kazuya Ishida



Bhagwat Prasad

Anand Desai

Other Collaborators



Dept. of Radiology: Jeanne Link, David Mankoff, Todd Richards, Janet Eary, Satoshi Minoshima, Ken Maravilla, Mark Muzi, Steve Shoner, Shirely Rene, David Lewis, Jean Lee and the PET suite team

Dept. of Medicine: Ann Collier and her team; Scott Lee and his team

Dept. of Anesthesiology: Karen Domino, Matthew Pennington

Dept. of Pharmaceutics: Bhagwat Prasad, Edward Kelly, Carol Collins, Joanne Wang

Kidney Research Institute: Jonathan Himmelfarb

Univ. of Greifswald: Stefan Oswald and team

Children's Mercy Hospitals: Steven Leeder and team

Aarhus University Hospital, Aarhus, Denmark, Dr. Gormsen

Acknowledgement: NIH P01DA032507, MH63641, P50 HD44404, RR 00166, HD47892, AG031485, RC1NS068904, BioIVT, Solvo Biotech., UWRAPT funded by Genentech, Merck, Biogen, Gilead, BMS, Takeda, Pfizer, Amgen

