

# **Clinical Significance and Regulatory Framework for the Evaluation of Organic Anion Transporting Polypeptide 1B-Based Drug-Drug Interactions**

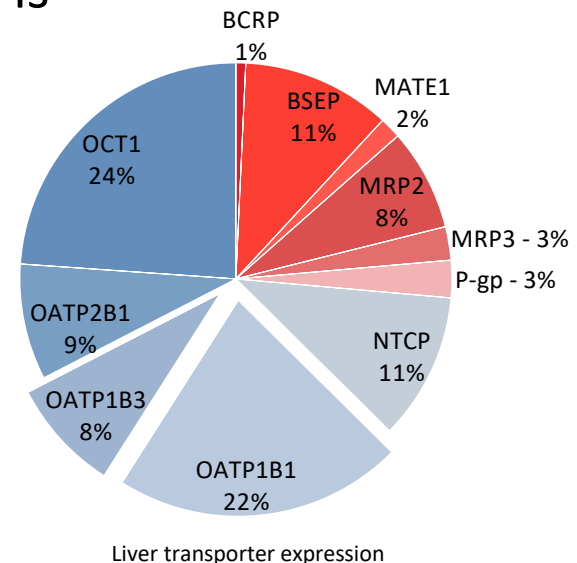
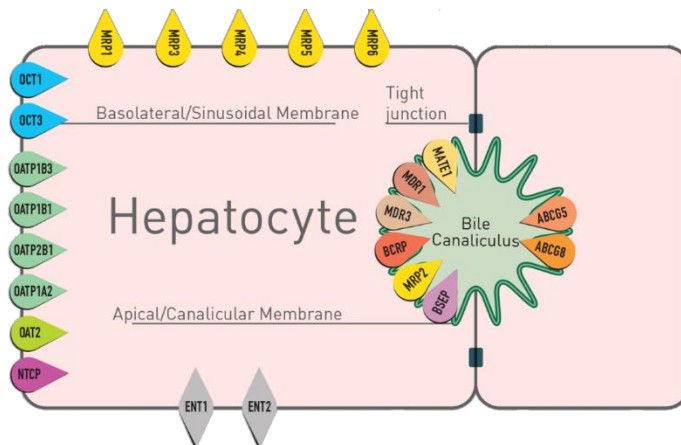
*Savannah McFeely, PhD*

*Research Scientist, UW Drug Interaction Solutions*

*5 September 2019*

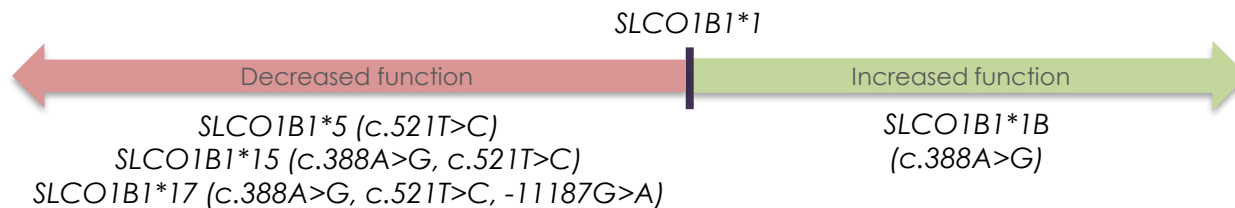
# OATP1B1/1B3

- The OATPs belong to the solute-carrier (SLC) family of transporters
  - Assumed to transport compounds based on concentration gradient or ion exchange
- OATP1B1 and 1B3 are uptake transporters exclusively expressed on sinusoidal membrane of hepatocytes
  - OATP1B1 and OATP1B3 share 80% amino acid identity
- Among the liver transporters, OATP1B1 shows the second highest expression (22%) and OATP1B3 expression is approximately one-third of that (8%)



# OATP1B1/1B3 Polymorphisms

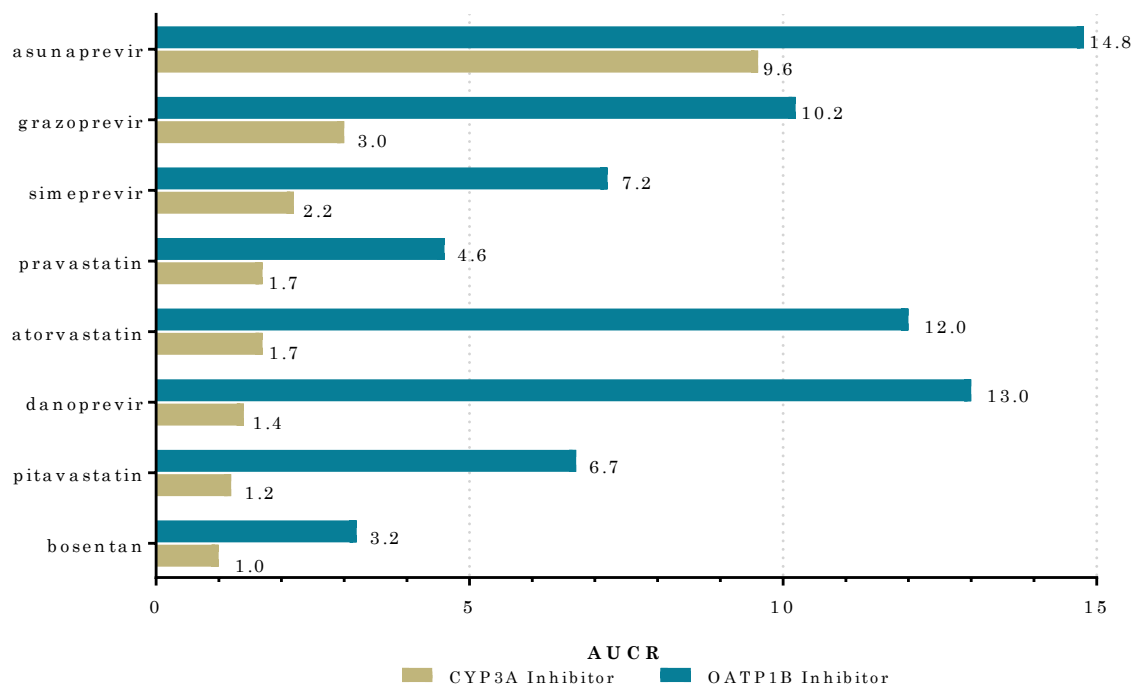
- Genetic variation in both *SLCO1B1* and *SLCO1B3* have been shown to affect function and therefore substrate exposure.
- 21 different *SLCO1B1* variant alleles have been identified to date with varying effects on transport efficiency relative to the wild type (*SLCO1B1*\*1)



- Variants of *SLCO1B3* are currently not as well characterized and while many have been identified, clinical effects are mostly unknown.
  - Decreased function *in vitro*: 334T>G, 699G>A, 1564G>T, -5035G>A

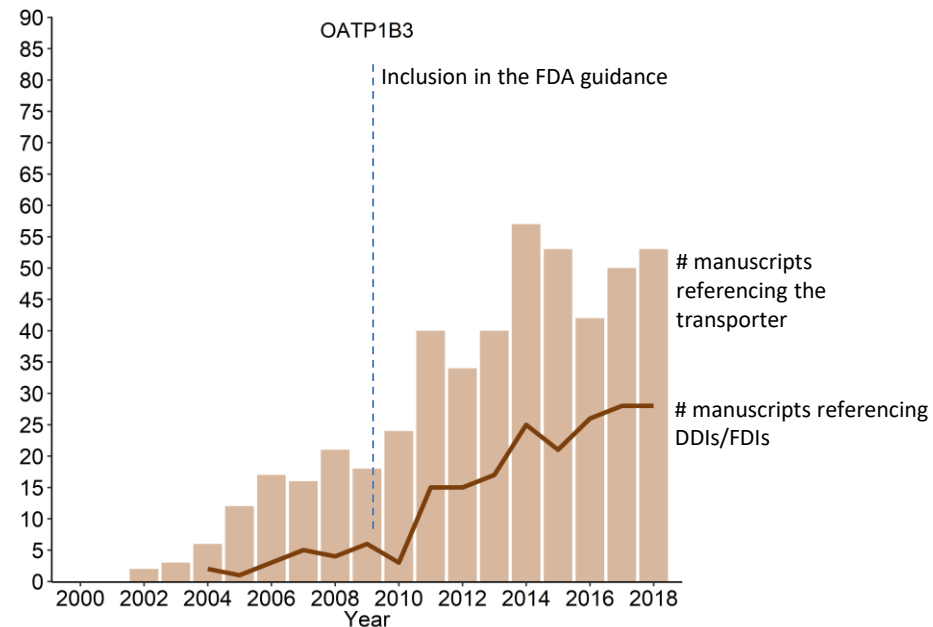
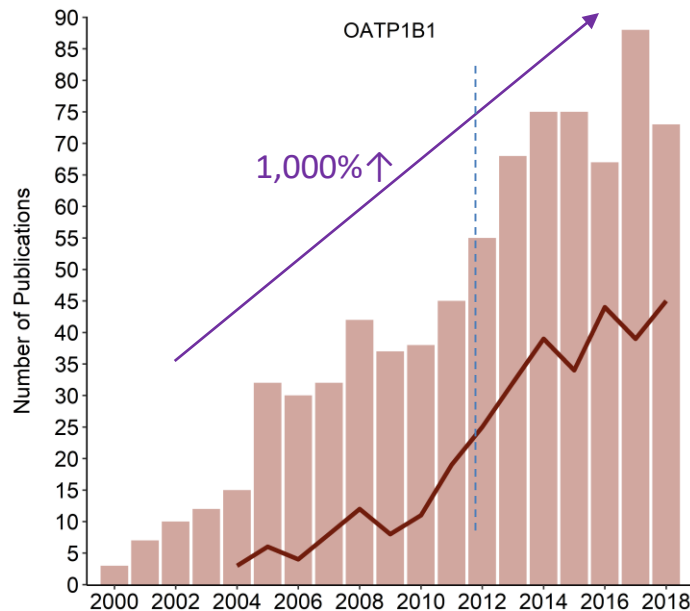
# Inhibition Causes Significant Changes in Exposure

- For many compounds, hepatic uptake is a rate-determining step and the effect of inhibition can meet or exceed that observed with CYP inhibition
  - Magnitude of change in exposure is also, on average, much higher than observed with other transporters



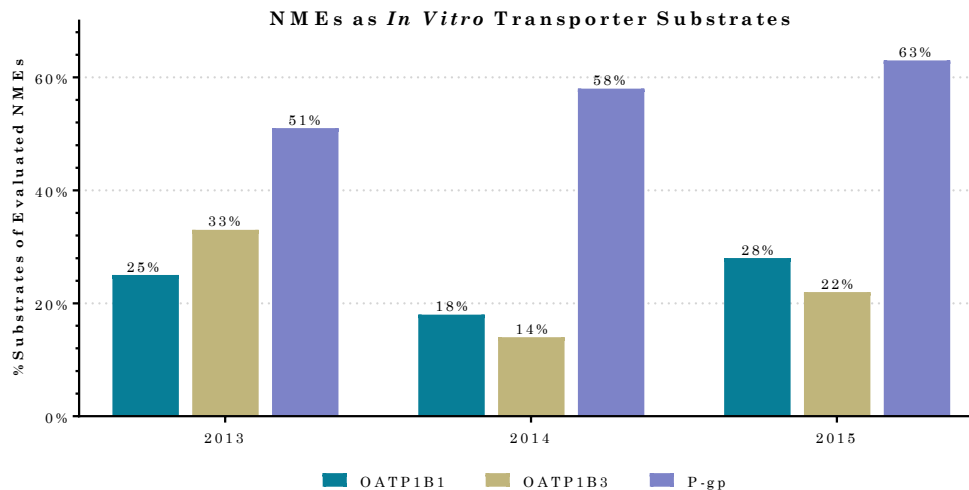
# OATP1B1/1B3 Research and Publications

- Since the transporters were identified in the early 2000s, the number of publications on the structure/function has steadily increased
- Recommended for evaluation during drug development in 2012
  - The number of reported drug-drug interactions (DDIs) and food-drug interactions (FDIs) continues to increase



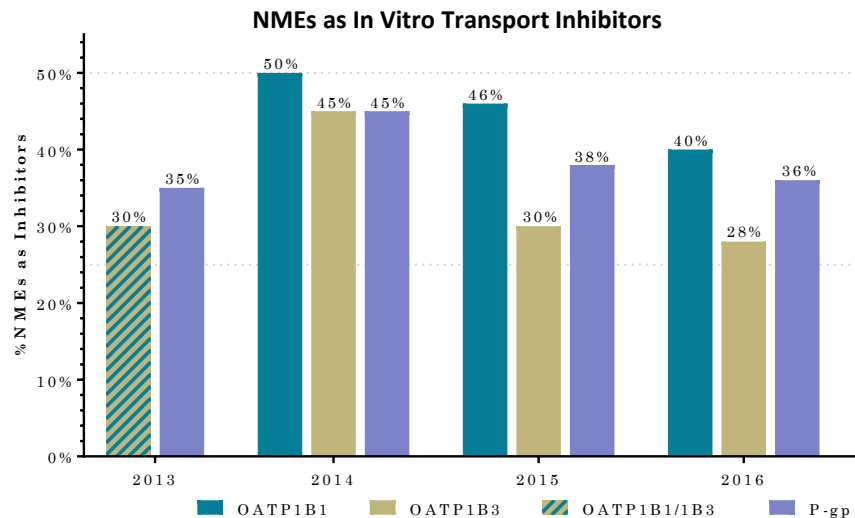
# OATPs in recent NDAs

- The 2012 revision to the FDA drug interaction guidance added six transporters, including OATP1B1/1B3, to be considered in the evaluation NMEs
- Based on *in vitro* data, less than 10 new drugs in the last four years are OATP1B1/1B3 substrates
  - P-gp is the most common (>40), followed by BCRP
- Overall, fewer drugs were tested as substrates of OATPs compared to P-gp



# OATPs in recent NDAs

- In the last four years, OATP1B1 is the transporter most commonly inhibited by NMEs (44 drugs) *in vitro*
  - Followed by P-gp (37 drugs) and OATP1B3 (33 drugs)



- When evaluated *in vivo*, only 10% significantly increased OATP1B1/1B3 substrate exposure

# Regulatory Guidance on Transporter Assessment

---

- *In Vitro*
  - As Substrate:  $\geq 25\%$  of  $CL_{\text{total}}$  is hepatic/biliary; site of action in the liver
    - Uptake Ratio  $\geq 2$ , decreases with known inhibitor by  $\geq 50\%$
  - As Inhibitor: all new compounds must be evaluated
    - R-value  $\geq$ ~~1.25~~ 1.1
- *In Vivo*
  - Positive *in vitro* result(s)
    - Change in AUC  $\geq 1.25$ -fold
- Recommended index drugs
  - Substrates: pitavastatin, pravastatin, rosuvastatin
  - Inhibitors: cyclosporine, single dose rifampin



# OATP1B1/1B3 Marker Compounds

- Recommended marker substrates and inhibitors are similar between agencies, but limited
  - Inhibitors: cyclosporine, single dose rifampin

*“Results from most transporter inhibition studies are not easily extrapolated to other drugs, because most inhibitors are not specific for a single transporter”*

- Substrates: pitavastatin, pravastatin, rosuvastatin

*“Several drugs are substrates of more than one transporter. For example, rosuvastatin is a substrate for BCRP and OATP.”*

Despite the increase in research on OATP1B1/1B3 since these transporters were included in the 2012 guidance, little has been updated regarding their evaluation

It has been well established in recent years that OATP1B1/1B3 are clinically relevant transporters for drug-drug interactions and should be considered during development, yet the current regulatory guidance offers a limited choice of selective substrates.

**By analyzing clinical and preclinical literature data, it is hypothesized that more sensitive and selective substrates and inhibitors of OATP1B1/1B3 can be identified, which can, in turn, be used to evaluate and improve the translatability of *in vitro* data to *in vivo* prediction.**

---

- *Aims of the evaluation:*

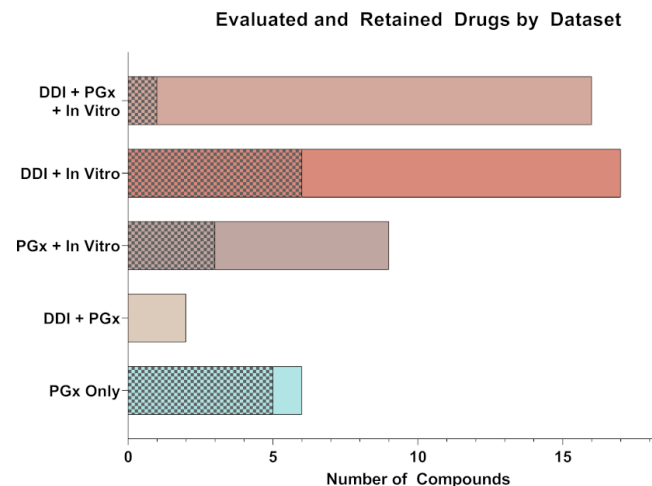
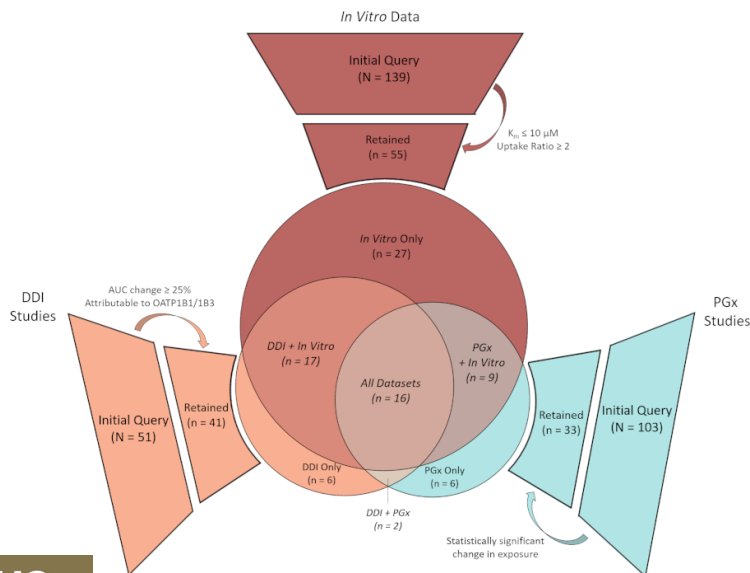
- Identify potential *in vivo* substrates of OATP1B1/1B3 and evaluate the identified compounds for clinical relevance using a novel indexing system
- Evaluate the sources of variability in the *in vitro* evaluation of OATP1B/1B3 inhibitors and the effect on clinical interaction predictions
- Identify potential inhibitors of OATP1B1/1B3 and evaluate the identified compounds for clinical relevance

# SUBSTRATE IDENTIFICATION

McFeely SJ et al. Identification and Evaluation of Clinical Substrates of Organic Anion Transporting Polypeptides 1B1 and 1B3. Clin Transl Sci. 2019 Jul 1;12(4):379–87.

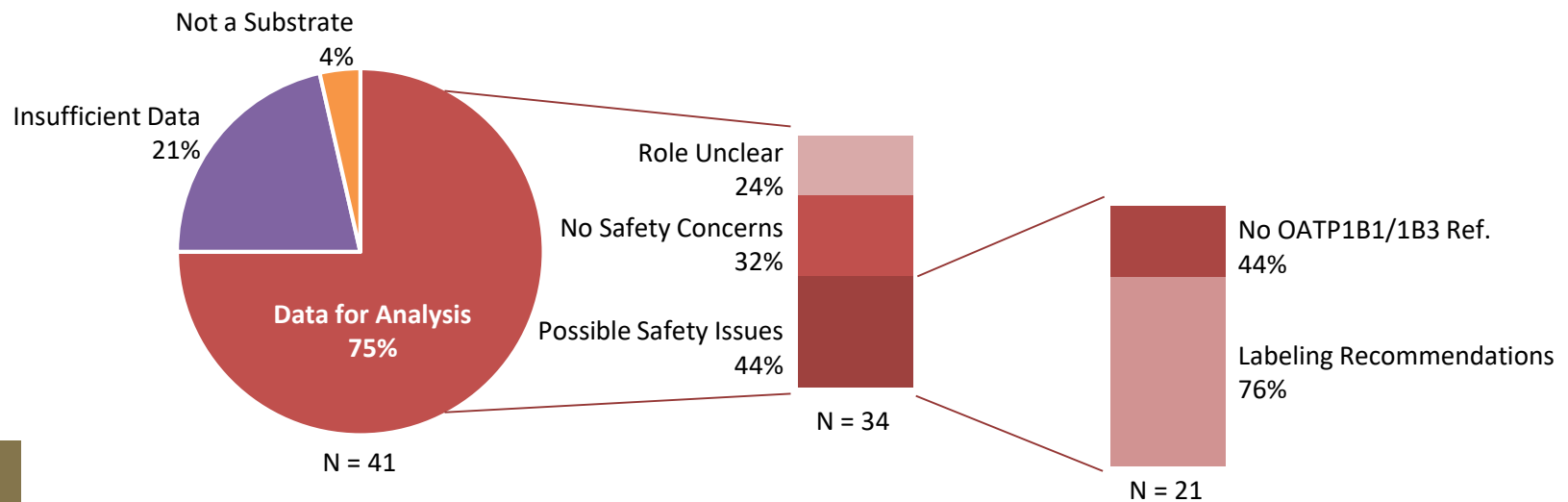
# Substrate Identification

- Queries of the *in vitro*, clinical DDI, and pharmacogenetic modules of the UW DIDB were completed to identify potential clinical substrates of OATP1B1/1B3
  - 53% of identified *in vitro* substrates did not have corresponding clinical data and were unable to be evaluated further.
  - 26% of substrates (22/83) had *in vitro* and either clinical DDI or PGx data
  - 19% of substrates (16/83) had data from all three sources



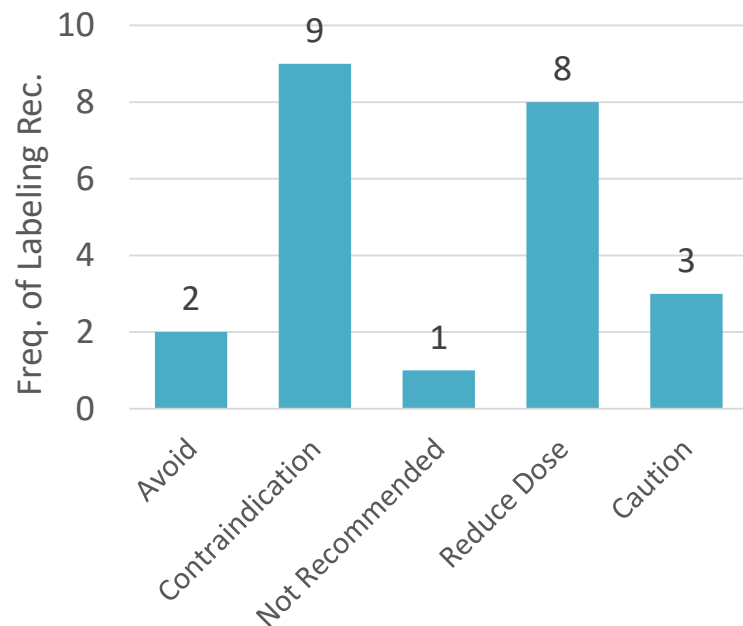
# Clinical Substrates of OATP1B1/1B3

- Of the 41 drugs identified as potential substrates, 34 (83%) had sufficient data to support a clinically significant role of OATP1B1/1B3
  - 21 show possible significant safety issues associated with OATP1B1/1B3 inhibition
  - 6 did not have sufficient data to determine the clinical impact of inhibition
  - 1 not a substrate of OATP1B1/1B3
- 16/21 identified substrates (76%) have labeling recommendations regarding OATP1B1/1B3 inhibition.



# Labeling Recommendations for Identified Substrates

- 16 of the 22 identified substrates (72%) have statements in the labeling regarding OATP1B1/1B3 inhibition.
  - 23 specific statements
  - Includes language towards “OATP inhibitors” (5) and specific inhibitors (11)
- 5 drugs (24%) do not currently have recommendations regarding OATPs



Identified substrates with no OATP1B1/1B3 labeling recommendations		
Drug	AUC Ratio	Possible Reason for Lack of Recommendation
caspofungin	1.6 (RIF)	caution recommended with CsA
danoprevir	15.6 (CsA)	not approved in US/Europe
docetaxel	1.6 (CsA)	reduce dose with strong CYP3A inhibitor (2.4-fold, keto)
lovastatin	5.0 (CsA)	avoid GEM or CsA (CYP3A)
SN-38	2.1 (PGx)	active metabolite of irinotecan

# Probe Index

An index was developed to quantitatively and objectively evaluate substrates for utility as an OATP1B1/1B3 probe substrate.

Primary  
Evaluation  
Categories

TOTAL SCORE	15	(top of each category + all positive criteria)
Sensitivity to OATP1B1/1B3 inhibition <sup>a</sup>	0	No PGX data or clinical studies with a specific inhibitor for OATP1B1/1B3 -or- AUC Ratio < 1.25
	1	$1.25 \leq \text{AUCR} < 2$
	2	$2 \leq \text{AUCR} < 3.5$
	3	$3.5 \leq \text{AUCR} < 5$
	4	$5 \leq \text{AUCR} < 7.5$
	5	$7.5 \leq \text{AUCR} < 10$
	6	$\text{AUCR} \geq 10$
Specificity <sup>b</sup>	0	Sensitive substrate for at least 2 metabolic enzymes or transporters (AUCR $\geq 5$ for each pathway) <sup>c,d</sup>
	1	Moderate sensitive substrate for at least 2 metabolic enzymes or transporters ( $2 \leq \text{AUCR} < 5$ for each pathway) <sup>c,d</sup>
	2	Sensitive substrate of one metabolic enzyme or transporter (AUCR $\geq 5$ )
	3	Weak substrate for at least 2 metabolic enzymes or transporters (AUCR < 2 for each pathway) <sup>c,d</sup>
	4	Moderate sensitive substrate of one metabolic enzyme or transporter ( $2 \leq \text{AUCR} < 5$ )
	5	Weak substrate of one metabolic enzyme or transporter (AUCR < 2)
	6	Only OATP1B1/1B3 contributes to the disposition of the compound
Safety Profile	-2	Unfavorable safety profile for a single dose (narrow therapeutic range or expected significant side effects) or clinical safety has not been fully evaluated at this time
	1	Can be administered as a single, low dose with a low risk of adverse events in a healthy population or is well tolerated over a wide dose range, no concerns administering to a healthy population

## Additional Criteria:

Additional  
Criteria

Positives	1	PGx studies completed showing an impact of <i>SLCO1B1</i> or <i>1B3</i> variants
	0.5	Microdosing validated
	0.5	Published and validated PBPK model
Negatives	-2	Only available as a combination therapy
	-0.5	Non-linear pharmacokinetics
	-0.5	Half-life longer than 24 h
	-0.5	Very low bioavailability ( $F < 5\%$ )

# Probe Index

- Six drugs are proposed as potential clinical marker substrates
  - High sensitivity towards OATP1B1/1B3 inhibition
  - Low or manageable contribution of other metabolism/transport
  - Favorable clinical safety profile

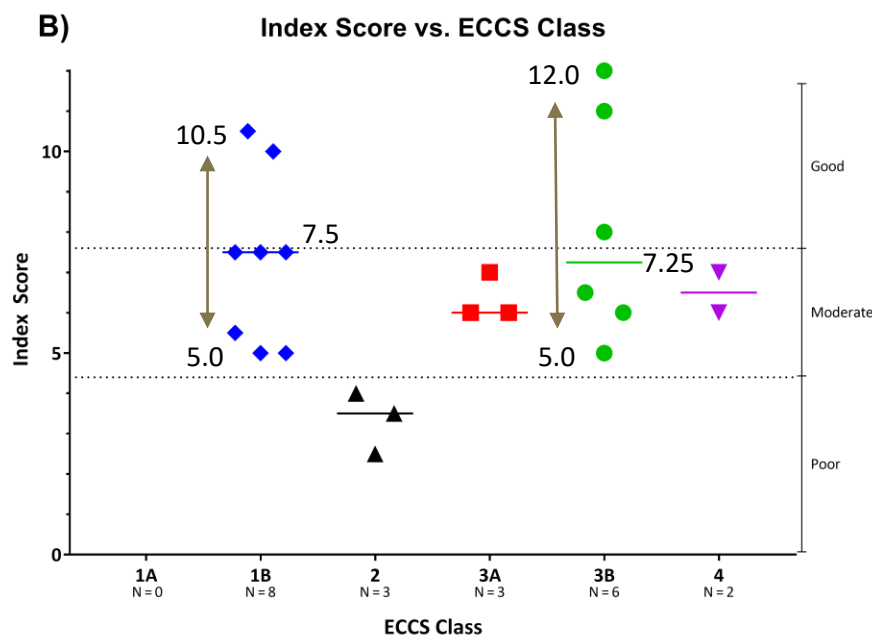
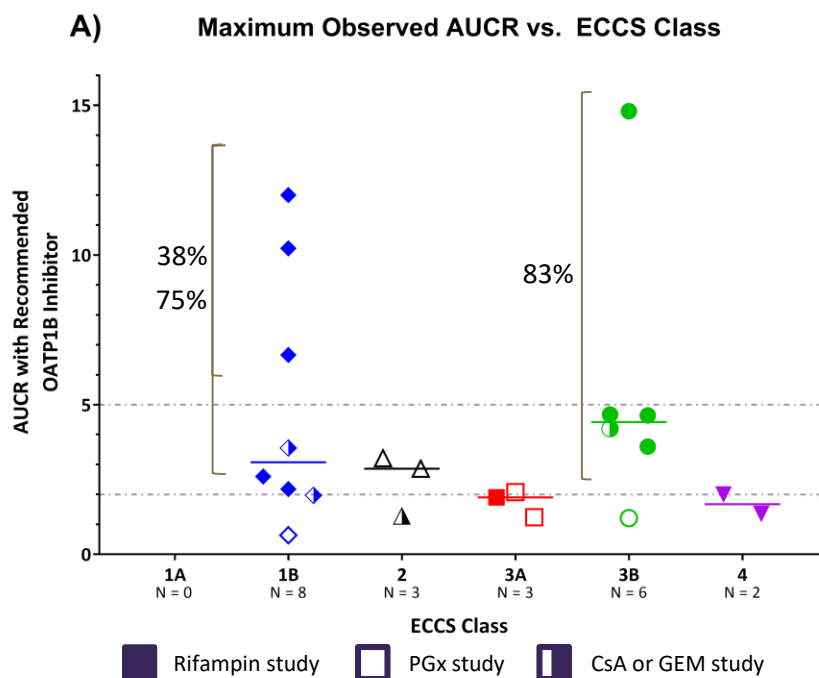
Drug	Rank	Index Score	ECCS Classification	Therapeutic Area	Highest Reported AUC Ratio	Highest Observed PGX Effect	Other Metabolism / Transport
pravastatin*	1	12.0	3B	statin	4.64	3.81	BCRP/OATP2B1/ P-gp
rosuvastatin*	2	11.0	3B	statin	4.67	2.18	CYP2C9 BCRP/OATP2B1/ P-gp
pitavastatin*	3	10.5	1B	statin	6.67	3.85	BCRP/OATP2B1/ P-gp
atorvastatin*	4	10.0	1B	statin	12.0	2.51	CYP3A BCRP/P-gp
eluxadoline	5	8.0	3B	GI agent	4.20 (CsA)	2.01	N/A
letermovir	5	8.0	--	antiviral	2.10 (CsA)	1.40	N/A

\*FDA/ITC Recommended Substrate



# Comparison to ECCS

- The ECCS evaluates drugs based on a combination of permeability, ionization state, molecular weight, and the separation of metabolic and transport rate- determining steps
  - The 1B and 3B classes should be the most promising OATP1B1/1B3 markers



# Substrate Summary

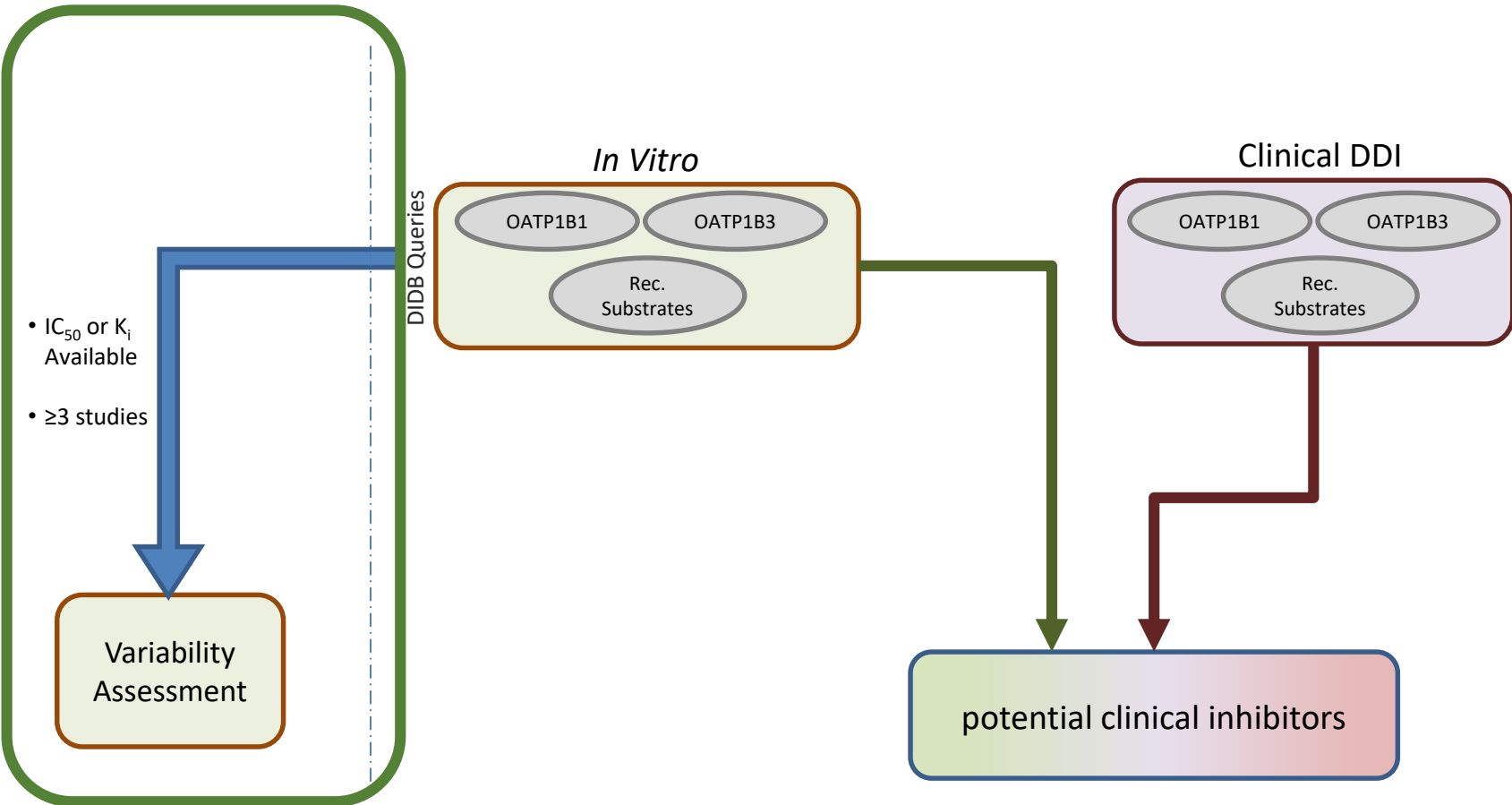
---

- 34 drugs were identified as clinical substrates of OATP1B1/1B3
  - Of these, 6 were identified using a novel ranking system as potential marker compounds
- A thorough understanding of the clinical disposition of these drugs allows for use of a fit-for-purpose marker
  - Isolate the contribution of OATP1B1/1B3 using a selective compound
    - Ex: pravastatin, pitavastatin, eluxadoline
  - Determine a “worst-case scenario” effect if the NME is an inhibitor of multiple pathways
    - CYP3A/P-gp/OATP1B: atorvastatin
- The current regulatory approach to *in vitro* substrate data has limitations
  - Uptake ratios are highly variable and currently do not have established acceptance or reporting criteria

# INHIBITOR IDENTIFICATION AND VARIABILITY

McFeely SJ, et al. Variability in *In Vitro* OATP1B1/1B3 Inhibition Data: Impact of Incubation Conditions on Variability and Subsequent Drug Interaction Predictions. Clin Transl Sci. 2019 [Epub ahead of print].

# Compound Identification



University of Washington Drug Interaction Database (DIB®, [www.druginteractioninfo.org](http://www.druginteractioninfo.org))

# *In Vitro* Variability

---

- 128 studies evaluated from 44 publications
  - Required to have  $\geq 3$  studies for retention
  - OATP1B1
    - $IC_{50}$  values: 21 substrate/inhibitor pairs
    - $K_i$  values: 7 substrate/inhibitor pairs
  - OATP1B3
    - $IC_{50}$  values: 2 substrate/inhibitor pairs
- Inhibitors: rifampin (27%), cyclosporine (25%), gemfibrozil (18%)
- Substrates: estradiol-17- $\beta$ -gluc (62%), atorvastatin (15%)
- Cell type: HEK293 (79%)

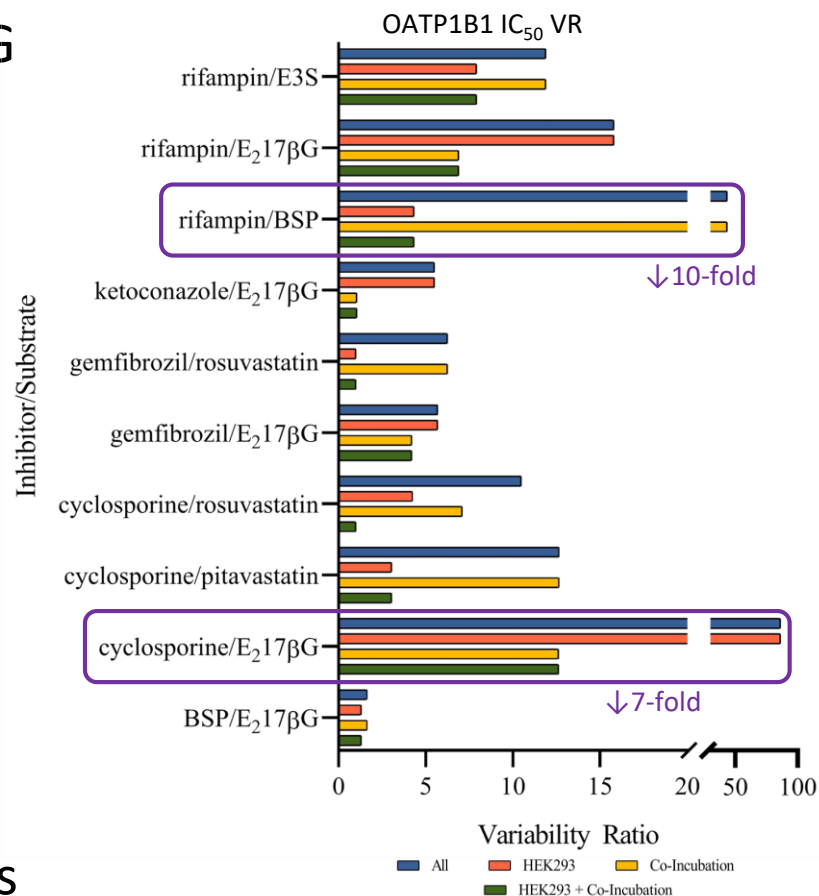
Variability ratios (highest  $IC_{50}$  or  $K_i$  relative to the lowest) were calculated for each pair

R-values were calculated from each inhibitor constant

# IC<sub>50</sub> and K<sub>i</sub> Variability

- The VR for the entire dataset = 12.4
  - VR<sub>IC50,OATP1B1</sub> = cyclosporine/E<sub>2</sub>-17β-G (86.4, n = 11)
  - VR<sub>Ki,OATP1B1</sub> = gemfibrozil/ E<sub>2</sub>-17β-G (7.2, n = 3)
  - VR<sub>IC50,OATP1B3</sub> = rifampin/ E<sub>2</sub>-17β-G (58.2, n = 7)

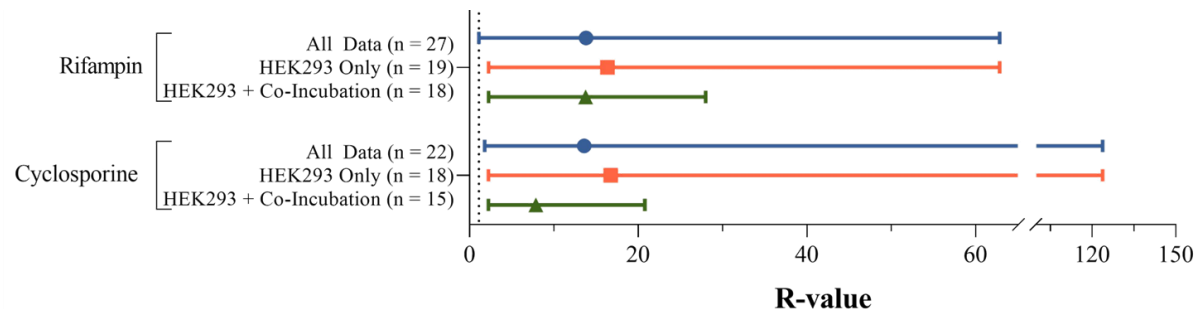
- Accounting for cell type and co-incubation reduced dataset variability (VR = 5.23)
- Substrate also contributed to variability
  - Highest VR for non-clinical substrates
    - cyclosporine/E<sub>2</sub>-17β-G (86.4, n = 11)
    - cyclosporine/pitavastatin (12.7, n = 4)



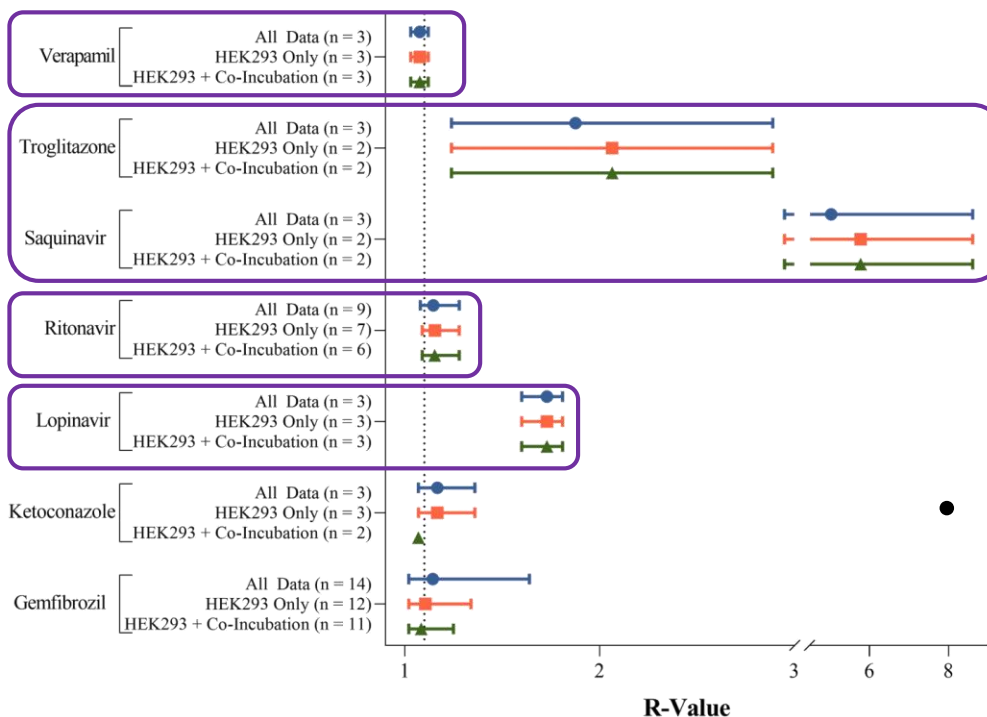
Reduced by controlling for cell type and co-incubation

# R-Value Variability

- Despite changes in VR when accounting for incubation conditions, resulting R-values did not show a significant shift relative to the FDA cut-off value of 1.1
- The recommended index inhibitors rifampin and cyclosporine had all R-values  $\geq 1.1$ 
  - Maximum fold-change was reduced when accounting for incubation conditions
    - Rifampin: 12.8  $\rightarrow$  5.7
    - Cyclosporine: 51.1  $\rightarrow$  8.6
- In contrast, only 5/14 (36%) of the R-values calculated for gemfibrozil met the FDA cut-off
  - Likely due to not accounting for inhibitory metabolites



# R-Value Variability



- Remaining drugs had mixed effects of incubation conditions on R-value
  - 4 showed all values  $\geq 1.1$  regardless of conditions
  - 2 resulted in R-values above and below the cutoff for all datasets
  - Ketoconazole did not have any R-values  $\geq 1.1$  for the most uniform dataset
- Very few drugs had clinical data available
  - Trend towards less variability in significance for strong inhibitors

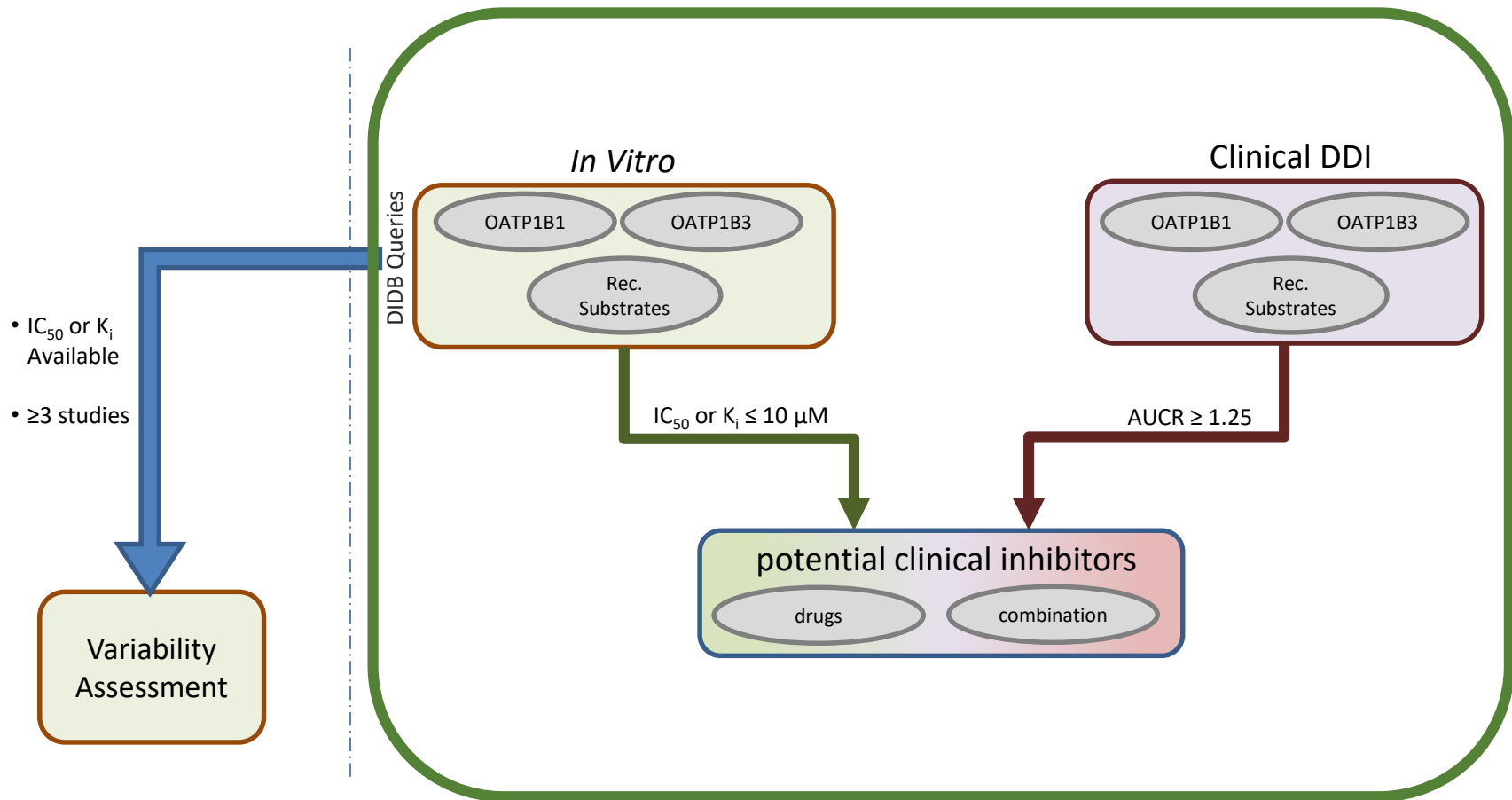


# Conclusions

---

- Two aspects of study design – cell type and preincubation – significantly contribute to *in vitro* variability
  - FDA recommends a 30-min preincubation as of the 2017 guidance
  - Over 80% of experiments performed in the last 5 years were completed in HEK293 cells
  - Substrate was also found to have an effect
- *In vitro* variability does not appear to have an effect on clinical predictions for the inhibitors evaluated
  - Weaker inhibitors may show predictions above and below the cut-off value
- Despite the broad range of values found in this work, the overall variability is lower than what has been observed for P-gp
  - P-gp showed over 700-fold variability for a single inhibitor/substrate pair

# Clinical Inhibitor Identification



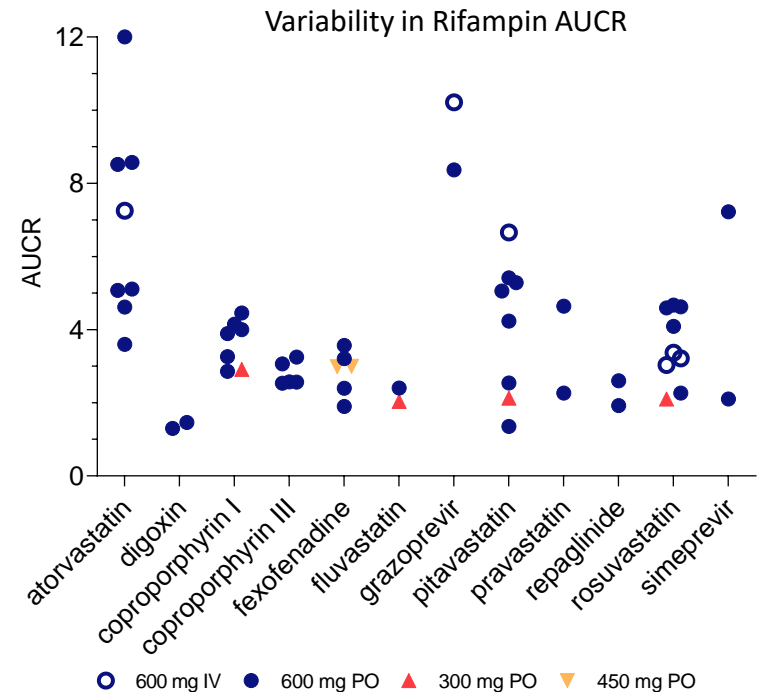
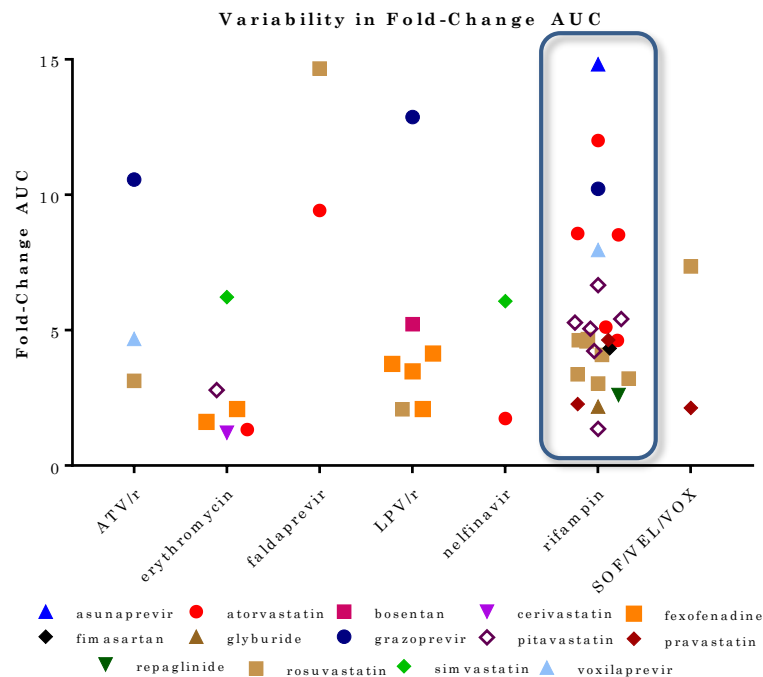
# Rifampin as a Marker Inhibitor

---

- Despite the long-standing use of rifampin, many aspects of its disposition and use as an index inhibitor have not been fully evaluated
  - Reproducibility and variability
  - Time-dependent inhibition
    - FDA now recommends a 30 min pre-incubation with inhibitor for *in vitro* evaluation
  - Induction / inhibition balance
- Areas for future research [regulatory perspective]
  - Impact of lower doses
  - Route / timing of administration
  - Use of PBPK modeling
- Clinical Use
  - Populations / regions where data is most relevant
  - What is known about the impact of RIF on co-meds

# Rifampin as a Marker Inhibitor

- In vivo*, there is high variability observed in the AUCR for a given inhibitor-substrate pair
  - pitavastatin – rifampin, 5-fold variation
  - atorvastatin – rifampin, 2.6-fold



# Rifampin as a Marker Inhibitor

---

- Currently, 89% of studies use a 600 mg dose of rifampin with 76% using a single, oral dose (68% overall)
  - Limited data for other doses
  - Alternate doses are almost exclusively multiple dose studies
- Static predictions for doses ranging from 300 mg - 900 mg show little difference for the sensitive substrate pravastatin (2.53 – 2.67)
  - Likely due to plasma concentrations >> lowest reported  $K_{i,OATP1B1}$
- Lower doses of rifampin could likely be used in inhibition studies, reducing risk to patients while still providing maximal inhibition

# Inhibitor Summary

---

- Over 60% of *in vitro* OATP1B1/1B3 inhibitors met the regulatory criteria for further clinical evaluation
  - Clinical data is limited – less than 40% of these compounds have study data available
- Using clinical data for identified sensitive substrates, 13 drugs and 16 combination treatments were identified as inhibitors of OATP1B1/1B3
  - Majority of interactions are weak (47%, AUCR < 2)
  - 14% of identified interactions have an AUCR ≥ 5
- No novel clinical index inhibitors were identified in this analysis, but these findings further support the utility of cyclosporine and rifampin as worst-case-scenario and targeted inhibitors, respectively
  - Despite the frequent use, many aspects of rifampin study design have not been fully evaluated
  - There is a limited understanding of the underlying causes of variability in AUCR for specific interactions

# Conclusions

---

- Thorough analysis of the clinical data identified 12 marker compounds for OATP1B1/1B3
  - Includes drugs from multiple therapeutic areas
  - 9/12 have labeling recommendations regarding OATP1B inhibition
- 13 clinical inhibitors have been identified from studies with known marker compounds
  - Most interactions result in AUCR < 2
  - A high number of potent *in vitro* inhibitors, yet clinical data are limited
  - Data supports the regulatory use of rifampin and cyclosporine
- Translating *in vitro* transport data to *in vivo* effects is inherently difficult
  - Uptake ratios are not currently well defined
  - High variability in IC<sub>50</sub>/K<sub>i</sub> values could contribute to poor predictions of clinical effect
  - Contribution of other metabolic and transport pathways confound clinical interpretation

# Acknowledgments

---

- University of Washington Drug Interaction Solutions
  - Isabelle Ragueneau-Majlessi
  - Tasha Ritchie
  - Rene Levy
  - Jingjing Yu
- Eva Gil Berglund
- Anna Nordmark

**Savannah McFeely, PhD**  
**sjkerr@uw.edu**  
**206.616.9751**