

A BioIVT Company

PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE
FROM DISCOVERY THROUGH CLINICAL SUPPORT

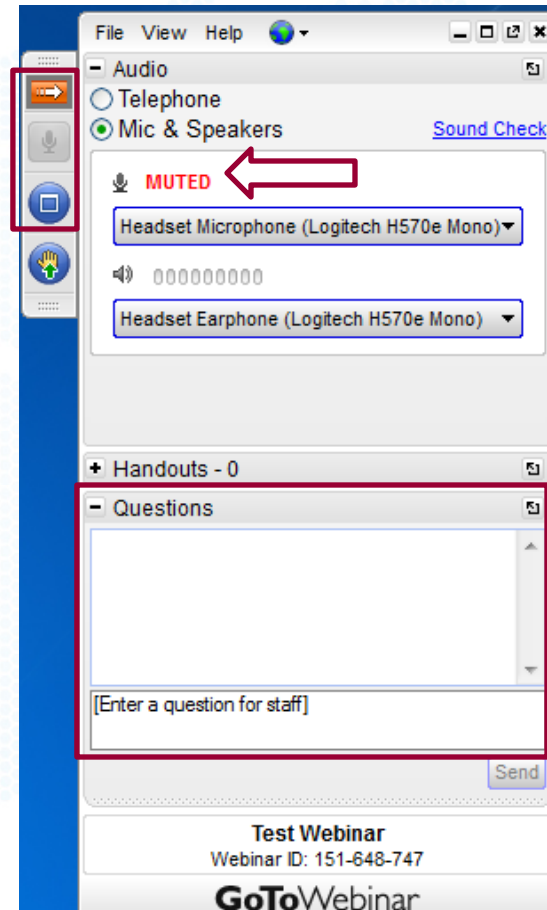
Evaluation of Ketoconazole and Its Alternative Clinical CYP3A4/5 Inhibitors as Inhibitors of Drug Transporters: The In Vitro Effects of Ketoconazole, Ritonavir, Clarithromycin and Itraconazole on 13 Clinically- Relevant Drug Transporters

Lydia Vermeer, Ph.D.
Senior Scientist, Drug Transport
lvermeer@xenotechllc.com



Before we begin...

- User console features
- Audience audio is muted
- Post questions to our staff anytime via “Questions” message pane
- Slides and recorded session available soon. Watch your e-mail for both.
- A short survey will appear at the conclusion of the webinar, we value your feedback!



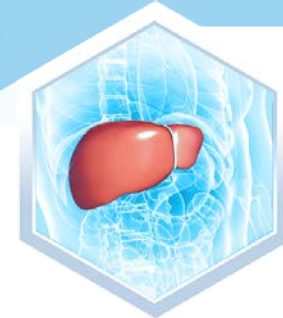
A BioIVT Company

PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE
FROM DISCOVERY THROUGH CLINICAL SUPPORT

Evaluation of Ketoconazole and Its Alternative Clinical CYP3A4/5 Inhibitors as Inhibitors of Drug Transporters: The In Vitro Effects of Ketoconazole, Ritonavir, Clarithromycin and Itraconazole on 13 Clinically- Relevant Drug Transporters

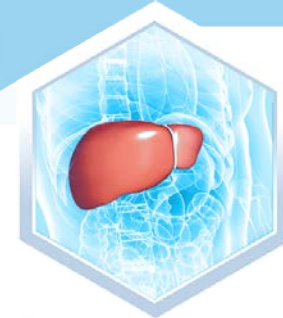
Lydia Vermeer, Ph.D.
Senior Scientist, Drug Transport
lvermeer@xenotechllc.com





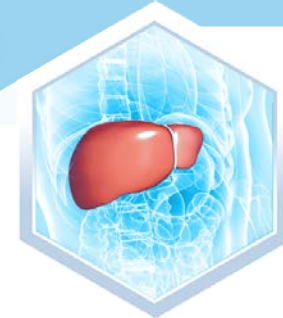
Drug Transport Technology at XenoTech

Transporter Category	Transporter	Experimental system	
SLC	OAT1, 3	HEK	
	OCT1, 2	HEK	
	OATP1B1, 1B3,	HEK	
ABC	MDR1 (human)	MDCK	Vesicle
		Caco-2 (BCRP also present)	
	MRP2-4	Vesicle	
	BCRP (human)	MDCK	Vesicle
	BSEP	Vesicle	



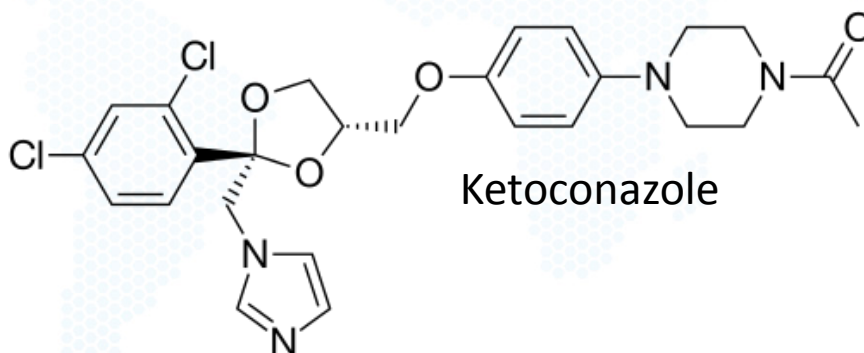
Drug Transport Technology at XenoTech

	FDA (2012)	EMA (2013)
Recommended transporters	P-gp, BCRP OATP1B1, OATP1B3, OAT1, OAT3, OCT2	P-gp, BCRP, BSEP OATP1B1, OATP1B3, OAT1, OAT3, OCT2
Transporters to consider	BSEP, MATE, MRP2 Others as necessary	OCT1, MATE1, MATE2 Others as necessary
Substrate evaluation	Yes (P-gp, BCRP) Others based on clearance/elimination routes	
Inhibitor evaluation	Yes, all of the recommended transporters (EMA and PDMA: in vitro data prior to Phase III trials)	

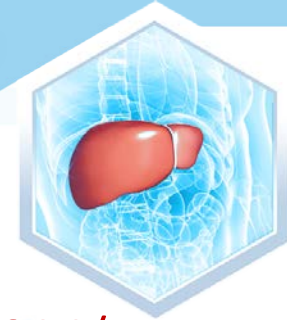


Ketoconazole

- Orally available
- Synthetic, broad spectrum, antifungal agent (imidazole)
- Approved in 1982 by FDA for use in fungal infections
- Known substrate and inhibitor of CYP3A4/5



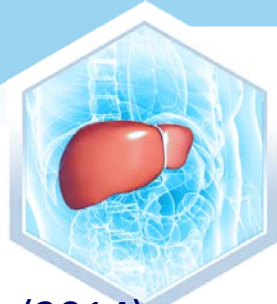
<https://www.nlm.nih.gov/>



Statement of the problem

High dose ketoconazole was previously the gold-standard strong CYP3A4/5 inhibitor used in clinical DDI studies

- Ketoconazole was essentially banned by FDA from clinical use in 2013
 - Typically asymptomatic, reversible liver function test abnormalities
 - 1984 estimate by Van Tyle: DILI in 0.1 to 1.0% of patients. No association with dosage, but increase with duration (e.g., months)
 - Later estimates: 1) 134.1 per 100,000 person-months; 2) 4.9 cases per 10,000 patients; 3) 3.6 to 4.2% (Reviewed by Greenblatt and Greenblatt, 2014)
- The FDA specifically recommended clarithromycin or itraconazole as alternatives for DDI studies, but noted other drugs may be used
- Ritonavir has been suggested as an alternative CYP3A4/5 inhibitor by some (Greenblatt et al., 2014 and 2015)
 - Excluded by Ke et al., due to non-specific CYP inhibition and induction



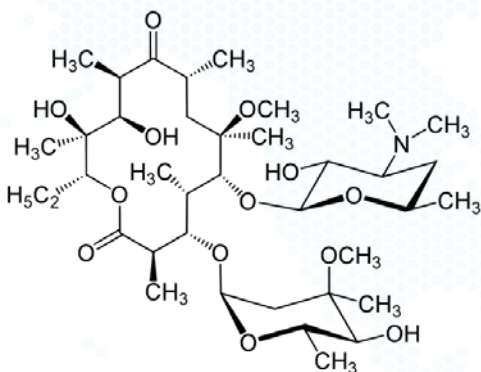
Statement of the problem (continued)

- 19 strong CYP3A inhibitors systematically evaluated by Ke and colleagues (2014). Only itraconazole and clarithromycin were deemed acceptable. Others excluded because:
 - 1) Drug not approved in U.S.
 - 2) Known to be a non-specific CYP inhibitor
 - 3) Significant safety issues
 - 4) Used exclusively in combination with ritonavir
 - 5) Are only moderate CYP3A4/5 inhibitors
- Some transporter data, but no previous comprehensive investigation of these alternative CYP3A4/5 inhibitors as transporter inhibitors
 - Ketoconazole IC_{50} values vary 37-fold for P-gp inhibition, 6.4-fold for OATP1B1 inhibition
 - Ritonavir: 85-fold variation for P-gp, 117-fold for OATP1B1 inhibition

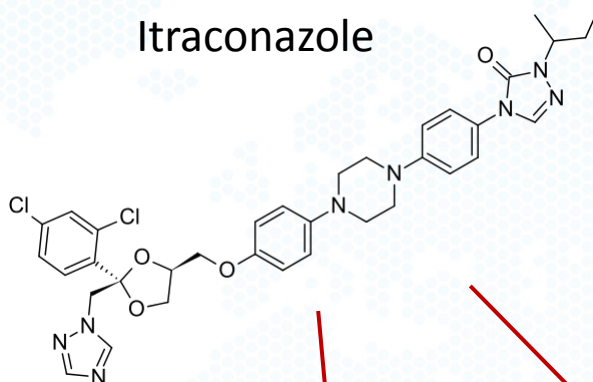
Goals of the study

Define the inhibition profile of the following compounds:

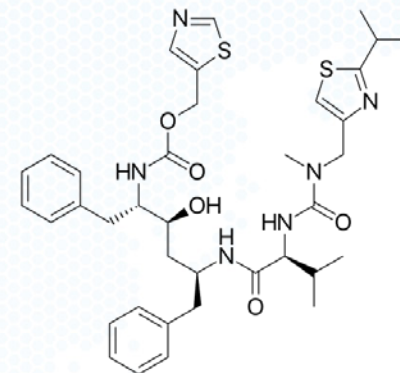
Clarithromycin



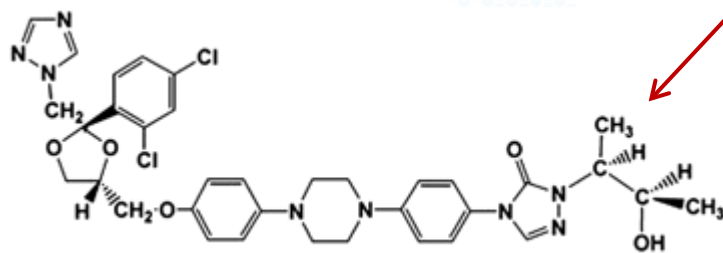
Itraconazole



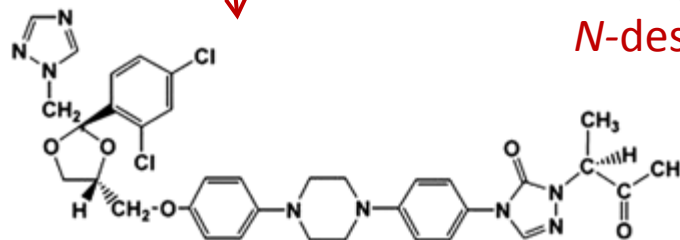
Ritonavir



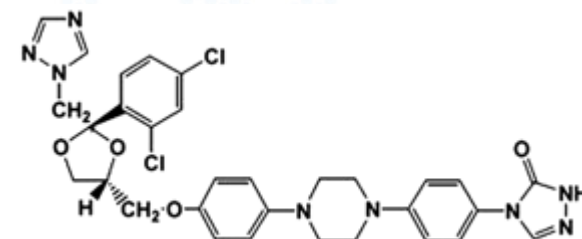
Hydroxyitraconazole

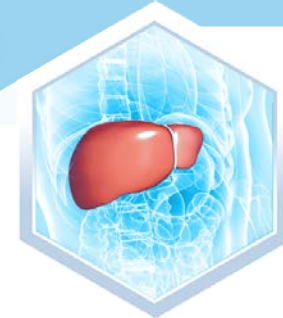


Ketoitraconazole



N-desalkylitraconazole



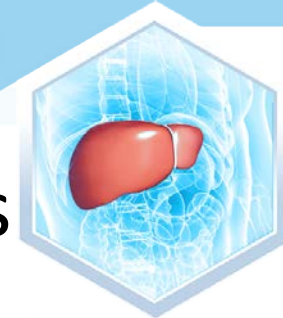


Goals of this study

To define the inhibition profile towards the following transporters:

OATP1B1	MATE2-K
OATP1B3	P-gp
OAT1	BCRP
OAT3	MRP2
OCT1	MRP3
OCT2	BSEP
MATE1	

To allow a more informed choice of a strong clinical CYP3A4/5 inhibitor for clinical DDI studies with a drug candidate known to be a substrate of one or more of these transporters: reduce confounding DDI results



Publication details / acknowledgements

1521-009X/44/3/1-7\$25.00

DRUG METABOLISM AND DISPOSITION

Copyright © 2016 by The American Society for Pharmacology and Experimental Therapeutics

<http://dx.doi.org/10.1124/dmd.115.067744>

Drug Metab Dispos 44:1-7, March 2016

Evaluation of Ketoconazole and Its Alternative Clinical CYP3A4/5 Inhibitors as Inhibitors of Drug Transporters: The In Vitro Effects of Ketoconazole, Ritonavir, Clarithromycin, and Itraconazole on 13 Clinically-Relevant Drug Transporters^[S]

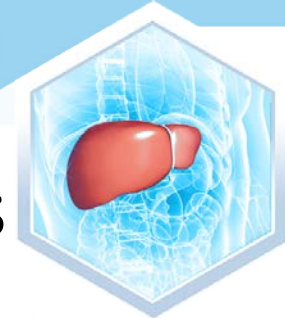
Lydia M. M. Vermeer,¹ Caleb D. Isringhausen,¹ Brian W. Ogilvie, and David B. Buckley

XenoTech, Kansas City, Kansas

Received October 8, 2015; accepted December 11, 2015

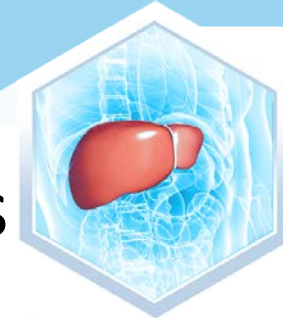
Drug Metab Dispos; Published ahead of print December 14, 2015;

doi:10.1124/dmd.115.067744



Methods – test systems and substrates

- HEK-293 assays:
 - Transfected (immortalized cell line with stable transfection) and control (wild type) cells were utilized (OATP1B1, OATP1B3, OAT1, OAT3, OCT1, and OCT2)
 - Transient transfection for MATE1 and MATE2-K
 - Seeded and cultured for 24 hours and supplemented with 2 mM sodium butyrate (OATP1B1, OAT1B3, MATE1, and MATE2-K) or fed with fresh supplemented DMEM (OAT1, OAT3, OCT1, and OCT2)
- Vesicle assays:
 - Membrane vesicles expressing P-gp, BCRP, MRP2, MRP3, and BSEP
 - ATP-dependent test system



Methods – test systems and substrates

HEK-293 assays

Membrane vesicle assays

Transporter	Substrate
OATP1B1	$[^3\text{H}]$ -Estradiol-17 β -glucuronide (50 nM)
OATP1B3	
OAT1	$[^3\text{H}]$ - <i>p</i> -Aminohippurate (1 μM)
OAT3	$[^3\text{H}]$ -Estrone 3-sulfate (50 nM)
OCT1	$[^3\text{H}]$ -Tetraethyl ammonium bromide (5 μM)
OCT2	$[^{14}\text{C}]$ -Metformin (10 μM)
MATE1	
MATE2-K	

Transporter	Substrate
P-gp	<i>N</i> -Methylquinidine (0.5 μM)
BCRP	$[^3\text{H}]$ -Estrone-3-sulfate (1 μM)
MRP2	$[^3\text{H}]$ -Estradiol -17 β -glucuronide (50 nM)
MRP3	
BSEP	$[^3\text{H}]$ -Taurocholic acid (0.4 μM)

- Inhibitors were pre-incubated with HEK-293 cells or vesicles for 15 min prior to addition of substrates to minimize effects of time-dependency and non-specific binding
- $[\text{Substrate}] = \sim 1\text{-}10\%$ of K_m for each assay (such that $\text{IC}_{50} \approx K_i$)

Methods – inhibitors in HEK-293 cells

CYP3A4 inhibitor	OATP1B1	OATP1B3	OAT1	OAT3	OCT1	OCT2	MATE1	MATE2-K
Ketoconazole	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.01 - 2 μ M	0.01 - 2 μ M
Itraconazole	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M
Hydroxy-itraconazole	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.001 - 0.3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M
Keto-itraconazole	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M
N-deskalkyl itraconazole	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M
Clarithromycin	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M
Ritonavir	0.03 - 10 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.1 - 20 μ M	0.1 - 20 μ M

- Concentrations based on clinical PK data for each inhibitor in commonly used dosing regimens in clinical CYP3A4/5 DDI studies, up to 10-fold higher than the average $C_{\max,ss}$
- Where concentrations are lower, an initial range-finding IC_{50} experiment was performed, and concentrations adjusted lower, as needed (values in red)



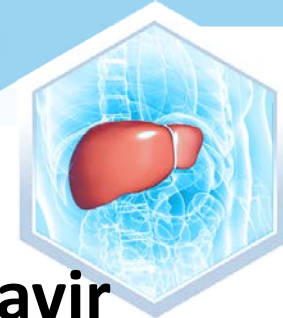
Methods – inhibitors in membrane vesicles

CYP3A4 inhibitor	P-gp	BCRP	MRP2	MRP3	BSEP
Ketoconazole	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M	0.1 - 20 μ M
Itraconazole	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M	0.03 - 10 μ M
Hydroxyitraconazole	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M
Ketoitraconazole	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M	0.01 - 3 μ M
N-deskalkyl itraconazole	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M	0.001 - 0.2 μ M
Clarithromycin	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M	0.3 - 50 μ M
Ritonavir	0.01 - 3 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.3 - 30 μ M	0.3 - 30 μ M

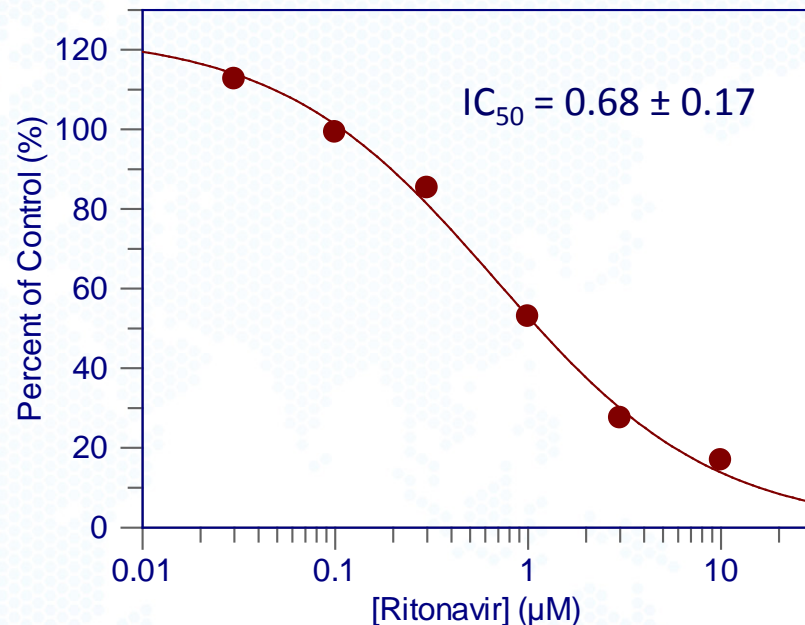
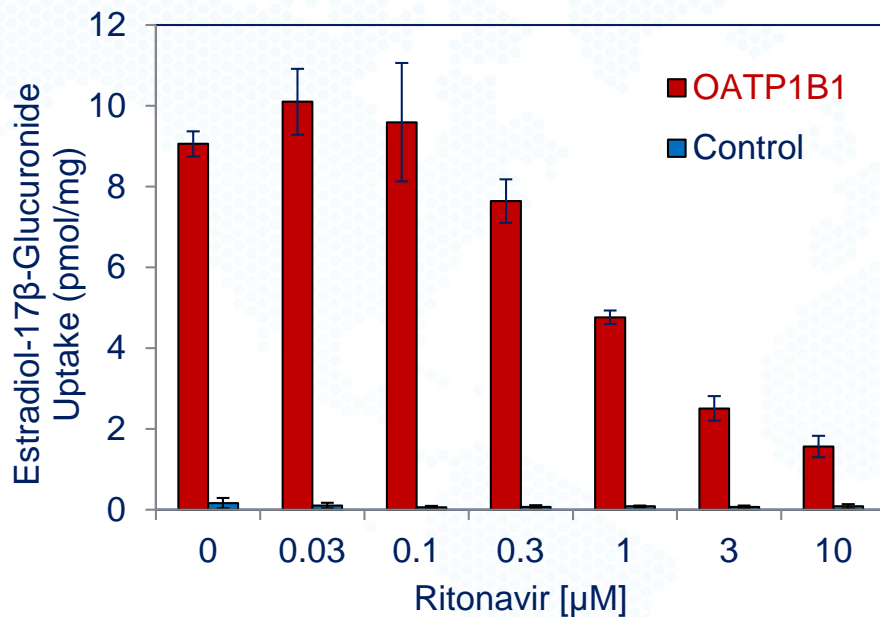
- Prototypical inhibitors used as positive controls for all assays

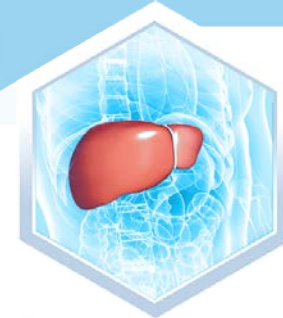
Methods – prototypical inhibitors

Transporter	Inhibitor
OATP1B1	Rifampin (10 μ M)
OATP1B3	
OAT1	Probenecid (100 μ M)
OAT3	
OCT1	Quinidine (300 μ M)
OCT2	
MATE1	Cimetidine (1000 μ M)
MATE2-K	
P-gp	Verapamil (60 μ M)
BCRP	Ko143 (1 μ M)
MRP2	Benzbromarone (100 μ M)
MRP3	
BSEP	Cyclosporine (20 μ M)

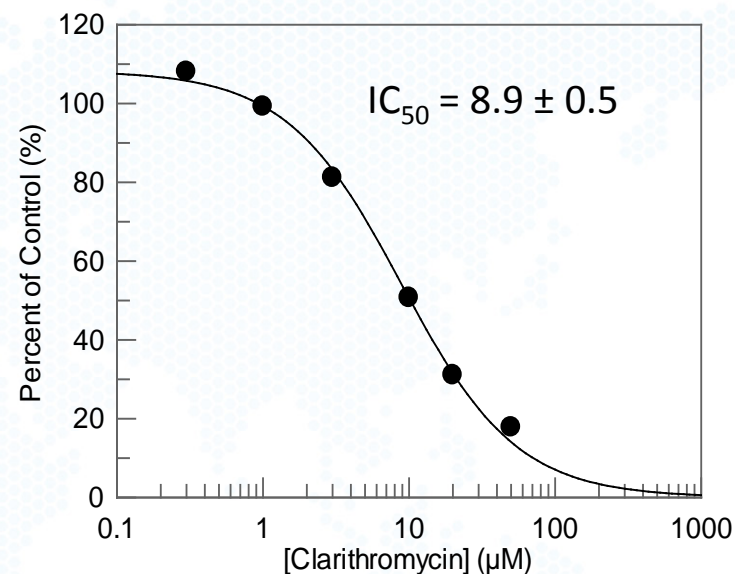
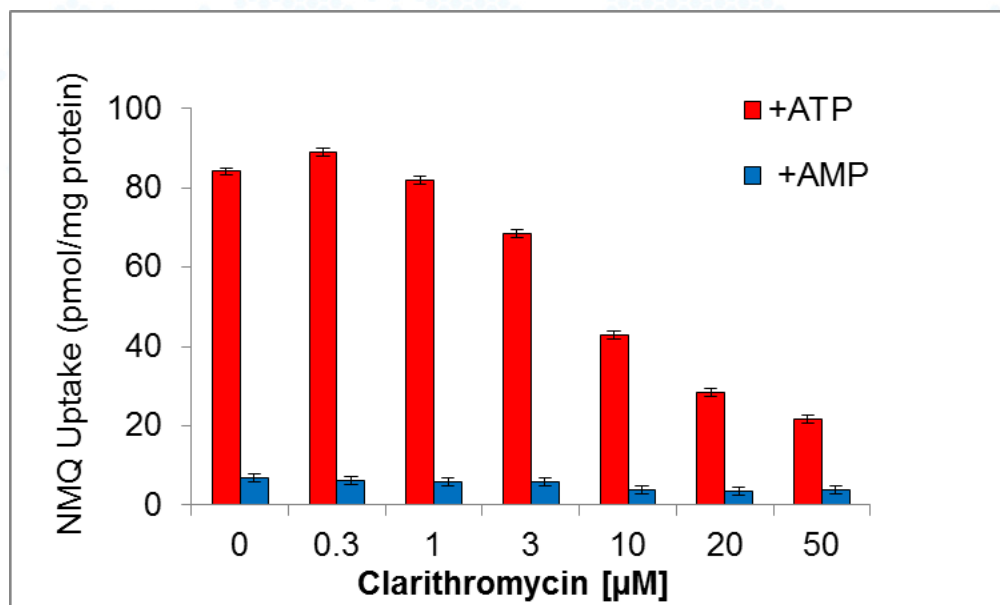


Results – Example: Inhibition of OATP1B1 by ritonavir





Results – Example: Inhibition of P-gp by Clarithromycin



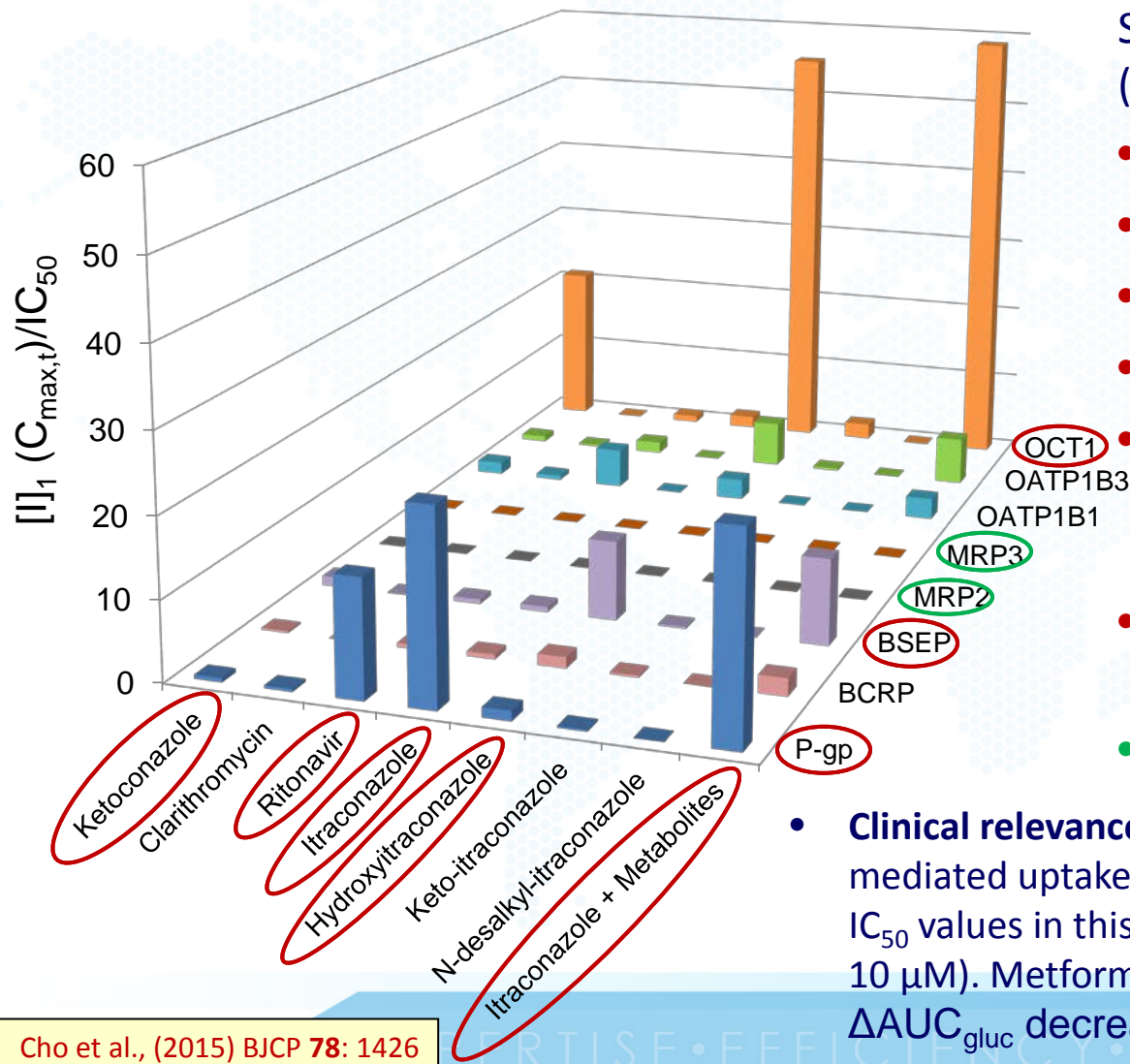
Results – IC50 values (μM)

Transporter	Ketoconazole	<u>Itraconazole</u>	Hydroxy- itraconazole	Keto- itraconazole	<u>N-Desalkyl- itraconazole</u>	<u>Clarithromycin</u>	Ritonavir
OATP1B1	1.8 ± 0.2	>10	0.23 ± 0.03	0.29 ± 0.04	>0.2	5.3 ± 1.3	0.68 ± 0.17
OATP1B3	3.9 ± 0.6	>10	0.10 ± 0.01	0.088 ± 0.035	>0.2	14 ± 2	2.3 ± 0.4
OAT1	5.7 ± 0.5	>10	>3	>3	>0.2	>50	17 ± 3
OAT3	0.86 ± 0.68	>10	2.0 ± 0.3	>3	>0.2	>50	>30
OCT1	0.13 ± 0.03	0.74 ± 0.24	0.01 ± 0.00	0.04 ± 0.01	>0.2	>50	4.1 ± 0.6
OCT2	0.89 ± 0.35	>10	>3	>3	>0.2	>50	>30
MATE1	0.37 ± 0.03	>10	0.84 ± 0.21	1.1 ± 0.2	>0.2	>50	1.2 ± 0.2
<u>MATE2-K</u>	>2	>10	>3	>3	>0.2	>50	15 ± 2
P-gp	5.6 ± 0.4	0.048 ± 0.04	0.49 ± 0.14	0.12 ± 0.12	0.26 ± 0.05	8.9 ± 0.5	0.24 ± 0.02
BCRP	12 ± 9	1.9 ± 0.3	0.44 ± 0.03	0.10 ± 0.01	>0.2	>50	6.6 ± 0.5
<u>MRP2</u>	>20	>10	>3	>3	>0.2	>50	>30
<u>MRP3</u>	>20	>10	>3	>3	>0.2	>50	>30
BSEP	2.4 ± 0.6	1.8 ± 0.0	1.2 ± 0.8	0.11 ± 0.02	>0.2	59 ± 8	6.1 ± 0.9

Compounds in **yellow** inhibited the fewest transporters
Transporters in **green** were inhibited by the fewest compounds



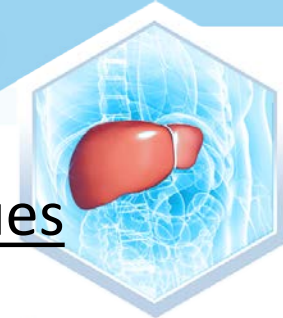
DDI predictions for hepatic uptake and efflux transport based on the FDA's basic model



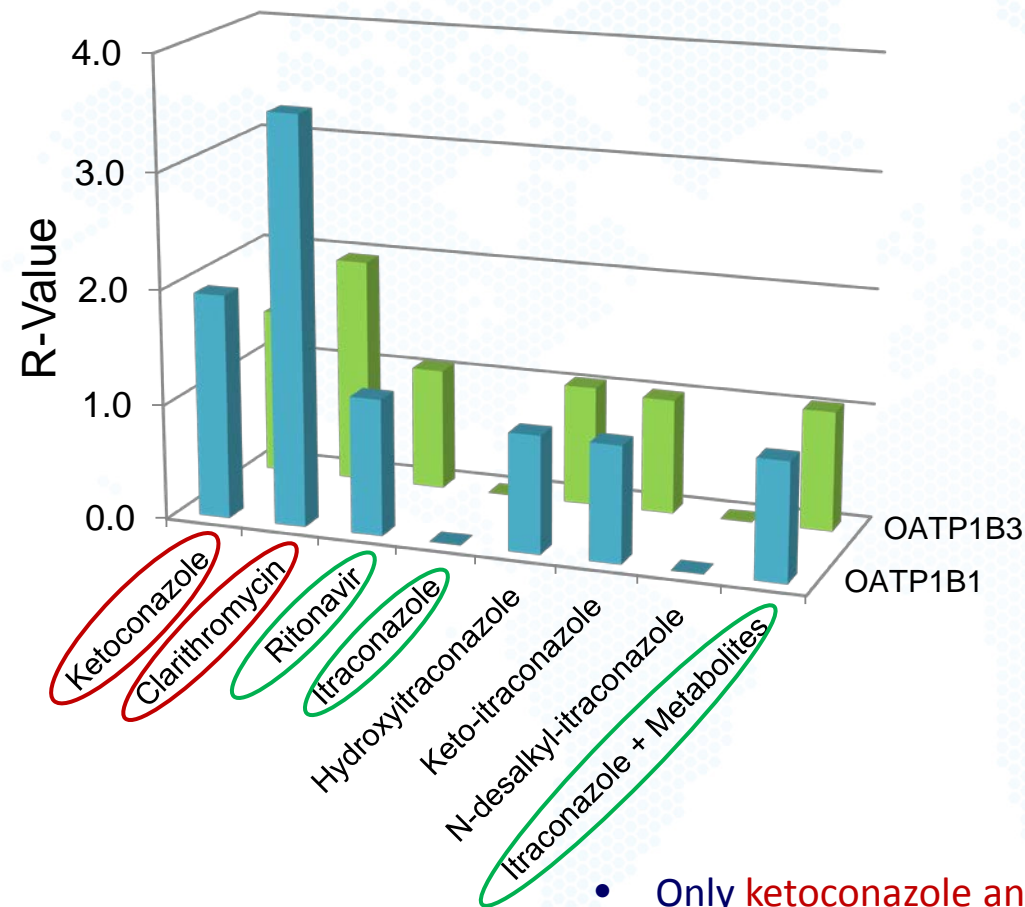
Some very high $[I]_1/IC_{50}$ values (>100-fold the 0.1 cut-off):

- Ketoconazole for OCT1 (~200x)
- Ritonavir for P-gp (~150x)
- Itraconazole for P-gp (~240x)
- OH-Itraconazole for OCT1 (550x)
- Sum of itraconazole + metabolites >250-fold cut-off for P-gp, BSEP and OCT1
- No inhibitors with all values <0.1
- MRPs not affected by any

- **Clinical relevance?** Verapamil inhibits OCT1-mediated uptake of metformin ($IC_{50}=13 \mu M$; Most IC_{50} values in this study much less than $10 \mu M$). Metformin plasma PK not altered, but ΔAUC_{gluc} decreased 240%.



DDI predictions for OATP-mediated hepatic uptake: R-values

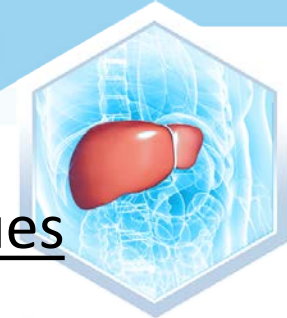


$$R \text{ value} = 1 + (f_u * \frac{I_{in,max}}{IC_{50}})$$

Where:

- f_u = fraction unbound of the inhibitor
- $I_{in,max}$ = estimated maximum inhibitor concentration at the inlet to the liver
- $I_{in,max} = C_{max} + (k_a * Dose * \frac{F_a F_g}{Q_h})$
- k_a = absorption rate constant of the inhibitor (assumed 0.1)
- $F_a F_g$ = fraction of the dose of inhibitor absorbed (assumed 1)
- C_{max} = maximum systemic plasma concentration of inhibitor

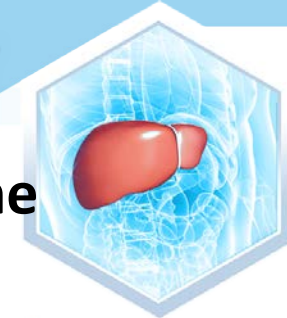
- Only **ketoconazole** and **clarithromycin** predicted to affect OATP1B1 and 1B3 using the R-value method with cut-off ≥ 1.25 .
- Neither **ritonavir** nor **itraconazole** had values > 1.25



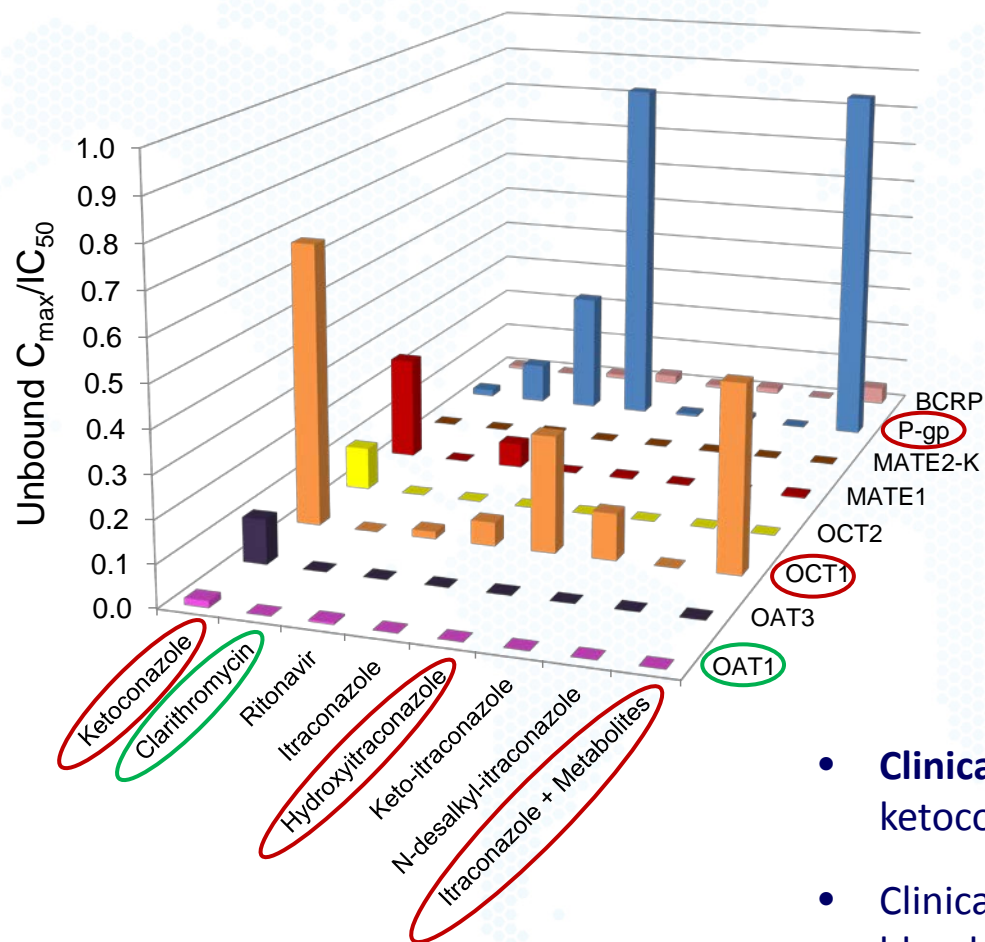
DDI predictions for OATP-mediated hepatic uptake: R-values

- Clinical relevance: Ketoconazole ↑ bosentan AUC by 122%
- Clarithromycin and:
 - Bosentan (273% ↑ AUC)
 - Glyburide (33% ↑ AUC)
 - Pravastatin (111% ↑ AUC)
- OATP inhibition explains only part of these interactions, with CYP3A4 inhibition also playing a significant role

University of Washington DIBD (2015)

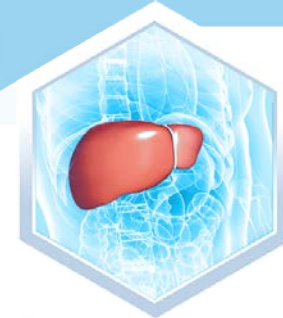


DDI predictions for renal or BBB transporters based on the FDA's basic model

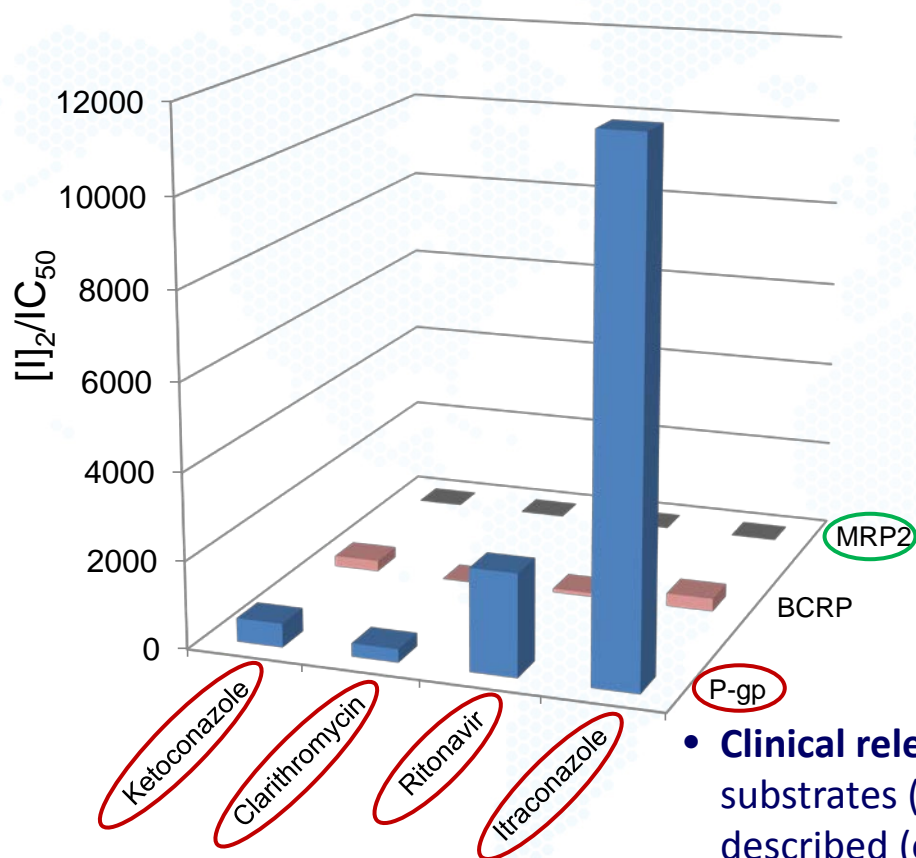


- With the unbound C_{\max}/IC_{50} value cut-off >0.1 all transporters except **OAT1** were predicted to be affected.
- All values were <10 -fold higher than the cut-off
- P-gp and OCT1 were most affected
- **Ketoconazole and OH-itraconazole** had the greatest impacts
- **Clarithromycin** did not have any values >0.1

- **Clinical relevance?** Only MATE1 inhibition by ketoconazole mentioned as possibly relevant
- Clinically relevant inhibition of drug efflux at the blood-brain-barrier is unlikely to occur



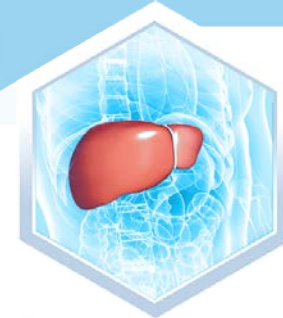
DDI predictions for intestinal efflux transporters based on the FDA's basic model



- With the $[I]_2/IC_{50}$ value cut-off >10, **MRP2** was not predicted to be affected by any inhibitors.
- For P-gp, $[I]_2/IC_{50}$ values were:
~50-, 30-, 230- and 1000-fold the cut-off for **ketoconazole**, **clarithromycin**, **ritonavir** and **itraconazole**, respectively
- For BCRP $[I]_2/IC_{50}$ values were:
25-, 8- and 30-fold the cut-off for **ketoconazole**, **ritonavir** and **itraconazole**, respectively
- **Clarithromycin** was not predicted to inhibit intestinal BCRP

• **Clinical relevance?** Interactions between all inhibitors and P-gp substrates (with minimal CYP3A4 contribution) are well described (e.g., fexofenadine, quinidine, dabigatran, digoxin - up to 2.5 fold AUC increase).

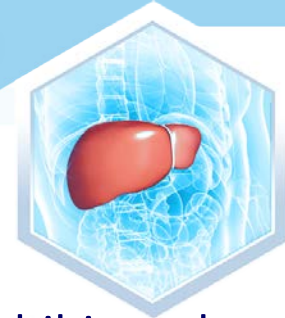
• BCRP mentioned in DDI between rosuvastatin and ritonavir



Potential limitations of the study

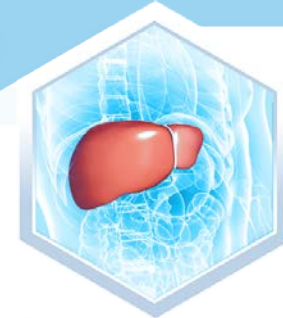
- The potential for non-specific binding of the inhibitors in the test systems was not directly measured
- If measured, should the free inhibitor concentration be evaluated with or without test system present and should pre-incubations be employed?
- Because of these questions, a consistent design was used throughout with a pre-incubation with inhibitor
- These studies were designed to use similar methodologies reported in recent publications (Brouwer et al., 2013; Izumi et al., 2013; Izumi et al., 2015; Shitara et al., 2013; Taub et al., 2011; Zamek-Gliszczynski et al., 2013)
- Designed to *minimize* effects of time-dependency and non-specific binding

Brouwer et al. (2013) CPT **94**: 95; Izumi et al. (2013) DMD **41**: 1859; Izumi et al. (2015) DMD **43**: 235; Shitara et al. (2013) JPS **102**: 3427; Taub et al. (2011) DMD **39**: 2093; Zamek-Gliszczynski et al. (2013) CPT **94**: 67.



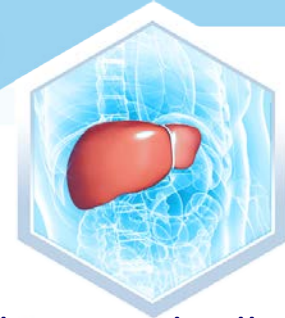
Summary – Part 1

- Of the alternatives to ketoconazole as a strong clinical CYP3A4/5 inhibitor, the following conclusions can be made:
 - MRP2 and MRP3 were not significantly inhibited by any at clinically relevant concentrations
 - MATE2-K was only inhibited by ritonavir
 - Itraconazole, *N*-desalkylitraconazole and clarithromycin inhibited the fewest transporters
- Based on the FDA's basic hepatic model $[I]_1/IC_{50}$:
 - **OCT1** was the most potently inhibited transporter, with values >250-fold the cut-off of 0.1 for **itraconazole + metabolites**
 - **P-gp** was most potently inhibited by **ritonavir**
 - No inhibitors had $[I]_1/IC_{50} < 0.1$ for all transporters, except MRPs



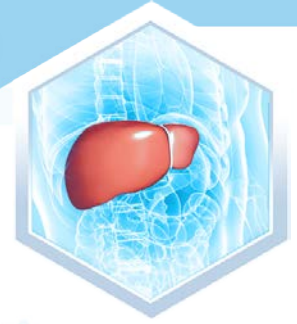
Summary – Part 2

- Based on the FDA's R-value hepatic model for OATP1B1 and 1B3:
 - **Ritonavir** and **itraconazole** have values <1.25
 - **Clarithromycin** most potently inhibits OATP1B1 (R value = 3.5)
- Based on the FDA's basic renal and BBB model (unbound C_{\max}/IC_{50}):
 - **OAT1** is predicted to be unaffected by all inhibitors
 - None of the unbound C_{\max}/IC_{50} values were >10-fold the cut-off
 - **Hydroxy-itraconazole** had the greatest impacts
 - **Clarithromycin** did not have any unbound C_{\max}/IC_{50} values >0.1
- Based on the FDA's basic intestinal model ($[I]_2/IC_{50}$):
 - **MRP2** is predicted to be unaffected by all inhibitors
 - **P-gp** and **BCRP** were predicted to be significantly affected by the inhibitors with the exception of **clarithromycin** for **BCRP**



Conclusions

- None of the alternatives to ketoconazole provided a clean inhibition profile towards all 13 drug transporters evaluated
- The alternatives to ketoconazole each have unique transporter inhibition profiles
- MRP2 and MRP3 were not inhibited by any alternative inhibitors
- Ritonavir and itraconazole may be the best alternatives for CYP3A4/5 substrates that are transported by OATP1B1 and 1B3
- CYP3A4/5 substrates that are transported by OAT1 may not be affected by any of these alternative inhibitors
- Clarithromycin may be the best choice for substrates of renal transporters
- For substrates of intestinal P-gp or BCRP, clarithromycin may be the best choice (although P-gp is still predicted to be affected)
- The best choice for a strong clinical CYP3A4/5 inhibitor depends on the unique transporter substrate profile of the drug candidate



Acknowledgments

- Andrea Wolff
- XenoTech's Analytical Services group

Upcoming Show

AAPS/FDA/ITC Joint Workshop on

Drug Transporters in ADME:
From the Bench to the Bedside

April 18-20, 2016

Renaissance Baltimore Harborplace Hotel



AAPS ITC Drug Transporter Workshop

April 18-20, 2016 – Baltimore, MD

Booth # 4

Dr. Lydia Vermeer & Lindsey Harston

Senior Scientist, Drug Transport
lvermeer@xenotechllc.com



Technical Account Manager
lhaston@xenotechllc.com

AAPS Drug Transporter Workshop Poster Presentation

Poster # M1032
Dr. Lydia Vermeer

Evaluation of Ketoconazole and Its Alternative Clinical CYP3A4/5 Inhibitors as Inhibitors of Drug Transporters: The In Vitro Effects of Ketoconazole, Ritonavir, Clarithromycin and Itraconazole on 13 Clinically-Relevant Drug Transporters

Tuesday, April 19th
5:15 – 6:30pm

Drug Transporter Services at XenoTech

- Over 50 transporters assays available (Tokai & Kansas site)
- Variety of models available
 - Cell lines
 - Vesicles
 - Transfected cells
 - Oocytes
 - Hepatocytes
 - Animal models
- GLP & non-GLP study options



Questions