

Ki determination in vitro and its impact on modelling and prediction of DDI

Yuichi Sugiyama

**Sugiyama Laboratory, RIKEN Innovation Center,
RIKEN, Research Cluster for Innovation,**

8th Meet the Experts Transporter Conference in Budapest

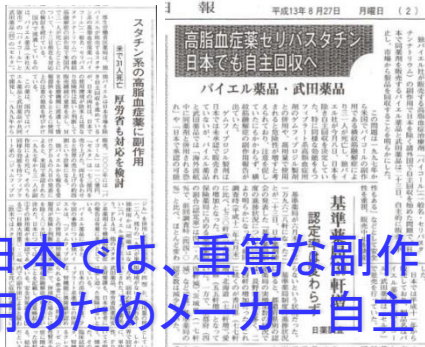


**April 26-27, 2018
Continental Hotel Budapest**

Contents

- 1) Introduction;**
Rate-determining process (focusing on the liver)
(Uptake, efflux, elimination, metabolism)
- 2) PBPK model based analysis of OATPs mediated drug-drug interaction (Top down + Bottom-up)
- 3) Other collaborations
Integrated model of rifampicin as a perpetrator
Time-dependent inhibition
Substrate-dependent inhibition

Drug-interaction between Cerivastatin and Gemfibrozil/CsA



日本では、重篤な副作用のためメーカー自主回収へ

薬事日報平成13年8月27日号

薬事日報平成13年8月15日号

WORLD HEALTH ORGANIZATION



ORGANISATION MONDIALE DE LA SANTE

QSM/MC/IEA.102

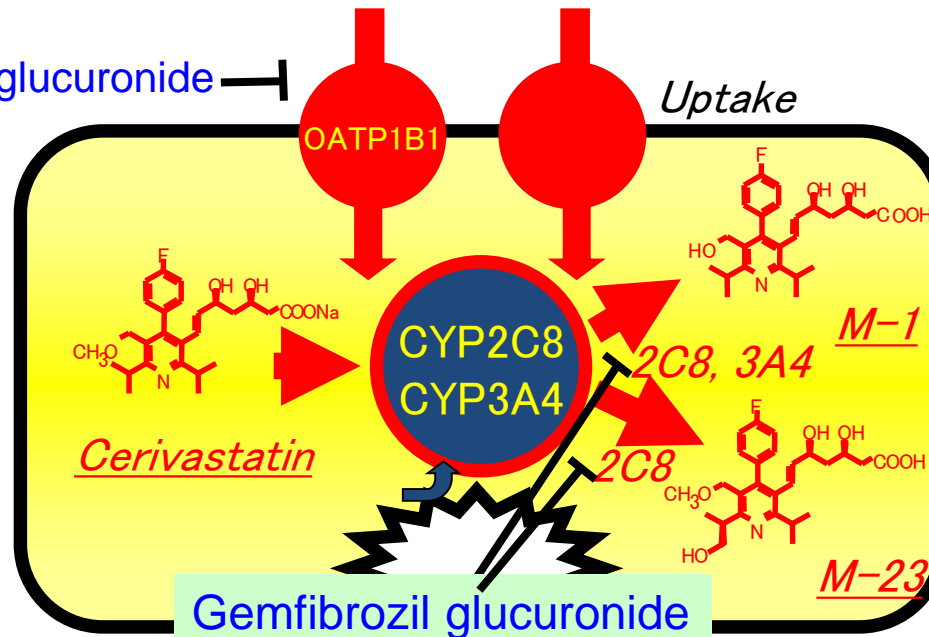
09 August 2001

Voluntary withdrawal of Cerivastatin – Reports of Rhabdomyolysis

Baycol (cerivastatin) was initially approved in the U.S. in 1997. It belongs to a group of cholesterol lowering drugs referred to as “statins”. While all statins could potentially cause this dangerous muscle reaction, rhabdomyolysis appears more frequent with cerivastatin, especially when used in high doses, in the elderly or, when taken along with gemfibrozil, another cholesterol lowering drug. In this connection it may be noted that Bayer has withdrawn all

- **52 patients died (US 31).**
- **Among 31 patients, 12 were given also gemfibrozil.**

Gemfibrozil glucuronide



Shitara, Y. et al.
J Pharmacol Exp Ther,
304(2): 610-6 (2003)

Shitara, Y. et al.
J Pharmacol Exp Ther,
311(1): 228-36 (2004)

Shitara, Y. and Sugiyama Y.
Pharmacol Ther,
112(1): 71-105 (2006)

Gemfibrozil glucuronide functions as **a dual inhibitor** of OATP1B1 and CYP2C8

Examples of substrates for uptake/efflux transporters and enzymes (1)

Substrates	Uptake transporter	Metabolic enzymes	Efflux transporter
Anti-Hyperlipidemic drugs (statins)			
atorvastatin	OATPs	CYP3A4	-
cerivastatin	OATPs	CYP2C8, 3A4	-
fluvastatin	OATPs	CYP2C9	-
pravastatin	OATPs	-	MRP2
rosuvastatin, pitavastatin	OATPs	-	BCRP
Anti-hypertension or -cardiovascular disease			
bosentan	OATPs	CYP3A4, 2C9	-
torasemide	OATPs	CYP2C9	-
telmisartan	OATP1B3	UGTs	-
valsartan	OATPs	-	MRP2
Anti-cancer drug			
docetaxel	OATP1B3	CYP3A4	-

Examples of substrates for uptake/efflux transporters and enzymes (2)

Substrates	Uptake transporter	Metabolic enzymes	Efflux transporter
Anti-diabetes			
repaglinide	OATPs	CYP2C8, 3A4	-
nateglinide, glibenclamide	OATPs	CYP2C9, 3A4	
Anti-HCV			
simeprevir, grazoprevir	OATP1B1	CYP3A4	-
asunaprevir, danoprevir, paritaprevir	OATPs	CYP3A4	Pgp
Others			
Montelukast	OATP2B1	CYP2C8, 2C9, 3A4	-
maraviroc	OATP1B1	CYP3A4	Pgp
fexofenadine	OATPs	-	Pgp

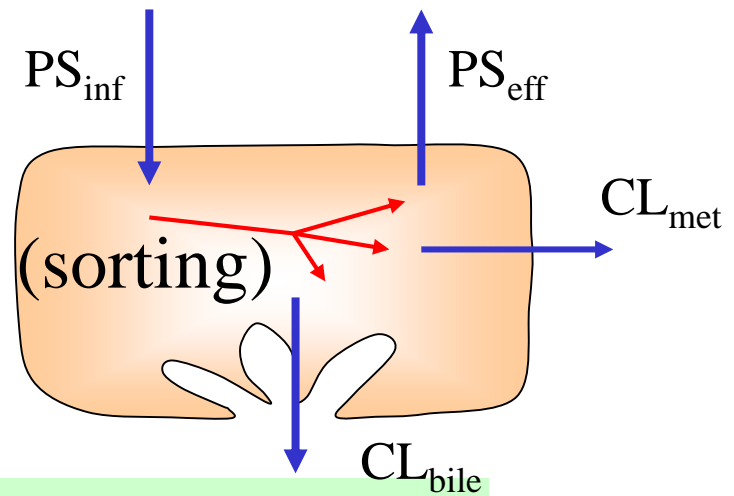
Extended clearance concept

$$CL_{int,all} = PS_{inf} \times \frac{CL_{bile} + CL_{met}}{PS_{eff} + CL_{bile} + CL_{met}}$$

R_{dif} values

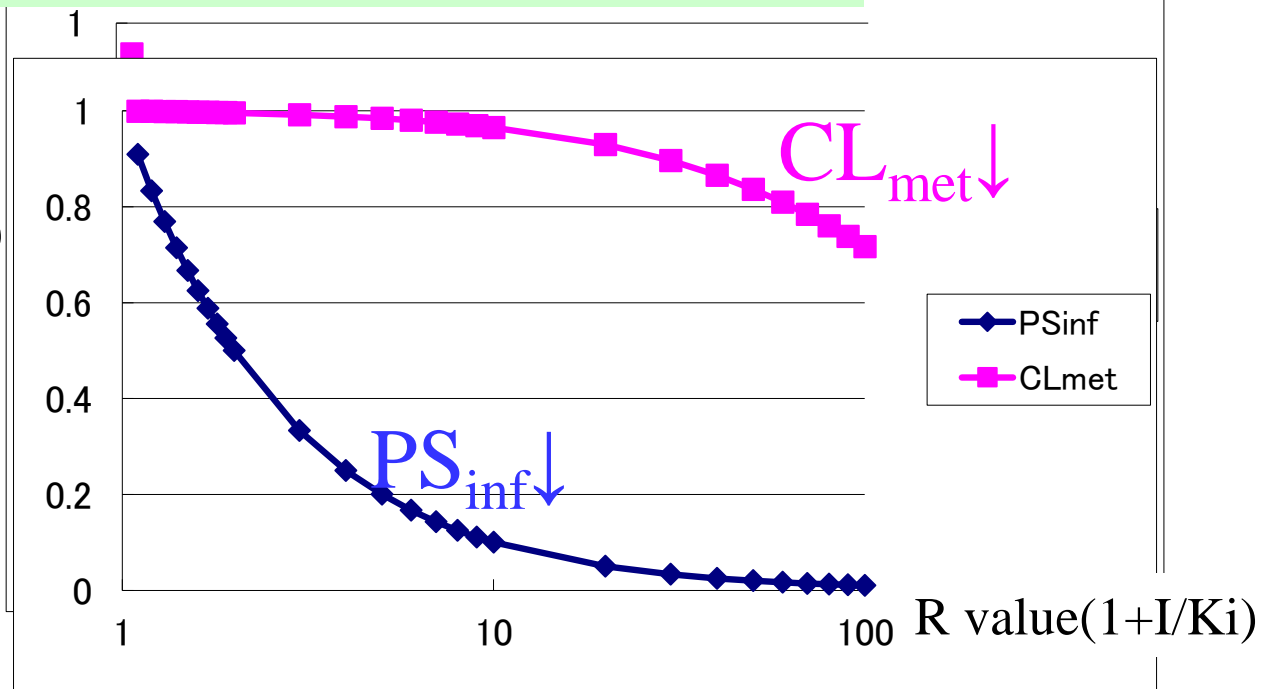
$$= PS_{inf,dif} / PS_{inf,act}$$

(β value)



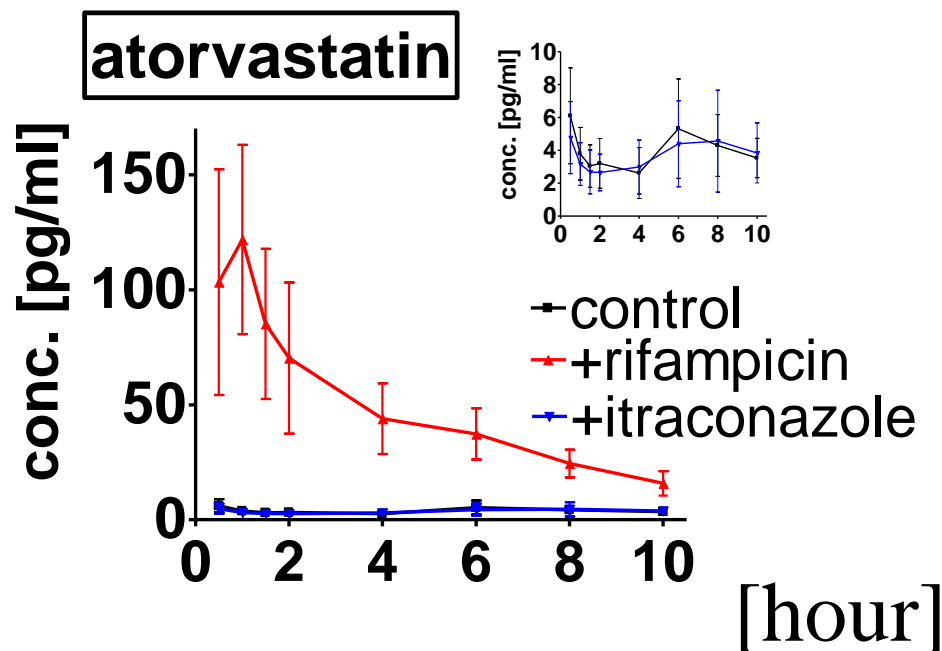
$PS_{inf}=100, PS_{eff}=2, CL_{bile}+CL_{met}=500$ (Case-1, limited by uptake) netabolism)

% of $CL_{int,all}(cont.)$
% $CL_{int,all}(cont.)$
(1/AUC)
(1/AUC)



Impact of the function of each pathway on the overall intrinsic clearance

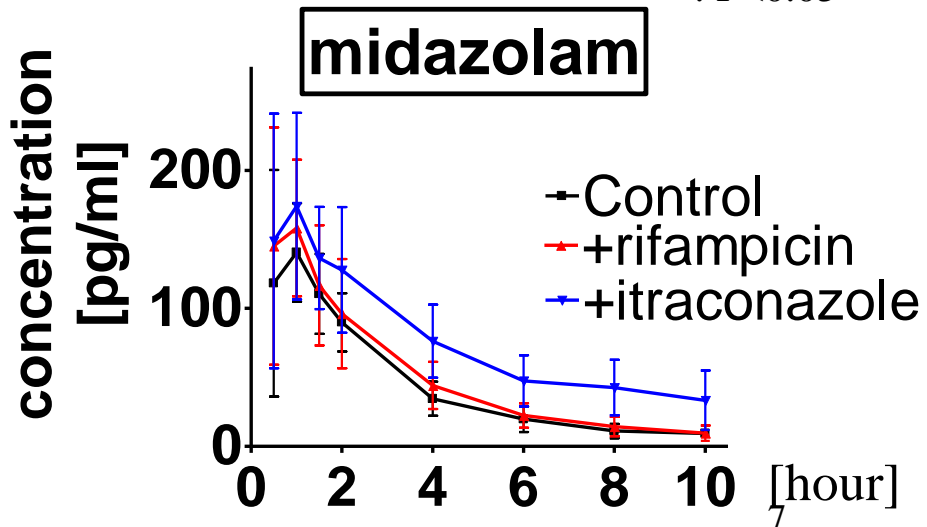
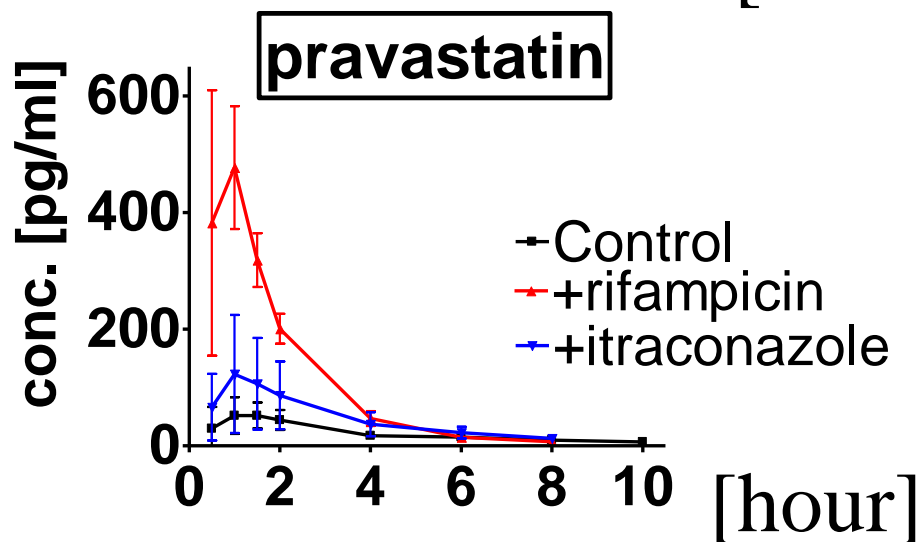
Plasma concentrations of atorvastatin was greatly increased by rifampicin, but not by itraconazole (Maeda K. et al., Clin Pharmacol Ther 90:575-581 (2011)).



AUC ₀₋₁₀ [pg*hr/ml]	ATV	PRV (AUC ₀₋₈)	MDZ
Cont.	38.5 ± 17.5	195 ± 78.7	434 ± 122
+RIF	439*** ± 134	949*** ± 179	471 ± 168
+ITZ	36.0 ± 19.2	386 ± 254	755* ± 276

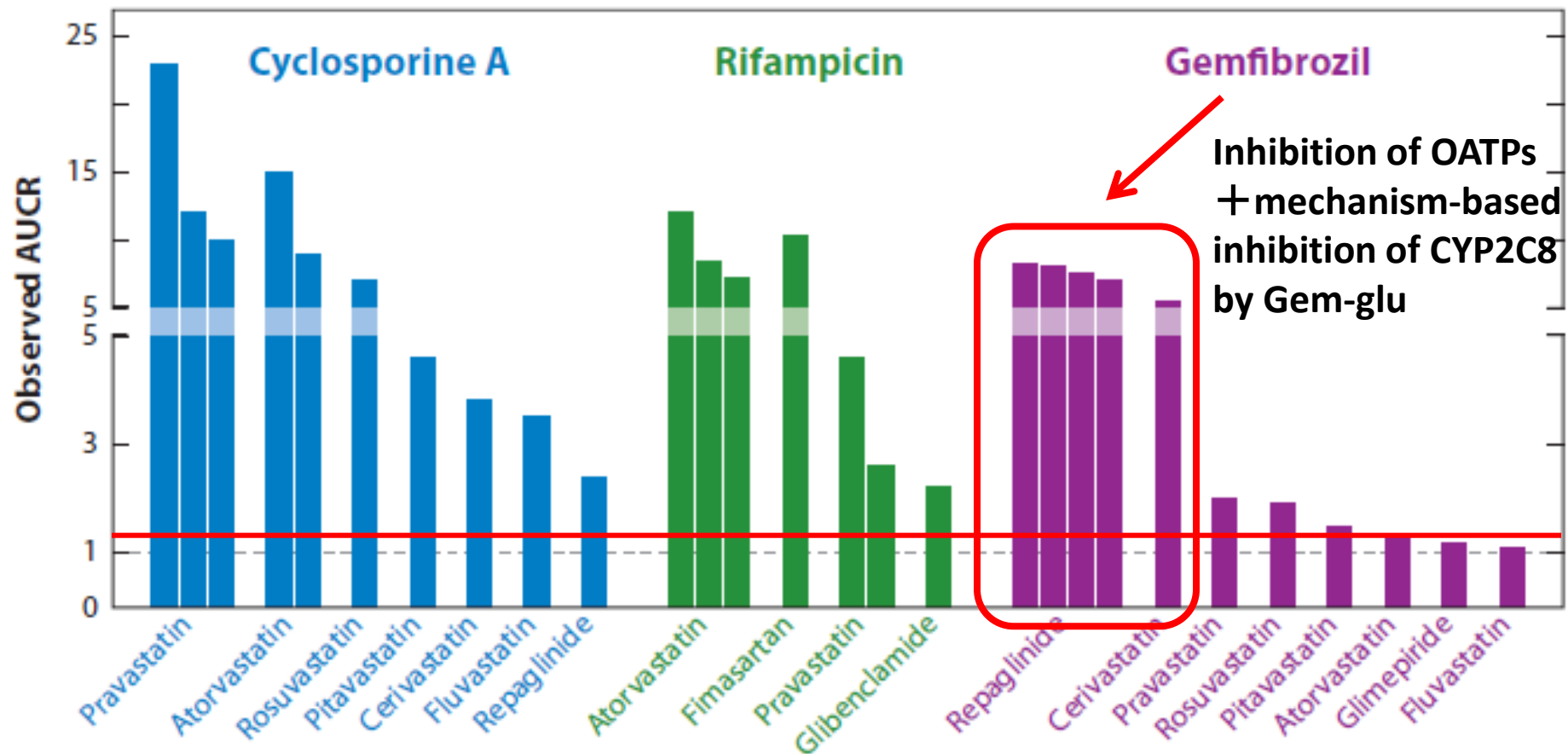
***: P<0.0005

*: P<0.05



*Doses of each substrates are 33μg

OATP1B1/1B3 - mediated DDIs



Atorvastatin, pravastatin, rosuvastatin exhibited relatively higher AUCR, while fluvastatin, repaglinide, and glibenclamide do lower AUCR.

The latter compounds are lipophilic and have higher R_{dif} and lower fo_{atp} values.

Separation of active uptake and passive uptake

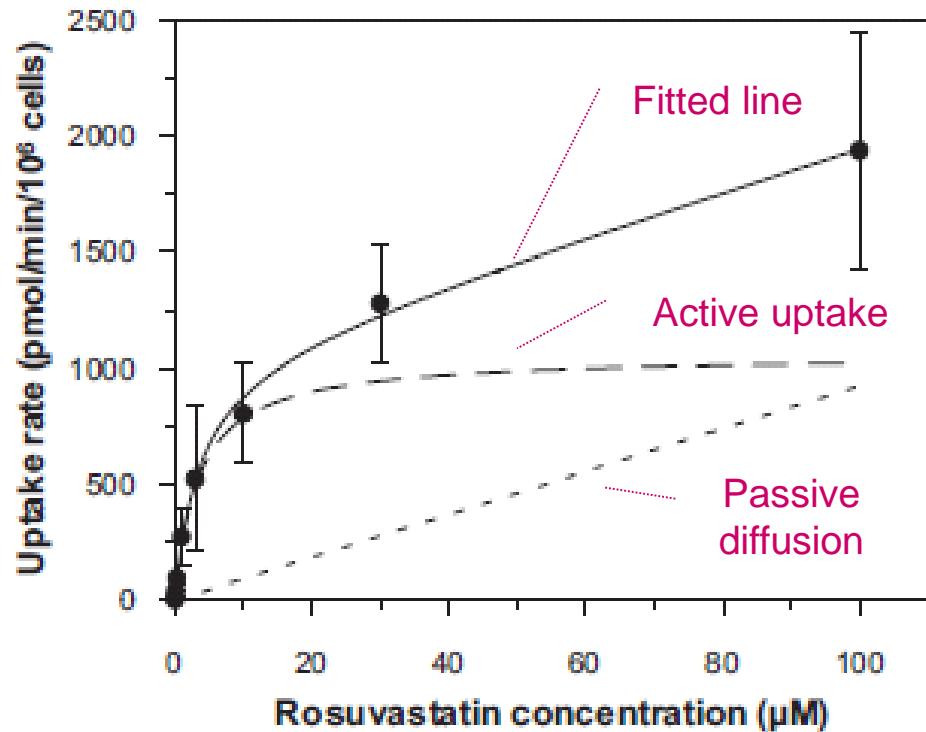
Yabe Y et al., *DMD* (2011) 39: 1808-14.

$$v = \frac{V_{\max} \times S}{K_m + S} + PS_{dif} \times S$$

R_{dif} values

$$= PS_{inf,dif} / PS_{inf,act}$$

$$= PS_{dif} / (V_{\max} / K_m)$$



Effect of fraction transported (foatp) by the affected transporter(OATP1Bs)

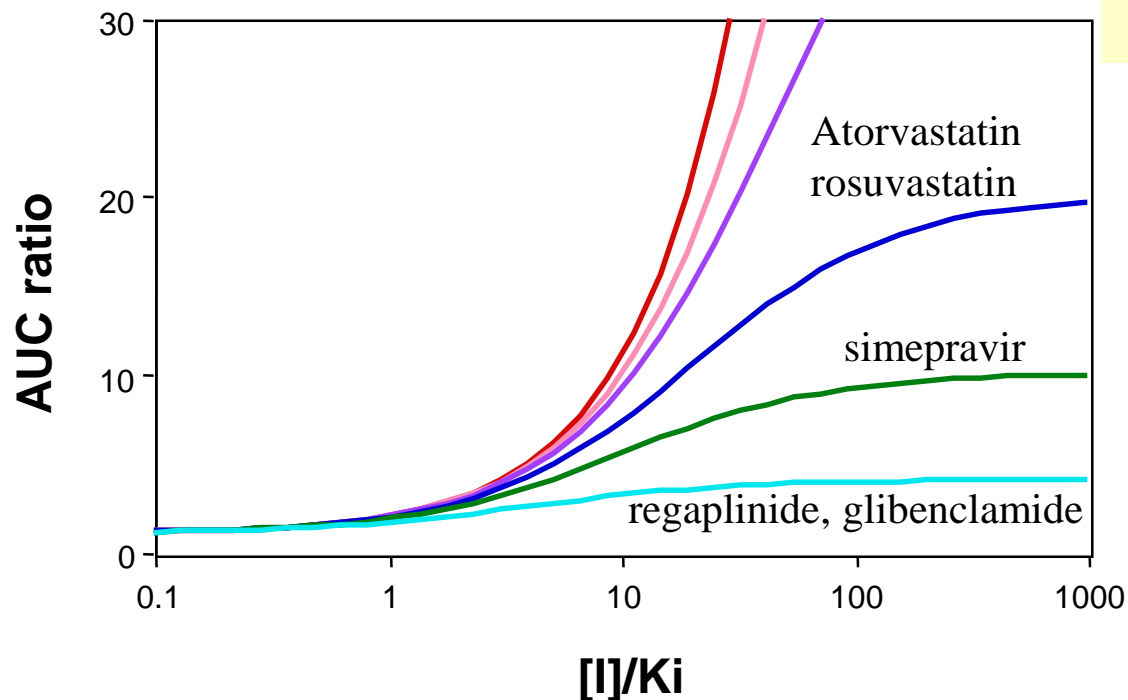
$$\frac{\text{AUC (+inhibitor)}}{\text{AUC (control)}} = \frac{1}{\frac{\text{fm}}{1+[I]/K_i} + (1 - \text{fm})}$$

$\text{fm} \Rightarrow \text{foatp}$

Lipophilicity of the substrate \uparrow

Then

the contribution of passive
diffusion increases, and foatp \downarrow



- foatp = 1
- foatp = 0.99
- foatp = 0.98
- foatp = 0.95
- foatp = 0.9
- foatp = 0.75

How to obtain the in vivo β value and R_{dif} values (=PS_{inf,dif}/PS_{inf,act}) from these clinical studies ?

- 1) R_{dif} values; Effect of single dose of rifampicin;
Atrovastatin AUCR=12, R_{dif} is known as 0.024 (in vitro; determined in my lab).
Then, from AUCR for other substrates together with the estimates of foatp(Rifampicin sensitive OATPs) value, **in vivo R_{dif} value of OATPs mediated uptake is estimated**

atorvastatin	AUCR=12.0	$R_{dif} = 0.024$
simeprevir	AUCR = 7.3	$R_{dif} = 0.07-0.10$
bosentan	AUCR=3.2	$R_{dif} = 0.3-0.4$
repaglinide	AUCR =2.0	$R_{dif} = 0.7-1.1$

- 2) β values; Effect of iv itraconazole(iv);
Midazolam AUCR=4.0 , fm(3A) is known as 0.9 (in vitro; determined in my lab)
Then, from AUCR for other substrates together with their fm values, in vivo β value of dual substrates of OATP and CYP3A4 is estimated (shown in the next slide)

(Summary)

$$\beta = \frac{CL_{bile} + CL_{met}}{PS_{eff} + CL_{bile} + CL_{met}}$$

ATV, Bosentan : Close to Case-1 (β high 0.8~1)

CER, REPG, SIMP: (β intermediate 0.3-0.7)

DAR (β low < 0.2)

From the clinical DDI studies so far done and also from the I/Ki value of ITZ (CYP3A4 inhibitor), we can estimate the β values of each compound. With high β value of substrate, the inhibition of metabolism and biliary excretion does not affect the blood AUC value of unchanged drug (uptake-limited case)

Yoshikado T et al., A Clinical Cassette Dosing Study for Evaluating the Contribution of Hepatic OATPs and CYP3A to Drug-Drug Interactions. Pharm Res. 34(8):1570-1583 (2017)

Contents

- 1) Introduction;
Rate-determining process (focusing on the liver)
(Uptake, efflux, elimination, metabolism)
- 2) **PBPK model based analysis of OATPs mediated drug-drug interaction (Top down + Bottom-up)**
- 3) Other collaborations
Integrated model of rifampicin as a perpetrator
Time-dependent inhibition
Substrate-dependent inhibition

Quantitative analyses of hepatic OATP-mediated interactions between statins and inhibitors using PBPK modeling with a parameter-optimization method.

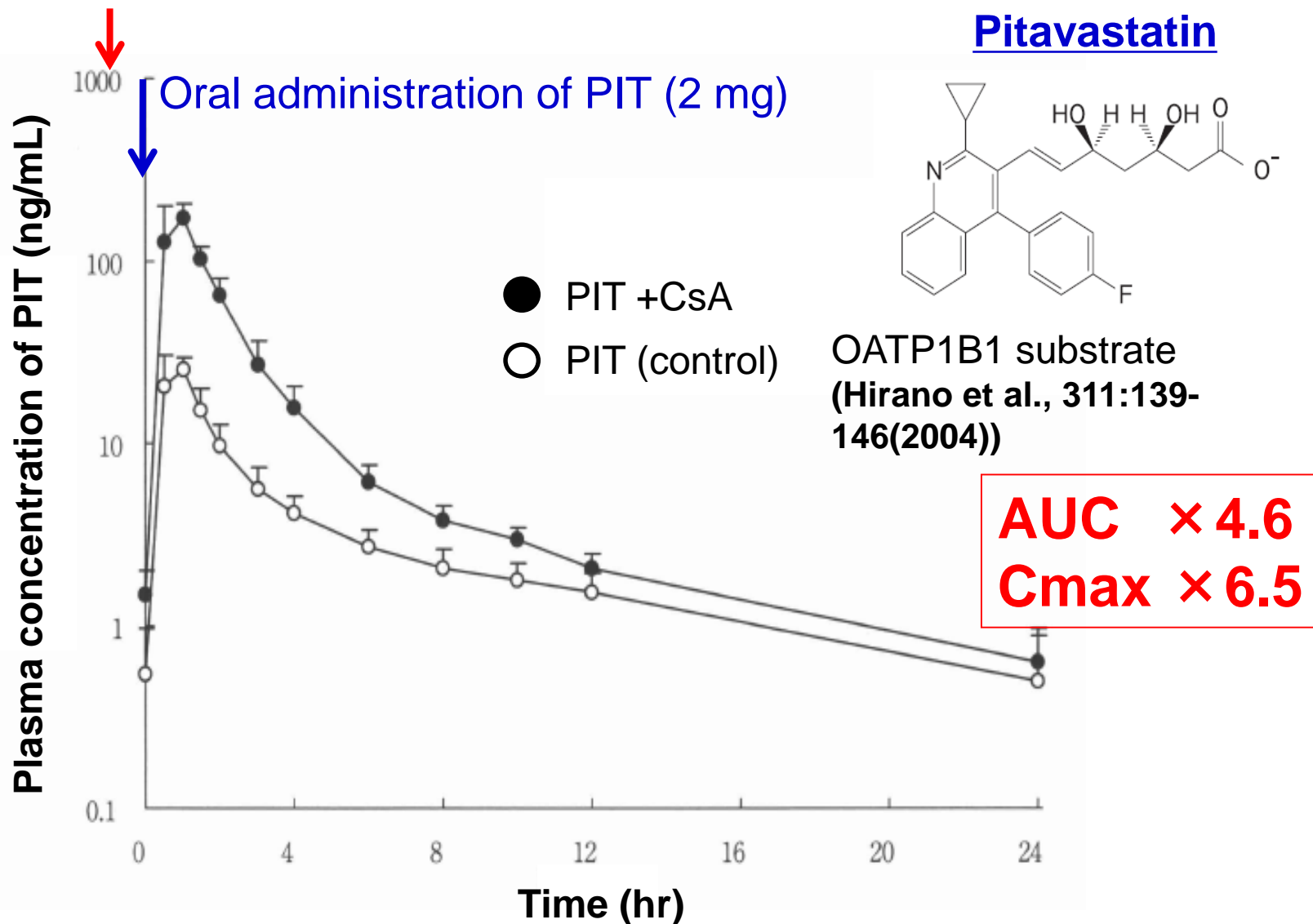
Yoshikado T, Yoshida K, Kotani N, Nakada T, Asaumi R, Toshimoto K, Maeda K, Kusuhashi H, Sugiyama Y.

Clin Pharmacol Ther. 100 (5):513-523 (2016)

Elevation of plasma pitavastatin (PIT) concentration by cyclosporine A (CsA)

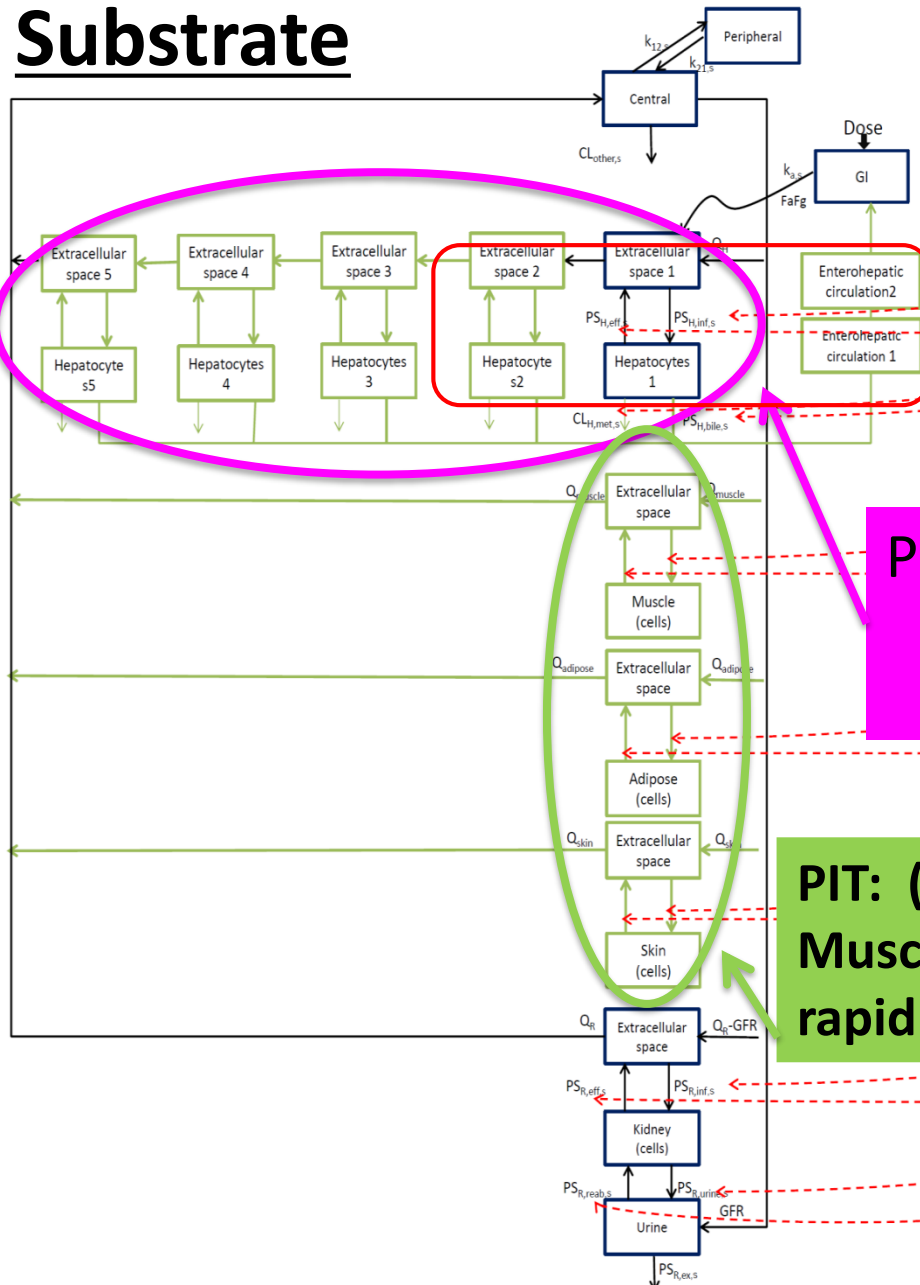
Prediction by dynamic model (PBPK model)

Oral administration of CsA (average 131 mg) 1 hr before PIT

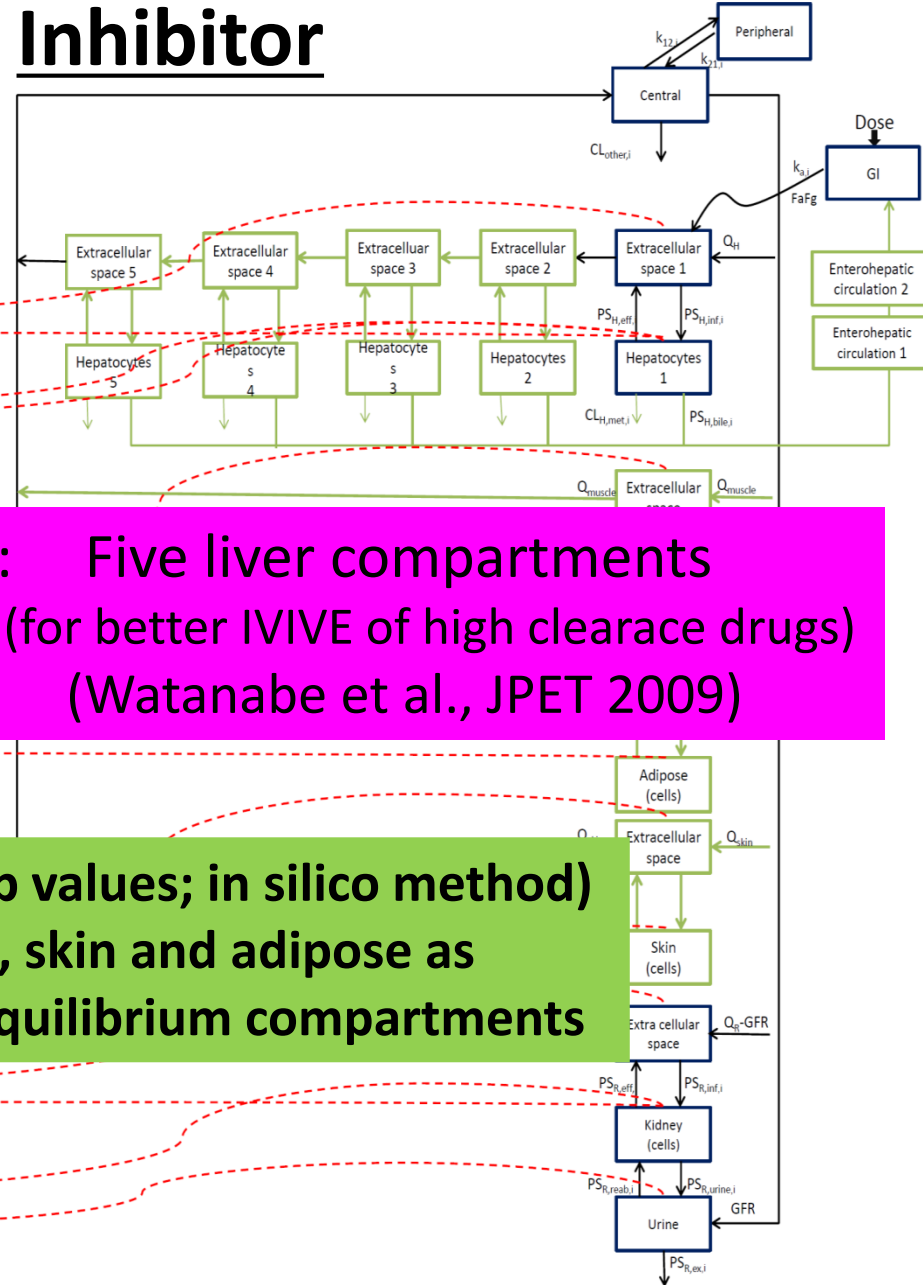


PBPK model for DDI analyses (PIT, CsA)

Substrate

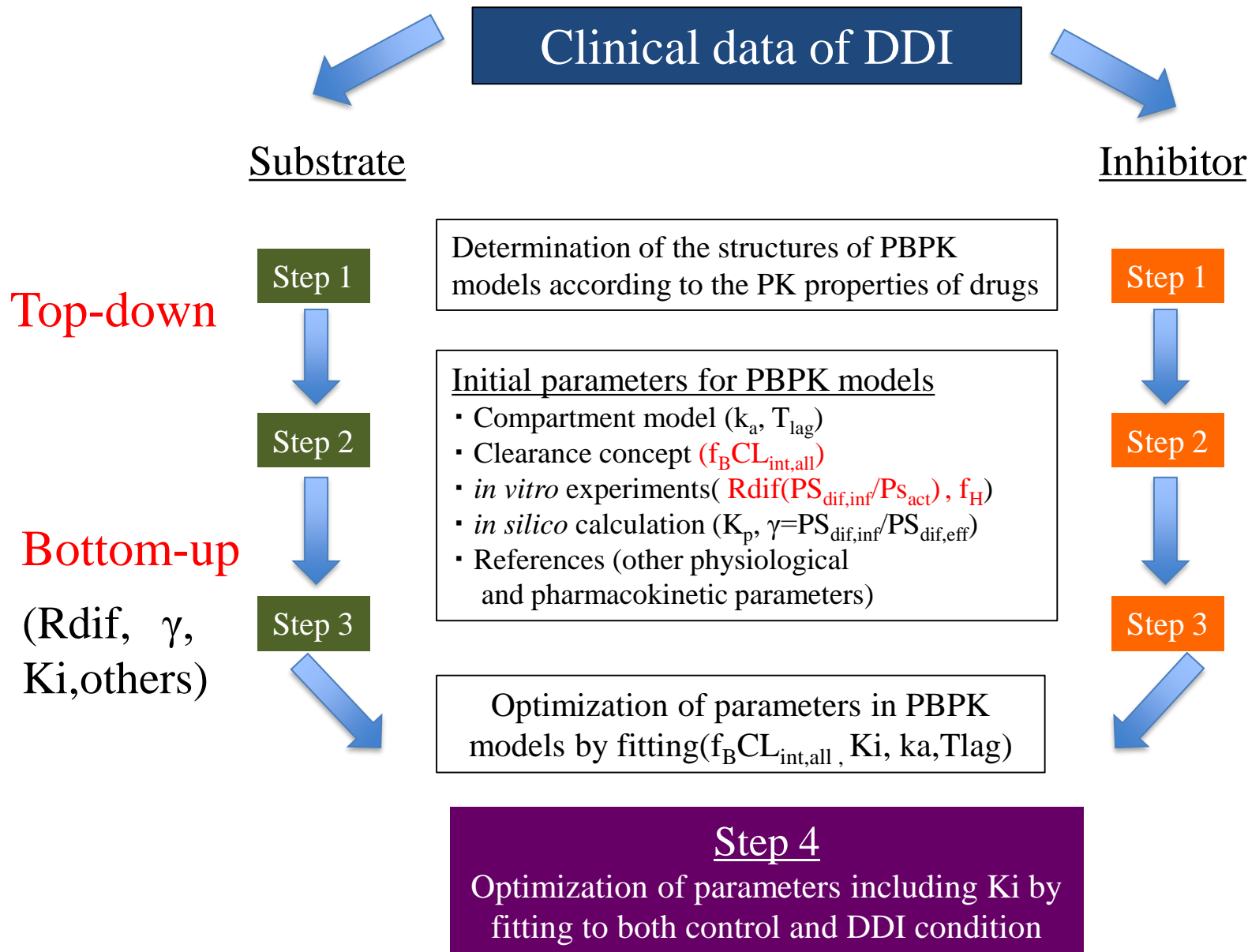


Inhibitor



PIT: Five liver compartments
(for better IVIVE of high clearance drugs)
(Watanabe et al., JPET 2009)

PIT: (K_p values; in silico method)
Muscle, skin and adipose as
rapid equilibrium compartments

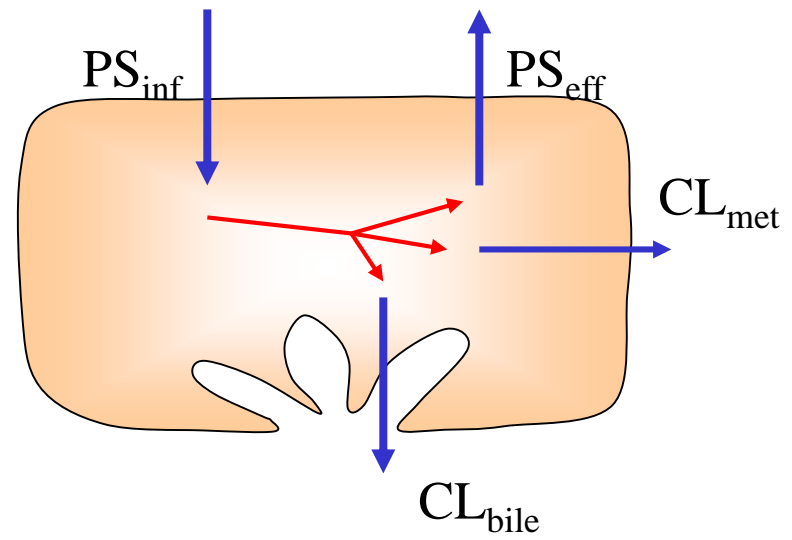


Scheme of the workflow of parameter optimization in the PBPK models to describe hepatic OATP-mediated DDIs.

Extended clearance concept

$$CL_{int,all} = PS_{inf} \times \frac{CL_{bile} + CL_{met}}{PS_{eff} + CL_{bile} + CL_{met}}$$

β value



Where $PS_{inf} = PS_{inf,act} + PS_{inf,dif}$

PS_{eff} assumed to be $PS_{eff,dif}$

$$R_{dif} = PS_{inf,dif} / PS_{inf,act}$$

$$K_{p,uu} = \frac{PS_{inf,act} + PS_{inf,dif}}{PS_{eff,dif}}$$

$$\gamma = PS_{inf,dif} / PS_{eff,dif}, \quad fbile = CL_{bile} / (CL_{met} + CL_{bile})$$

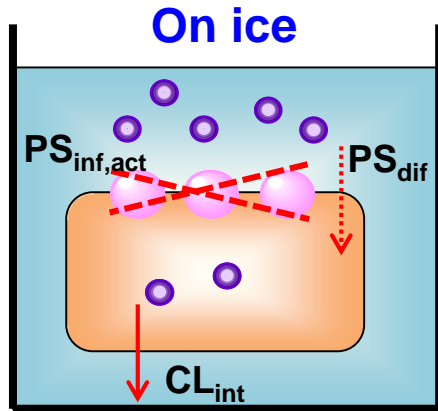
Parameters for elementary steps

$$PS_{inf,act}, PS_{inf,dif}, PS_{eff,dif}, CL_{bile}, CL_{met}$$

Hybrid parameters for describing hepatobiliary elimination steps

$$CL_{int,all}, \beta, K_{p,uu}, R_{dif}, fbile$$

How to determine $K_{p,uu(true)}$ from in vitro experiment?



$$C/M \text{ ratio}(\text{on ice}) = \left[\frac{C_{cell}}{C_{medium}} \right]_{\text{on ice}} = \frac{PS_{dif,inf}}{f_T \cdot PS_{dif,eff}}$$

Assumption: membrane potential is completely lost under on ice condition (recently demonstrated experimentally)

$$PS_{dif,inf} = PS_{dif,eff} \quad (\gamma_{\text{on ice}} = \frac{PS_{dif,inf}}{PS_{dif,eff}} = 1)$$

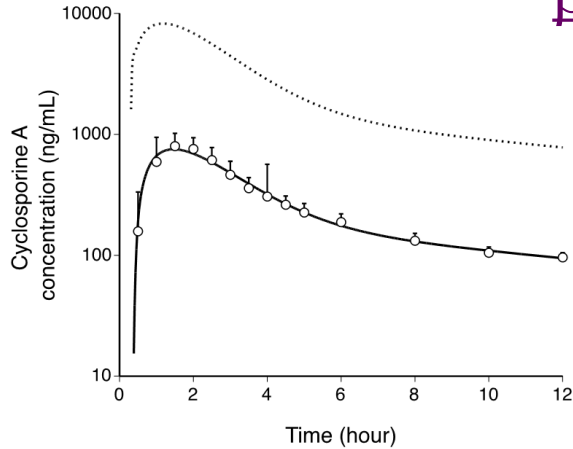
$$\frac{C/M \text{ ratio}(37^\circ\text{C})}{C/M \text{ ratio}(\text{on ice})} = \frac{PS_{act,inf} + PS_{dif,inf}}{f_T \cdot (PS_{dif,eff} + CL_{int,met})} \cdot f_T = K_{p,uu(true)}$$

Comparison of methods for estimating unbound intracellular-to-medium concentration ratios in rat and human hepatocytes using statins

Yoshikado et al., Drug Metab Dispos. 45(7):779-789 (2017).

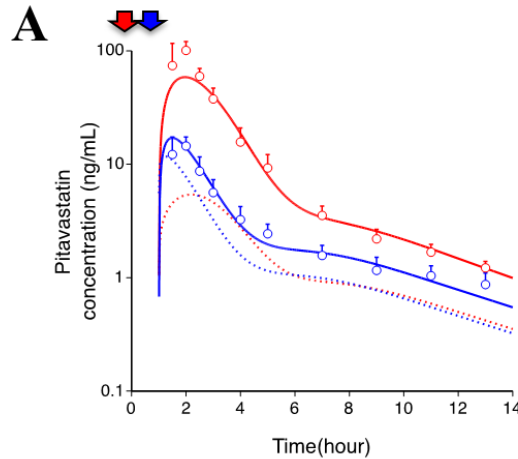
Simulated time course of plasma conc. of cyclosporine A and pitavastatin/fluvastatin

Cyclosporine A

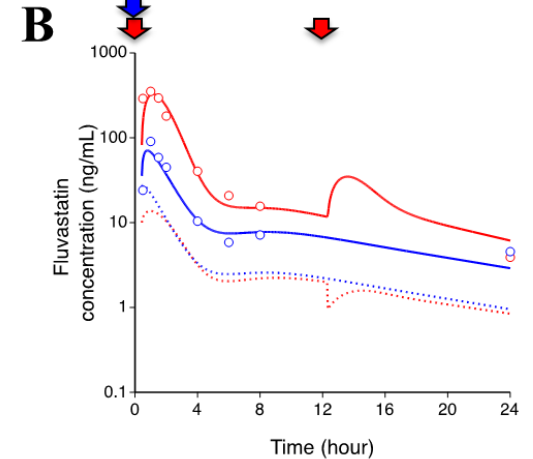


$\beta=0.8$

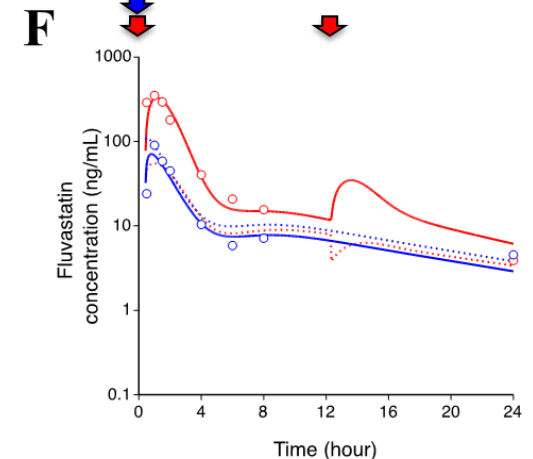
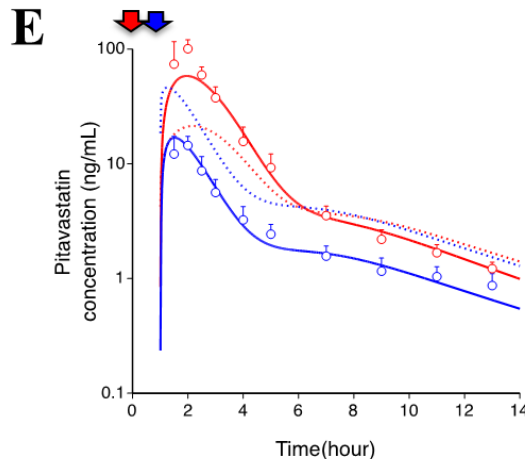
Pitavastatin (EHC model)



Fluvastatin (EHC model)



$\beta=0.2$



Plasma conc..

— control
— +CsA

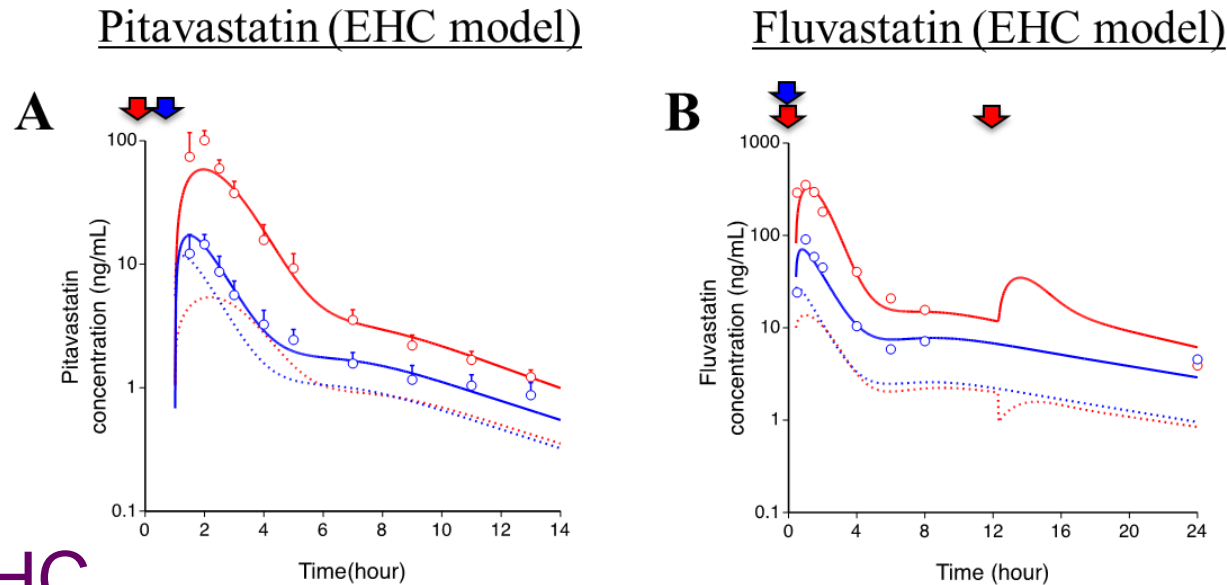
Hepatic conc..

--- control
--- +CsA

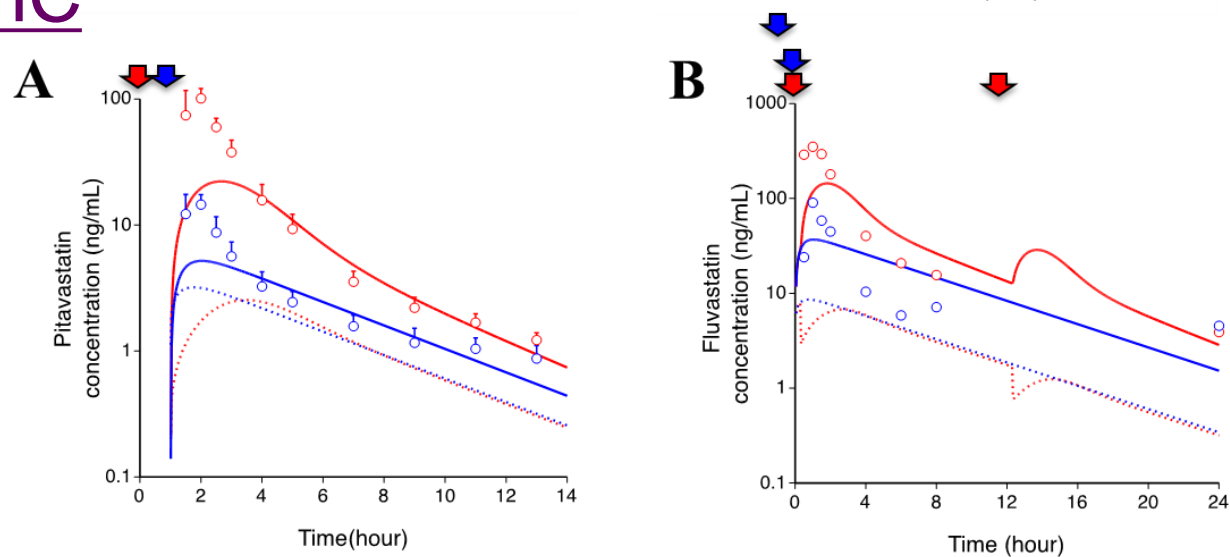
Impact of the consideration of enterohepatic circulation on the simulated time course of plasma conc. of pitavastatin/fluvastatin

(EHC model gave much better fit to the data)

EHC



Non-EHC



- PBPK models were constructed for pitavastatin, fluvastatin and pravastatin as substrates and cyclosporin A (CsA) and rifampicin (RIF) as inhibitors, where enterohepatic circulations (EHC) of statins were incorporated. Without EHC, good fitting was not obtained for either substrate.

(In vitro measured R_{dif} values, γ values were well incorporated into this modeling)

- Similar *in vivo* inhibition constant (K_i) values of each inhibitor against OATPs were obtained, regardless of the substrates.

CsA: 0.012 μ M (pitavastatin) 0.010 μ M (fluvastatin)

Rifampicin; 0.23 μ M (pitavastatin) 0.19 μ M (pravastatin)

- Estimated K_i values of CsA were comparable to reported *in vitro* values with the preincubation of CsA, while those of RIF were 3-5 folds smaller than reported *in vitro* values .

CsA(+preincubate) ; 0.014–0.080 μ M, Trans-inhibition mechanism

Rifampicin; 0.65–1.1 μ M,

(Mechanism of lower K_i value in vivo of Rif. compared with in vitro are not known yet.)

Conclusion

- * Standardized protocol of top-down analyses of complex DDI (where transporters and enzymes are inhibited) based on PBPK modeling were established. The *in vivo* K_i values were obtained, leading to the prediction of complex DDI of other substrates
- * We need some *in vitro* measured parameters such as R_{dif} value, K_{puu} value, γ values (index of the asymmetry of passive diffusion via basolateral membrane) in this model based analyses
- * An important parameter, β which determines rate-determining process of drugs is set to different values (0.2, 0.5, 0.8) in the model analyses, and the outcome of analyses are not so much different as far as the plasma-concentration time profiles are analyzed.
- * However, this β values should affect the hepatic concentration time profiles. Therefore, this should be estimated in near future from the *in vitro* experiments
(isolated and sandwich cultured hepatocytes, HLMs)

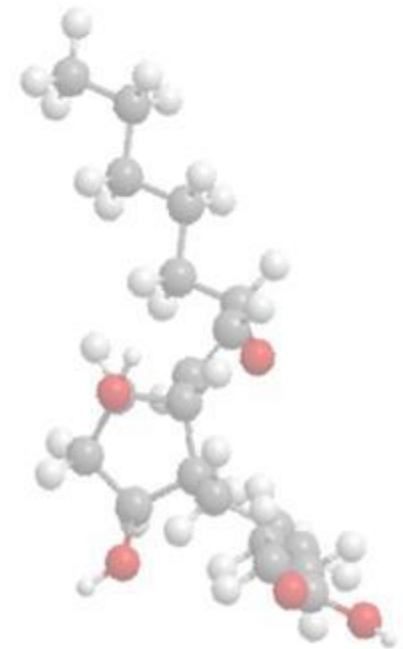
Contents

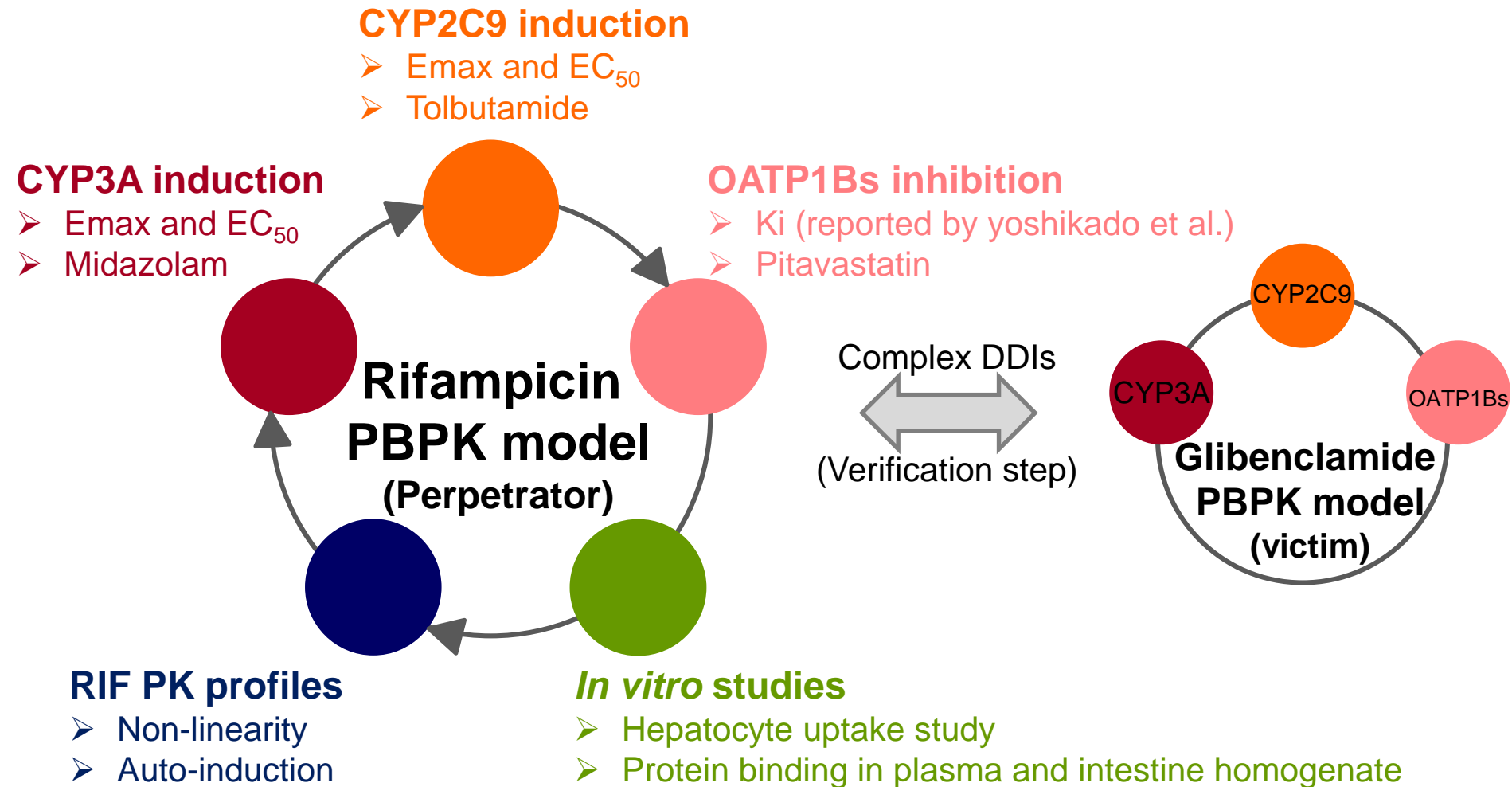
- 1) Introduction;
Rate-determining process (focusing on the liver)
(Uptake, efflux, elimination, metabolism)
- 2) PBPK model based analysis of OATPs mediated drug-drug interaction (Top down + Bottom-up)
- 3) Other collaborations
Integrated model of rifampicin as a perpetrator
Time-dependent inhibition
Substrate-dependent inhibition

PBPK model of rifampicin i
hepatic uptake and auto-in
various types of DDIs such
and/or OATPs inhibitions

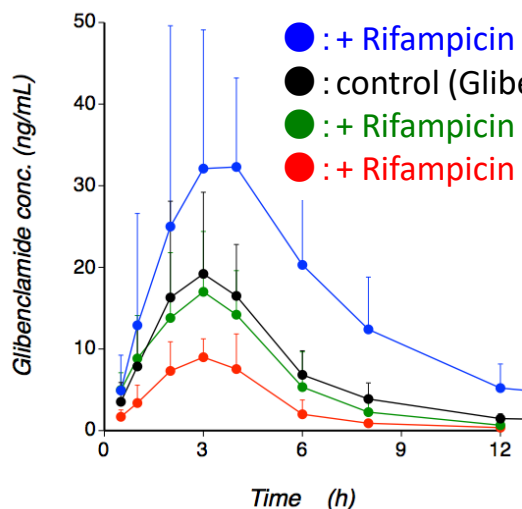
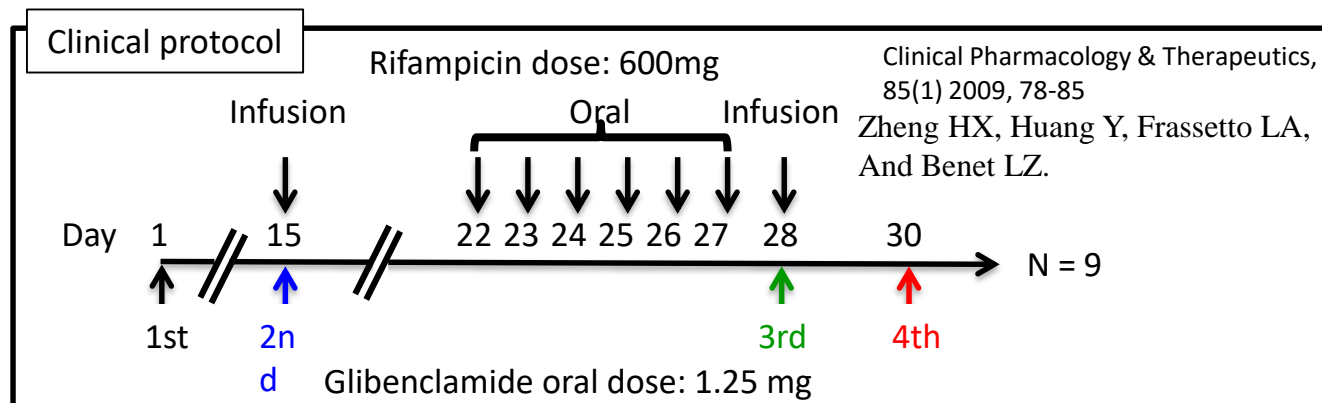
Asaumi R, Toshimoto K, Tobe Y, Hashizume K,
Comprehensive PBPK Model of Rifampicin for
Drug-Drug Interactions: CYP3A/2C9 Induction
CPT Pharmacometrics Syst Pharmacol. (2018) 7

CPT Pharmacometrics Syst Pharmacol. (2018) 7, 186–196

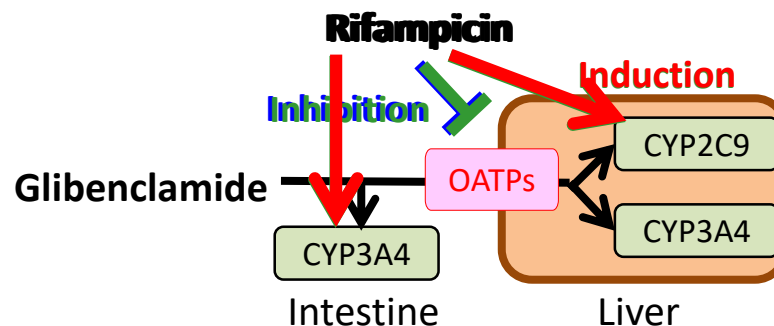




Complex rifampicin DDI effects on glibenclamide pharmacokinetics



Collaboration with Asaumi (Ono Pharm)



Further studies recently done on a collaboration base on the OATP-mediated DDIs

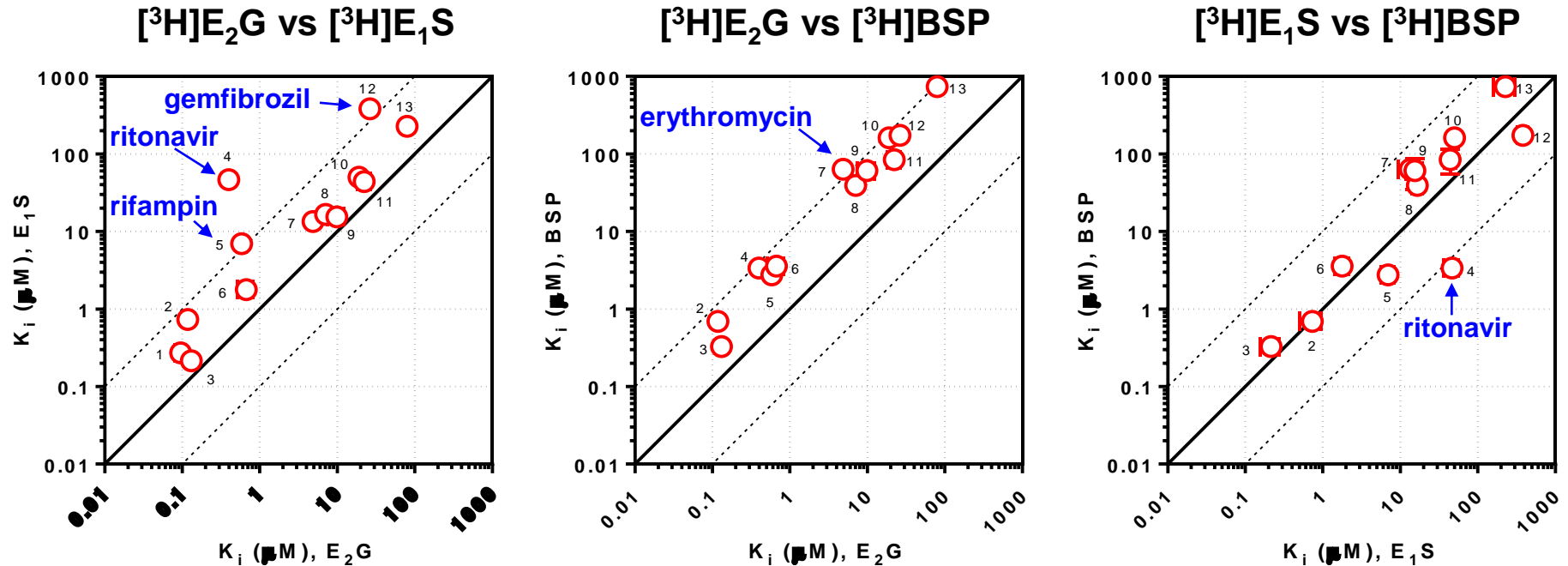
A) Time-dependent inhibition of OATPs by cyclosporine A

Shitara Y et al., (Sanofi) *Pharmacol Therap.* S0163-7258: 30066-9 (2017)

B) Substrate-dependent inhibition. Collaboration with Eisai

- 1) **Izumi S, Nozaki Y**, Komori T, Maeda K, Takenaka O, Kusano K, Yoshimura T, Kusuhara H and Sugiyama Y. Substrate-dependent inhibition of organic anion transporting polypeptide 1B1: comparative analysis with prototypical probe substrates estradiol-17beta-glucuronide, estrone-3-sulfate, and sulfobromophthalein. *Drug Metab Dispos* 41:1859-1866 (2013).
- 2) **Izumi S, Nozaki Y**, Maeda K, Komori T, Takenaka O, Kusuhara H and Sugiyama Y. Investigation of the Impact of Substrate Selection on In Vitro Organic Anion Transporting Polypeptide 1B1 Inhibition Profiles for the Prediction of Drug-Drug Interactions. *Drug Metab Dispos* 43:235-247 (2015).
- 3) **Izumi S, Nozaki Y**, Komori T, Takenaka O, Maeda K, Kusuhara H, Sugiyama Y. Investigation of Fluorescein Derivatives as Substrates of Organic Anion Transporting Polypeptide (OATP) 1B1 To Develop Sensitive Fluorescence-Based OATP1B1 Inhibition Assays. *Mol Pharm.*13: 438-48 (2016)
- 4) **Izumi S, Nozaki Y**, Komori T, Takenaka O, Maeda K, Kusuhara H, and Sugiyama Y. Comparison of the Predictability of Human Hepatic Clearance for Organic Anion Transporting Polypeptide (OATP) Substrate Drugs Between Different In Vitro-In Vivo Extrapolation Approaches *J.Pharm.Sci., in press*

Comparison of K_i values for OATP1B1 between prototypical probe substrates

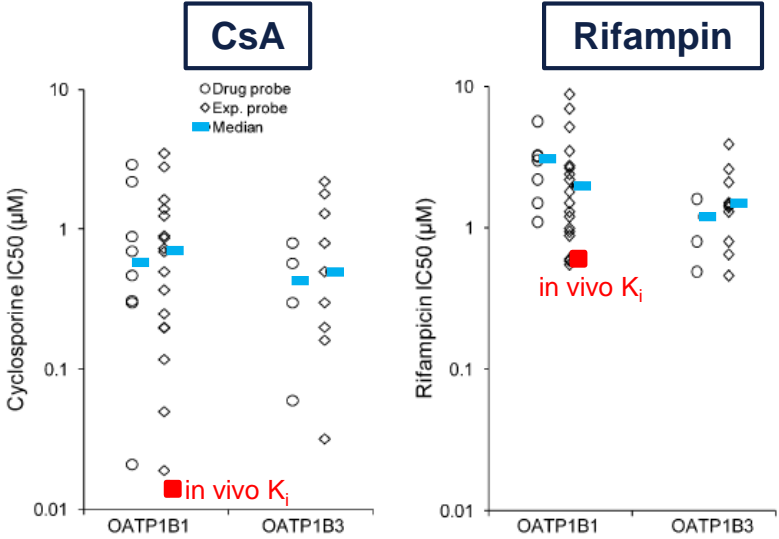


1, E₁S; 2, CsA; 3, BSP; 4, Ritonavir; 5, Rifampin; 6, Tacrolimus; 7, Erythromycin; 8, E₂G; 9, Ketoconazole; 10, TCA; 11, Verapamil; 12, Gemfibrozil; 13, Probenecid.

Izumi et al. (2013) *Drug Metab Dispos* 41: 1859-1866

- Some inhibitors (eg, ritonavir, gemfibrozil, and rifampin) showed >10-fold (117-fold for ritonavir) variations in the K_i values, depending on the substrates used.
- Of the 3 substrates, [³H]E₂G provided the lowest K_i values for all inhibitors examined.
- [³H]E₂G can be used as a sensitive probe substrate of OATP1B1, which could offer conservative K_i values and help mitigate the risk of false-negative DDI prediction.

Comparison of IC₅₀ (or K_i) values of CsA and rifampicin obtained in vitro and in vivo (Collaboration with Izumi, Nozaki et al)



IC₅₀ values of CsA, rifampin for OATP1B1 and OATP1B3

In vitro IC₅₀ values were higher than in vivo K_i values obtained from PBPK model, particularly for CsA.

Li et al. (2014) Clin Pharmacokinet 53: 659-678

Transporters	Substrates	Pre-incubation time (min)	K _i or IC ₅₀ of CsA		Ratio
			Without CsA	With CsA	
			(µM)	(µM)	
OATP1B1 ^a	Atorvastatin	60	0.47 ± 0.34	0.021 ± 0.004	22.4
	[³ H]E ₂ G		0.0458 ± 0.0041	0.0139 ± 0.0066	3.3
	[³ H]E ₁ S		0.134 ± 0.017	0.0264 ± 0.0085	5.1
OATP1B1 ^b	[³ H]BSP	60	0.252 ± 0.057	0.0799 ± 0.0273	3.2
	Pitavastatin		0.0985 ± 0.0360	0.0252 ± 0.0038	3.9
	Atorvastatin		0.0986 ± 0.0250	0.0229 ± 0.0033	4.3
OATP1B1 ^c	[³ H]E ₂ G	30	0.198 ± 0.069	0.019 ± 0.007	10.4
OATP1B3 ^c	[³ H]E ₂ G		0.162 ± 0.062	0.032 ± 0.003	2.6

In vitro IC₅₀ values of CsA for OATP1B1 and 1B3 obtained w/ or w/o CsA pre-incubation.

K_i (or IC₅₀) = 0.01 ~ 0.08 µM

- a) Amundsen et al. (2010) DMD 38: 1499-1504
- b) Izumi et al. (2015) DMD 43: 235-247
- c) Gertz et al. (2913) Pharm Res 30: 761-780

Experimentally obtained K_i value of CsA after pre-incubation with CsA was comparable to the in vivo K_i value.

Dialogue and Debate session

"The unfolded story of Long-Lasting OATP Transporter Inhibition"



Aleksandra Galetin, PhD,
Centre for Applied Pharmacokinetic Research,
University of Manchester, UK

Title: Translation of prolonged OATP1B1 inhibition *in vitro*
to clinical DDI risk assessment



Yuichi Sugiyama, PhD,
Head of Sugiyama Laboratory
RIKEN, Yokohama, Japan

Title: Preincubation Time-Dependent and Long-Lasting
Inhibition of Organic Anion Transporting Polypeptides
(OATPs)

Moderators

Wei Yue, Ph.D.
Assistant Professor of
Pharmaceutical Sciences
The University of Oklahoma
Health Sciences Center



Dan Bow, Ph.D.
Senior Principal Research
Scientist at AbbVie



Preincubation-dependent and long-lasting inhibition of organic anion transporting polypeptide (OATP) and its impact on drug-drug interactions



Yoshihisa Shitara^a, Yuichi Sugiyama^{b,*}

^a Pharmacokinetics, Dynamics and Metabolism, Translational Medicine and Early Development, R&D Operations, Sanofi, Tokyo, Japan

^b Sugiyama Laboratory, RIKEN Innovation Center, RIKEN, Yokohama, Japan

ARTICLE INFO

Available online 27 February 2017

Keywords:

OATP1B1

Drug-drug interactions

Hepatic uptake

Time-dependent inhibition

Long-lasting inhibition

Physiologically based pharmacokinetic model

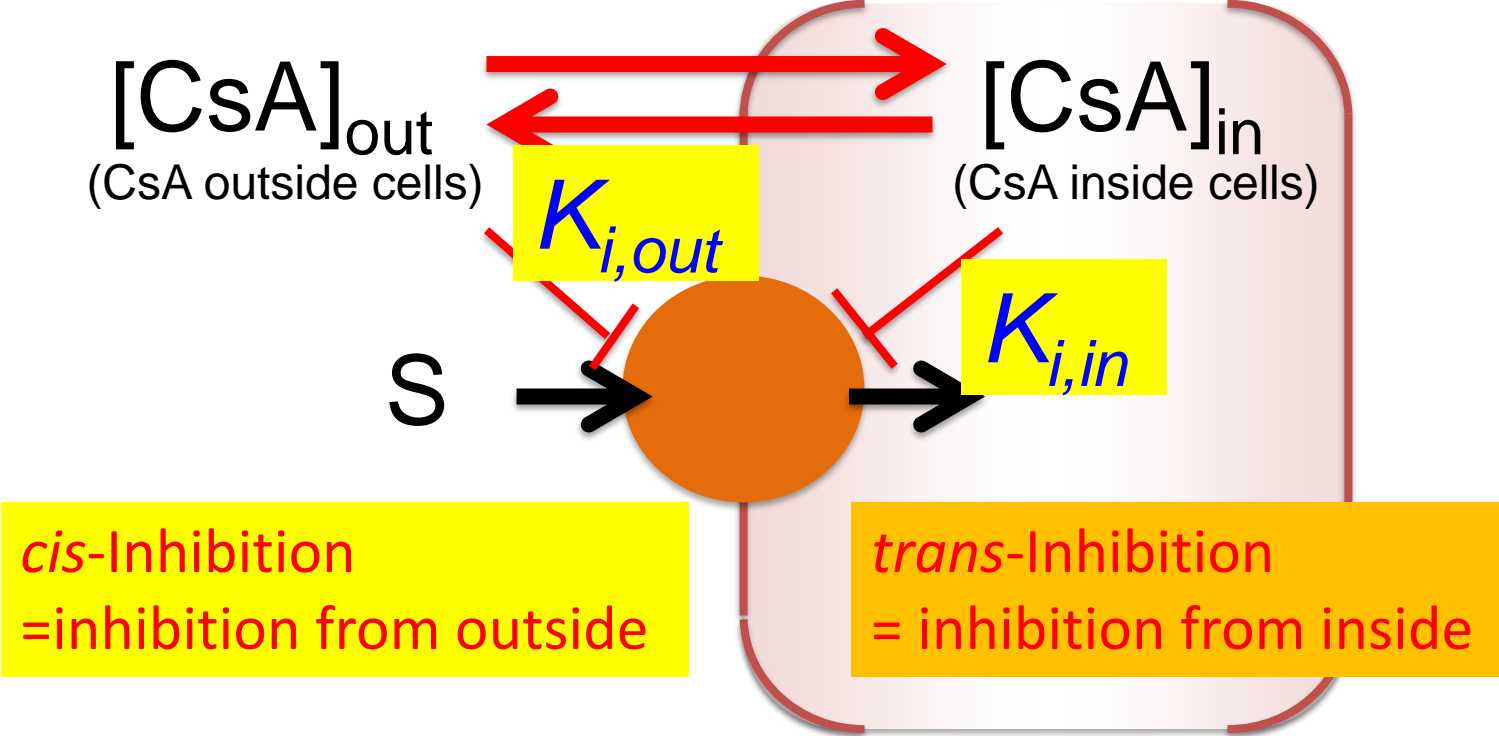
Modeling & simulation

ABSTRACT

Preincubation with cyclosporin A (CsA), a potent inhibitor of organic anion transporting polypeptide 1B1 (OATP1B1) and OATP1B3, enhanced its inhibitory effects on these transporters *in vitro*. A similar effect was observed upon preincubation with some other inhibitors. Removing these from the incubation media did not readily reverse the inhibition on OATP1B1 and OATP1B3. This preincubation-dependent long-lasting inhibition appeared to be related to CsA concentration in the cells in addition to that in the incubation media. Thus, we hypothesized that CsA inhibits OATP1B1 and OATP1B3 from inside (*trans*-inhibition) as well as outside (*cis*-inhibition) the cells and constructed the *cis*- and *trans*-inhibition model. The enhanced inhibitory effect of CsA on OATP1B1 observed after preincubation was quantitatively described using $K_{i,OUT}$ and $K_{i,IN}$ as inhibition constants for *cis*- and *trans*-inhibitions, respectively. In addition, a long-lasting inhibition was also described by this model. Additional factors taken into consideration when simulating *in vivo* pharmacokinetic alterations by CsA are potential inhibition by AM1, a major metabolite of CsA, which has been reported to inhibit OATP1B1 and OATP1B3. Based on the physiologically based pharmacokinetic model incorporating *trans*- and *cis*-inhibition of OATP1B1 by CsA, the simulation showed that OATP1B1-mediated drug-drug interaction with CsA was suggested to be time-dependent also *in vivo* although further clinical studies are required for confirmation.

© 2017 Elsevier Inc. All rights reserved.

Cis- and trans-inhibition of CsA on OATP1B1



Uptake clearance of OATP1B1 substrates in the liver when co-administered with CsA:

$$CL_{\text{uptake}}(+I) = \frac{V_{\text{max}}}{K_m \cdot (1 + f_{u,B} \cdot I_{EH} / K_{i,out}) + S}$$

$(1 + f_{u,H} \cdot I_H / K_{i,in})$

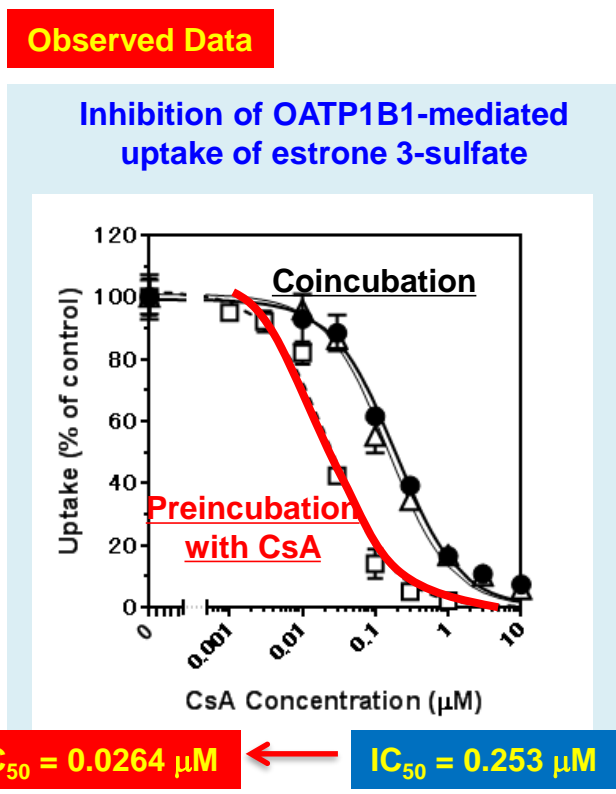
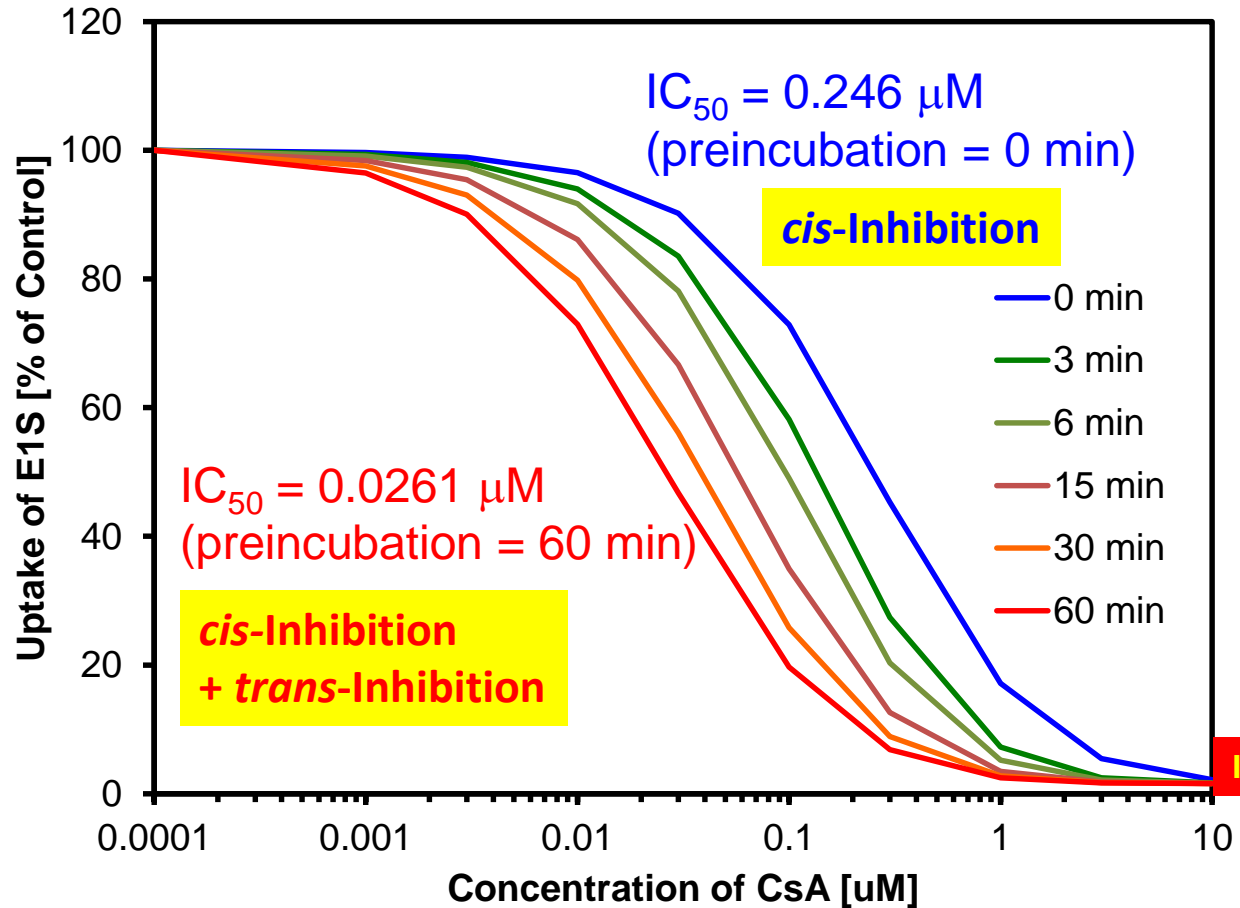
trans-Inhibition (non-competitive)

$(1 + f_{u,B} \cdot I_{EH} / K_{i,out})$

cis-Inhibition (competitive)

* I_H , I_{EH} : inhibitor (CsA) concentrations in the liver and at the extracellular space of the liver, respectively

Simulation analysis of time-dependent enhancement effect of inhibition of OATP1B1 by CsA



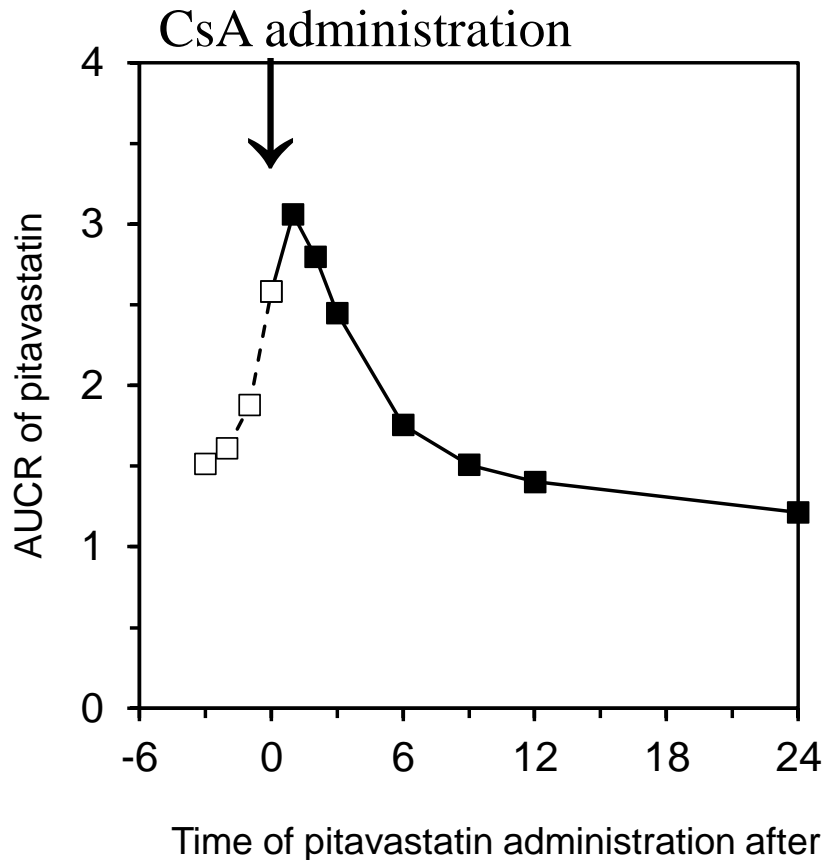
Izumi S et al. (2015)
Drug Metab Dispos 43, 235

$K_{i,out}$: 0.296 (μ M), $K_{i,in}$: 0.0119 (μ M): fitted values

$$1/(1 + I_{out}/IC_{50,app}) = 1/[(1 + I_{out}/K_{i,out}) \times (1 + I_{u,in,unbound}/K_{i,in})]$$

:with respect to time $I_{u,in} \uparrow, K_{i,in} \rightarrow \Rightarrow IC_{50,app} \downarrow$

When pitavastatin is administered one hour after the CsA administration, AUCR of pitavastatin exhibited the maximum value



Estimation of pitavastatin AUCR by coadministration with CsA for different time interval between two drugs.

AUCR values of pitavastatin by coadministration with CsA with several intervals were estimated from the PBPK modeling considering trans-inhibition mechanism.

What is the best translational approach until the mechanism is fully elucidated?

- 1) Benefit in using 'shifted' IC50 in the PBPK models as conservative approach?
- 2) More mechanism based PBPK modeling; transfer of inhibitor to intracellular sites \Rightarrow estimation of on/off rate of inhibitors \Rightarrow estimation of intracellular kinetics of inhibitors

COI

I am a scientific advisory board member of the following companies.

- 1) SimCYP
- 2) SEKISUI Medical

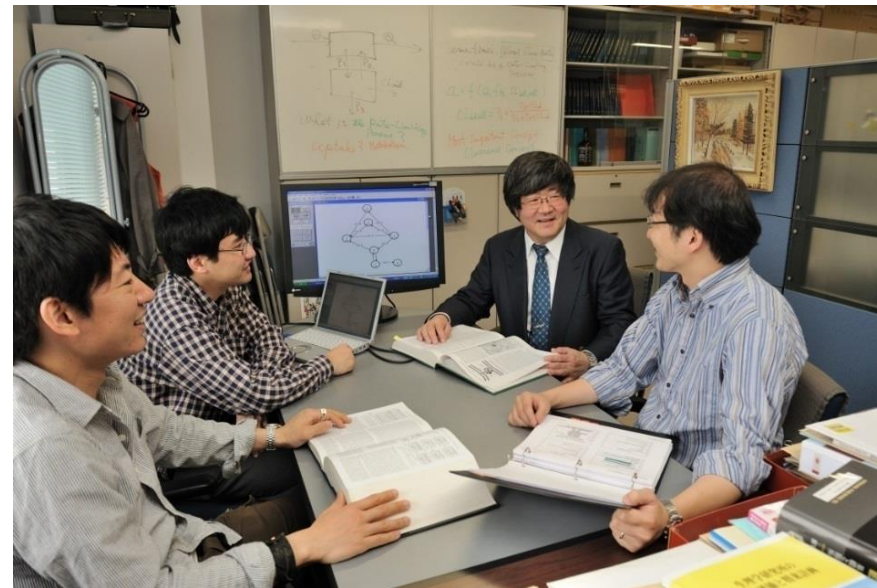
I am and have been a consultant of and collaborating with 30 domestic and global pharmaceutical industries



RIKEN Acknowledgement

Sugiyama Lab Main Members

Takashi Yoshikado (wet +dry, IVIVE DDI)
 Kota Toshimoto (dry, VCT (Virtual clinical study))
 Kim, Soo-Jin: (DDI, PGx, IVIVE)
 Lee, Kyeong Ryoong (Daewoong) (DDI, IVIVE)
 Kim, Sae Hee (Daewoong) (DDI, IVIVE)
 Atsuko Tomaru (IVIVE, Bioanalysis with LC/MS/MS)
 Aya Kiriake (IVIVE, transporter expression systems)
 Kiyoe Morita (IVIVE, isolated hepatocyte (suspension, plated))
 In-Soo Yoon (DDI, isolated hepatocytes (suspension, plated))



Kazuya Maeda, Hiroyuki Kusuvara (Univ of Tokyo)

Azusa Futatsugi (DDI, PGx)
 Keiko Ishigame (IVIVE, Bioanalysis)
 Hisako Motokawa (IVIVE, Bioanalysis)

Secretaries

Sachie Sato
 Chinami Yamada
 Rie Abe
 Aya Miyamoto

Acknowledgement

The members of PKPD seminar

PKPD seminar held once a month (18:30 – 22:00)

Headed by Yuichi Sugiyama

