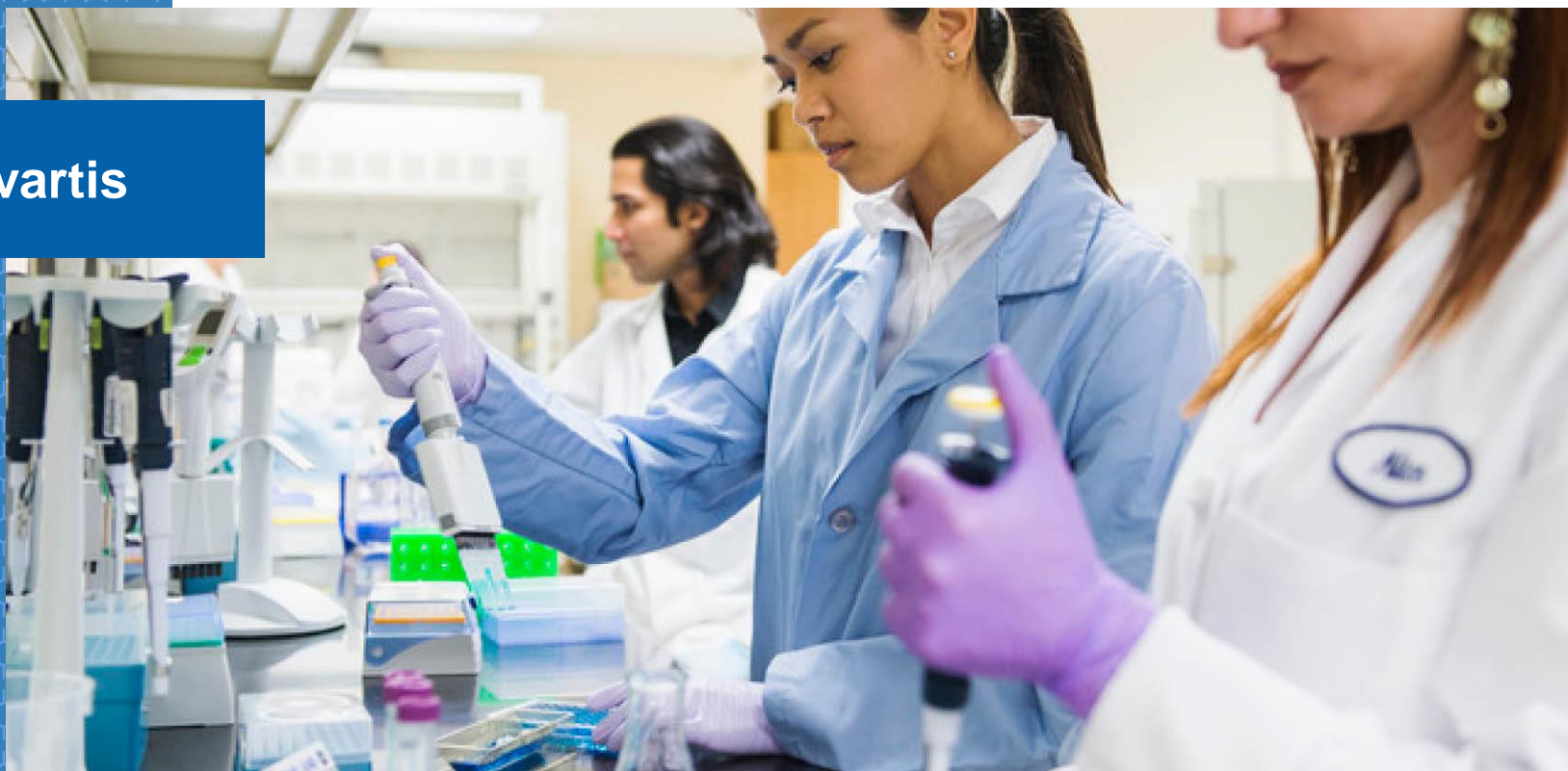


Novartis



Drug disposition classification systems: A comparative review of BDDCS, ECCS and ECCCS

Birk Poller, Gian Camenisch - Novartis

SOLVO - Meet The Experts Transporter Conference

April 26, 2018



Drug disposition classification systems

BCS

	High Solubility	Low Solubility
High Permeability	Class 1 Metabolism	Class 2 Metabolism
Low Permeability	Class 3 Renal and/or Biliary Elimination of Unchanged Drug	Class 4 Renal and/or Biliary Elimination of Unchanged Drug

Amidon et al, 1995, *Pharm Res*;12:413-20

BDDCS

	High Solubility	Low Solubility
Extensive Metabolism	Class 1 High Solubility Extensive Metabolism (Rapid Dissolution and $\geq 70\%$ Metabolism for Biowaiver)	Class 2 Low Solubility Extensive Metabolism
Poor Metabolism	Class 3 High Solubility Poor Metabolism	Class 4 Low Solubility Poor Metabolism

Wu and Benet, 2005, *Pharm Res*;22:11-23

ECCCS

ECC class 1	ECC class 2
<p>ECCCS: $PS_{app} \leq CL_{int}$ BDDCS: Solubility \times Dose/250 mL</p> <p>ECCCS: $PS_{app} > 3 \times CL_{int}$ BDDCS: Metabolism $> 70\%$</p> <p>Hepatic elimination primary via metabolism</p> <p>Transporter effects minimal</p>	<p>ECCCS: $PS_{app} \gg CL_{int}$ BDDCS: Solubility $<$ Dose/250 mL</p> <p>Hepatic elimination via metabolism and possibly biliary secretion of unchanged drug</p> <p>Efflux transporter effects can occur</p>
<p>ECCCS: $PS_{app} < 3 \times CL_{int}$ BDDCS: Metabolism $< 30\%$</p> <p>Renal and/or biliary elimination of unchanged drug</p> <p>Uptake transporter effects predominant</p>	<p>Biliary and/or renal elimination of unchanged drug</p> <p>Uptake and efflux transporter effects may be important</p>
ECC class 3	ECC class 4

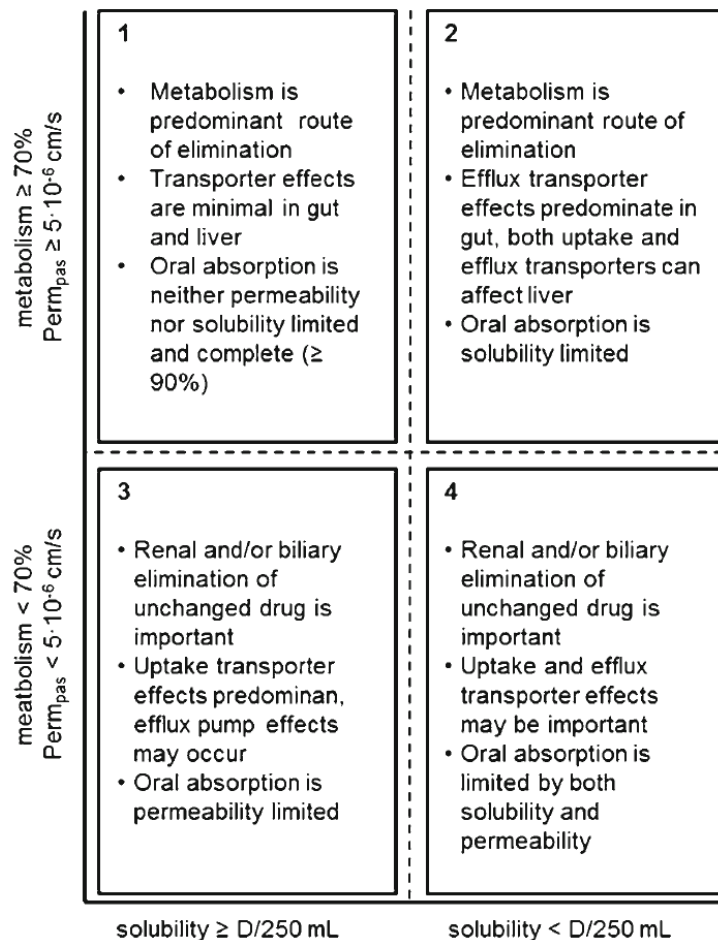
Camenisch et al, 2015, *ADMET&DMPK*;1:1-14
Camenisch, 2016, *Pharm Res*;33:2583-93

ECCS

High Permeable	<p>Class 1</p> <p>MW ≤ 400 Class 1A Metabolism</p> <p>MW > 400 Class 1B Hepatic uptake</p>	Class 2 Metabolism
Low Permeable	<p>Class 3</p> <p>MW ≤ 400 Class 3A Renal</p> <p>MW > 400 Class 3B Hepatic uptake (or) Renal</p>	Class 4 Renal
	<u>Acids/Zwitter</u>	<u>Bases/Neutrals</u>

Varma et al, 2015, *Pharm Res*;32:3785-802

Biopharmaceutics Drug Disposition Classification System (**BDDCS**)



Classification based on **human *in vivo* metabolism (or passive permeability)** and **soluble dose**

→ *Rather applicable in late drug development phases*

→ *Provides information about involvement of potential transport processes in absorption and elimination*

→ *Observation based classification system*

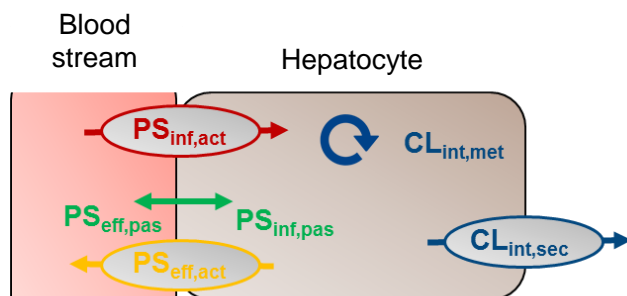
Extended Clearance Classification System (ECCS)

$\text{Perm}_{\text{pas}} \geq 5 \cdot 10^{-6} \text{ cm/s}$	1A <ul style="list-style-type: none"> • Clearance determined by metabolism • Eliminated as metabolites ($\geq 70\%$) • Absorption is not permeability-limited 	1B <ul style="list-style-type: none"> • Clearance determined by (active) hepatic uptake • Eliminated as metabolites ($\geq 70\%$) • Absorption is not permeability-limited 	2 <ul style="list-style-type: none"> • Clearance determined by metabolism • Eliminated as metabolites ($\geq 70\%$) • Absorption is not permeability-limited
	3A <ul style="list-style-type: none"> • Clearance determined by renal • Eliminated as parent in urine ($\geq 70\%$) • Permeability-limited absorption 	3B <ul style="list-style-type: none"> • Clearance determined by hepatic uptake or renal • Eliminated as parent in bile or urine ($\geq 70\%$) • Permeability-limited absorption 	4 <ul style="list-style-type: none"> • Clearance determined by renal • Eliminated as parent in urine ($\geq 70\%$) • Absorption is permeability-limited
	MW ≤ 400 acids/zwitter	MW > 400	bases/neutrals

Classification based on *in vitro* permeability and physicochemical properties (MW, charge)

- Applicable in early drug development phases
- Allows to identify the rate-limiting clearance processes (absorption, distribution and elimination model)
- Observation based classification system (based on the extended clearance concept)

Extended Clearance Concept Classification System (**EC3CS** = **EC3S**)



$$CL_{h,int} = \frac{PS_{inf,act} + PS_{inf,pas}}{PS_{eff,act} + PS_{eff,pas} + CL_{int}} \times CL_{int}$$

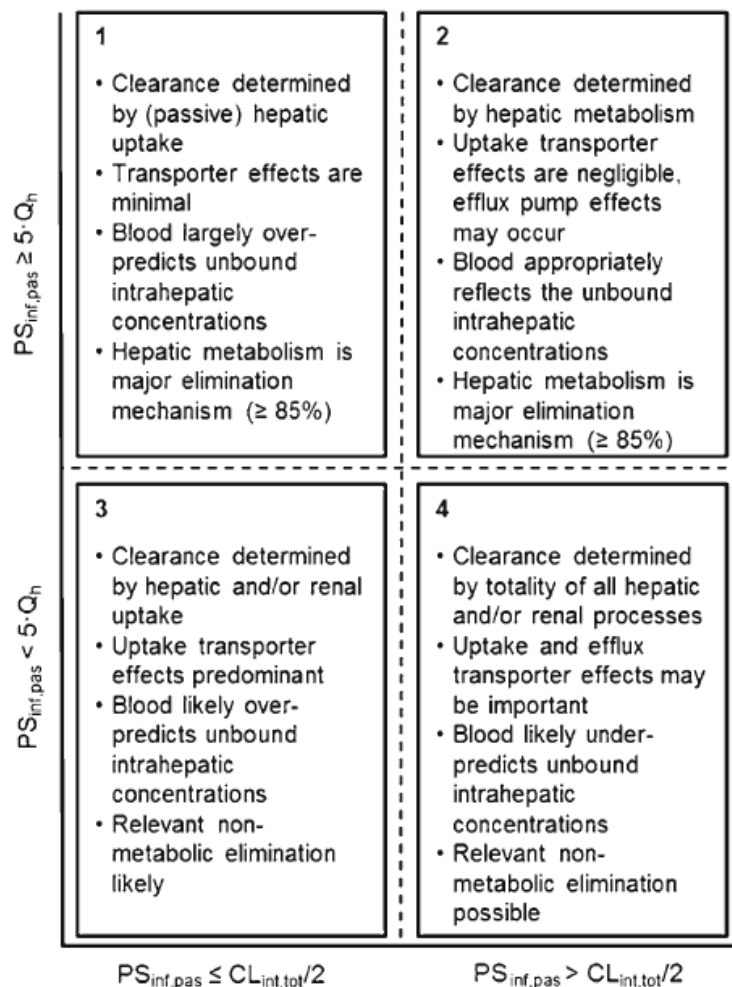
	$PS_{inf,pas} < CL_{int}$	$PS_{inf,pas} \geq CL_{int}$
$PS_{inf,pas} \geq 3-5 Q_h$	EC3S class 1 $PS_{inf,pas}$	EC3S class 2 CL_{int}
$PS_{inf,pas} < 3-5 Q_h$	PS_{inf} EC3S class 3	$\frac{PS_{inf} \times CL_{int}}{PS_{inf,pas}}$ EC3S class 4

in vitro input parameters

$PS_{inf,pas}$	Hepatic uptake / MDCK permeability
$PS_{inf,act}$	Hepatic uptake
$CL_{int,met}$	Liver microsomes / Hepatocytes / S9
$CL_{int,sec}$	Sandwich-cultured hepatocytes
$PS_{eff,act}$	= $PS_{inf,pas}$

Camenisch, 2016, *Pharm Res*;33:2583-93;
 Shitara et al, 2005, *Annu Rev Pharmacol Toxicol*;45:689-723
 Sirianni and Pang, 1997, *J Pharmacokinet Biopharm*;25:449-70

Extended Clearance Concept Classification System (**ECCCS** = **EC3S**)



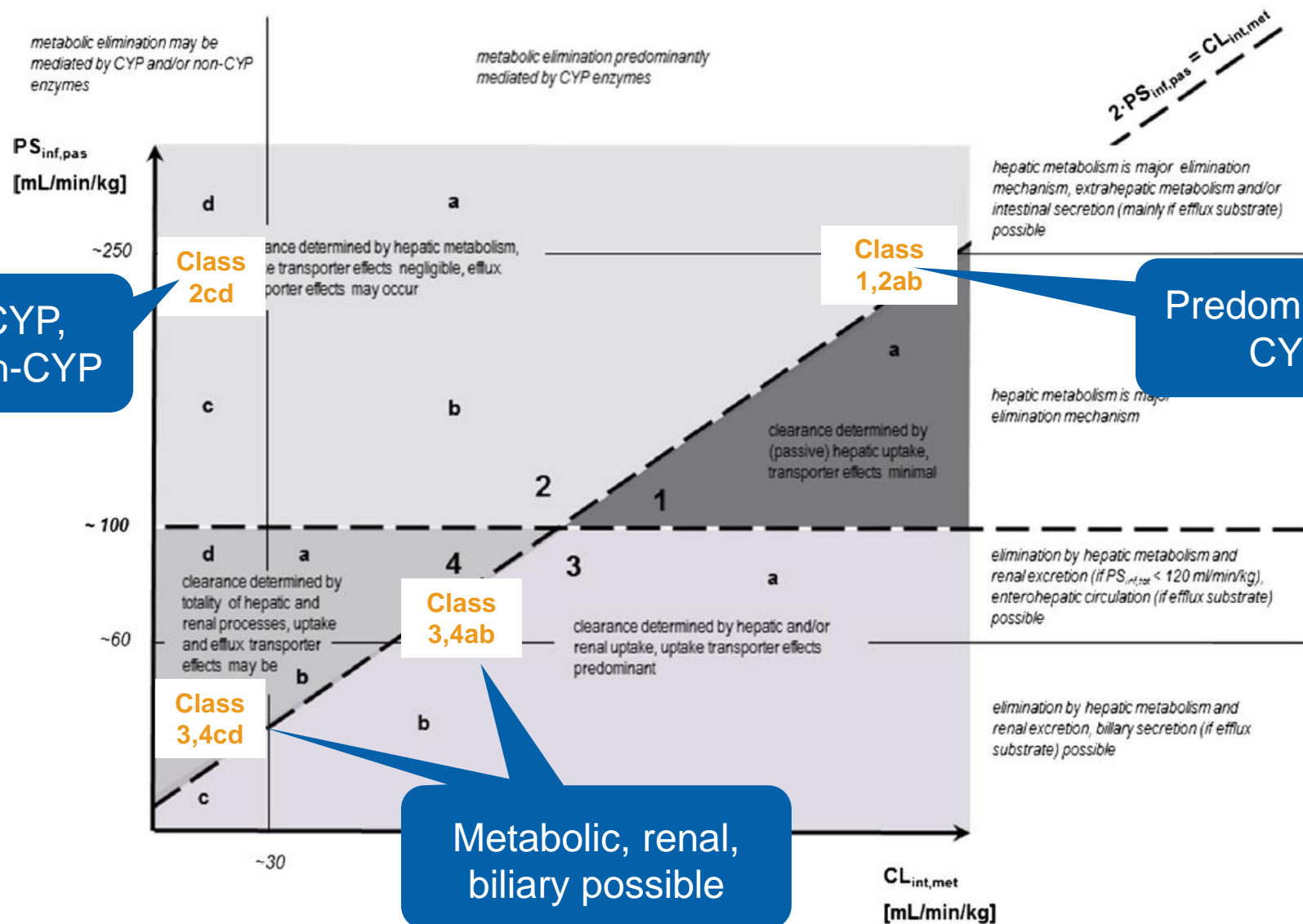
Classification based on *in vitro* permeability and *in vitro* metabolic and biliary clearance data

→ Allows to identify the rate-limiting clearance processes (absorption, distribution and elimination model)

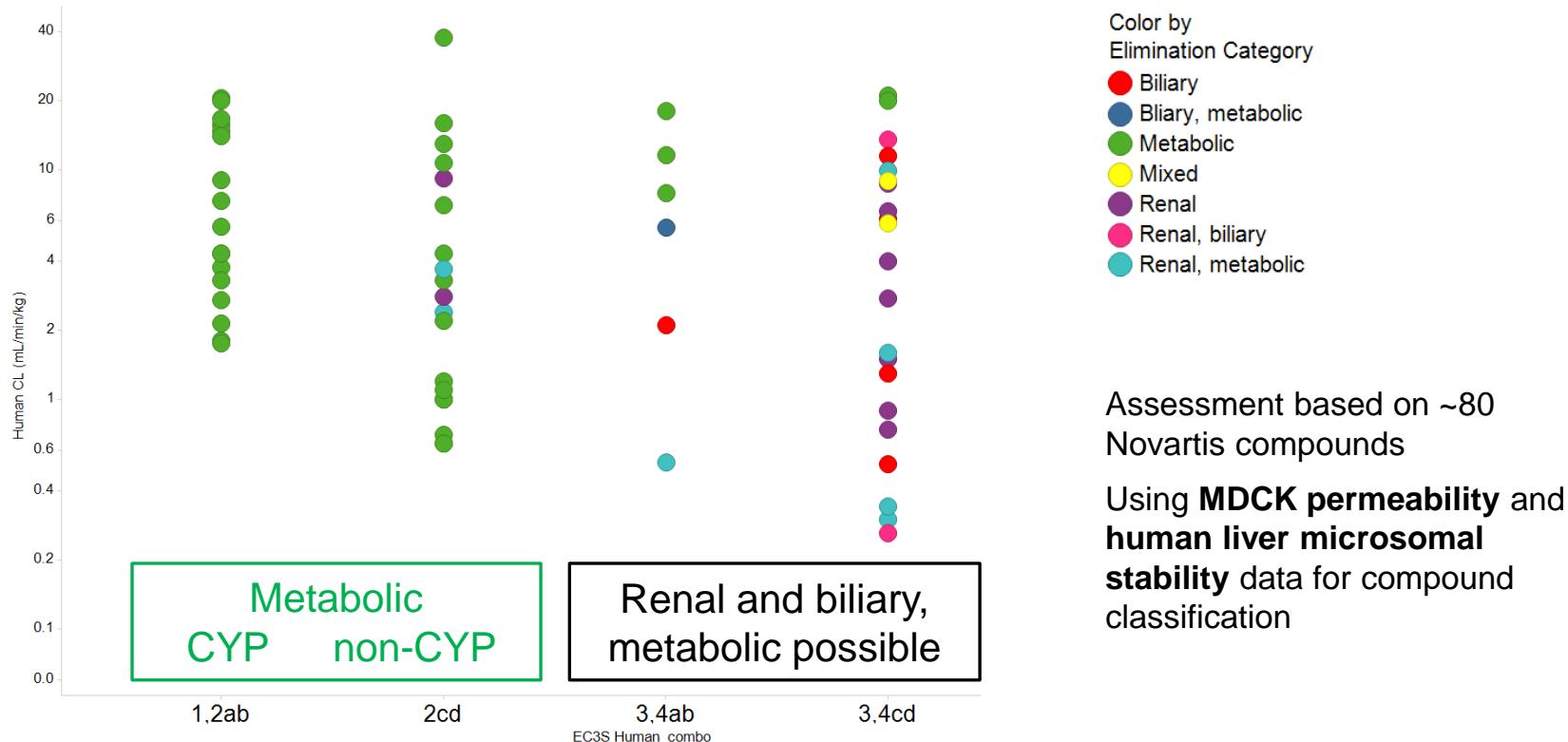
→ Rate-limiting step of hepatic elimination

→ **Model-based drug absorption, distribution and elimination drug classification system**

EC3S – Elimination mechanisms



EC3S – Elimination mechanisms

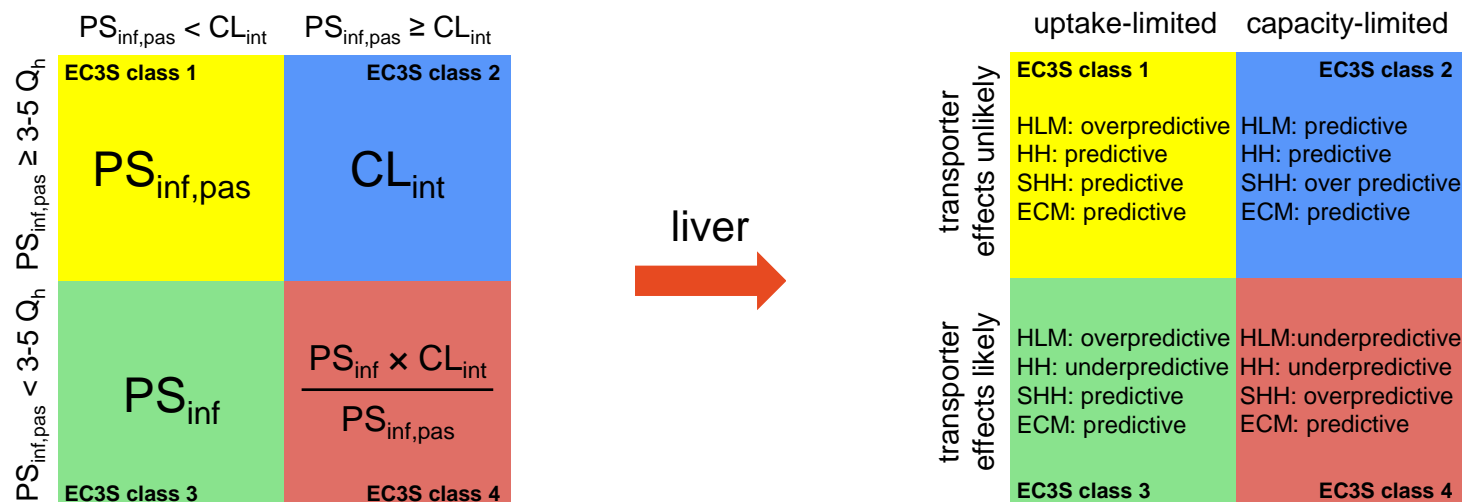


- Metabolic elimination generally well predicted (MDCK-LE $P_{app} > 5 \cdot 10^{-6}$ cm/s)
- EC3S provides information for CYP (Class 1,2ab) vs non-CYP (class 3,4ab)

→ Prediction of major elimination mechanisms in early development phase

EC3S – Hepatic clearance IVIVE

Rate-determining process



Hypothesis: knowing the rate-limiting process of hepatic elimination will facilitate selection of the most predictive clearance prediction tool

EC3S – Hepatic clearance IVIVE

Mechanism:	In vitro assay:
sinusoidal influx/efflux	suspended hepatocytes (SHH)
metabolism	liver microsomes (HLM), hepatocytes (HH)
biliary secretion	sandwich-cultured hepatocytes (SCH)
Extended Clearance Model (ECM)	HLM, SHH, SCH
plasma protein binding	ultrafiltration, ultracentrifugation or or equilibrium-dialysis

$$CL_{int,all} = CL_{met,u}$$

HH, HLM



$$CL_h = Q_h \cdot E_h = \frac{Q_h \cdot f_{u,b} \cdot CL_{int,all}}{Q_h + f_{u,b} \cdot CL_{int,all}}$$



$$CL_{int,all} = \frac{PS_{inf} \times (CL_{sec,u} + CL_{met,u})}{PS_{inf,pas} + (CL_{sec,u} + CL_{met,u})}$$

ECM

$$CL_{int,all} = PS_{inf}$$

SHH



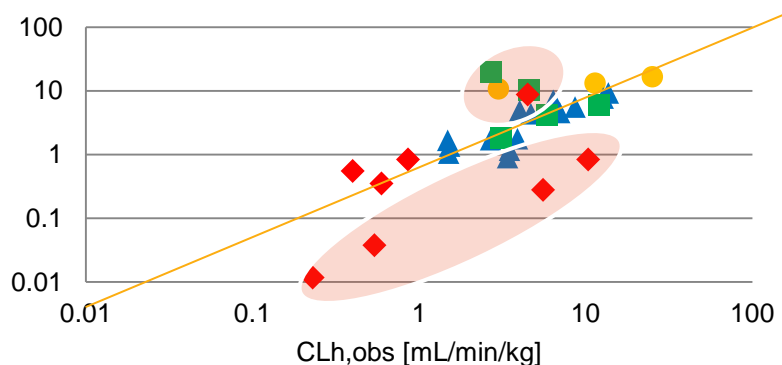
$$CL_{int,all} = \frac{PS_{inf} \times CL_{met,u}}{PS_{inf,pas} + CL_{met,u}}$$

ECM (-)

→ **Expectation: different outcomes depending on rate-limiting clearance mechanism (EC3S class-dependent)**

EC3S – Hepatic clearance IVIVE

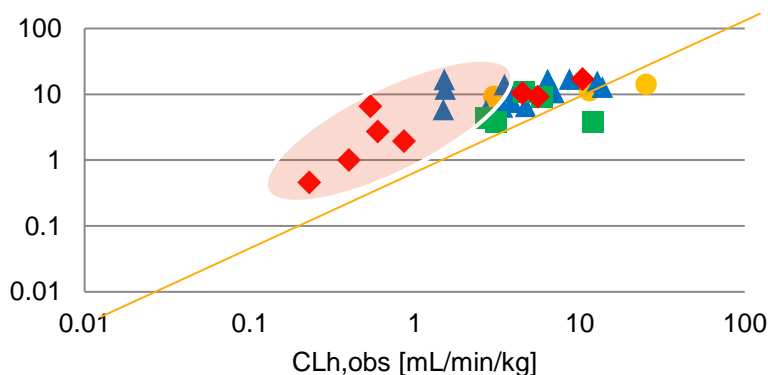
HLM:



$$CL_{int,all} = CL_{met}$$

- Class 2 generally well predicted
- Often under-predictive for class 4
- Tendency for being over-predictive for class 1 and class 3

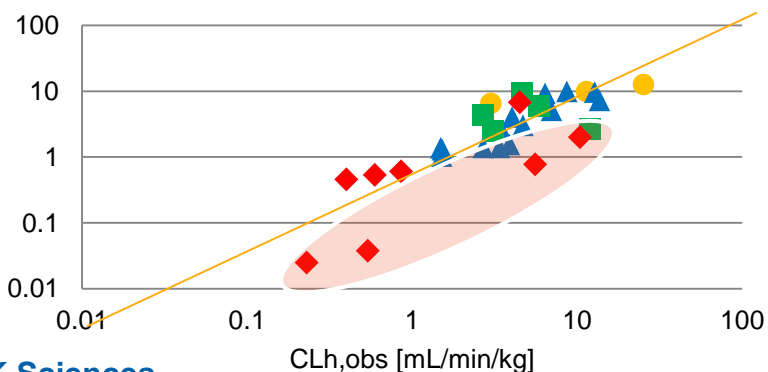
SHH



$$CL_{int,all} = PS_{inf}$$

- Class 1 and class 3 generally well predicted
- Over-predictive for some class 2 and class 4 cpds

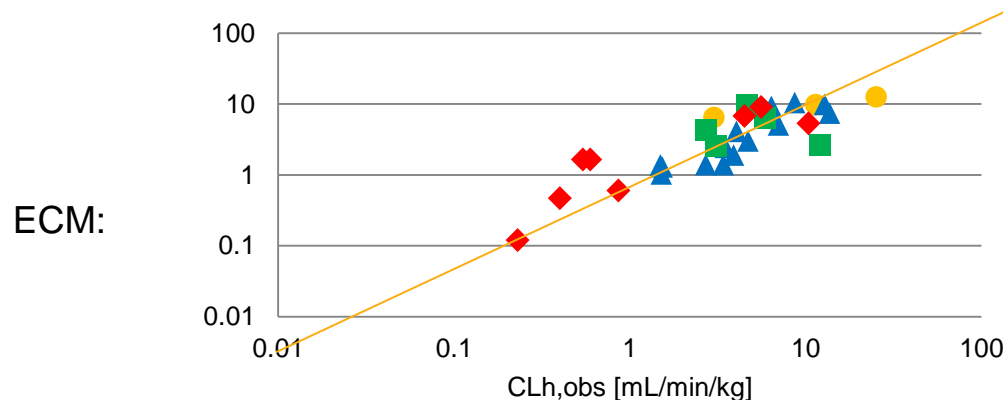
ECM (-):



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

- Class 1, class 3 and class 2 generally well predicted
- Under-predictive for some class 4 cpds

EC3S – Hepatic clearance IVIVE



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

→ Predictive for all EC3S classes

IVIVE recommendations:

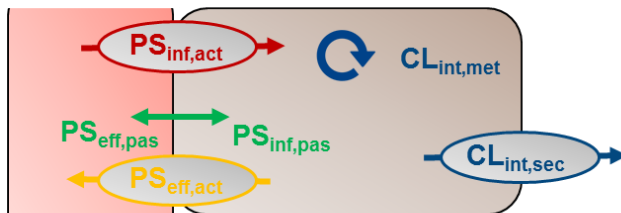
- **HH** is the method of choice for IVIVE of EC3S **class 1** cpds
- **HLM** or **HH** is the method of choice for IVIVE of EC3S **class 2** cpds
- **SHH** is recommended for EC3S **class 3** cpds (**HH** is the best alternative)
- **ECM** is needed for EC3S **class 4** cpds (no real alternative available)

EC3S – Total Clearance IVIVE

Estimation of fractional hepatic elimination

Total drug clearance (CL_{tot})

= hepatic drug clearance (CL_h) + renal drug clearance (CL_{ren})



Extended Clearance Model

no appropriate renal
in vitro model available

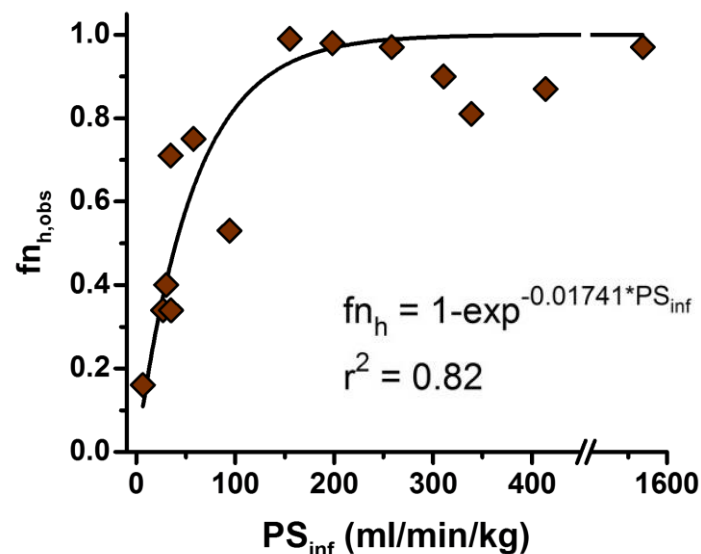
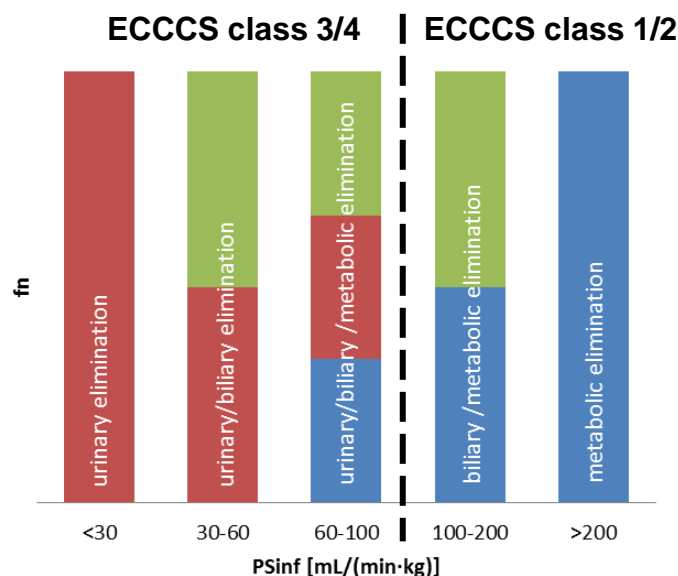
→ Is it possible to estimate relative contributions of hepatic (fn_h) and non-hepatic elimination pathways?

$$CL_{tot} = \frac{CL_h}{fn_h}$$

EC3S – Total Clearance IVIVE

Estimation of fractional hepatic elimination

Observation: hepatic uptake permeability correlates with elimination pathway



$$1 = fn_{ren} + \underbrace{fn_{sec} + fn_{met}}_{fn_h}$$

Camenisch et al, 2015, ADMET&DMPK; 3:1-14
Riede et al, 2016, Eur J Pharm Sci;86:96-102.

EC3S – Total Clearance IVIVE

Estimation of fractional hepatic elimination

ECCCS class 1/2:

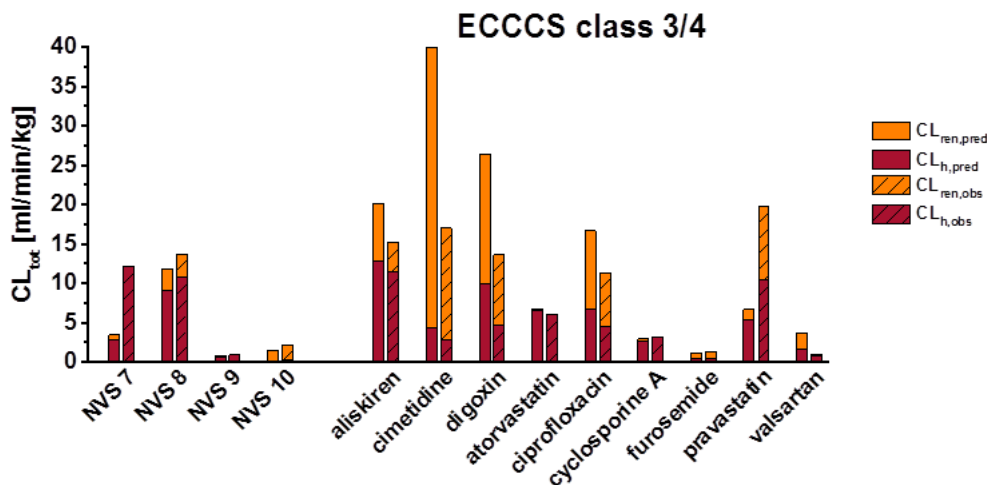
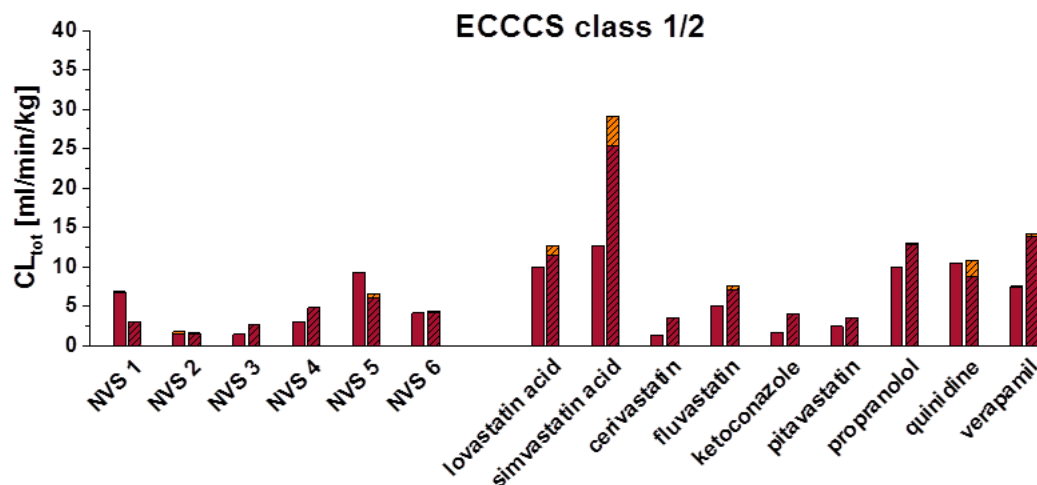
hepatic drug elimination

$$CL_{tot} = \frac{CL_h}{f_{n_h}}$$

ECCCS class 3/4:

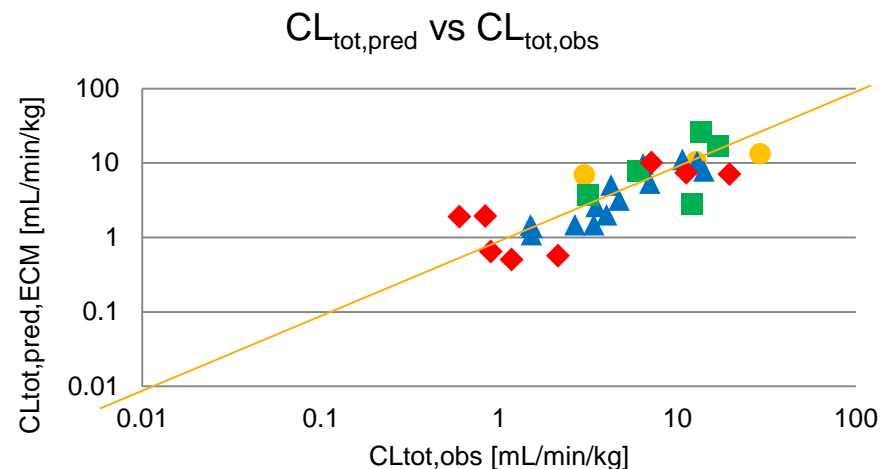
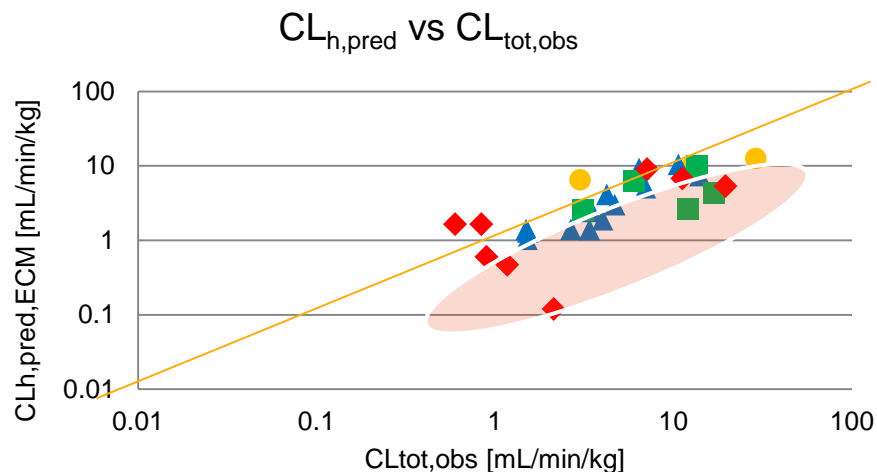
hepatic and **renal** drug elimination

→ Accurate prediction of total drug clearance independent of elimination pathways



EC3S – Total Clearance IVIVE

Estimation of fractional hepatic elimination



Prediction of total human clearance

- 1) $f_{n,h}$ estimated from $PS_{inf,pas}$
- 2) CL_{tot} calculated with : $CL_{tot,pred} = CL_{h,pred} / f_{n,h,pred}$

(assuming absence of other elimination routes)

Applications for drug classification

Guidance for the drug development process

- What is the recommendation with regards to metabolism investigations (*in vivo* or *in vitro*) ?
- What is the most appropriate clearance prediction tool for IVIVE?
- Are there opportunities to waive any animal studies (e.g. bile-duct cannulation studies)?
- What is the potential leverage with regards to in silico PK work?
- Is it recommended to synthesize a radiolabel in an early development phase?
- Which DDI follow-up studies (cpd as perpetrator vs victim) are recommended?

Applications for drug classification

	BDDCS	ECCS	EC3S
IVIVE		<ul style="list-style-type: none"> – Varma et al, Pharm Res (2015) 	<ul style="list-style-type: none"> – Umehara and Camenisch, Pharm Res (2012) – Camenisch and Umehara, Biopharm Drug Dispo (2012) – Riede et al, Eur J Pharm Sci (2016)
Elimination Mechanism	<ul style="list-style-type: none"> – Hosey et al, The AAPS Journal (2016) 	<ul style="list-style-type: none"> – Varma et al, Pharm Res (2015) – El Kattan et al, Pharm Res (2016) 	<ul style="list-style-type: none"> – Riede et al, Eur J Pharm Sci (2016)
DDI	<ul style="list-style-type: none"> – Shugarts and Benet, Phar Res (2009) 	<ul style="list-style-type: none"> – El Kattan et al, Pharm Res (2016) 	<ul style="list-style-type: none"> – Kunze et al, Drug Metab Pers Ther (2015)
$K_{p_{uu}}$			<ul style="list-style-type: none"> – Riede et al, Drug Metab Dispos (2017)
Food effect	<ul style="list-style-type: none"> – Custodio et al, Adv Drug Deliv Rev (2008) – Himbach et al, The AAPS Journal (2012) 		

Summary

- All compound classification systems provide information on drug disposition and the interplay between metabolic enzymes and transporters.
- ECCs and EC3S use *in vitro* data only. BDDCS requires information of a clinical dose and may therefore be positioned at a later stage in the drug development process.
- EC3S provides directly enables quantitative estimates of hepatic clearance and disposition processes given the required *in vitro* parameters are generated.
- All three classification systems may facilitate the compound class-dependent drug development process by guiding the selection of the most appropriate *in vitro* and *in vivo* studies

Acknowledgments

- Gian Camenisch
- Dallas Bednarczyk
- Sujal Deshmukh
- Bernard Faller
- Imad Hanna
- Anett Kunze
- Julia Riede
- Patrick Schweigler
- Kenichi Umehara

Thank you for your attention

