

# Timing and application of transporter studies in drug discovery & development

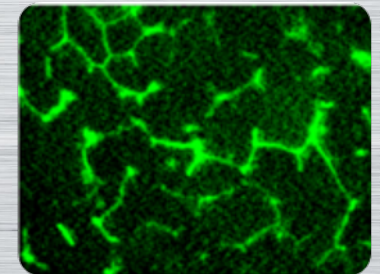
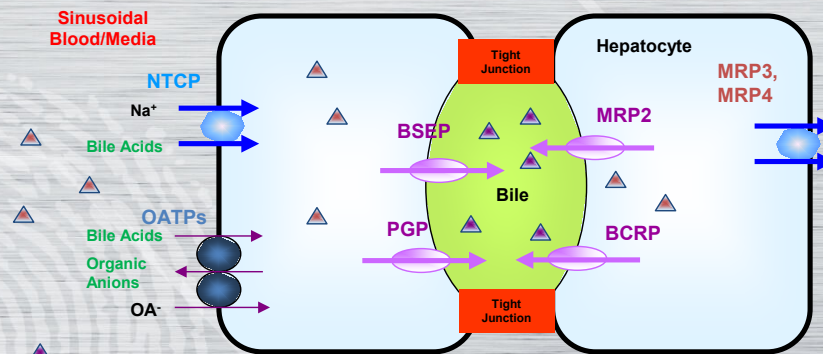
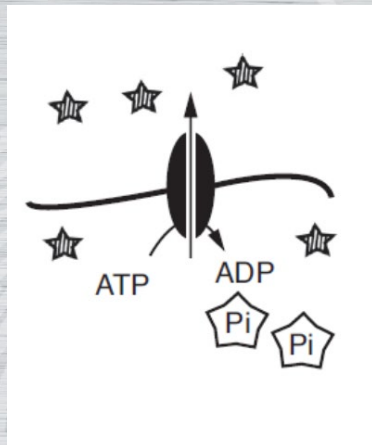
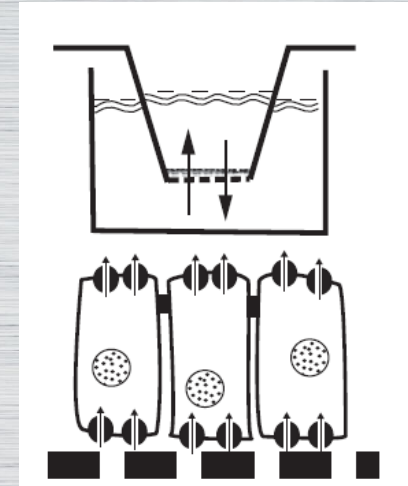
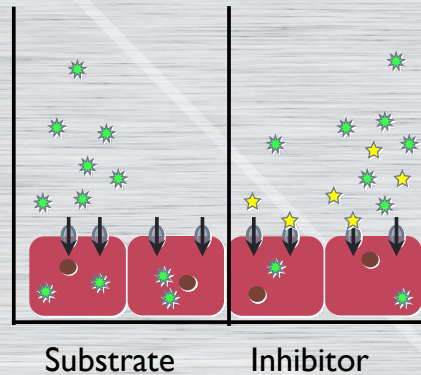
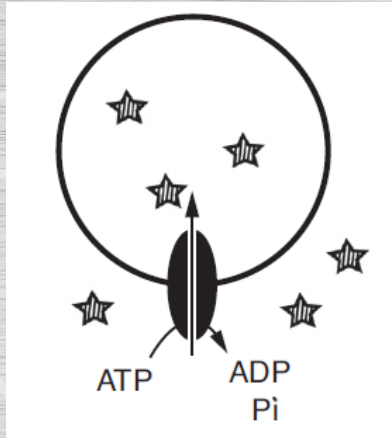
**Joe Zolnerciks, Ph.D.**

VP, Business Development

SOLVO Biotechnology | A Charles River Company  
Seattle, WA



# Multiple methods available for studying transporters





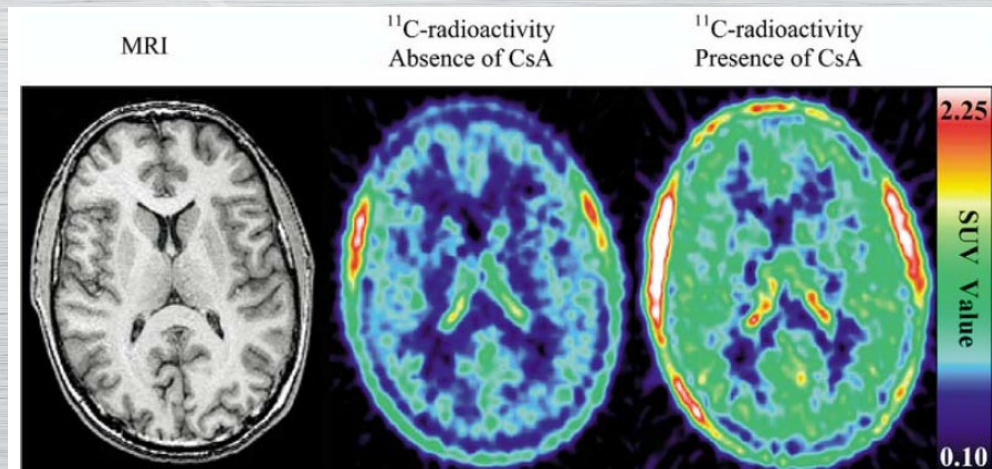
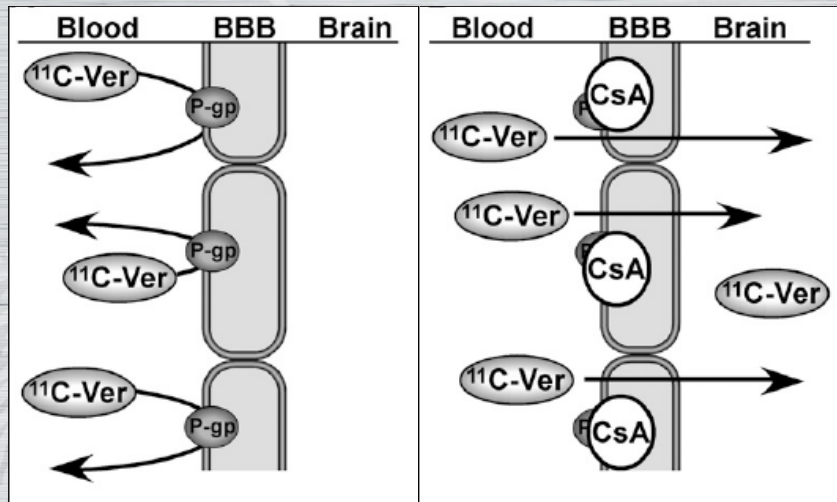
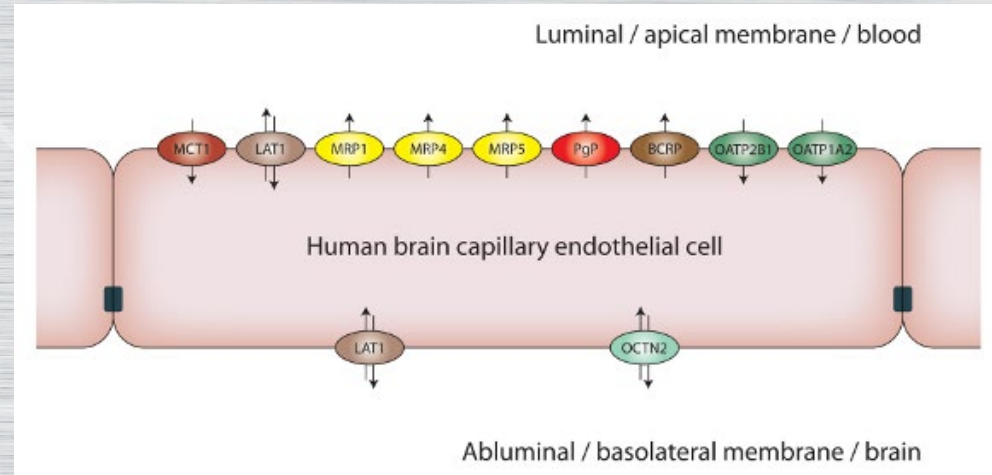
Why

~~When~~ should transporters be studied?

# Transporters can limit tissue distribution

## CNS Drugs

- P-glycoprotein (P-gp/MDR1;ABCB1)
  - Efflux transporter expressed on luminal side of blood-brain barrier
  - Prevents entry to CNS for transported substrates
  - PET imaging of  $^{11}\text{C}$ -labeled verapamil (P-gp substrate) shows marked increase in brain accumulation following co-administration with cyclosporine A (P-gp inhibitor)

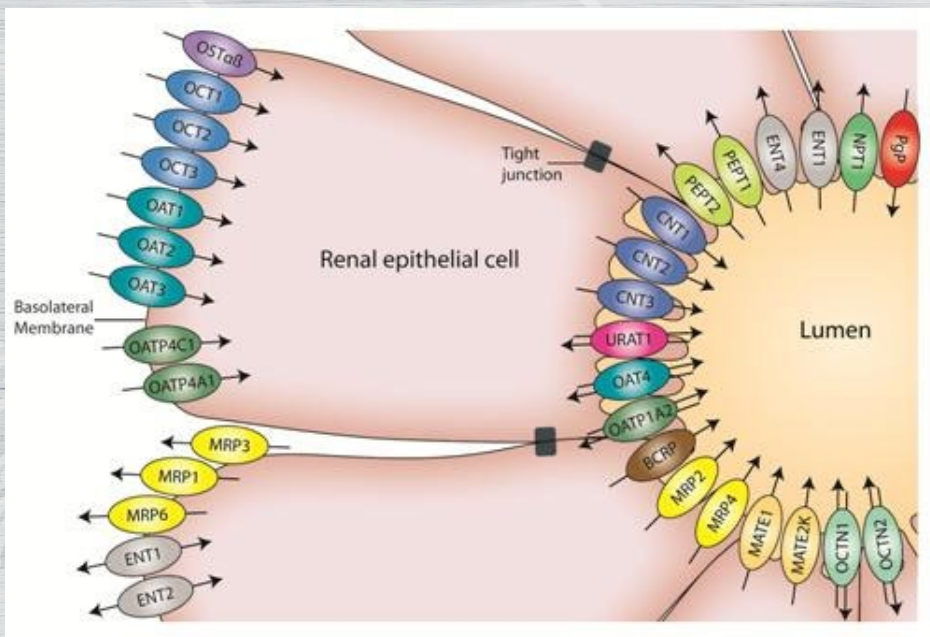
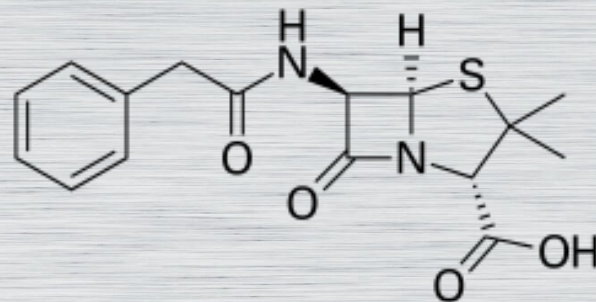




# Transporters can mediate drug clearance

## Benzylpenicillin (Penicillin G)

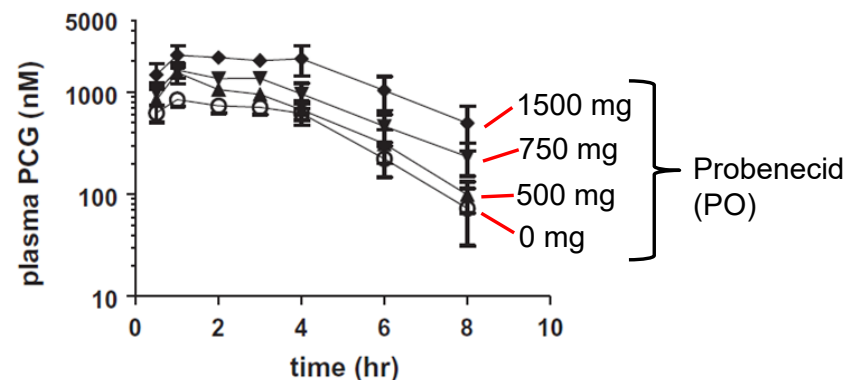
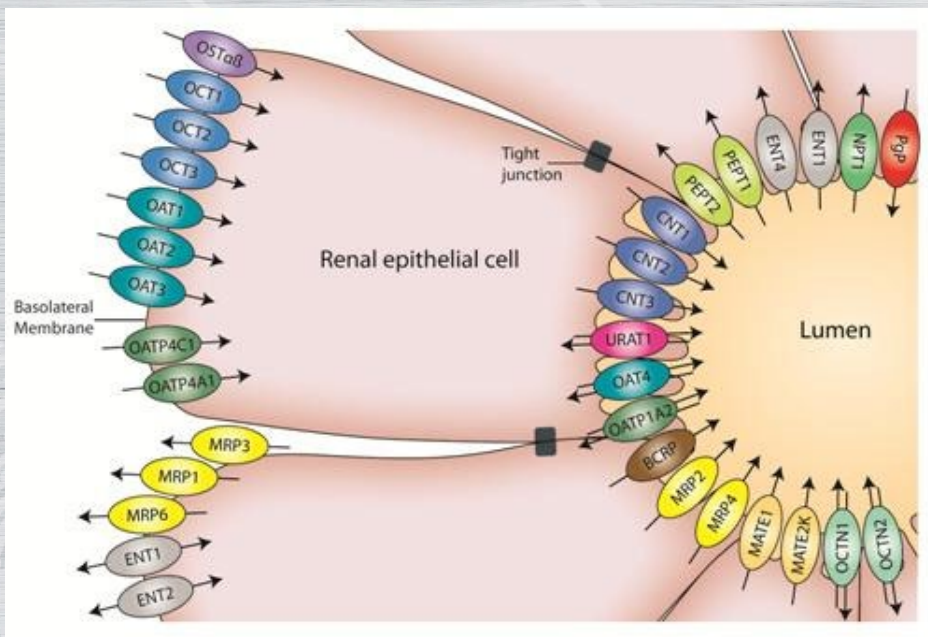
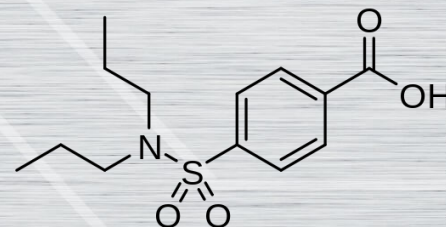
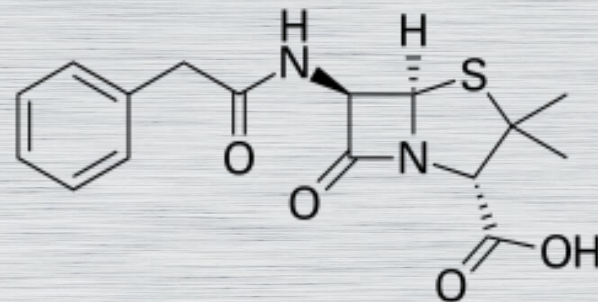
- Discovered in 1929
- Rapidly cleared via renal elimination (half life ~30 minutes)
- Supply so limited during early use, drug was re-isolated from urine of patients



# Transporters can mediate drug clearance

## Benzylpenicillin (Penicillin G) + Probenecid

- 3.3-fold increase in benzylpenicillin (PCG) AUC at highest dose of probenecid
- Renal clearance equal to GFR at highest dose (no active secretion of PCG)



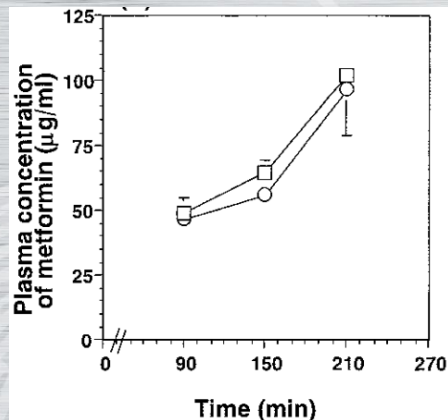


# Transporters can mediate tissue distribution

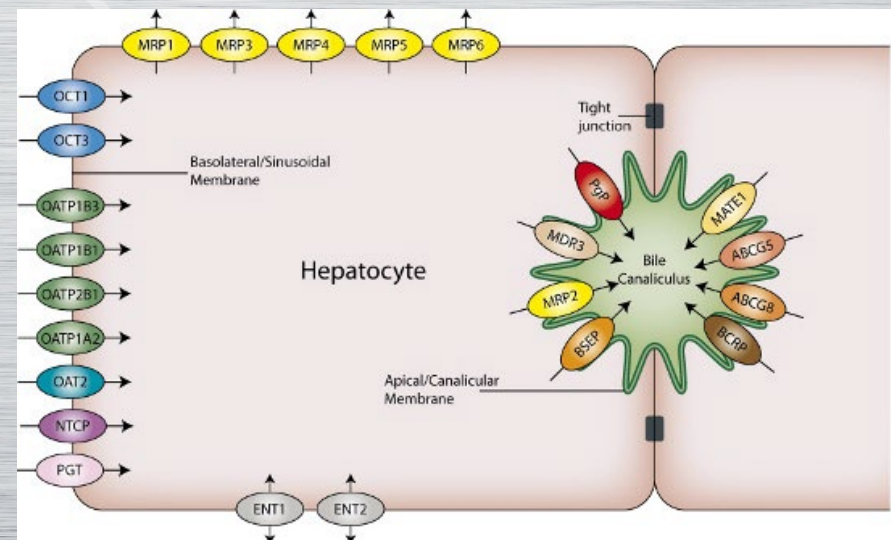
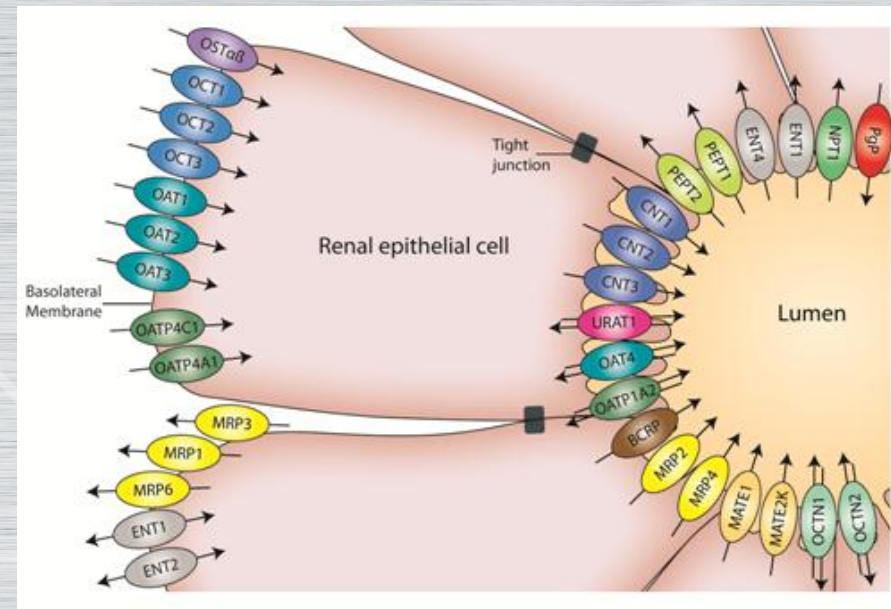
## Metformin

- First line therapy for type 2 diabetes
- Site of action- liver
- Primarily excreted in urine (active secretion by OCTs & MATEs)

	C <sub>plasma</sub>	Liver
	$\mu\text{g/ml}$	C <sub>liver</sub> *
	$\mu\text{g/g}$	$\mu\text{g/g}$
Wild-type	102 ± 11	417 ± 178
Oct(-/-)	97.7 ± 15.1	49.3 ± 10.4



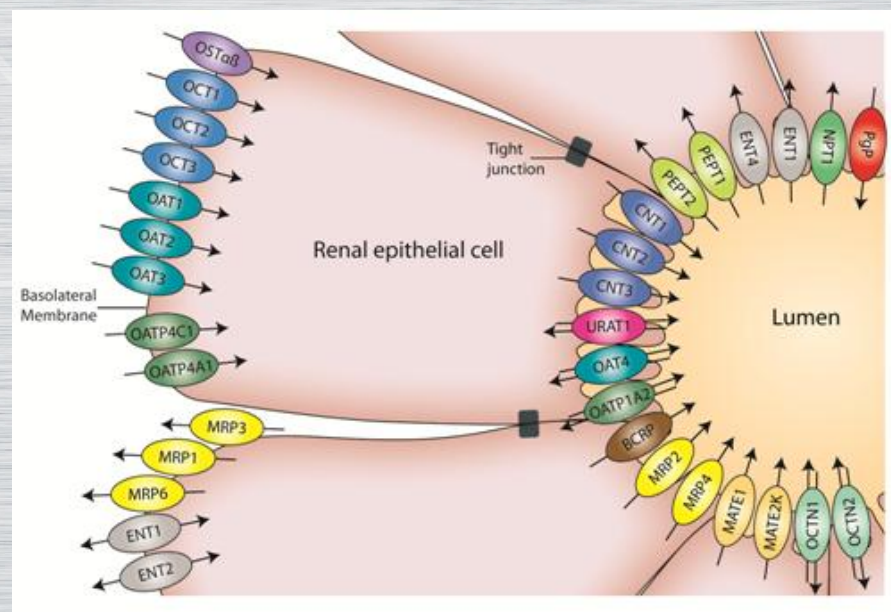
- OCT1(-/-) mice show no change in metformin plasma AUC
- ~90 % reduction in liver distribution



# Transporters as targets

## Transporters as targets

- Transporters are being harnessed as drug targets
  - Sodium-glucose cotransporter 2 (SGLT2; SLC5A2)
    - Canagliflozin (Invokana)
    - Dapagliflozin (Farxiga)
    - Empagliflozin (Jardiance)
  - Uric acid transporter 1 (URAT1; SLC22A12)
    - Lesinurad (Zurampic)
- Transporters are being utilized to improve absorption/distribution profile
  - Peptide transporter 1 (PEPT1; SLC15A1)
    - Valacyclovir
    - valganciclovir





# A changing regulatory environment

## Guidance for Industry

**2006:** FDA Draft Guidance (P-gp)

**2007:** Formation of International Transporter Consortium (ITC)

## REVIEWS

**2010:** ITC Transporter White Paper (P-gp, BCRP, OATPIBI, OATPIB3, OCT2, OAT1, OAT3)

**2010:** EMA Draft Guidance (P-gp, BCRP, OATPIBI, OATPIB3, OCT2, OAT1, OAT3, BSEP, OCT1)

**2012:** Revised FDA Draft Guidance (P-gp, BCRP, OATPIBI, OATPIB3, OCT2, OAT1, OAT3, BSEP, MATEs, MRPs)

**2013:** Seven ITC Whitepapers Published

**2013:** Final EMA Guidance (more detailed)

**2014:** PMDA Guidance published

**2017:** EMA Guidance Concept Paper Released

**2017:** Revised FDA Draft Guidance Released

**Drug Interaction Studies —  
Study Design, Data Analysis, and  
Implications for Dosing and Labeling**

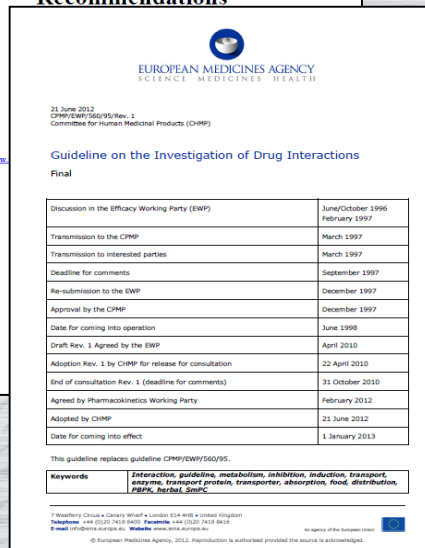
This guide  
Comments and suggestions  
publication in the FDA  
guidance. Submit to  
Drug Administration  
should be identified  
the Federal Register  
For questions regarding  
or (CBER) Toxicology

Memorandum  
developed

The International  
Abstract/Memo  
and efficacy  
including which  
which in vitro  
In addition, which  
should be considered  
International  
intended to have  
transporter in  
development  
transporter in  
questions (for  
understanding  
drug molecules



## Guidance for Industry Drug Interaction Studies — Study Design, Data Analysis, Implications for Dosing, and Labeling Recommendations



# EMA Guidance Update

Update to Guideline on the investigation of drug interactions

Concept paper released **Q1 2017**

Proposed changes include:

- Recommendations for *in vitro* studies on
  - Transport
  - Time-dependent inhibition
  - Induction
- Update of transporter list (for inhibition screening)
- Update on cut-offs for transporter inhibition
- Transport as rate-limiting step
- Clarifications of guideline text
- Discussion of DDI study requirement with contraceptive steroids



# FDA Draft Guidance Revision

Revision to Guidance for Industry Drug Interaction Studies – Study Design, Data Analysis, Implications for Dosing, and Labeling Recommendations

Released **October 2017**

Changes include:

- Preparation of two separate guidance documents
  - *In vitro*
  - *In vivo*
- Update of transporter list (for inhibition screening)
- Update on cut-offs for transporter inhibition
- Additional recommendations for *in vitro* experiments

---

## In Vitro Metabolism- and Transporter- Mediated Drug-Drug Interaction Studies Guidance for Industry

### DRAFT GUIDANCE

This guidance document is being distributed for comment purposes only.

Comments and suggestions regarding this draft document should be submitted within 90 days of publication in the *Federal Register* of the notice announcing the availability of the draft guidance. Submit electronic comments to <https://www.regulations.gov>. Submit written comments to the Division of Dockets Management (HFA-305), Food and Drug Administration, 5630 Fishers Lane, rm. 1061, Rockville, MD 20852. All comments should be identified with the docket number listed in the notice of availability that publishes in the *Federal Register*.

For questions regarding this draft document, contact (CDER) Office of Clinical Pharmacology, Guidance and Policy Team at [CDER\\_OCP\\_GPT@fda.hhs.gov](mailto:CDER_OCP_GPT@fda.hhs.gov).

U.S. Department of Health and Human Services  
Food and Drug Administration  
Center for Drug Evaluation and Research (CDER)

October 2017  
Clinical Pharmacology

---

100417

# Types of transporter study

## Inhibition studies

- “Perpetrator” studies
- What is the risk of your NME causing a clinically significant change in disposition of co-administered drug(s)
- Same basic panel required for all NMEs

## Substrate studies

- “Victim” studies
- What is the risk of co-administered drug(s) changing the disposition of your NME
- Transporter panel based upon important clearance pathways of NME





# Regulatory Requirements

		INHIBITION STUDIES		SUBSTRATE STUDIES	
		EMA	FDA	EMA	FDA
EFFLUX	Transporter				
	P-gp	yes	yes	consider	yes
	BCRP	yes	yes	consider	yes
	BSEP	prefer	no	consider	no
UPTAKE	MRPs	no	no	consider	no
	OAT1	yes	yes	consider	≥25 % of elimination is active renal
	OAT3	yes	yes	consider	
	OATP1B1	yes	yes	≥25 % of elimination hepatic	≥25 % of elimination hepatic or biliary
	OATP1B3	yes	yes		
	OCT1	consider	no	consider	no
	OCT2	yes	yes	consider	≥25 % of elimination is active renal
	MATE1	consider	yes	consider	
	MATE2-K	consider	yes	consider	

# NDA Approvals - Transporter Information

Period	Percentage of Total NDA Approvals *							
	Pgp	BCRP	OATPs	OATs	OCTs	MATES	BSEP	MRPs
<b>2006-2011 (n= 183)**</b>	86 %	15 %	18 %	8 %	15 %	-	-	-
<b>2012 (n=30)</b>	73 %	37 %	40 %	23 %	30 %	3 %	3 %	13 %
<b>2013 (n=21)</b>	95 %	38 %	52 %	24 %	33 %	10 %	14 %	14 %
<b>2014 (n=25)</b>	96 %	68 %	72 %	64 %	72 %	16 %	20 %	44 %
<b>2015 (n=32)</b>	84 %	66 %	75 %	59 %	69 %	28 %	34 %	38 %
<b>2016 (n=14)</b>	93 %	93 %	86 %	71 %	79 %	36 %	43 %	43 %
<b>2017 (n=35)</b>	86 %	77 %	77 %	74 %	69 %	20 %	34 %	20 %
<b>to Q3 2018 (n=36)</b>	89 %	89 %	86 %	83 %	83 %	53 %	42 %	22 %

\* Total approvals adjusted for 2013 onwards to exclude “Biologicals” (MABs and Peptide therapeutics), imaging agents and topical products.

\*\* Tweedie et al, CPT, 94(1), 113-25, 2013



# NDA Approvals - Transporter Information

- Predominantly in vitro data
  - Predominantly inhibition (perpetrator) studies reported
  - All transporter families significantly represented
- Total number of clinical investigations broadly similar
  - 19, 35, 28, 11, 29 and 30 each year for 2013- Q3 2018, respectively
- Range of transporters tested clinically has broadened markedly
  - Consequence of increased range of transporters investigated in vitro
- Occasional post-marketing commitments
- Negative and positive findings referred to in labels
- Trend to combine metabolism and transporter interactions
  - Increasing use of PBPK modelling approaches

When should transporters be studied?



# When should transporters be studied?

- Regulatory considerations
- Patient population
  - Therapeutic indication and likely co-meds
- Physicochemical properties of compound
- Distribution/site of action
- Safety assessment

# Regulatory considerations - EMA Viewpoint

## Substrate studies

*In vitro* information supporting the prediction of the effects by other medicinal products on the pharmacokinetics of the investigational drug should preferably be available before introducing the investigational product to patients (phase II) and is generally required before starting phase III.

## Inhibition studies

*In vitro* data on the effects of the investigational parent drug on the pharmacokinetics of other drugs should preferably be available before starting phase II studies unless all concomitant drug treatments at risk of being clinically relevantly affected can be avoided in these studies. The *in vitro* information should be available before starting phase III.

- Preferably available before starting phase II, required before phase III

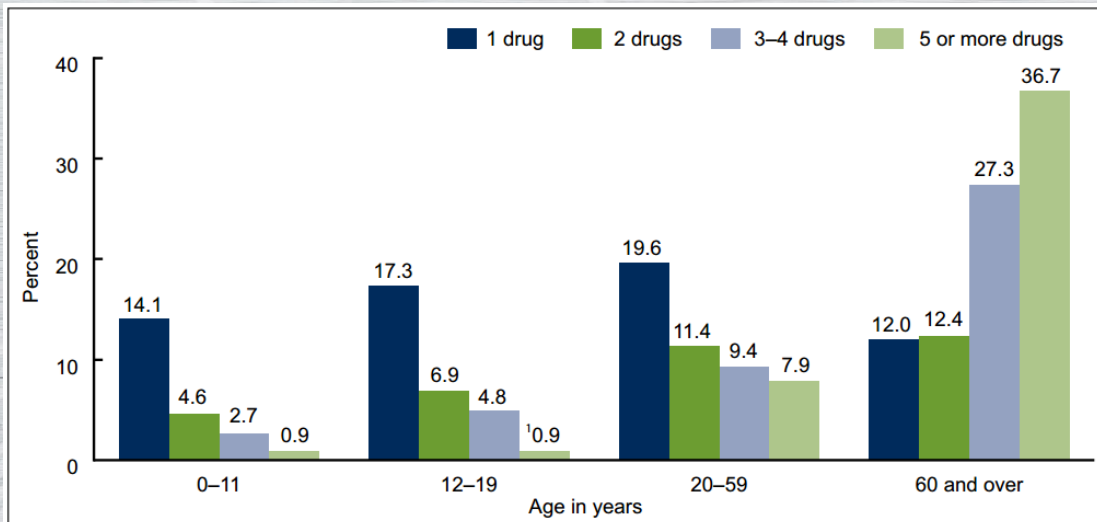
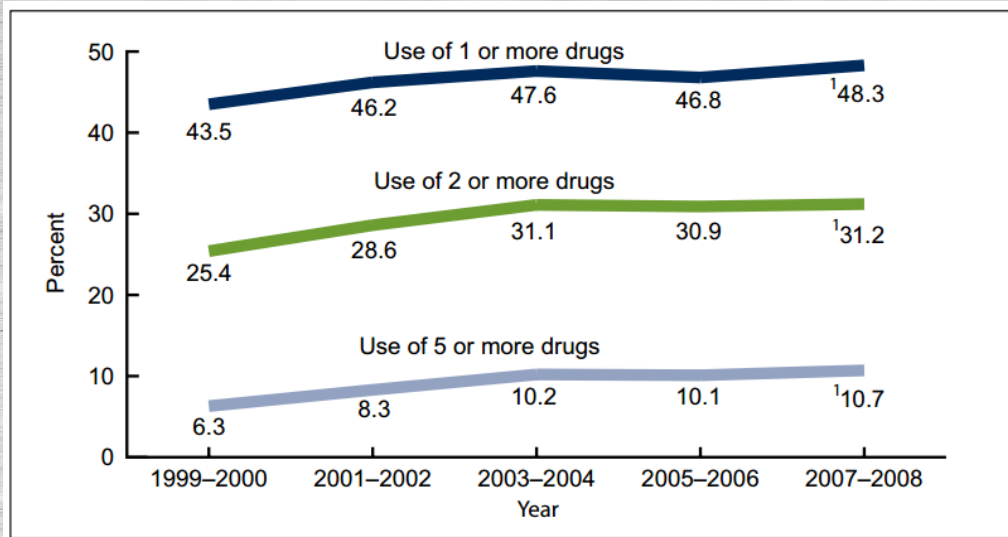


# Regulatory considerations - FDA Viewpoint

The timing of the in vitro evaluation of each transporter may vary depending on the therapeutic indications of the investigational drug. For example, if the intended population is likely to use statins, the sponsor should examine the potential of the investigational drug to interact with OATP1B1/1B3 before clinical studies in patients. If in vitro experiments indicate a low potential for an interaction between the transporter and investigational drug, subjects taking statins may be included in clinical studies to better represent the intended patient population.

- Timing of in vitro studies depends on therapeutic indication and patient population
- Can be before clinical studies

# Patient population - Polypharmacy

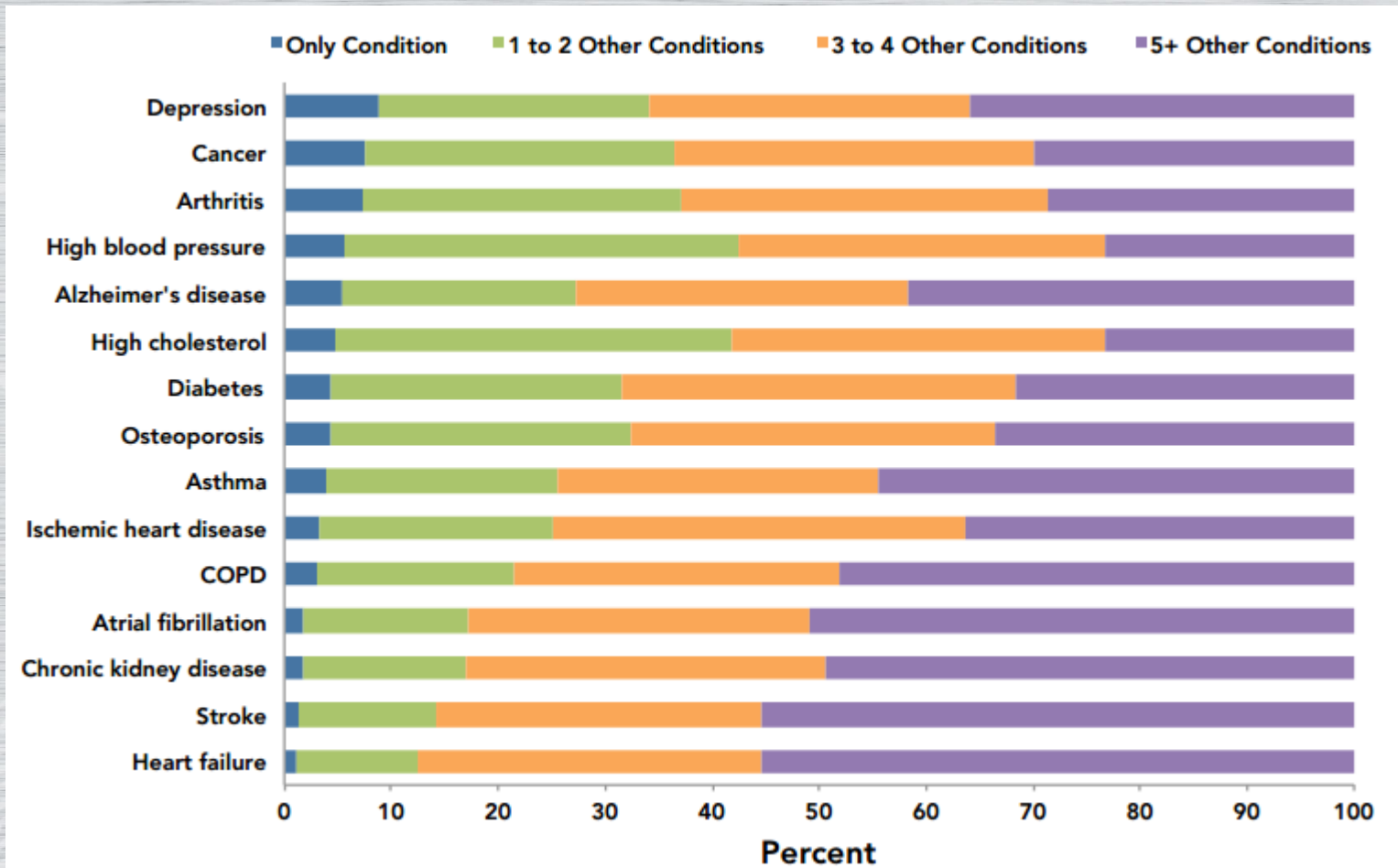


<sup>1</sup>Estimate is unstable; the relative standard error is greater than 30%.

- Increasing prevalence of polypharmacy
- Increased risk of DDI and ADR
- 'Prescribing cascade'



# Patient population – Co-morbidities



# Patient population – Co-morbidities

## FIVE MOST PREVALENT DYADS

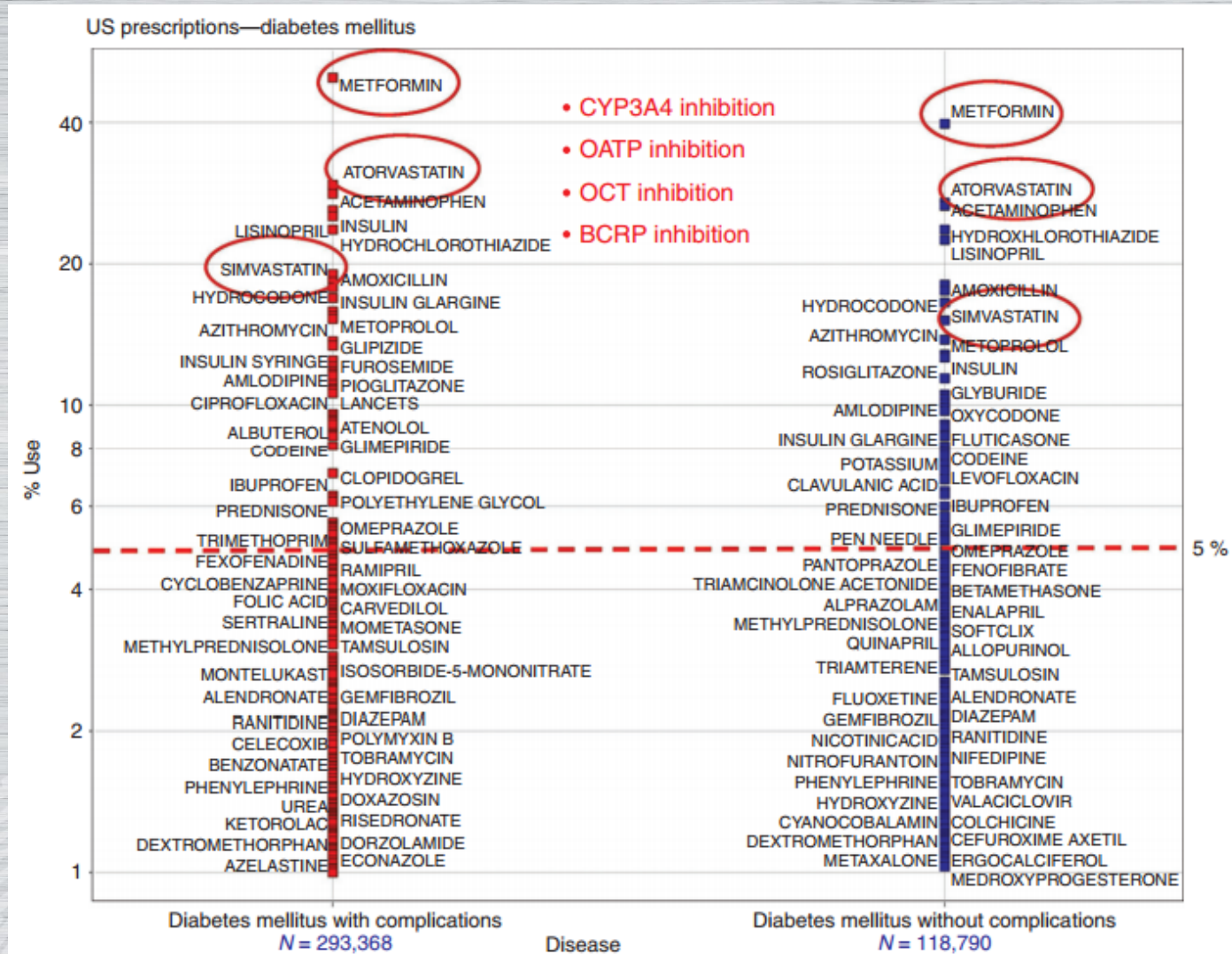
Dyads	Prevalence (%)
High cholesterol and High blood pressure	52.9
High cholesterol and Ischemic heart disease	36.2
High cholesterol and Diabetes	32.3
High cholesterol and Arthritis	31.1
Ischemic heart disease and High blood pressure	29.6

## FIVE MOST PREVALENT TRIADS

Triads	Prevalence (%)
High cholesterol and High blood pressure and Ischemic heart disease	33.7
High cholesterol and High blood pressure and Diabetes	29.9
High cholesterol and High blood pressure and Arthritis	25.7
High cholesterol and Diabetes and Ischemic heart disease	21.5
High cholesterol and Ischemic heart disease and Arthritis	19.3



# Patient population – Co-meds



# Victim Drugs - Therapeutic Index

Victim DDI for NTI drugs	Pgp	BCRP	OATP	OAT	OCT
Severe clinical risk (toxicity severe in nature, options for management include contraindication, dose adjustment, directed monitoring)	Digoxin	Topotecan	Repaglinide	Methotrexate	Dofetilide Pilsicainide Procainamide
Moderate clinical risk (potential consequences of toxicity are less severe in nature and/or risk is monitorable before serious sequelae occur)		Atorvastatin Fluvastatin Rosuvastatin Simvastatin	Atorvastatin Pitavastatin Pravastatin Rosuvastatin Simvastatin Glyburide		Metformin



# Victim Drugs - Therapeutic Index

Therapeutic window	Wide	<b>1</b> <b>No impact of DDIs</b> No dedicated studies	<b>2</b> <b>Low impact of DDIs</b> Focussed DDI evaluation; may defer to later phases of clinical development
	Narrow	<b>3</b> <b>Low/medium impact of DDIs</b> Focussed DDI evaluation; may defer to later phases of clinical development; restriction of use may be required	<b>4</b> <b>High impact of DDIs</b> Comprehensive early DDI evaluation; seek NCE without DDI liability if feasible
		Low	High
		Frequency of co-medication use	

# Physicochemical Properties – BCS and BDDCS

## Biopharmaceutical Drug Disposition Classification System (BDDCS)

## Biopharmaceutics Classification System (BCS)

		High Solubility	Low Solubility
Permeability	High	<b><u>Class 1</u></b> High Solubility High Permeability Rapid Dissolution	<b><u>Class 2</u></b> Low Solubility High Permeability
	Low	<b><u>Class 3</u></b> High Solubility Low Permeability	<b><u>Class 4</u></b> Low Solubility Low Permeability

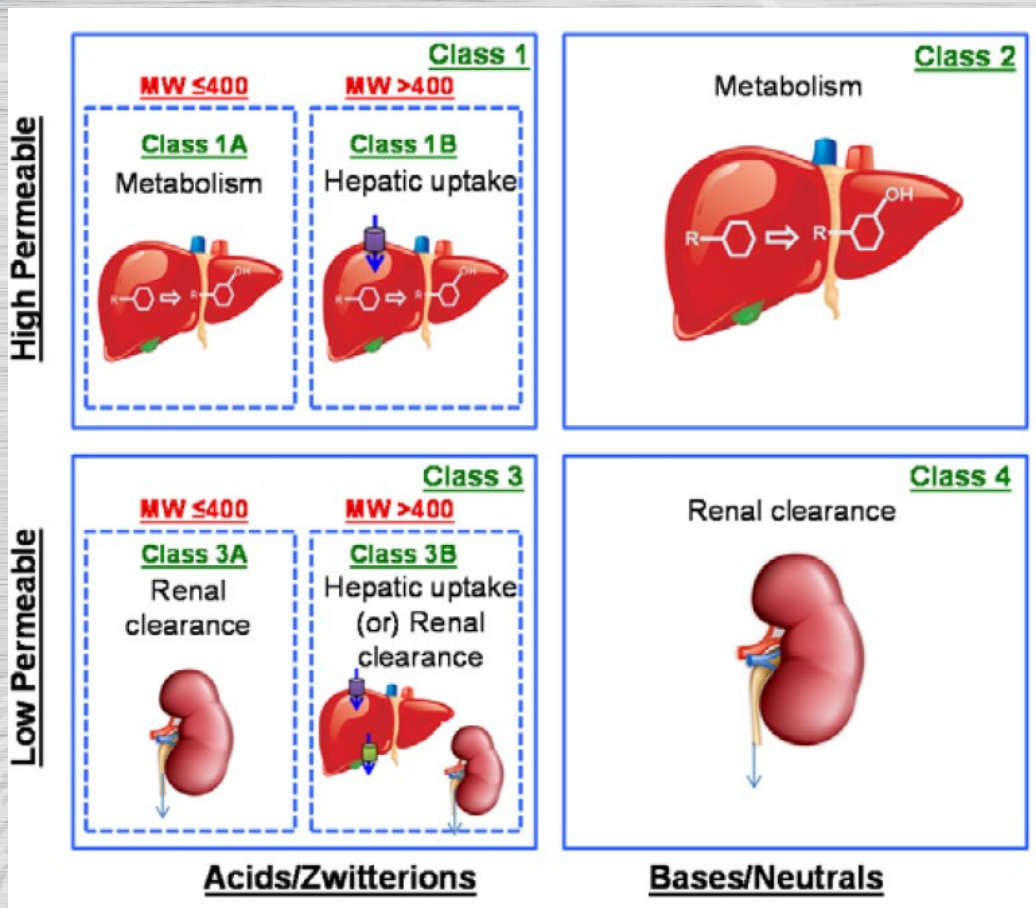
		Oral Dosing: Transporter Effects By BDDCS Class	
		High Solubility	Low Solubility
Metabolism	Extensive	<b><u>Class 1</u></b> Transporter effects minimal	<b><u>Class 2</u></b> Efflux transporter effects predominate in the gut, while absorptive and efflux transporter effects occur in the liver
	Poor	<b><u>Class 3</u></b> Absorptive transporters effects predominate (but may be modulated by efflux transporters)	<b><u>Class 4</u></b> Absorptive and efflux transporters effects could be important

Fig. 5. Transporter effects, following oral dosing, by BDDCS class.



# Physicochemical Properties – ECCS

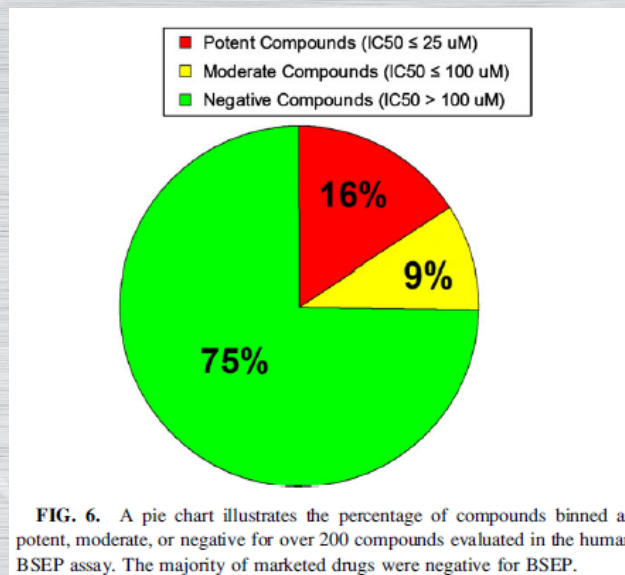
## Extended Clearance Classification System (ECCS)



- Prediction of clearance rate-determining step using physicochemical properties (permeability, molecular weight, ionization)

# Safety considerations - cholestasis

- Ranking cpds based on BSEP  $IC_{50}$ , 25  $\mu M$  cutoff
- 79 % of cpds with BSEP  $IC_{50} < 25 \mu M$  associated with DILI
- $C_{ss} / BSEP IC_{50} > 0.1$  gave 95 % correlation with DILI incidence
- Recommendations:
  - BSEP-VT screen for potent inhibitors
  - Confirmatory transporter assays (MRP2, MRP3 and MRP4) may be helpful



**TABLE 2**  
Number of Compounds With Evidence of Liver Injury/Total Number of Compounds Fitting Column and Row Criteria (%)

Transporter Assay	$C_{ss} / IC_{50}$ Ratio < 0.01	$C_{ss} / IC_{50}$ Ratio < 0.1	$C_{ss} / IC_{50}$ Ratio $\geq 0.1$
BSEP	18/44 (41%)	34/70 (49%)	36/38 (95%)
MRP2	6/9 (67%)	9/13 (69%)	1/1 (100%)
MRP3	7/11 (64%)	17/23 (74%)	5/6 (83%)
MRP4	10/23 (53%)	26/39 (67%)	14/17 (82%)

*Notes.* The closer exposure values in humans approach *in vitro* potency values in the transporter assays, the stronger the association with liver injury. Conversely, as the exposure values fall further below the *in vitro* potency values, the weaker the association with liver injury.



# Safety considerations - cholestasis

## Transporter Panel Flow Scheme for Hazard Identification

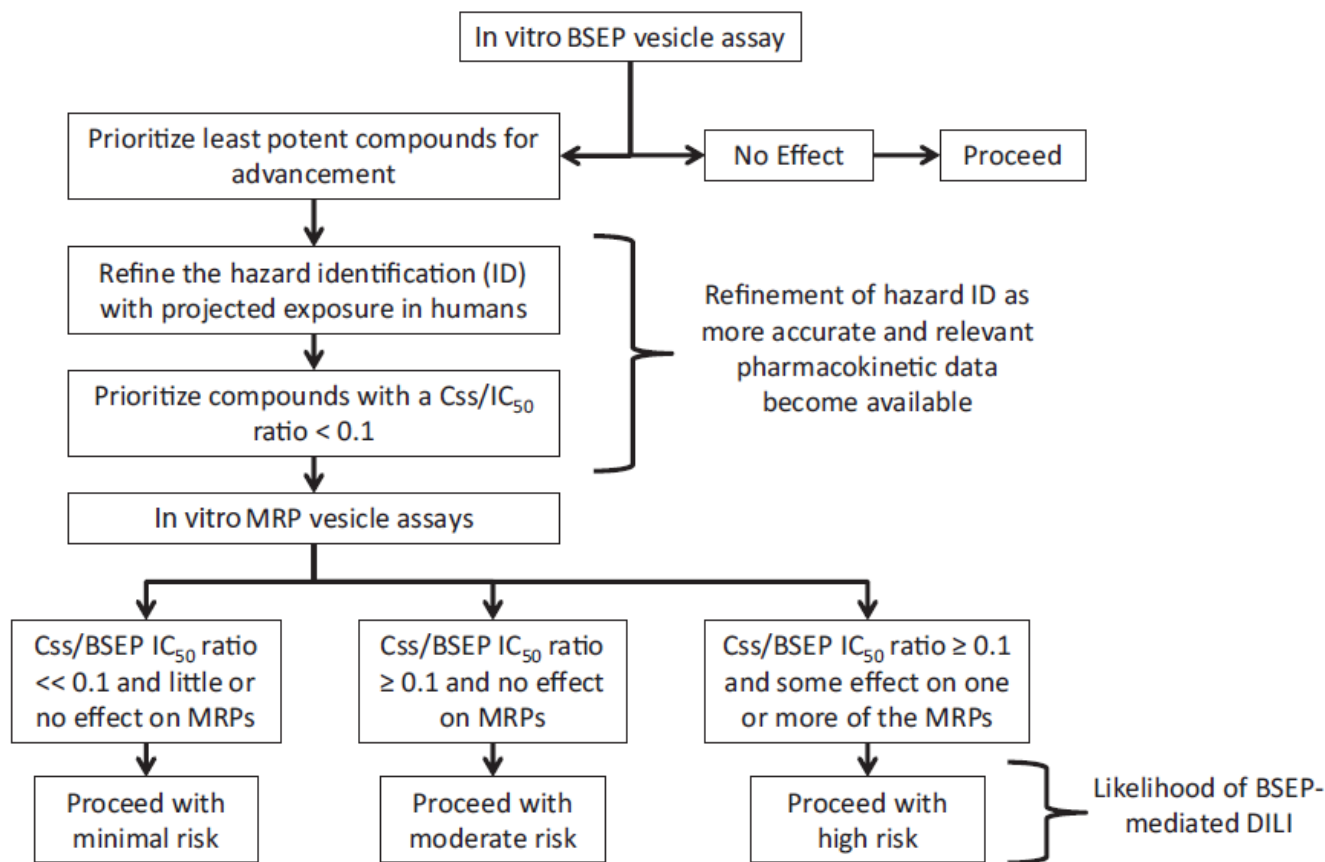


FIG. 7. Flow scheme for deploying a transporter panel during early therapeutic compound development. Abbreviations: BSEP, bile salt export pump; DILI, Drug-induced liver injury; MRP, multidrug resistance-associated protein.

# Safety considerations - cholestasis



SOT | Society of  
Toxicology  
www.toxsci.oxfordjournals.org

TOXICOLOGICAL SCIENCES, 167(2), 2019, 458–467

doi: 10.1093/toxsci/kfy253

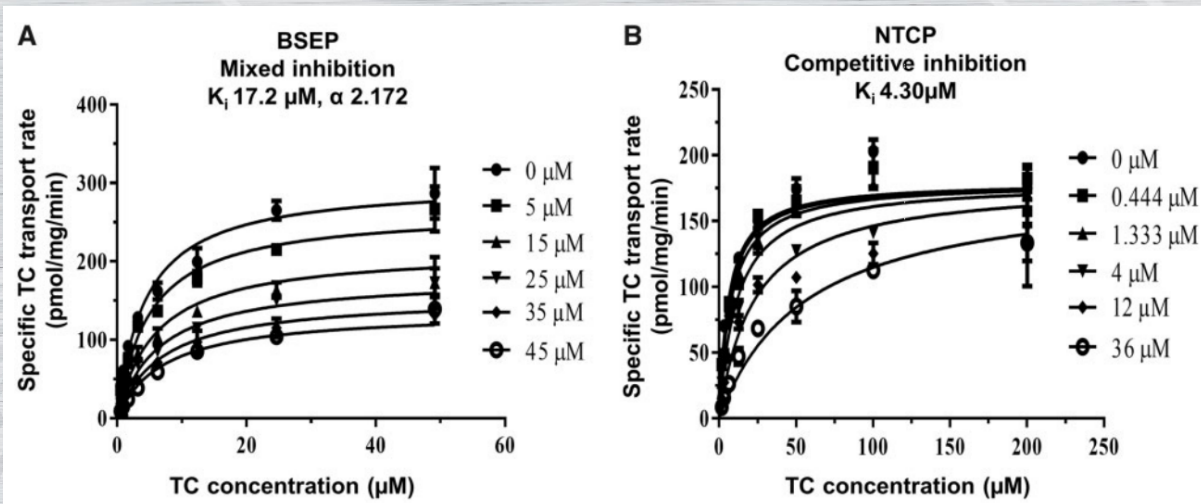
Advance Access Publication Date: October 5, 2018

Research Article

## Quantitative Systems Toxicology Analysis of *In Vitro* Mechanistic Assays Reveals Importance of Bile Acid Accumulation and Mitochondrial Dysfunction in TAK-875-Induced Liver Injury

Diane M. Longo,<sup>\*,1</sup> Jeffrey L. Woodhead,<sup>\*</sup> Paul Walker,<sup>†</sup> Krisztina Herédi-Szabó,<sup>‡</sup> Károly Mogyorósi,<sup>‡</sup> Francis S. Wolenski,<sup>§</sup> Yvonne P. Dragan,<sup>§</sup> Merrie Mosedale,<sup>¶,||</sup> Scott Q. Siler,<sup>\*</sup> Paul B. Watkins,<sup>\*,¶,||</sup> and Brett A. Howell<sup>\*</sup>

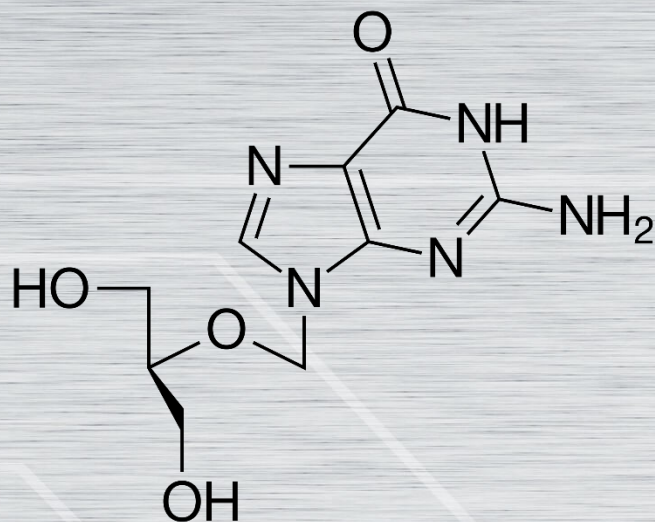
- TAK-875 withdrawn in Ph.III due to liver toxicity
- Used *in vitro* mechanistic data together with PBPK modeling (DILIsym)
- Mode of transporter inhibition found to be significant for predicting tox



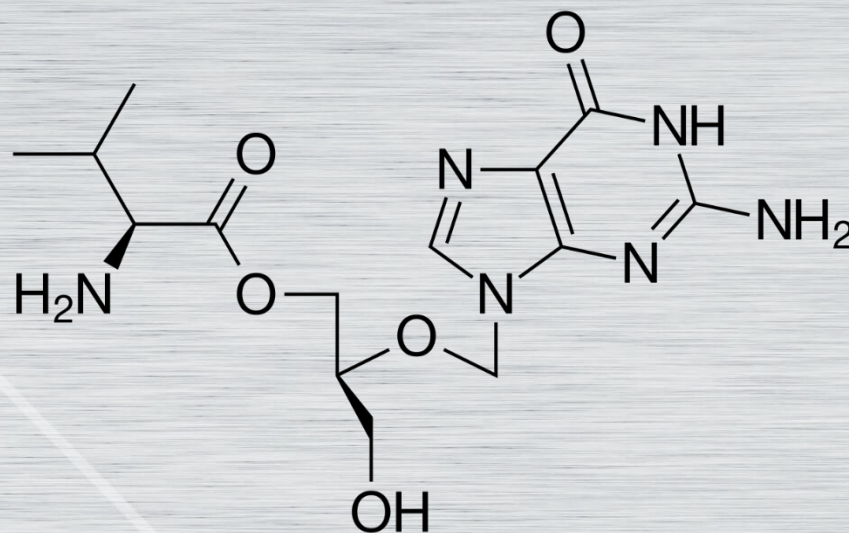
Longo, Toxicol Sci 2019



# Distribution – Prodrugs



Ganciclovir  
~ 9% bioavailability

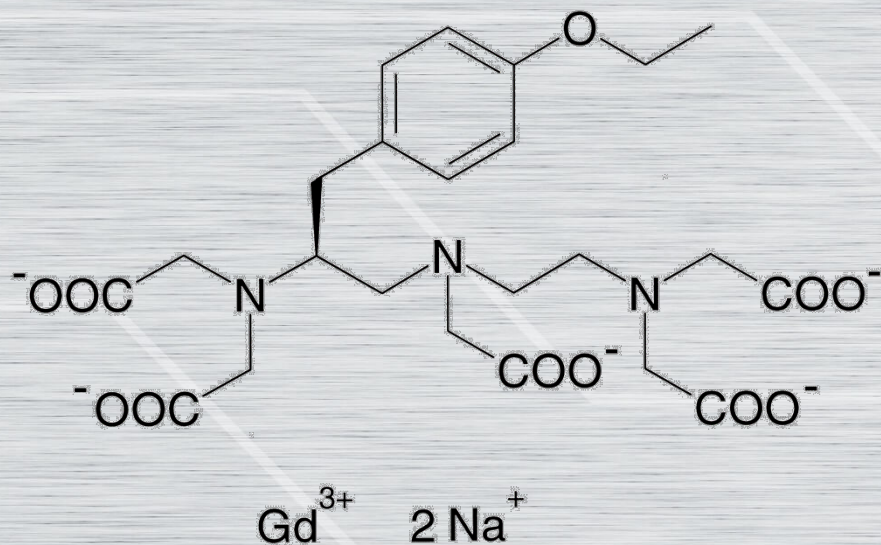


Valganciclovir  
~ 60% bioavailability

- Addition of valine moiety makes Valganciclovir a substrate of the PEPT1 uptake transporter in the gut, increasing bioavailability following oral administration

# Distribution – transporter targeting

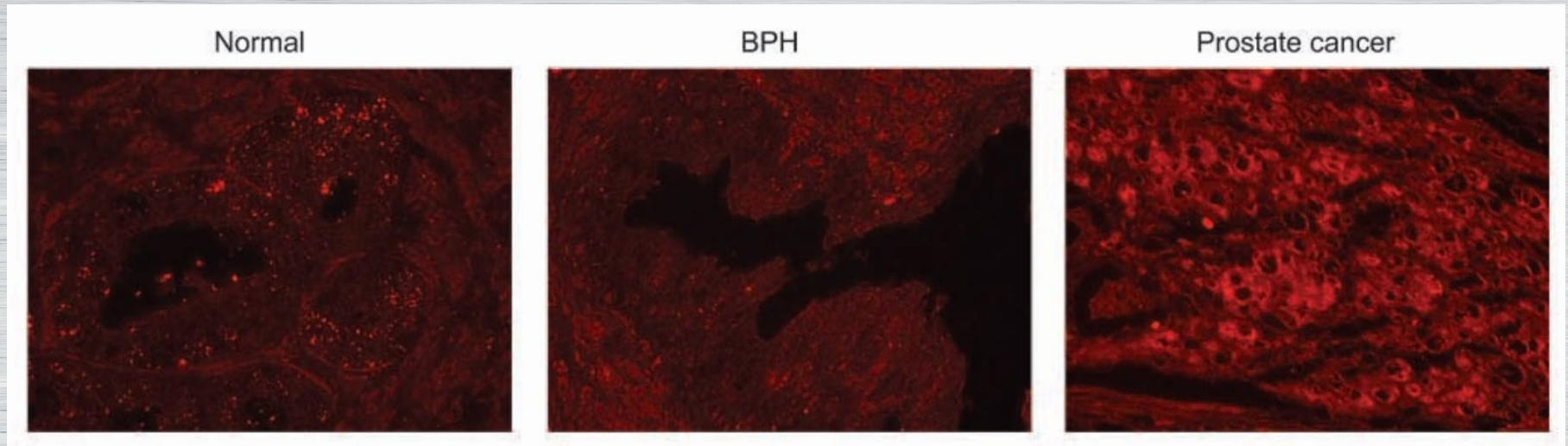
## Gadoxetate (Eovist)



- MRI contrast agent
- Used for imaging hepatocellular carcinoma and liver metastases
- Substrate for OATPIB1 and OATPIB3



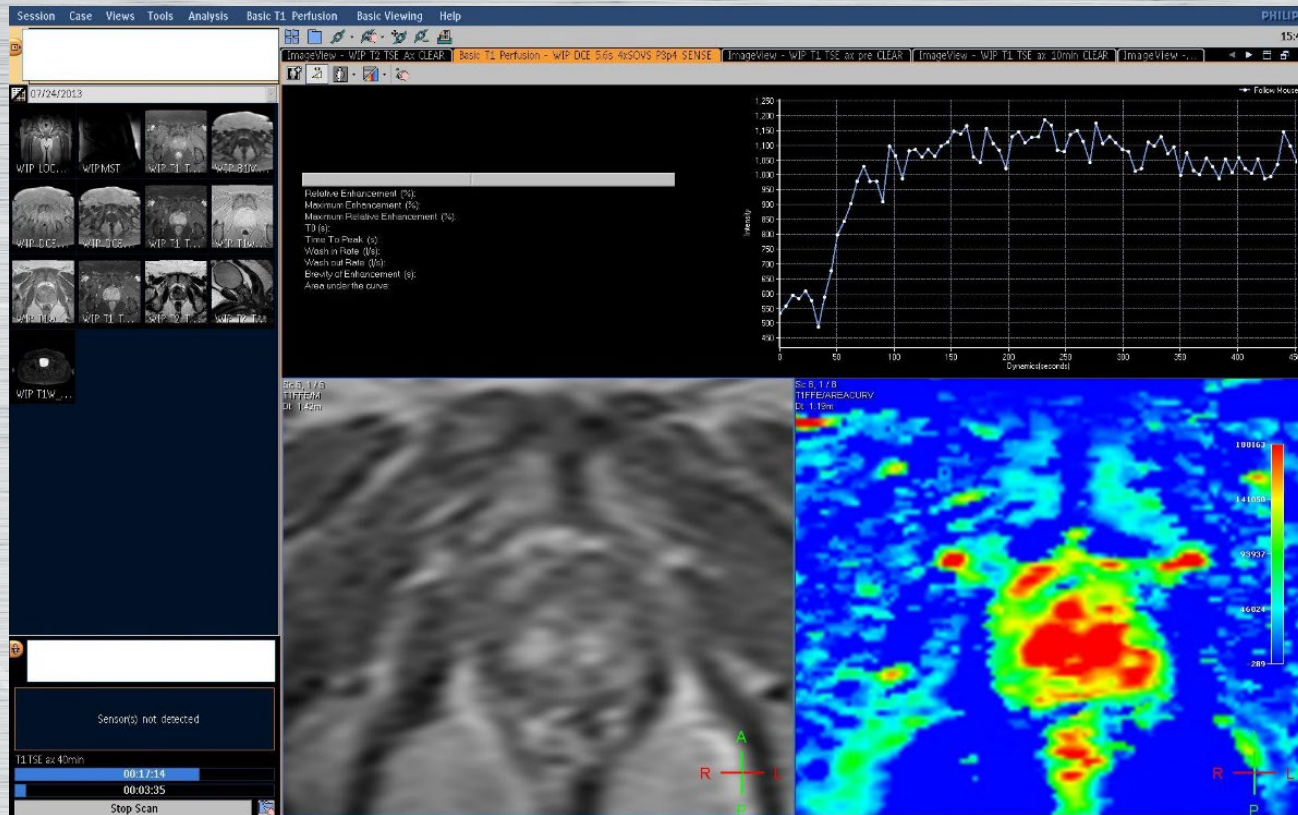
# Distribution – transporter targeting



- OATP1B3 expression is normally liver specific
- OATP1B3 expression is upregulated in prostate cancer



# Distribution – transporter targeting



- Gadoxetate (MRI contrast agent and OATP substrate)
- Accumulation of gadoxetate in prostate cancer



# Development Stage - Preclinical

- Screening for (hepato)tox
  - BSEP, (& MRP2-4, NTCP)
  - Incorporating *in vitro* data into PBPK models
- Transporter as a target
  - Substrate/inhibition assessment – ie. SGLT2 (eg. Canagliflozin/Invokana) or URATI (eg. Lesinurad/Zurampic)
- Route of administration
  - Oral - MDRI & BCRP substrate assessment
- Target biology (tissue/site of action)
  - Central nervous system? MDRI & BCRP
  - Organ targeting (ie. liver via OATPs)
- Patient population/co-meds
  - Statins? OATPs
  - Metformin? OCTs & MATEs

# Development Stage - Preclinical

- Physicochemical properties
  - BDDCS
  - ECCS
- Patient population/co-meds
  - Statins? OATPs
  - Metformin? OCTs & MATEs
- Narrow therapeutic index
  - ie. Oncology drugs
  - Substrate assessment to determine extent of DDI liability as victim



# Development Stage – Phases I & II

## Phase I

- Will have animal PK data (indication of extent of metabolism, renal/hepatic elimination)
- Little/no human PK/PD data
- Adding to transporter data information on substrate studies (test compound as victim of DDI)

## Phase II

- Will have human PK data (indication of extent of metabolism, renal/hepatic elimination)
- Additional transporter data required - Substrate studies (test compound as victim of DDI)
- Complete information on transporter DDI typically required for End-of-Ph.II meeting

# Summary

- Early transporter screening utilizing vesicle/cell-based assays useful to address specific questions
  - Tissue targeting
  - Mechanism of action
  - BSEP (& MRP) inhibition and hepatotoxicity
- Prioritization of inhibition studies for patient population with high rates of co-morbidities/co-medications, or likely transporter-interacting co-meds
- Prioritization of substrate studies for patient population with high rates of co-morbidities/co-medications, narrow therapeutic index
- Some transporter inhibition data may be required pre-IND