

In Vitro DDI Drug Transporter Studies: **Efflux and Uptake** **Transporters**

Andrew G. Taylor, Ph.D.

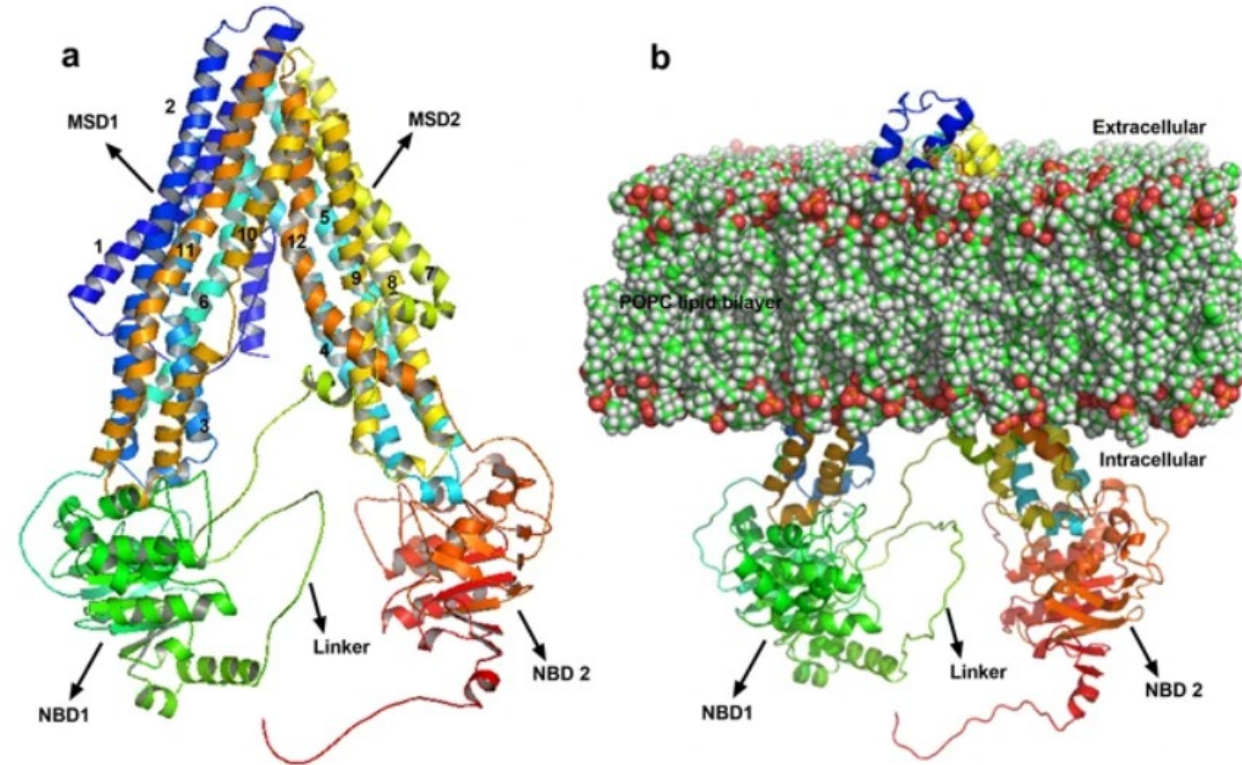
Manager, Technical Support for Services

XenoTech



What are Drug Transporters?

- Proteins bound to membranes that prevent or assist in movement of drugs in and out of cells
- Located throughout the body
 - Intestine
 - Liver
 - Kidney
 - Blood-brain barrier
- Two superfamilies
 - ATP-binding cassette (ABC) Transporters (Efflux)
 - P-gp (MDR1), BCRP, BSEP, MRPs
 - Solute carrier (SLC) Transporters (Uptake)
 - OATP1B1/3, OCT1/2, OAT1/3, MATE1/2K

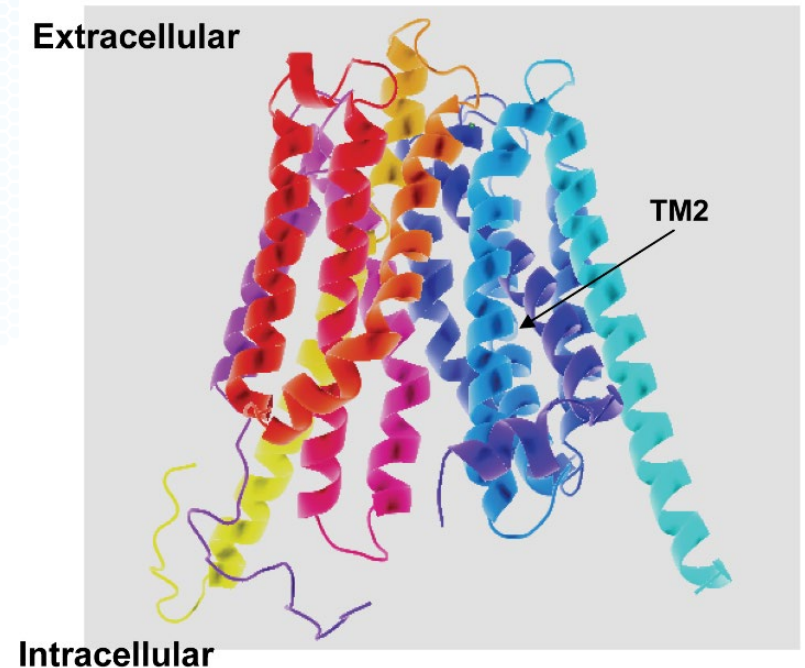


Syed, S.B., Arya, H., Fu, I. *et al.* Targeting P-glycoprotein: Investigation of piperine analogs for overcoming drug resistance in cancer. *Sci Rep* **7**, 7972 (2017). <https://doi.org/10.1038/s41598-017-08062-2>

Why are Transporters Important?

The AD&E in ADME

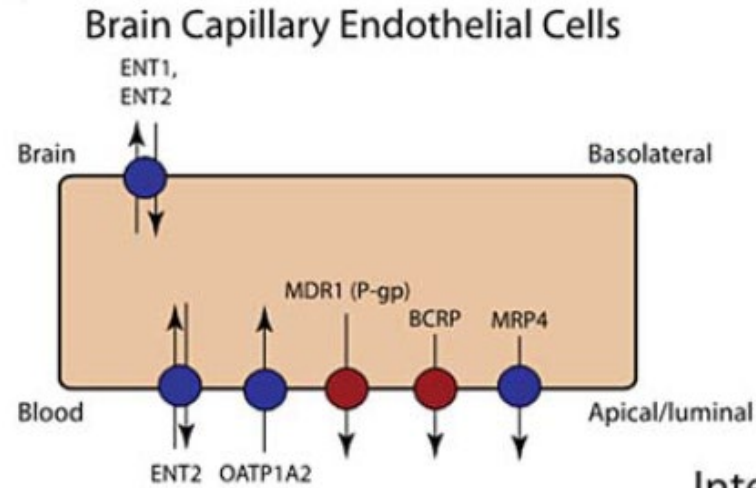
- Drug absorption, distribution, tissue-specific drug targeting, and elimination
- Drug-drug interactions
 - Clearance of transporter substrates (Victims) can be impacted by transporter inhibitors or inducers (Perpetrators)
 - Toxicity or loss of efficacy
- Real world example - Statins
 - Hepatic uptake transporter (OATPs) substrates: taken up in the liver, reduce cholesterol
 - Cyclosporine inhibits OATPs: up to 10-fold increase in statin exposure
 - Toxic side effect: rhabdomyolysis (skeletal muscles break down, cells released into bloodstream, can lead to kidney failure and possibly death)



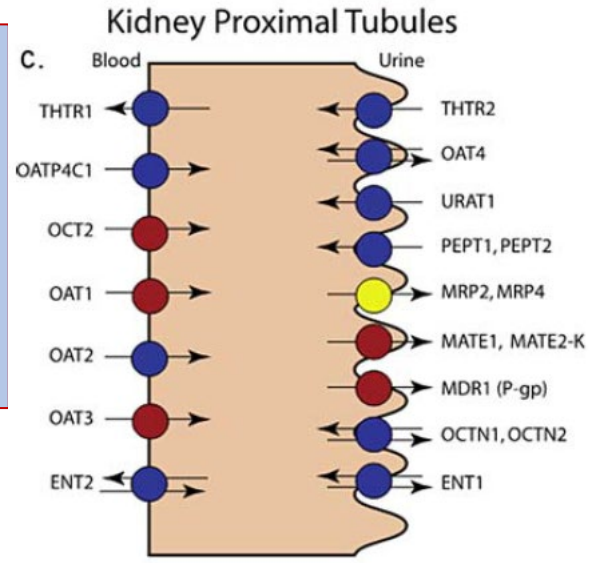
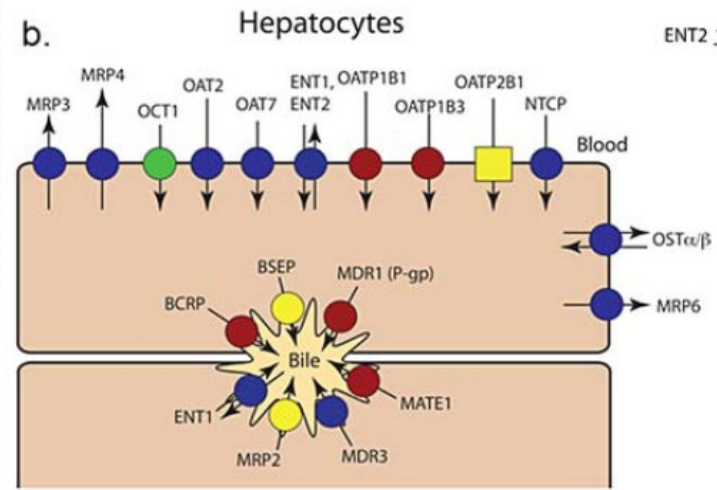
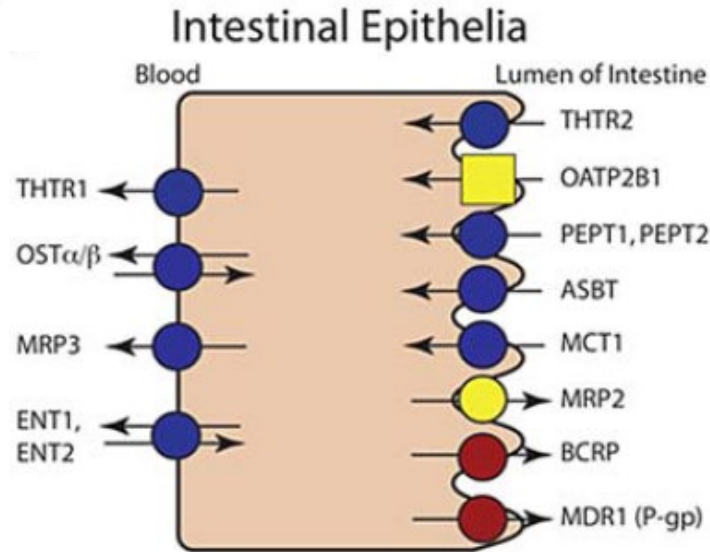
Li N, Hong W, Huang H, Lu H, Lin G, et al. (2013) Correction: Identification of Amino Acids Essential for Estrone-3-Sulfate Transport within Transmembrane Domain 2 of Organic Anion Transporting Polypeptide 1B1. PLOS ONE 8(5): 10.1371/annotation/35943d4c-3455-4818-976a-be73a2e07fff.

Transporter Expression and Function

Liver (hepatocytes) and kidney (proximal tubules): uptake transporters (SLC) remove compounds from the blood and efflux (ABC) or SLC transporters pump them into the bile or urine



Intestine and BBB: efflux (ABC) transporters predominate (particularly P-gp & BCRP) regulating absorption or entry into the brain



Figures apted from Zamek-Gliszczynski et al., Clin Pharm Ther 2018

Regulatory Guidance on Transporters

	FDA (2020)	EMA (2013)	PMDA (2014)
Recommended transporters	P-gp, BCRP OATP1B1, OATP1B3, OAT1, OAT3, OCT2, MATE1, MATE2-K	P-gp, BCRP, BSEP OATP1B1, OATP1B3, OAT1, OAT3, OCT2	P-gp, BCRP OATP1B1, OATP1B3, OAT1, OAT3, OCT2, MATE1, MATE-2K
Transporters to consider	BSEP, MRP2, OCT1 Others as necessary	OCT1, MATE1, MATE-2K Others as necessary	Others as necessary
Substrate evaluation	Yes (P-gp, BCRP) Others based on clearance/elimination routes and knowledge of drugs in same therapeutic class EMA: Two systems for P-gp, 4 conc (100-fold range)		
Inhibitor evaluation	Yes, all of the recommended transporters (EMA and PMDA: in vitro data prior to Phase III trials) EMA: Two inhibitors as positive controls		
Induction	FDA and EMA: In vivo studies Yes, for CYP3A4, PXR, CAR inducers (P-gp)		
Metabolites	≥ 25% of parent drug AUC		

General Transporter Study Design: **Inhibition**

- Solubility: Typically DMSO, 0.1% v/v (avoid 1% or more)
- Preliminary experiments
 - Cytotoxicity, stability in media, non-specific binding to plates
- Inhibition
 - Probe substrate concentration $\ll K_m$ ($IC_{50} \approx K_i$)
 - Short incubation time (in linear range with acceptable fold uptake)
 - Highest [TA] based on EMA/FDA guidance
 - Minimal design: 2 [TA] + no solvent and solvent control
 - Definitive design: 7 [TA] + no solvent and solvent control to determine IC_{50}



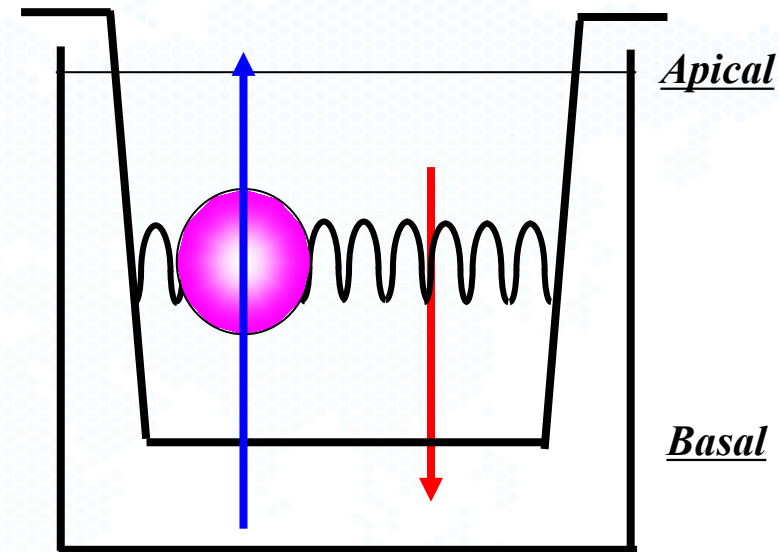
General Transporter Study Design: **Substrate**



- Substrate
 - [TA] pharmacologically relevant (too low, then BLQ; too high, then saturate)
 - Transwell
 - Minimal design: 1 [TA] , 3 times, \pm inhibitor
 - Definitive design: 4 [TA] , 3 times, \pm inhibitor and control cells at 1 [TA]
 - Uptake
 - Minimal design: 1 [TA] , 2 times, \pm inhibitor, transporter and control cells
 - Definitive design: 4 [TA] , 3 times, transporter and control cells, \pm inhibitor at 1 [TA]

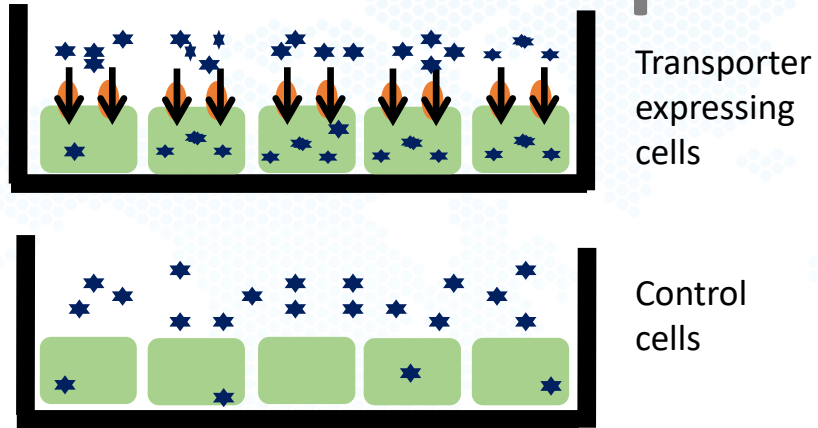
Efflux Transporter: Transwell Assays

- **Test system:** polarized cells grown on transwell plates
- P-gp and BCRP pump in basal to apical direction. Restrict permeation in apical to basal direction.
- **Substrate:** measure bidirectional permeability of test article across cells
- **Inhibition:** measure effect of test article on bidirectional permeability of probe substrate
- P-gp test systems: Caco-2 and MDCKII-MDR1
- BCRP test system: MDCKII-BCRP



Transporter	Probe substrate	Positive control inhibitors
P-gp	Digoxin	Valspodar, Verapamil
BCRP	Prazosin	Ko143, Lopinavir

SLC Transporter Uptake Assays

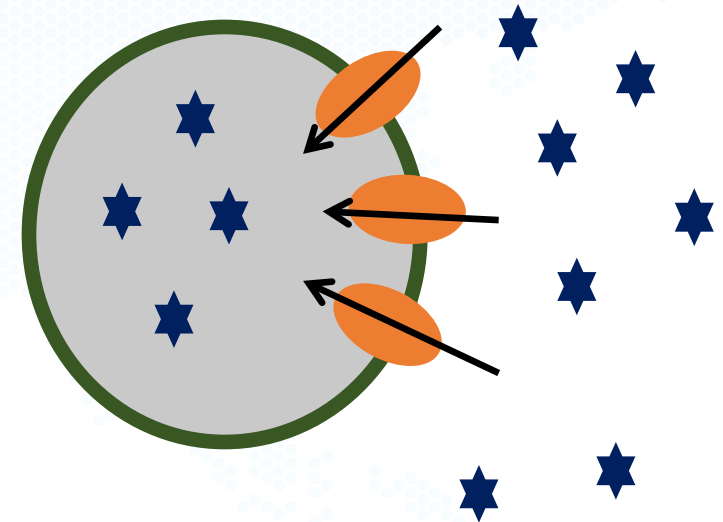


- **Test system:** transfected HEK293 cells grown in cell culture plates
- **Substrate:** measure the accumulation of the test compound in transfected and control cells
- **Inhibition:** measure the effect of the test article on the accumulation of a probe substrate

Transporter	Probe substrate	Positive control inhibitors
OATP1B1/ OATP1B3	³ H-Estradiol-17β-glucuronide	Rifampin, Cyclosporine
OAT1	³ H-p-Aminohippurate	Probenecid, Novobiocin
OAT3	³ H-Estrone-3-sulfate	Probenecid, Ibuprofen
OCT2	¹⁴ C-Metformin	Quinidine, Cimetidine
OCT1	¹⁴ C-Tetraethylammonium bromide	Quinidine, Verapamil
MATE1/2K	¹⁴ C-Metformin	Cimetidine, Pyrimethamine

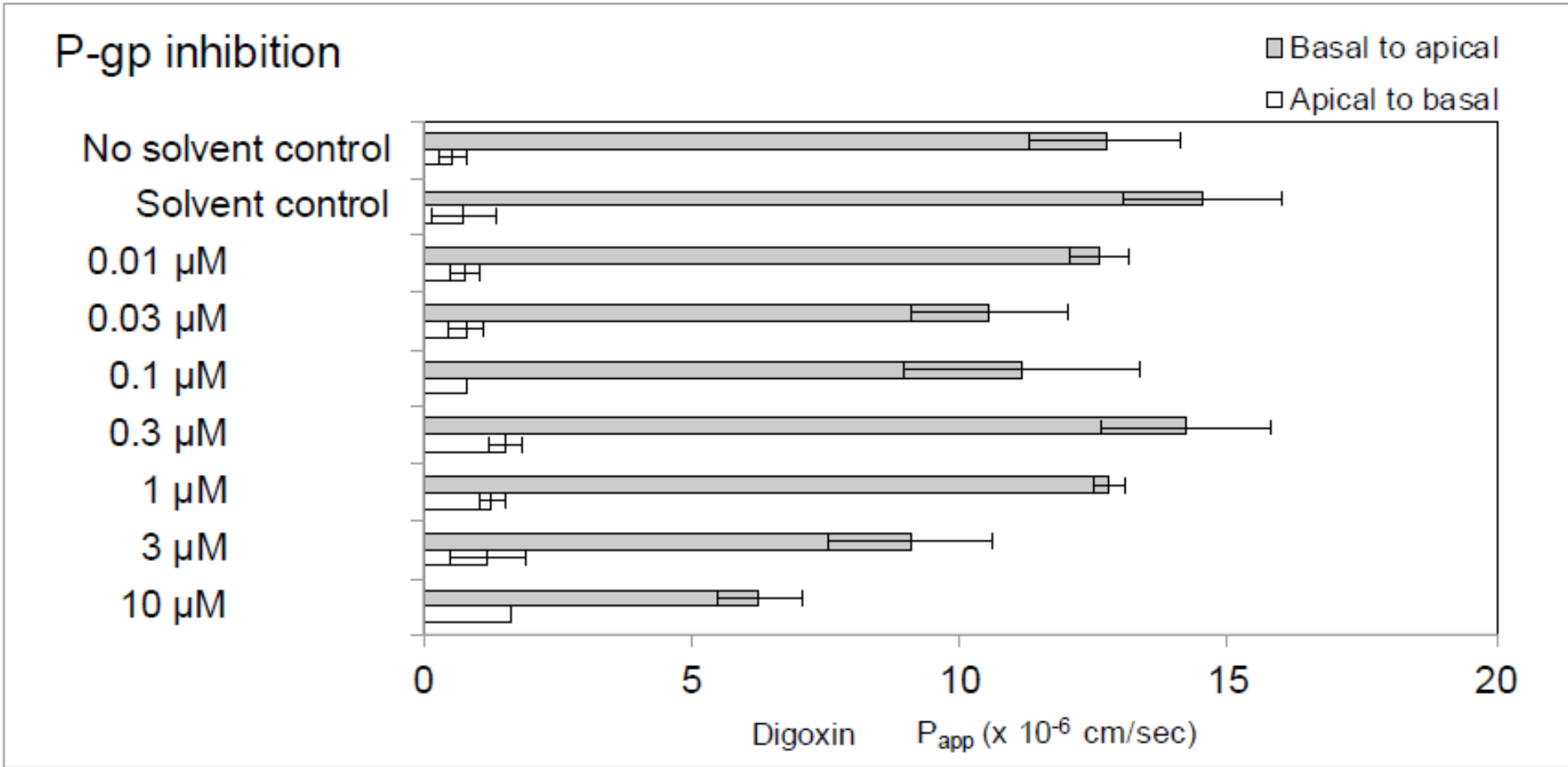
BSEP and MRP2 (Vesicle assays)

- **Test system:** Inverted plasma membrane vesicles, from cells over expressing transporter
- **Substrate:** measure accumulation of test article in transfected vesicles \pm ATP
- **Inhibition:** measure effect of test article on accumulation of probe substrate \pm ATP
- Used for other efflux transporters (e.g., P-gp, BCRP, MRP1, MRP3, MRP4)

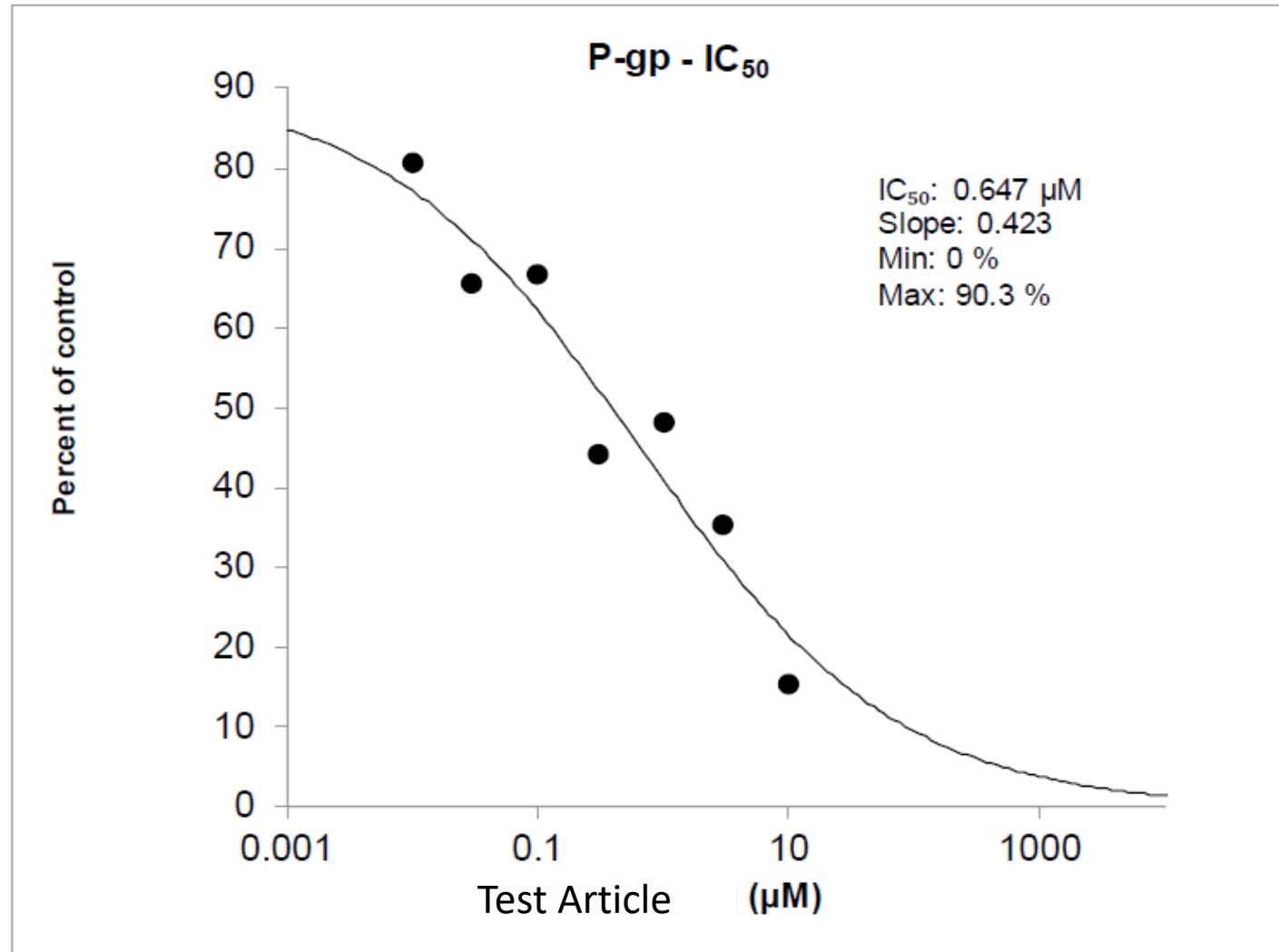


Transporter	Probe substrate	Positive control inhibitors
BSEP	³ H-Taurocholate	Cyclosporine, Glyburide
MRP2	³ H-Estradiol-17 β -glucuronide	Benzbromarone

Transporter Results Example



Transporter Results Example





Transporters Available

Description of construct	Parental Cell line	Gene	Substrates	Kinetic parameters	Kinetic parameters for inhibitors
MDR1	LLC-PK1	ABCB1	1 µM Digoxin	Km: >100 µM	IC50 Verapamil : 1.6 µM IC50 Cyclosporin A: 0.94µM
BCRP	LLC-PK1	ABCG2	10 nM Prazosin	KM: 19.6 µM	IC50 Ko143: 0.0344 µM
	Vesicle		1 µM Estrone sulfate	KM: 6.77 µM (14.2)	IC50 Ko143: 0.0186 µM (0.026)
BSEP	Vesicle	ABCB11	2 µM Taurocholic acid	KM: 19.1 µM	IC50 Cyclosporin A: 6.68 µM
MRP2	Vesicle	ABCC2	0.1 µM Leukotriene C ₄	KM: 5.79 µM (0.24-5.9)	IC50 BSP: 10.0 µM IC50 Benzbromarone: 11.1 µM (7.3-8.3)
OATP1B1	HEK	SLCO1B1	50 nM Estradiol glucuronide	KM: 3.94 µM (3.7-8.3)	IC50 Rifampicin: 0.243 µM(0.94)
OATP1B3	HEK	SLCO1B3	50 nM Estradiol glucuronide	KM: 26.5 µM (5-25)	IC50 Cyclosporin A: 0.215 µM IC50 Rifampicin: 0.388µM (1.5)
			20 nM CCK-8	KM: 18.6 µM	IC50 Cyclosporin A: 0.386 µM IC50 Rifampicin: 1.06 µM
OATP2B1	HEK	SLCO2B1	50 nM Estrone sulfate	KM: 14.5 µM (7-21)	IC50 Cyclosporin A: 0.612 µM IC50 Rifampicin: 20.6µM (90)
OAT1	HEK	SLC22A6	1000 nM p-Aminopipurate	KM: 11.9 µM (4-22)	IC50 Cyclosporin A: 1.05 µM IC50 Probenecid: 4.70 µM (6.3-12.5)
	S ₂		1000 nM p-Aminopipurate	KM: 24.6 µM (4-22)	IC50 Indomethacin: 1.95 µM IC50 Probenecid: 11.7 µM (6.3-12.5)
OAT2	S ₂	SLC22A7	50 nM PGF2α	KM: 0.137 µM (0.4-0.8)	IC50 Mefenamic acid: 11.7 µM (21.7)
OAT3	HEK	SLC22A8	50 nM Estrone sulfate	KM: 22.8 µM (2-7)	IC50 Probenecid: 2.70 µM (4-30)
	S ₂		50 nM Estrone sulfate	KM: 1.59 µM (2-7)	IC50 Probenecid: 5.33 µM (4-30)
OAT4	HEK	SLC22A11	50 nM Estrone sulfate	KM: 15.8 µM (1-10)	IC50 BSP: 0.377 µM (<5)
	S ₂		50 nM Estrone sulfate	KM: 7.42 µM (1-10)	IC50 Probenecid: 81.8 µM (25.4)
OAT7	HEK	SLC22A9	50 nM Estrone sulfate	KM: 12.9 µM (8.7-40)	IC50 BSP: 0.121 µM
OCT1	HEK	SLC22A1	5 µM Tetraethylammonium bromide	KM: 576 µM (160-1600)	IC50 Quinidine: 4.57 µM (5.4-23.4)
	S ₂		10 µM Metformin	KM: 899 µM	IC50 Verapamil: 0.272 µM IC50 Quinidine: 13.8 µM
OCT2	HEK	SLC22A2	5 µM Tetraethylammonium bromide	KM: 453 µM (160-1600)	IC50 Quinidine: 17.9 µM (5.4-23.4)
			10 µM Metformin	KM: 3600 µM (990)	IC50 Quinidine: 88.6 µM (1 mM 5-20%)
OCT3	HEK	SLC22A3	5 µM Tetraethylammonium bromide	KM: 723 µM	IC50 Quinidine: 93.7 µM
			10 µM Metformin	KM: 12500 µM (990)	IC50 Verapamil : 192 µM IC50 Quinidine: 104 µM
OCTN1	S ₂	SLC22A3	5 µM Tetraethylammonium bromide	KM: 35.1 µM	IC50 Verapamil : 298 µM IC50 Quinidine : 12.7 µM
			5 µM Histamine	KM: 26.1 µM(180-420)	IC50 Verapamil: 11.6 µM (24)
OCTN2	S ₂	SLC22A4	5 µM Tetraethylammonium bromide	KM: 315 µM (199-1280)	IC50 Verapamil: 12.3 µM (8.4)
NTPC1	HEK	SLC10A1	30 nM Carnitine	KM: 4.73 µM (2-66)	IC50 Verapamil: 5.19 µM (500 µM 2.4%)
NTCP2 (ASBT)	HEK	SLC10A2	50 nM Taurocholate	KM: 10.2 µM (6.2-10)	IC50 Cyclosporin A: 1.20 µM (1)
	HEK	SLC10A2	200 nM Taurocholate	KM: 15.8 µM (9.4-20)	IC50 Chenodeoxycholic acid: 3.48 µM (Ki 3.3)
NPT1	HEK	SLC17A1	50 nM Estradiol glucuronide	KM: 6.88 µM	IC50 Indomethacin: 2.17 µM IC50 Probenecid: 69.1 µM
PEPT1	HEK	SLC15A1	100 nM Glycylsarcosine	KM: 556 µM (290-3130)	IC50 Cephalixin: 1460 µM (5200-13700)
PEPT2	HEK	SLC15A2	100 nM Glycylsarcosine	KM: 59.4 µM (74)	IC50 Cephalixin: 124 µM (10mM 0%)
URAT1	HEK	SLC22A12	50 µM Uric acid	KM: 1050 µM (371)	IC50 Benzbromarone: 0.0334 µM (0.3)
MATE1	HEK	SLC47A1	10 µM Metformin	KM: 73.4 µM (220)	IC50 Probenecid: 80.0 µM IC50 Cimetidine: 1.32 µM (<20)
			5 µM Tetraethylammonium bromide	KM: 29.9 µM	IC50 Cimetidine : 1.10 µM IC50 Quinidine : 1.40 µM
MATE2-K	HEK (Transient)	SLC47A2	10 µM Metformin	KM: 252 µM	IC50 Cimetidine: 4.95 µM IC50 Cimetidine : 36.6 µM
			5 µM Tetraethylammonium bromide	KM: 109 µM	IC50 Cimetidine : 36.6 µM IC50 Quinidine : 16.3 µM
rUrat1	HEK	Slc22a12	Uric acid	KM: 677 µM	IC50 Quinidine : 16.3 µM IC50 Benzbromarone : 1.88 µM
OSTα/β	HEK	SLC51	Taurocholic acid	KM: 7960 µM	IC50 Indomethacin : 262 µM

P-gp (MDR1)	LLC-PK1, MDCKII, Caco-2	✓			
MRP1		✓			
MRP2		✓	OATP1B1		
MRP3		✓			
MRP4		✓			
BCRP	LLC-PK1, MDCKII, Caco-2	✓			
BSEP		✓			
OATP1A2				✓	
OATP1B1	HEK293		MRP2	✓	✓
OATP1B3	HEK293			✓	✓
OATP2B1	HEK293			✓	
OAT1	HEK293, S ₂			✓	
OAT2	S ₂			✓	
OAT3	HEK293, S ₂			✓	
OAT4	HEK293, S ₂			✓	
OAT7	HEK293			✓	
OCT1	HEK293, S ₂			✓	
OCT2	HEK293, S ₂			✓	
OCT3	HEK293, S ₂			✓	
OCTN1	S ₂			✓	
OCTN2	S ₂			✓	
OST α/β	HEK293			✓	
PEPT1	HEK293			✓	
PEPT2	HEK293			✓	
URAT1	HEK293			✓	
NTCP1	HEK293			✓	✓
NTCP2 (ASBT)	HEK293			✓	
MATE1	HEK293			✓	
MATE2-K	HEK293 (TEA & Metformin)			✓	
NPT1	HEK293			✓	
ENT1	MDCKII			✓	
ENT2	MDCKII			✓	
ENT4	MDCKII			✓	
CNT1	MDCKII			✓	
CNT2	MDCKII			✓	
CNT3	MDCKII			✓	
GLUT1				✓	
GLUT2				✓	
Mouse Mdr1a	LLC-PK1				KO animal
Mouse Mdr1b	LLC-PK1				KO animal
Rat Mdr1a	LLC-PK1				KO animal
Rat Mdr1b	LLC-PK1				
Rat Urat1	HEK293			✓	
Rat Oatp4/Oatp1b2				✓	
Dog Mdr1	LLC-PK1				

XT Products (Transporters)

Primary Human Hepatocytes

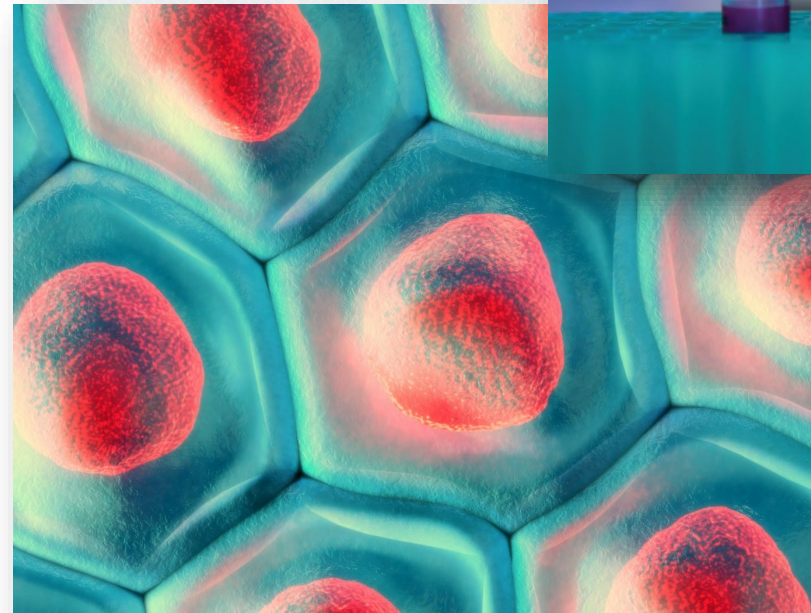
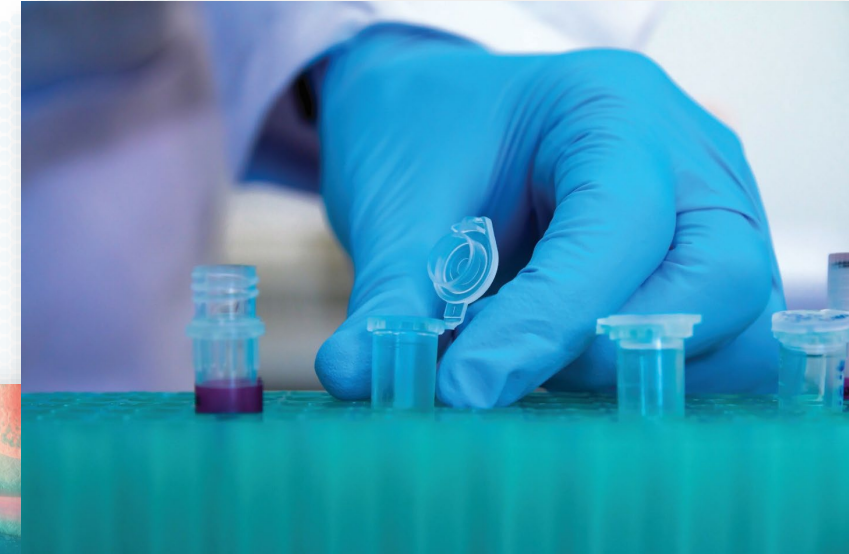
- Suspension (uptake)
 - Some lots with uptake characterization
 - Individual donor and Pooled CryostaX
- Plated
 - CryostaX (plated uptake)
 - No individual lots are characterized, but no reason not to believe that they do not maintain uptake activity.
 - No Efflux characterized products

Primary Animal Hepatocytes

- No lot characterization, but have R&D data to show that our cell preparation protocols maintain uptake activity (rat and mouse)
 - Suspension
 - Plated

Support Reagents

- PHH
 - OptiTHAW
 - OptiPLATE
 - OptiCULTURE
 - OptiINCUBATE



Thank you for watching!

Questions? Get in touch through the **Contact Us** tab on our website

Please contact your regional account manager if you are interested in a placing a contracted drug transporter study or have interest in high-quality test systems for your assays

