

TRANSPORTER INDUCED PROTEIN BINDING SHIFT (TIPBS): IMPACT ON OATP1B1/1B3 MEDIATED DRUG TRANSPORT AND DRUG-DRUG INTERACTIONS

Xuexiang Zhang, Jason Baik, Mirza Jahic, Wenjie Jiang and Yong Huang*



* Correspondence email: yhuang@optiviabio.com

ABSTRACT

BACKGROUND: We recently proposed a Transporter-Induced Protein Binding Shift (TIPBS) hypothesis to describe the effects of serum proteins on transporter-mediated drug transport. This work substantiates our previous findings and theory by demonstrating drug-dependent discrepancies between predicted and measured OATP1B1/1B3 mediated transport and inhibition in human serum.

METHODS: OATP1B1 and OATP1B3 mediated transport of highly protein-bound substrates, and rifampicin OATP1B1 IC₅₀s were measured in protein-free HBSS and 100% human serum, using CHO cells stably expressing the transporters. Serum unbound fraction (fu) was used to calculate substrate transport or inhibitor IC₅₀s in serum from the constants measured in HBSS. The predicted values of OATP mediated transport and its inhibition in human serum using fu adjustment method were contrasted to that measured from assays conducted in human serum.

RESULTS: The fu adjustment method generally under-estimated substrate uptake and inhibitor potency in serum. The extent of under-estimation appeared to be compound and transporter dependent. For examples, actual OATP1B1 mediated uptake of atorvastatin, pitavastatin, simvastatin and E17bG in serum is significantly higher than that of predicted values based on measured uptake in HBSS and fu; while such difference was not observed with fluvastatin and lovastatin. In terms of rifampicin inhibition of OATP1B1 mediated transport of these substrates, measured IC₅₀s in serum were generally significantly lower than that of predicted values based on measured rifampicin IC₅₀s in HBSS and fu, however, the extent of discrepancies was substrate dependent. On various substrates and inhibitors with different fu indicated that the extent of underestimation were drug AND transporter dependent, possibly due to difference in drug binding affinities to serum proteins and transporters as predicted by our TIPBS models.

CONCLUSION: Our work further confirms previously reported experimental observations and theoretical modeling that suggest drug transport and inhibition in serum may not be described by the simple unbound drug model, raising the question on whether/when is appropriate to use the conventional fu adjustment method for predicting in vivo drug clearance and DDIs.

INTRODUCTION

Transporters play a crucial role in drug clearance and drug-drug interactions (DDI). Currently *in vitro* cell-based assays are conducted to predict potential clinical DDI. Typically, in these assays transporter kinetic parameters (V_{max}, K_m, and K_i/IC₅₀) are measured in protein free buffers (e.g. HBSS without albumin). However, the presence of proteins in serum may affect significantly the free concentration of drugs. Thus conventional "equilibrium model" is applied by applying drug unbound fraction at equilibrium (fu), in predicting *in vivo* drug clearance and DDIs. This approach assumes that drugs are in binding equilibrium with serum proteins and only the free fraction (fu) is transported or has inhibitor effect.

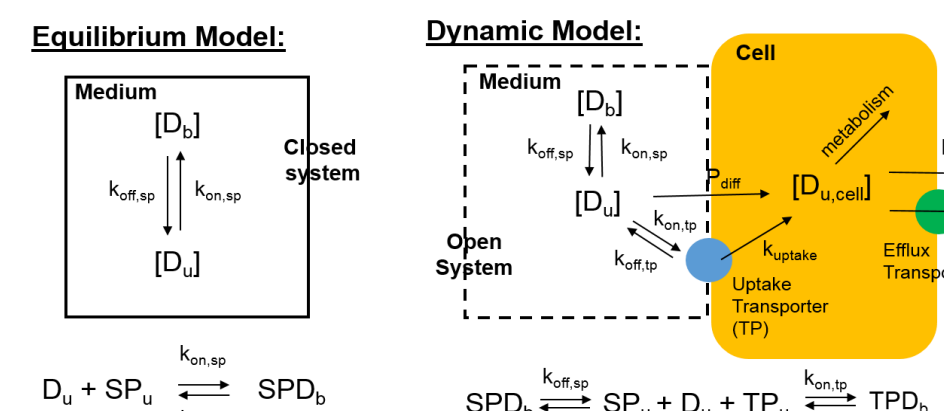


Figure 1. Schematic illustration of the conventional rapid equilibrium model (left) whereas [D_u] = [D_{total}] × fu, whereas $f_u = \frac{1}{1 + [A]/K_d}$, and our dynamic TIPBS model (right) in which [D_u] at cell surface may not be determined by the fu determined under equilibrium conditions

We have previously shown for certain highly protein bound transporter substrates and inhibitors, the apparent transport rate or inhibition potency measured in 100% human serum were significantly higher than those calculated based on equilibrium fu and the intrinsic transporter IC₅₀s assessed in HBSS [1]. We further hypothesized that such discrepancies may arise from transporter induced protein binding shift (TIPBS) for compounds which exhibit higher binding affinities toward transporters than serum binding proteins; thus the transporters may pull compounds away from binding proteins then clear them out, creating a "sink" condition under which binding equilibrium may not be reached (Open system) [2]. Here we aim to further interrogate experimentally TIPBS effects on OATP1B1/1B3 mediated transport of highly protein bound statins and its inhibition by Rifampicin. The extent of shift in active uptake is correlated with estimated compound binding affinities toward the transporter and serum albumin, to interrogate our mathematical modeling that suggests binding affinity difference is a major cause of TIPBS.

MATERIALS AND METHODS

CELL CULTURE: CHO cells stably expressing human OATP1B1 and OATP1B3 transporters were maintained in Dulbecco's modified Eagle's medium (DMEM) supplemented with 10% fetal bovine serum, 100 U/mL penicillin, 100 µg/mL streptomycin and 500 µg/mL G418 at 37 °C in a humidified atmosphere with 5% CO₂. Approximately 48 hr before assay, cells were seeded at 20K per well on Corning BioCoat™ 96-well flat bottom plates in medium without G418. CHO parental cells were also used as a control.

TRANSPORT ASSAY: Briefly, cells were washed and pre-incubated with HBSS for 15 min at 37°C. Radioisotope-labelled statins (atorvastatin, pitavastatin, lovastatin, simvastatin and fluvastatin) in HBSS or human serum were incubated with cells for 2 min based on the previously determined linear range for each statin. At the end of dosing, statins were removed and ice-cold PBS was used to wash the cells. The intracellular contents of statins were extracted by adding 50% acetonitrile in water and quantified by radiometry. Net transporter-mediated uptake of statins was calculated by subtracting uptake in the CHO-parental cells from uptake in the CHO-OATP1B1 and CHO-OATP1B3 cells. To determine the intrinsic K_m and V_{max} values for each statin as a substrate of OATP1B1 and OATP1B3, concentrations of up to 100 µM in HBSS were tested for each statin as a free monomer. To evaluate the IC₅₀ of rifampicin against the transport of each statin, a series concentrations of rifampicin (0.1-100 µM) was mixed with a fixed concentration of each statin tested at 100nM in HBSS and 10µM in serum (except for Atorvastatin tested at 5µM). The mixture was in either HBSS buffer or human serum. To provide better direct comparisons, transport assays in HBSS and human serum were conducted side-by-side on same plate. Actual concentrations of inhibitor Rifampicin was quantified by LC-MS/MS and used in calculation of IC₅₀ values by non-linear regression using GraphPad Prism.

DETERMINATION OF DRUG SERUM UNBOUND FRACTION F_u USING RAPID EQUILIBRIUM DIALYSIS: Rapid Equilibrium Dialysis (RED) kit (Thermo Scientific, Rockford, IL) was used to estimate the extent of statins and rifampicin binds to serum proteins. A series concentrations of drugs were prepared in 100% human serum and 200 µL was loaded the sample chamber. 350 µL PBS was added to the buffer chamber. The device was sealed and incubated at 37°C on an orbital shaker at approximately 250 rpm for 4 hr. 50 µL samples was taken each chamber and measured by either scintillation counter or LC-MS/MS. The fraction unbound was calculated by dividing the concentration in the buffer chamber by the concentration in corresponding serum sample chamber.

F_u ADJUSTMENT METHOD FOR PREDICTING SUBSTRATE TRANSPORT AND INHIBITOR POTENCY IN SERUM: To distinguish transport and inhibition constants in HBSS and serum, we hereby denote constants measured in protein-free HBSS as *intrinsic* as they reflect the "true" kinetic properties without influence from protein binding. The fu adjustment method as defined by the following equations are used to predict substrate transport and apparent inhibitor IC₅₀ in serum.

$$\text{Substrate: } V_{\text{serum}} = \frac{V_{\text{max}}}{K_m} \times C_{\text{total}} \times f_u, \text{ assuming } C_{\text{total}} \times f_u \ll K_m \quad \text{Inhibitor: } IC_{50, \text{serum}} = IC_{50} / f_u$$

F_{TIPBS} AND EXTENT OF TIPBS EFFECT: We define F_{TIPBS} as the ratio of measured and predicted transport rate or inhibition potency in serum. Larger F_{TIPBS} value suggests more profound shift in serum protein binding.

$$\text{Substrate: } F_{\text{TIPBS}} = V_{\text{serum, measured}} / V_{\text{serum, predicted}} \quad \text{Inhibitor: } F_{\text{TIPBS}} = IC_{50, \text{serum, predicted}} / IC_{50, \text{serum, measured}}$$

ESTIMATION OF COMPOUND DISSOCIATION CONSTANT K_d FOR TRANSPORTER AND ALBUMIN: For simplicity of calculation, it is assumed that albumin is the dominant serum binding protein and serum albumin concentration is 660µM. Albumin K_d is calculated from measured equilibrium fu. Transporter K_d is approximated by K_m assuming K_{off,tp} >> k_{uptake} (Fig. 1).

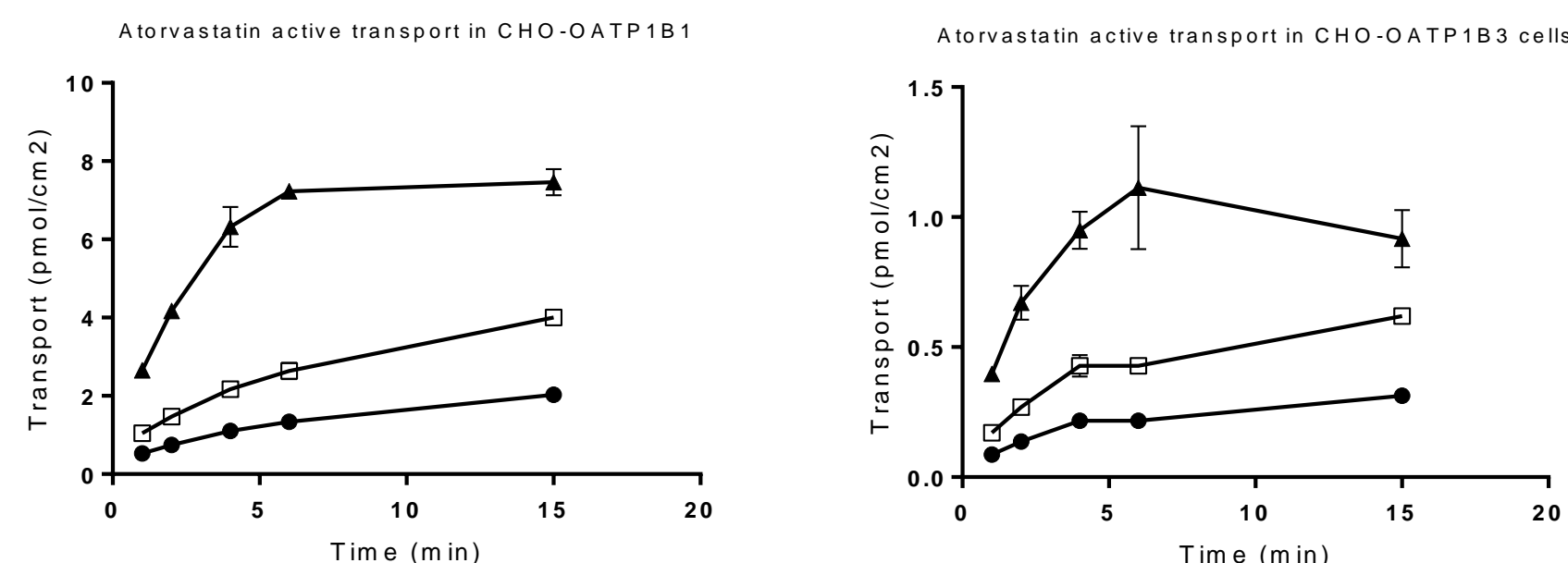


Figure 2. Time-dependent active transport of 100nM Atorvastatin in HBSS (solid circle) and 5µM Atorvastatin in 100% human serum (measured fu 3.95% mediated by OATP1B1 (left) and OATP1B3 (right)). Measured Atorvastatin active uptake (solid triangle) in serum is significantly higher than that of predicted (open square).

RESULTS, DISCUSSION AND CONCLUSIONS

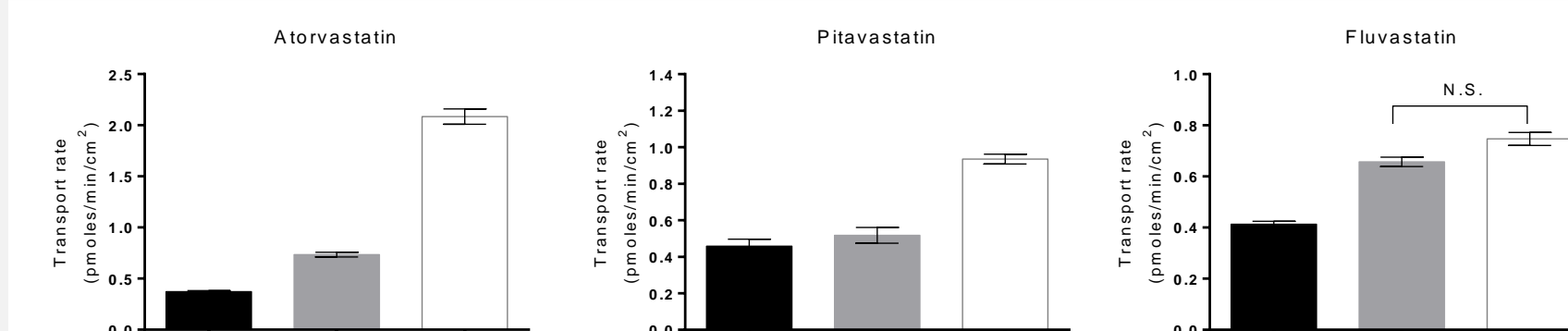


Figure 3. Representative measured active transport rate of 100nM statins in HBSS (black), predicted (grey) and actual (open) uptake rate of 5µM Atorvastatin, 10µM Pitavastatin and Fluvastatin, in 100% human serum

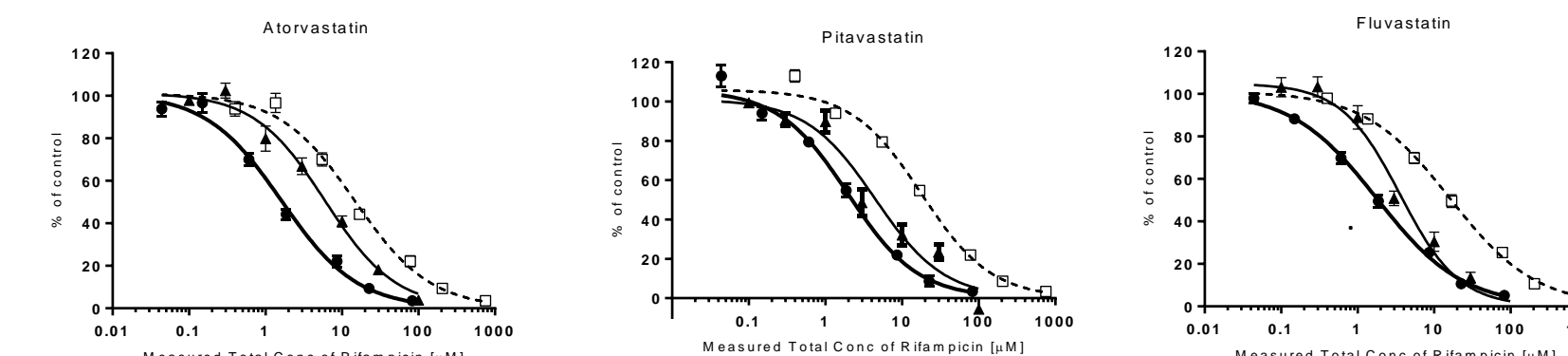


Figure 4. Representative dose-dependent inhibition of OATP1B1 mediated transport of different statins by Rifampicin. Solid circle: in HBSS, open square: predicted in serum, solid triangle: measured in serum

OATP1B1 Substrate	Meas'd f _u	Est. AB K _d (µM)	Meas'd OATP1B1 K _m (µM)	Affinity ratio (K _d /K _m)	Calc'd F _{TIPBS}	OATP1B1 Substrate	Meas'd HBSS IC ₅₀ (µM)	Meas'd HS IC ₅₀ (µM)	Pred'd HS IC ₅₀ (µM)	Calc'd F _{TIPBS}
Atorvastatin	3.95%	27.1	1.9	14.3	2.84	Atorvastatin	1.6	6.0	14.8	2.5
Pitavastatin	1.13%	7.5	1.7	4.5	1.81	Pitavastatin	2.0	4.4	18.0	4.1
Simvastatin	3.17%	21.6	6.0	3.6	1.60	Simvastatin	3.3	7.6	30.0	3.9
Lovastatin	5.92%	41.5	9.8	4.3	1.19 (ns)	Lovastatin	1.5	5.3	13.6	2.6
Fluvastatin	1.59%	10.7	10.6	1.0	1.14 (ns)	Fluvastatin	1.8	3.6	16.3	4.5
E17βG	2.62%	17.8	3.0	5.9	2.2	E17βG	0.6	1.7	5.5	3.2

Table 1. Summary of protein binding, active transport constants and TIPBS effect of various statins and reference OATP1B1 substrate E17βG (ns: no statistical difference)

Table 2. Summary of measured and predicted Rifampicin IC₅₀s of different substrates in HBSS and human serum (measured rifampicin fu 11%)

The presented *in vitro* data substantiate our previous experimental findings and TIPBS hypothesis, by demonstrating that both OATP1B1 and OATP1B3 mediated transport of highly protein bound substrates in human serum, and their inhibition by Rifampicin, can be under-estimated using the conventional fu adjustment approach, which can attribute to the reported compound-dependent scaling factors required in predicting *in vivo* hepatic clearance of these drugs and DDIs. Moreover, except for Lovastatin, there is a positive correlation between the albumin/OATP1B1 binding affinity difference and the extent of TIPBS effect (Table 1), supporting our previous hypothesis, based on mathematical simulations using a dynamic binding model [2], that high protein binding itself may not be sufficient to indicate whether a compound is subject to binding shift due to its interactions with transporters. It is intriguing that although the fu adjustment method led to underestimation of Rifampicin inhibition potencies in human serum for all OATP1B1 substrates, it is not clear when the extent of underestimation is substrate dependent.

Our current and previous experimental studies and theoretical modeling collectively suggest that high affinity binding to a transporter by its substrate or inhibitor may displace the compound's binding to serum binding proteins, hence, the conventional approach of using equilibrium unbound fraction fu may not be valid in predicting *in vivo* clearance and DDI for such compound. We therefore call for more research on this subject in order to better understand and predict drug transport in complex, protein-rich biological matrices.

REFERENCES

- M. Jahic, J. Baik, et. al., the 13th EU ISSX, P164, Glasgow, 2015
- J. Baik and Y. Huang, the 20th Intl. ISSX meeting, P272, Orlando, 2015