112

# Investigation of inhibitor preincubation condition on human OATP1B1, P-gp and BCRP transporter *in vitro* inhibitory potencies



Hayley Atkinson, Valentin Clements, Jennifer Meadows, Sam Outteridge, Sarah Pickering, Amy Radcliffe and Robert Elsby Cyprotex, No 24 Mereside, Alderley Park, Macclesfield, Cheshire, United Kingdom

# Background

Previously it has been demonstrated that preincubation with inhibitor enhances in vitro determined inhibitory potency versus OATP1B1 and that this phenomenon likely reflects a requirement of time for specific inhibitors to accumulate to a sufficient intracellular concentration in order to exert trans-inhibition from inside the cell, alongside the normal cis-inhibition from outside [1,2]. Furthermore the 2017 FDA In vitro metabolism and transporter mediated drug-drug interaction studies guidance for industry advises that a 30 minute preincubation with test compound is included in the experimental set-up for OATP1B1 and OATP1B3 inhibition assessment [3].

## Purpose

To evaluate whether the determined in vitro  $IC_{50}$  versus OATP1B1 following a 15 min preincubation step with inhibitor is comparable to that determined following a 30 min preincubation step with inhibitor [1], and determine whether inhibitor preincubation impacts on in vitro  $IC_{50}$  values determined in polarised cell monolayers versus P-gp and BCRP.

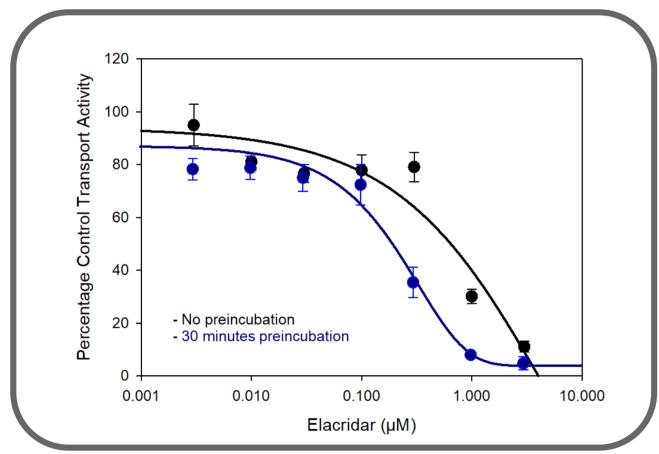


Figure 1: Comparison of in vitro determined  $IC_{50}$  values against P-gp-mediated [ $^{3}H$ ]digoxin ( $^{5}\mu$ M) transport following a 30 min preincubation without inhibitor (vehicle buffer) and a 30 min preincubation with elacridar

Table 1: Inhibitory potency (IC $_{50}$ ) determined against OATP1B1-mediated transport of [ $^{3}$ H]estradiol glucuronide (0.02 $\mu$ M) with inhibitor preincubation for 15 and 30 min or 15 min preincubation without inhibitor (vehicle buffer)

Compound		Fold change (none vs 15		
	No Preincubation (vehicle buffer only)	Preincubation 15 minutes	Preincubation 30 minutes	minutes preincubation)
Cyclosporin A	0.699 ± 0.461	0.372 ± 0.0899	0.354 ± 0.0235	1.78
Atorvastatin	0.473 ± 0.165	0.235 ± 0.0597	0.248 ± 0.0562	2.00

Table 2: Inhibitory potency (IC<sub>50</sub>) determined against BCRP-mediated transport of [<sup>3</sup>H]estrone 3-sulfate (1 μM) with inhibitor preincubation (30 min) or 30 min preincubation without inhibitor (vehicle buffer)

	BCRP IC		
Compound	No Pre-incubation (vehicle buffer only)	Preincubation	Fold change
Novobiocin	2.01 ± 0.722	2.06 ± 0.884	1.01
Fumitremorgin C	0.273 ± 0.0711	0.250 ± 0.0540	1.09

Table 3: Inhibitory potency (IC<sub>50</sub>) determined against P-gp-mediated transport of [ $^{3}$ H]digoxin (5  $\mu$ M) with inhibitor preincubation (30 min) or 30 min preincubation without inhibitor (vehicle buffer). (NS - not significant)

	P-gp IC50 (μM)			
Compound	No Pre-incubation (vehicle buffer only)	Preincubation	Fold change	
Ketoconazole	14.9 ± 5.20	8.83 ± 4.09	1.86 (NS)	
Cyclosporin A	1.60 ± 0.330	0.931 ± 0.0574	1.71 (NS)	
Verapamil	78.4 ± 15.0	54.7 ± 10.3	1.44 (NS)	
Elacridar	0.814 ± 0.0427	0.284 ± 0.0452	2.92 ( <i>p</i> ≤0.001)	

#### Methods

Inhibition of OATP1B1-mediated [3H]estradiol glucuronide transport

HEK293 Corning® TransportoCells<sup>TM</sup> overexpressing OATP1B1 and corresponding vector control cells were seeded onto 24-well poly-D-lysine coated plates at 3 x 10<sup>5</sup> cells per well and cultured for 24 hours. All incubations were carried out in uptake buffer (HBSS containing 10mM HEPES, pH 7.4) at 37<sup>o</sup>C.

Investigations compared in vitro inhibitory potential ( $IC_{50}$ ) determined following either a 15 min preincubation step with vehicle-containing (1 % v/v DMSO) buffer only, a 15 min preincubation with inhibitor or a 30 min preincubation with inhibitor. Following this, uptake of probe substrate [ $^3$ H]estradiol glucuronide (0.02  $\mu$ M) was determined over 2 min in the absence and presence of either cyclosporin A (0.01-10  $\mu$ M) or atorvastatin (0.01-30  $\mu$ M). Each set of incubation conditions utilised the same inhibitor solutions and were performed using triplicate wells per inhibitor concentration over 3 experimental occasions (n=9).

Uptake of [³H]estradiol glucuronide was terminated by washing the cells twice in ice cold uptake buffer, cells were lysed using M-PER reagent and the amount of radioactivity inside the cells determined by liquid scintillation counting. In parallel, protein concentration for each well at each experimental condition were determined using the BCA method. Uptake (pmol/mg) in control cells was subtracted from that determined in OATP1B1 overexpressing cells to give corrected uptake. Corrected uptake was converted to a percentage of vehicle control uptake, plotted against inhibitor concentration and the resulting curve fitted to determine inhibitory potency (IC<sub>50</sub>).

Inhibition of P-gp-mediated [3H]digoxin transport and BCRP-mediated [3H]estrone 3-sulfate transport

MDCK-MDR1 (for P-gp inhibition studies) and Caco-2 cells (for BCRP inhibition studies) were seeded onto Multiscreen<sup>™</sup> plates (Millipore, MA, USA) at a density of 3.4 x 10<sup>5</sup> and 1 x 10<sup>5</sup> cells/cm<sup>2</sup> and cultured for 4 and 20 days, respectively. All incubations were carried out in transport buffer (HBSS containing 25mM HEPES, 4.45mM glucose, pH 7.4) at 37°C.

Unidirectional (basolateral-to-apical) flux of probe substrate [ $^3$ H]digoxin (5  $\mu$ M), or [ $^3$ H]estrone sulfate (1  $\mu$ M), was assessed across polarised MDCK-MDR1, or Caco-2 cells, respectively. Incubations (90 min) were performed in the absence or presence of inhibitor, following a 30 min preincubation step with either inhibitor or vehicle-containing (1 % v/v DMSO) buffer alone. Inhibitors included cyclosporin A (0.01-10  $\mu$ M), elacridar (0.003-3  $\mu$ M), ketoconazole (0.1-100  $\mu$ M) or verapamil (0.1-100  $\mu$ M) for P-gp and novobiocin (0.1-100  $\mu$ M) and fumitremorgin C (0.01-10  $\mu$ M) for BCRP. Each set of incubation conditions utilised the same inhibitor solutions and were performed using triplicate wells per inhibitor concentration over 3 experimental occasions (n=9).

Vectorial flux of probe substrate was measured by monitoring its appearance on the apical side of the membrane by liquid scintillation counting and was used to determine apparent permeability ( $P_{app}$ ) in the absence and presence of inhibitor. The passive  $P_{app}$  of probe substrate (observed when P-gp/BCRP is completely inhibited) was subtracted from all  $P_{app}$  values to give a corrected transporter-mediated B-A  $P_{app}$ . This was subsequently converted to percentage vehicle control, plotted against inhibitor concentration and the resulting curve fitted to determine inhibitory potency ( $IC_{50}$ ).

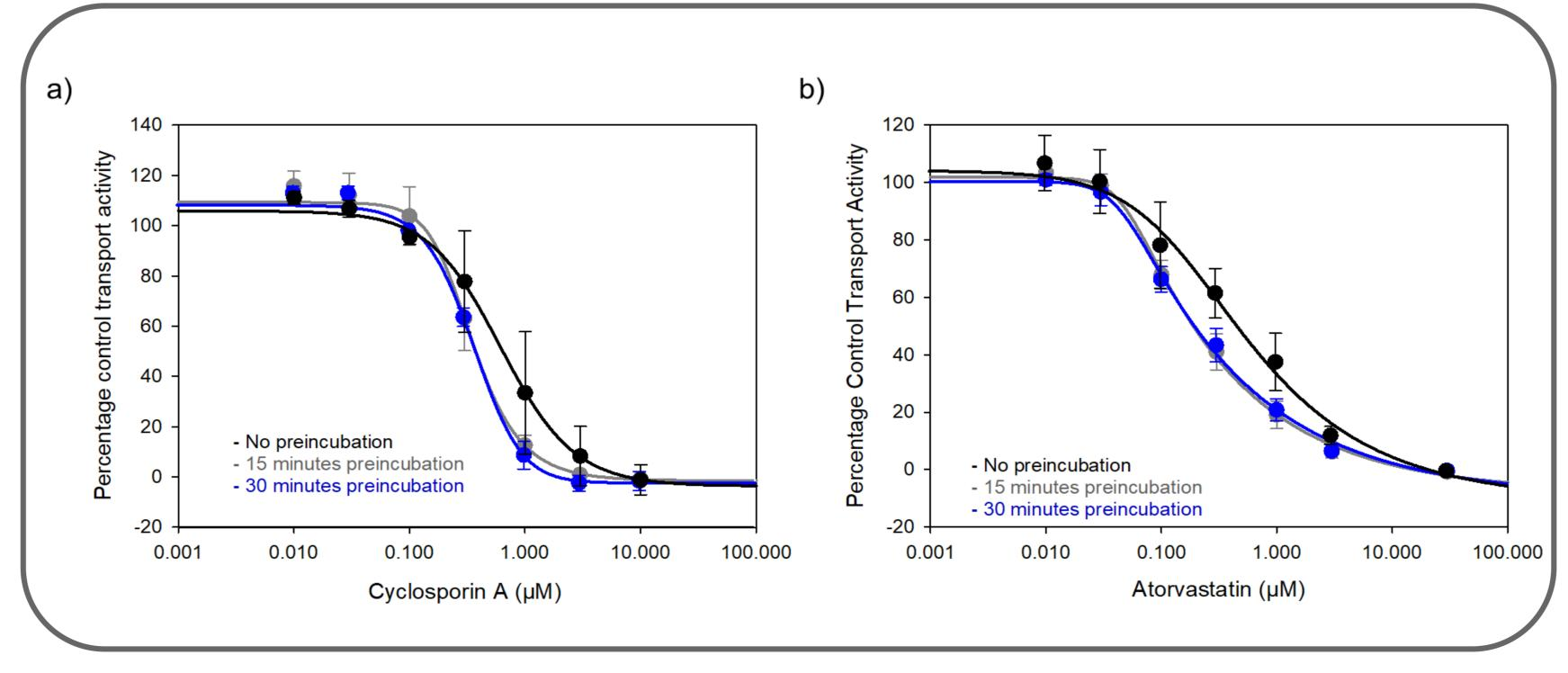


Figure 2: Comparison of in vitro determined IC<sub>50</sub> values against OATP1B1-mediated [βH]estradiol glucuronide (0.02μM) transport following a 15 min preincubation without inhibitor (vehicle buffer) and a 15 or 30 min preincubation with inhibitor. (a) concentration-dependent inhibition for cyclosporin A, (b) concentration dependent inhibition for atorvastatin

### Results

Cyclosporin A caused equipotent inhibition of OATP1B1-mediated transport following either a 15 min preincubation step with inhibitor (mean  $IC_{50}=0.372~\mu M$ ) or a 30 min preincubation step with inhibitor (mean  $IC_{50}=0.354~\mu M$ ; Table 1). A similar result was observed for atorvastatin which inhibited OATP1B1 with a mean  $IC_{50}$  value of 0.235  $\mu M$  or 0.248  $\mu M$ , determined following a 15 min or 30 min preincubation with inhibitor, respectively (Table 1). Preincubation with inhibitor for both 15 and 30 minutes enhanced the inhibitory potency of cyclosporin A and atorvastatin in vitro by decreasing their determined  $IC_{50}$  values 1.78 and 2.00-fold when compared to preincubation with vehicle buffer alone (Figure 2).

For BCRP inhibition, preincubation of Caco-2 cells with vehicle buffer alone resulted in mean  $IC_{50}$  values of 2.01  $\mu$ M and 0.273  $\mu$ M for novobiocin and fumitremorgin C, respectively. Furthermore, inclusion of inhibitor in the preincubation step had no impact on the determined  $IC_{50}$  values for novobiocin (2.06 $\mu$ M) and fumitremorgin C (0.250 $\mu$ M; Table 2).

Conversely for P-gp, whilst preincubation of MDCK-MDR1 cells with inhibitor resulted in only a small shift (decrease) in the determined mean IC $_{50}$  values for cyclosporin A (1.60  $\mu$ M to 0.931  $\mu$ M) and ketoconazole (14.9  $\mu$ M to 8.83  $\mu$ M) compared to preincubation with vehicle buffer alone, these changes were not statistically significant. Furthermore, no difference in IC $_{50}$  was observed for verapamil (Table 3). Only elacridar exhibited a statically significant 2.9-fold decrease in mean IC $_{50}$  (0.814  $\mu$ M to 0.284  $\mu$ M) following inhibitor preincubation (p ≤ 0.001; unpaired t-test assuming unequal variances; Table 3, Figure 1).

#### Conclusions

In agreement with our previous findings and with the literature we have shown that preincubation with inhibitor enhances the inhibitory potency determined against OATP1B1 in vitro [1,2].

Based on the OATP1B1 probe substrate and reference inhibitor combinations utilised in this study, our observations indicate that a 15 min preincubation step with inhibitor produces the same degree of inhibition ( $IC_{50}$ ) against OATP1B1 as the 30 min preincubation period suggested in the current FDA draft guidance (2017).

Furthermore, whilst an inhibitor preincubation step may possibly be deemed necessary for certain compounds when assessing inhibition of P-gp-mediated transport in MDCK-MDR1 cell monolayers, it does not appear to be a necessary requirement when studying inhibition of BCRP in Caco-2 cell monolayers based on the probe substrate and reference inhibitor combinations utilised in this study.

#### References

[1] Elsby R, Chidlaw S, Outteridge S, Sullivan R and Pickering S. (2016) Further investigation of the impact of inhibitor pre-incubation on human OATP1B1, OAT3, OCT2 and MATE1 transporter in vitro inhibitory potencies. The AAPS Journal (M1022), available from http://www.aapsj.org.

[2] Shitara Y, Sugiyama Y. (2017) Preincubation-dependent and long-lasting inhibition of organic anion transporting polypeptide (OATP) and its impact on drug-drug interactions. Pharmacology & Therapeutics 177: 67-80.

[3] Draft FDA Guidance for Industry – In vitro metabolism and transporter mediated drug-drug interaction studies, October 2017.