# EXAMPLE A BiolVT Company

# In Vitro System-Dependent Inhibition of Cytochrome P450 Enzymes (CYP), UDP-Glucuronosyl Transferases (UGT) and Transporters by Oligonucleotides Faraz Kazmi, Phyllis Yerino, Ellis Bixler, Chase McCoy and David B. Buckley

XenoTech, LLC, 16825 W 116th St, Lenexa, KS, USA

## Introduction

Oligonucleotides represent a growing class of biotherapeutics currently being developed for a variety of indications. They are often developed as antisense molecules (of approximately 20 nucleotides in length) designed to hybridize to a target mRNA molecule, preventing its translation to a protein. In many cases, these large polyanionic molecules are developed with phosphorothioate backbones to improve their biological stability. Phosphorothioate oligonucleotides are biopharmaceutics classification system (BCS) class III compounds, such that they have high aqueous solubility but low intestinal permeability, and are highly bound to plasma proteins (Geary, 2009; Yu *et al.*, 2013). Despite the increased pharmaceutical development of phosphorothioate oligonucleotides, there are limited drug-drug interaction (DDI) data available on this class of biotherapeutics.

We previously examined the cytochrome P450 (CYP) inhibition potential of oligonucleotides and reported that the phosphorothioate, but not phosphodiester, backboned molecules caused potent inhibition of CYP1A2 and CYP2C8 in human liver microsomes (HLM) but not in cryopreserved human hepatocytes (CHH) (Buckley *et al.*, 2009). In this study we expanded upon our previous work to include direct, time- (TDI) and metabolism-dependent

## Results (cont.)

There was evidence of time-dependent inhibition (TDI) of CYP enzymes in HLM in the absence of NADPHwith oligo #2 and #4 for CYP2B6, CYP2C8, CYP2C19, CYP2C9, CYP2D6 and CYP3A4/5. Little to no TDI was observed for UGT enzymes in HLM (data not shown). In contrast to HLM, as shown in **Table 3** and **Figure 3**, we observed little to no direct inhibition by any oligo in CHH (with the exception of oligo #3 with CYP2C19 and TDI by oligo #2 with CYP2C8), demonstrating system-dependent outcomes. There was moderate inhibition of CYP3A4/5 in CHH by oligo #2 after a 90 min pre-incubation; however, this inhibition was not concentration dependent and not considered significant despite yielding an IC<sub>50</sub> value (49  $\mu$ M).

The oligos were also tested as inhibitors of uptake and efflux transporters and the results are shown in **Table 4** and **Figure 4**. In summary, oligo #1 was found to inhibit OAT1 and OCT2; oligo #2 was found to inhibit OATP1B1, OATP1B3, OAT1, OAT3 and OCT2; oligo #3 was found to inhibit OAT1, OAT3 and OCT2; and oligo #4 was found to inhibit OATP1B3, OAT1 and OCT2. While phosphorothioate oligos overall inhibited more uptake transporters, phosphodiester oligos were inhibitory towards OAT1, OAT3 and OCT2. None of the oligos significantly inhibited the efflux transporters Pgp and BCRP.

#### **Table 4**. $IC_{50}$ values for uptake and efflux transporter inhibition with oligonucleotides

Transporter	Qubaturata	IC <sub>50</sub> (μΜ)									
	Substrate	Oligo #1	Oligo #2	Oligo #3	Oligo #4						
OATP1B1	Estradiol 17-β-glucuronide	>100	90	>100	>100						
OATP1B3	Estradiol 17-β-glucuronide	100	14	100	92						
OAT1	<i>p</i> -Aminohippuric acid	46	12	52	16						
OAT3	Estrone 3-sulfate	> 100	26	69	>100						
OCT1	Tetraethylammonium bromide	>100	>100	>100	-100						
OCT2	Metformin	17	14	27	29						
Pgp	Digoxin	>100	>100	>100	>100						
BCRP	Prazosin		> 100	> 100	>100						

### inhibition (MDI) of 7 CYP and 8 UGT enzymes in pooled HLM and pooled CHH. Additionally, inhibition of the transporters OATP1B1, OATP1B3, OAT1, OAT3, OCT1, OCT2, Pgp and BCRP was also examined.

Legend

## Materials & Methods

#### <u>Oligonucleotides</u>

Two ~20 base oligonucleotides with partial sequences towards glutathione peroxidase (5'-GCTCGTTCATCTGGGTGTAGT-3') and beta actin (5'-CCTCCTCTTTGTTCCCTTCT-3'), each with either a phosphodiester or phosphorothioate backbone (generic structures shown in **Figure 1**) were purchased from Integrated DNA Technologies (Coralville, IA).

#### <u>Chemicals</u>

Chenodeoxycholic acid (CDCA), digoxin, estradiol glucuronide, estrone sulfate, morphine, 1-naphthol, saccharic acid 1,4-lactone, oxazepam, *p*-aminohippuric acid, prazosin, propofol, testosterone and trifluoperazine were all purchased from Sigma-Aldrich (St. Louis, MO). Levomedetomidine was a gift from Orion Corporation (Espoo, Finland). [3H]-Estradiol glucuronide and [3H]-Estrone sulfate were purchased from Perkin Elmer (Waltham, MA). [14C]-Metformin was purchased from Moravek Biochemicals (Brea, CA). [14C]-Tetraethylammonium bromide and [3H]-*p*-aminohippuric acid were purchased from American Radio-labeled Chemicals (St. Louis, MO). d₅-Testosterone 17-O-glucuronide,



Oligonucleotide

d<sub>5</sub>-oxazepam-*N*-glucuronide and prochlorperazine glucuronide used as internal standards were prepared in-house. All other deuterated glucuronides were purchased from Toronto Research Chemicals (Toronto, Canada). The sources of all other reagents have been described previously (Parkinson *et al.*, 2011; Kazmi *et al.*, 2014).

#### Table 1. Experimental conditions for measurement of CYP and UGT activity for HLM and CHH studies

Enzyme	CYP or UGT activity	Substrate concentration (µM)	HLM (mg/mL)	CHH (million cells/mL)	Incubation time in HLM (min)	Incubation time in CHH (min)	Mass transition monitored (m/z)	Internal Standard
CYP1A2	Phenacetin O-deethylation	40	0.1				152 / 110	d₄-Acetaminophen
CYP2B6	Bupropion hydroxylation	50	0.1	0.5	5		256 / 238	d <sub>6</sub> -Hydroxybupropion
CYP2C8	Amodiaquine <i>N</i> -dealkylation	1.5 (7 for CHH)	0.0125				328 / 283	d <sub>5</sub> - <i>N</i> -Desethylamodiaquine
CYP2C9	Diclofenac 4'-hydroxylation	6				10	310 / 266	d <sub>4</sub> -4'-Hydroxydiclofenac
CYP2C19	S-Mephenytoin 4'-hydroxylation	40	]				233 / 190	d <sub>3</sub> -4'-Hydroxymephenytoin
CYP2D6	Dextromethorphan O-demethylation	7.5	]				258 / 157	d <sub>3</sub> -Dextrorphan
CYP3A4/5	Midazolam 1'-hydroxylation	4	0.1				342 / 324	d₄-1′-Hydroxymidazolam
UGT1A1	Estradiol 3-O-glucuronidation	9		NA	5		447 / 271	d <sub>5</sub> -Estradiol 3-O-glucuronide
UGT1A3	Chenodeoxycholic acid 24-O-glucuronidation	20			10		567 / 391	d <sub>5</sub> -Chenodeoxycholic acid 24-O-glucuronide
UGT1A4	Trifluoperazine glucuronidation	12			5		584 / 408	Prochlorperazine glucuronide
UGT1A6	1-Naphthol glucuronidation	1	0.0125			NA	319 / 143	d7-Naphthol glucuronide
UGT1A9	Propofol glucuronidation	20			5		353 / 177	d <sub>17</sub> -Propofol glucuronide
UGT2B7	Morphine 3-O-glucuronidation	400					462 / 286	d <sub>3</sub> -Morphine 3-O-glucuronide
UGT2B10	Levomedetomidine glucuronidation	7	0 1				377 / 201	d₄-1′-Hydroxymidazolam
UGT2B15	S-Oxazepam glucuronidation	50			10		463 / 286	d <sub>5</sub> -S-Oxazepam glucuronide
UGT2B17	Testosterone 17-O-glucuronidation	5	]				465 / 289	d <sub>5</sub> -Testosterone 17-O-glucuronide
UGT	4-Methylumbelliferone glucuronidation	70		0.5	NA	10	353 / 177	d <sub>5</sub> -7-Hydroxycoumarin glucuronid
NA: Not ap	plicable HLM: Human I	iver microsom	es	CHF	I: Cryoprese	erved huma	an hepatocytes	

#### **Table 2**. $IC_{50}$ values for CYP and UGT inhibition with oligonucleotides in HLM

	Substrate		IC <sub>50</sub> (μΜ)											
Enzyme			Oligo #1			Oligo #2			Oligo #3			Oligo #4		
		Direct	TDI	MDI	Direct	TDI	MDI	Direct	TDI	MDI	Direct	TDI	MDI	
CYP1A2	Phenacetin				0.8	0.8	1.8				4.2	3.4	7.3	
CYP2B6	Bupropion				15	15	13	>100	>100	>100	39	20	22	
CYP2C8	Amodiaquine		× 100	> 100	1.1	0.6	0.6				12	8.3	3.8	
CYP2C9	Diclofenac		>100	>100	97	36	47				>100	53	57	
CYP2C19	S-Mephenytoin				21	8.7	19	29	36		80	39	98	
CYP2D6	Dextromethorphan				81	41	>100		>100		>100	52	>100	
CYP3A4/5	Midazolam				>100	38	28					66	63	
UGT1A1	Estradiol	>100	I		4.5						5.4			
UGT1A3	Chenodeoxycholic acid				51						48			
UGT1A4	Trifluoperazine				>100						>100			
UGT1A6	1-Naphthol				52			>100			55			
UGT1A9	Propofol		N	D	19	Ν	٧D		ND		>42	ND		
UGT2B7	Morphine				>100						>100			
UGT2B10	Levomedetomidine				26						41			
UGT2B15	S-Oxazepam			n 1	38						45			
UGT2B17	Testosterone				7.2						18			

Figure 3.  $IC_{50}$  plots of CYP1A2, CYP2C8 and UGT inhibition by oligonucleotides in CHH



#### <u>Test system</u>

Pooled human liver microsomes (HLM, n = 200, mixed gender) and pooled cryopreserved human hepatocytes (CHH, n = 100, mixed gender) were prepared from non-transplantable livers and characterized at XenoTech, LLC (Lenexa, KS) as described previously (Pearce *et al.*, 1996; Parkinson *et al.*, 2004). HEK293 (Human embryonic kidney 293) cells expressing transporter (HEK293 cells transfected with vectors containing human transporter cDNA for OATP1B1, OATP1B3, OAT1, OAT3 and OCT2) and control cells (HEK293 cells transfected with only vectors) will be used. HEK293 cells were supplied from American Type Culture Collection (Manassas, VA) and transfected with the vector containing transporter cDNA by ADME & Tox. Research Institute, Sekisui Medical Co., Ltd. (Tokai, Japan); MDCKII cells transfected with BCRP were obtained from Netherlands Cancer Institute (Amsterdam, Netherlands); and Caco-2 cells were obtained from American Type Culture Collection (Manassas, VA).

#### In vitro CYP and UGT inhibition

The effects of oligonucleotides on CYP enzymes in HLM were evaluated in IC<sub>50</sub> experiments with and without a preincubation step (in the presence and absence of NADPH) as described previously (Parkinson *et al.*, 2011). For CHH, incubations were conducted in 100  $\mu$ l mixtures at 0.5 million cells/mL in KHB. Reactions were initiated with the addition of CHH to the oligos and incubations were conducted at 37°C with 95% humidity and 5/95% CO<sub>2</sub>/air for 0, 30 and 90 min, followed by addition of a CYP marker substrate (or the broad specific UGT substrate 4-methylumbelliferone; 4-MU) at a concentration approximately equal to its  $K_m$  (see **Table 1**). Reactions were terminated after 10 min by the addition of 100  $\mu$ l of acetonitrile containing the appropriate internal standard. Precipitated protein was removed by centrifugation (920 RCF for 10 min at 10°C) followed by LC/MS/MS analysis.

For UGT inhibition assays, oligos were incubated at 37°C in 150 µl incubation mixtures containing pooled HLMs (<0.1 mg/mL), Tris buffer (100 mM, pH 7.7), MgCl<sub>2</sub> (10 mM), EDTA (1 mM, pH 7.4), saccharic acid 1,4-lactone (0.1 mM), UDPGA (10 mM), and a UGT marker substrate at a concentration approximately equal to its  $K_m$ , at the final concentrations indicated in **Table 1**. Reactions were initiated by the addition of UDPGA and terminated after 5 or 10 min by the addition of 175 µl of acetonitrile containing the appropriate internal standard. Precipitated protein was removed by centrifugation (920 RCF for 10 min at 10°C) followed by LC/MS/MS analysis. TDI for UGT enzymes was assessed with a single concentration of each oligo (30 µM) with and without a 30 min preincubation step (in the presence and absence of UDPGA) with the methodology described above. IC<sub>50</sub> values were determined with GraFit (version 7.0.2; Erithricus Software Ltd., Surrey, UK).

#### In vitro uptake and efflux transporter inhibition

To evaluate inhibition of uptake transporters, HEK293 cells over-expressing various uptake transporters were plated in 24-well plates and grown in an incubator (37°C, 95% relative humidity, 5% CO<sub>2</sub>) for two days. On the day of the assay, cells were pre-incubated with oligos or solvent control for 15 min followed by 2 min incubation with a combination of probe substrate (a mixture of cold and radiolabeled compound) and oligo. 50 nM estradiol 17-β-glucuronide substrate was used for OATP1B1 and OATP1B3 assays (with approximate equimolar radiolabeled and cold substrate); 1 µM p-aminohippuric acid for OAT1 (1:4 ratio of radiolabeled to cold substrate); 50 nM estrone 3-sulfate for OAT3 (approximately equimolar radiolabeled and cold substrate); and 5 µM tetraethylammonium bromide and 10 µM metformin for OCT1 and OCT2 respectively (all radiolabeled). Uptake assays were terminated by aspiration of the incubation media followed by washing the cells three times, first a rinse with ice cold PBS containing 0.2% BSA then two rinses with ice cold PBS. For analysis, cells were extracted with 0.1N NaOH for scintillation counting on a Microbeta 2 instrument (Perkin Elmer, Waltham, MA). Evaluation of efflux transporter inhibition was conducted in Caco-2 (MDR1) or MDCK-II cells (transfected with BCRP) plated on 24-well trans-well plates. The bidirectional permeability of specific probe substrates (10 µM digoxin for Caco-2 cells and 1 µM prazosin for MDCK-II cells) was measured in the presence of oligonucleotides. Oligos were added to both apical and basolateral sides for 120 min. After the incubation, samples were stopped with 50:50 v/v methanol:water and analyzed by LC-MS/MS. IC<sub>50</sub> values were determined with GraFit (version 7.0.2; Erithricus Software Ltd., Surrey, UK).

**Figure 2.** IC<sub>50</sub> plots of CYP1A2, CYP2C8 and UGT1A1 inhibition by oligonucleotides in HLM



#### **Figure 4.** The inhibition of the uptake transporters OATP1B3, OAT1, OAT3 and OCT2 by oligonucleotides



## Conclusions

- In HLM, phosphorothioate but not phosphodiester oligonucleotides were found to be inhibitors of CYP and UGT enzymes.
- In CHH, inhibition of CYP and UGT enzymes was dramatically less potent than in HLM, suggesting test system-dependent inhibitory effects.
- Both phosphodiester and phosphothioate oligonucleotides were found to be inhibitors of uptake but not efflux transporters.

## Results

As shown in **Table 2**, **Figure 2** and **Table 3**, the results indicated that oligos #1 and #3 with the phosphodiester backbones had little to no inhibitory effect on all CYP and UGT enzymes in HLM and CHH with the exception of oligo #3 in HLM for CYP2C19. Conversely, the oligos with the phosphorothioate backbone caused direct inhibition of CYP and UGT enzymes. The rank order of direct inhibition in HLM for oligo #2 was CYP1A2 > CYP2C8 > UGT1A1 > UGT2B17 > CYP2B6 > UGT1A9 > CYP2C19 > UGT2B10 > UGT2B15 > UGT1A3 > UGT1A6 > CYP2D6 > CYP2C9. Similar results were observed for oligo #4: CYP1A2 > UGT1A1 > CYP2C8 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B17 > CYP2C8 > UGT2B10 > UGT2B10 > UGT1A1 > CYP2C8 > UGT2B17 > CYP2B6 > UGT2B15 > UGT1A3 > UGT1A6 > CYP2B6 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B17 > CYP2C8 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B10 > UGT2B17 > CYP2C8 > UGT2B10 > UG

Table 3. IC<sub>50</sub> values for CYP and UGT inhibition with oligonucleotides in CHH

	Substrate	-	IC <sub>50</sub> (μΜ)												
Fnzvme			Oligo #1			Oligo #2			Oligo #3			Oligo #4			
		Direct	MDI 30 min	MDI 90 min	Direct	MDI 30 min	MDI 90 min	Direct	MDI 30 min	MDI 90 min	Direct	MDI 30 min	MDI 90 min		
CYP1A2	Phenacetin		>100	0 >100	>100	>100	>100 >100	- >100 - 36		>100	>100	>100	>100		
CYP2B6	Bupropion									77					
CYP2C8	Amodiaquine					35	44								
CYP2C9	Diclofenac	>100				>100	>100		>100						
CYP2C19	S-Mephenytoin									>100					
CYP2D6	Dextromethorphan									- 100					
CYP3A4/5	Midazolam						49	>100							
UGT	4-Methylumbelliferone						>100	1							

 Overall this study has implications on the design of *in vitro* DDI studies for oligonucleotide based therapeutics based upon test system-dependent differences.

## References

- Buckley DB, Kazmi F, Yerino P, Ogilvie BW, and Parkinson A (2009) Inhibition of cytochrome P450 (CYP) Enzymes, CYP1A2 and CYP2C8, by oligonucleotides in human liver microsomes (HLM): A system-dependent outcome. *Drug Metab Rev* **41 (S3):**94-95.
- 2 Geary RS (2009) Antisense oligonucleotide pharmacokinetics and metabolism. *Expert Opin Drug Metab Toxicol* **5**:381-391.
- 3 Kazmi F, Haupt LJ, Horkman JR, Smith BD, Buckley DB, Wachter EA, and Singer JM (2014) *In vitro* inhibition of human liver cytochrome P450 (CYP) and UDP-glucuronosyltransferase (UGT) enzymes by rose bengal: system-dependent effects on inhibitory potential. *Xenobiotica* **44**:606-614.
- 4 Parkinson A, Kazmi F, Buckley DB, Yerino P, Paris BL, Holsapple J, Toren P, Otradovec SM, and Ogilvie BW (2011) An evaluation of the dilution method for identifying metabolism-dependent inhibitors of cytochrome p450 enzymes. *Drug Metab Dispos* **39**:1370-1387.
- 5 Parkinson A, Mudra DR, Johnson C, Dwyer A, and Carroll KM (2004) The effects of gender, age, ethnicity, and liver cirrhosis on cytochrome P450 enzyme activity in human liver microsomes and inducibility in cultured human hepatocytes. *Toxicol Appl Pharmacol* **199**:193-209.
- 6 Pearce RE, Rodrigues AD, Goldstein JA, and Parkinson A (1996) Identification of the human P450 enzymes involved in lansoprazole metabolism. *J Pharmacol Exp Ther* **277**:805-816.
- 7 Yu RZ, Grundy JS, and Geary RS (2013) Clinical pharmacokinetics of second generation antisense oligonucleotides. *Expert Opin Drug Metab Toxicol* **9:**169-182.