XENOTECH A BiolVT Company

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 inhibition (MDI) of 7 CYP and 8 UGT enzymes in pooled HLM and pooled CHH.

Materials & Methods

Oligonucleotides

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).

Chemicals

Chenodeoxycholic acid (CDCA), estradiol glucuronide, estrone sulfate, morphine, 1-naphthol, saccharic acid 1,4-lactone, oxazepam, p-aminohippuric acid, propofol, testosterone and trifluoperazine were all purchased from SigmaAldrich (St. Louis, MO). Levomedetomidine was a gift from Orion Corporation (Espoo, Finland). d₅-Testosterone 17-O-glucuronide, d₅--oxazepam-*N*-glucuronide and prochlorperazine glucuronide used as internal standards were prepared inhouse. 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).

| Legend | Oligonucleotide |
|----------|---------------------------------|
| Oligo #1 | Glutathione peroxidase phospho |
| Oligo #2 | Glutathione peroxidase phospho |
| Oligo #3 | Beta actin phosphodiester back |
| Oligo #4 | Beta actin phosphorothioate bac |



Test system

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 (Kansas City, KS) as described previously (Pearce et al., 1996; Parkinson et al., 2004).

In vitro CYP and UGT inhibition

The effects of oligonucleotides on CYP enzymes in HLM were evaluated in IC_{50} 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 µl 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% CO₂/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 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.

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), MgCl₂ (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. IC₅₀ values were determined with GraFit (version 7.0.2; Erithricus Software Ltd., Surrey, UK).

Results

As shown in **Table 2**, **Figure 2** and **Table 3**, the results indicated that phosphodiester-backbone oligonucleotides (#1 and #3) had little to no inhibitory effect on CYP and UGT enzymes in HLM and CHH with the exception of oligo #3 in HLM for CYP2C19. Conversely, the phosphorothioate-backbone oligonucleotides (#2 and #4) caused direct inhibition of CYP and UGT enzymes. The rank order of direct inhibition in HLM for oligo #2 and oligo #4 were similar. CYPs 1A2, 2B6, 2C8, and UGTs 1A1 and 2B17 were most highly inhibited by both phosphorothioatebackbone oligonucleotides.

Maciej Czerwinski, Faraz Kazmi, Phyllis Yerino, and David B. Buckley XenoTech, LLC, Kansas City, KS, USA

odiester backbone prothioate backbone oone ckbone

There was evidence of time-dependent inhibition (TDI) of CYP enzymes in HLM in the absence of NADPH with 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_{50} value (49 μ M).

| 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 | | | | 256 / 238 | d ₆ -Hydroxybupropion | |
| CYP2C8 | Amodiaquine <i>N</i> -dealkylation | 1.5 (7 for CHH) | 0.0125 | 0.5 | | | 328 / 283 | d ₅ - <i>N</i> -Desethylamodiaquine | |
| CYP2C9 | Diclofenac 4'-hydroxylation | 6 | | | 5 | 10 | 310 / 266 | d ₄ -4'-Hydroxydiclofenac | |
| CYP2C19 | S-Mephenytoin 4'-hydroxylation | 40 | | | | | 233 / 190 | d ₃ -4'-Hydroxymephenytoin | |
| CYP2D6 | Dextromethorphan O-demethylation | 7.5 | | | | | 258 / 157 | d ₃ -Dextrorphan | |
| CYP3A4/5 | Midazolam 1'-hydroxylation | 4 | 0.1 | | | | 342 / 324 | d₄-1′-Hydroxymidazolam | |
| UGT1A1 | Estradiol 3-O-glucuronidation | 9 | | | 5 | | 447 / 271 | d ₅ -Estradiol 3- <i>O</i> -glucuronide | |
| UGT1A3 | Chenodeoxycholic acid 24-O-glucuronidation | 20 | | | 10 | | 567 / 391 | d₅-Chenodeoxycholic acid 24-O-glucuronide | |
| UGT1A4 | Trifluoperazine glucuronidation | 12 | | | | | 584 / 408 | Prochlorperazine glucuronide | |
| UGT1A6 | 1-Naphthol glucuronidation | 1 | 0.0125 | NA | 5 | NA | 319 / 143 | d ₇ -Naphthol glucuronide | |
| UGT1A9 | Propofol glucuronidation | 20 | | | 5 | | 353 / 177 | d ₁₇ -Propofol glucuronide | |
| UGT2B7 | Morphine 3-O-glucuronidation | 400 | | | | | 462 / 286 | d ₃ -Morphine 3-O-glucuronide | |
| UGT2B10 | Levomedetomidine glucuronidation | 7 | 0.1 | | 10 | | 377 / 201 | d₄-1′-Hydroxymidazolam | |
| UGT2B15 | S-Oxazepam glucuronidation | 50 | 0.1 | | | | 463 / 286 | d₅-S-Oxazepam glucuronide | |
| UGT2B17 | Testosterone 17-O-glucuronidation | 5 | | | | | 465 / 289 | d ₅ -Testosterone 17-O-glucuronide | |
| UGT | 4-Methylumbelliferone glucuronidation | 70 | | 0.5 | NA | 10 | 353 / 177 | d ₅ -7-Hydroxycoumarin glucuronide | |
| NA: Not applicable HLM: Human liver microsomes CHH: Cryopreserved human hepatocytes | | | | | | | | | |

Table 2. IC₅₀ values for CYP and UGT inhibition with oligonucleotides in HLM Oligo #4 Oligo #3 TDI MDI Direct TDI MDI 4.2 3.4 7.3 22 39 20 >100 100 3.8 12 8.3 >100 >100 57 53 29 39 98 36 80 >100 52 >100 >100 66 63 5.4 48 >100 100 55 ND ND >42 >100 41 45 18

| | | ΙC ₅₀ (μΝ | | | | | | | | | |
|--------------|-----------------------|----------------------|----------|-------|----------|-----|------|---|--|--|--|
| Enzyme | Substrate | | Oligo #1 | 1 | Oligo #2 | | | | | | |
| | | Direct | TDI | MDI | Direct | TDI | MDI | D | | | |
| CYP1A2 | Phenacetin | | | | 0.8 | 0.8 | 1.8 | > | | | |
| CYP2B6 | Bupropion | 7 | | | 15 | 15 | 13 | | | | |
| CYP2C8 | Amodiaquine | 7 | >100 | > 100 | 1.1 | 0.6 | 0.6 | | | | |
| CYP2C9 | Diclofenac |] | | >100 | 97 | 36 | 47 | | | | |
| CYP2C19 | S-Mephenytoin | 7 | | | 21 | 8.7 | 19 | | | | |
| CYP2D6 | Dextromethorphan | 7 | | | 81 | 41 | >100 | | | | |
| CYP3A4/5 | Midazolam | 7 | | | >100 | 38 | 28 | | | | |
| UGT1A1 | Estradiol | >100 | | | 4.5 | | | | | | |
| UGT1A3 | Chenodeoxycholic acid |] | | | 51 | 1 | | | | | |
| UGT1A4 | Trifluoperazine |] | ND | | >100 | ND | | > | | | |
| UGT1A6 | 1-Naphthol |] | | | 52 | | | | | | |
| UGT1A9 | Propofol | 7 | | | 19 | | | | | | |
| UGT2B7 | Morphine |] | | | >100 | | | | | | |
| UGT2B10 | Levomedetomidine | | | | 26 | | | | | | |
| UGT2B15 | S-Oxazepam | | | | 38 | | | | | | |
| UGT2B17 | Testosterone |] | | | 7.2 | | | | | | |
| ND: Not dete | ermined | | | | | | | | | | |



Hepatocytes Are the Preferred Test System to Evaluate **Oligonucleotide-CYP Interactions** In Vitro

| Table 3. IC_{50} values for CYP and UGT inhibition with oligonucleotides in CHH | | | | | | | | | | | | | | |
|---|-----------------------|--------|---------------|---------------|--------|---------------|---------------|--------|--------------------|-------------------|--------|---------------|---------------|--|
| | IC ₅₀ (μM) | | | | | | | | | - | | | | |
| Enzyme | Substrate | | 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 | >100 | > 100 | > 100 | >100 | | >100 | | >100 | >100 | |
| CYP2B6 | Bupropion | | | | | >100 | >100 | | >100 >100 36 | 77 >10 >100 |] | | | |
| CYP2C8 | Amodiaquine | | >100 | | | 35 | 44 | | | | >100 | | | |
| CYP2C9 | Diclofenac | >100 | | | | | | 1 | | | | | | |
| CYP2C19 | S-Mephenytoin | | | | | | >100 | 36 | | | | | | |
| CYP2D6 | Dextromethorphan | | | | | >100 | | | | | | | | |
| CYP3A4/5 | Midazolam | | | | | | 49 |] >100 | | | | | | |
| UGT | 4-Methylumbelliferone | | | | | | >100 | | | | | | | |



Conclusions

- In HLM, phosphorothioate but not phosphodiester oligonucleotides were found to be inhibitors of CYP and UGT enzymes.
- system-dependent inhibitory effects.
- Overall this study indicates that hepatocytes, a test system physiologically more relevant than isolated microsomes, are the preferred model to evaluate oligonucleotide-CYP interactions in vitro.

References

- 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.
- system-dependent effects on inhibitory potential. Xenobiotica 44:606-614.
- p450 enzymes. Drug Metab Dispos 39:1370-1387.
- human hepatocytes. Toxicol Appl Pharmacol 199:193-209.
- 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.

In CHH, inhibition of CYP and UGT enzymes was dramatically less potent than in HLM, suggesting test

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

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:

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

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

6 Pearce RE, Rodrigues AD, Goldstein JA, and Parkinson A (1996) Identification of the human P450 enzymes