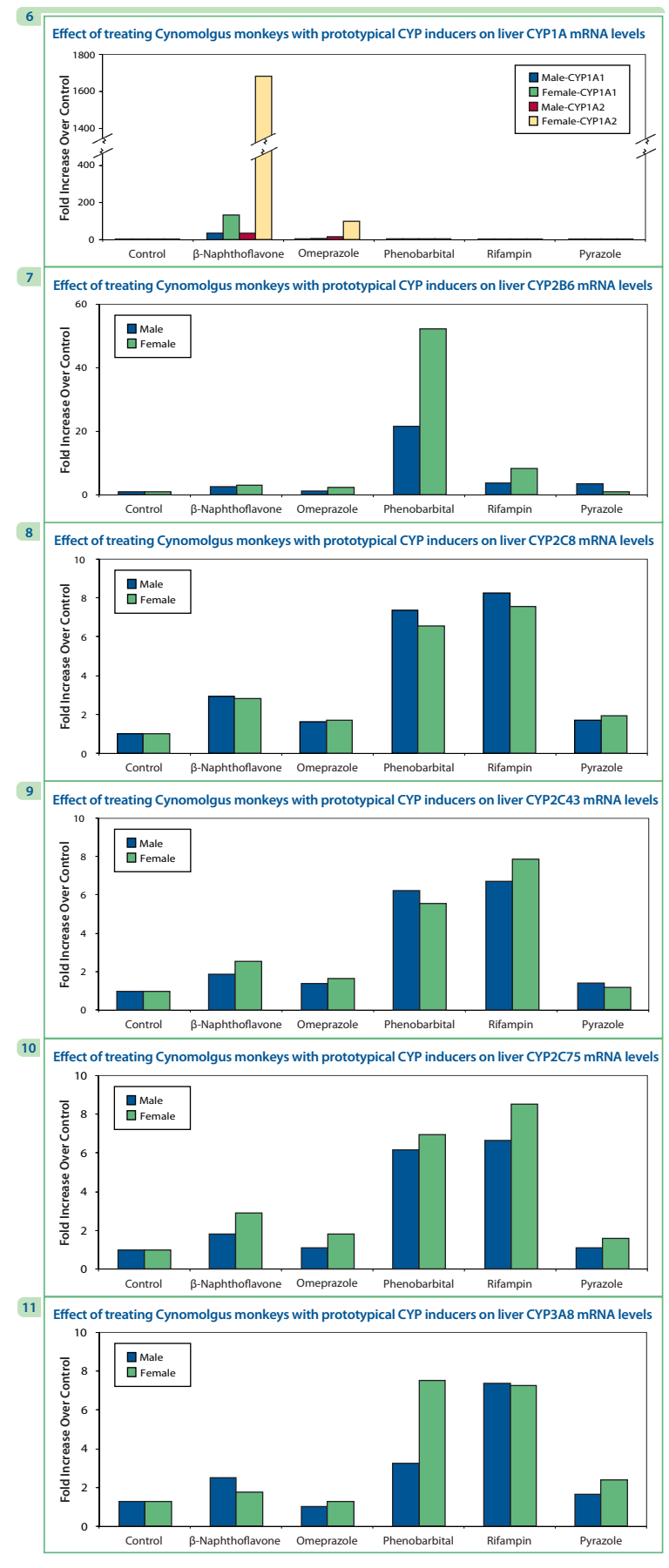


FIGURE 6-11



REFERENCES

Bullock P, Pearce R, Draper A, Podval J, Braxken W, Veltman J, Thomas P and Parkinson A. 1995. Induction of liver microsomal cytochrome P450 in cynomolgus monkeys. *Drug Metab Dispos* 23:736-748.

Burke MD, Thompson S, Weaver RJ, Wolf CR and Mayer RT. 1994. Cytochrome P450 specificities of alkoxyresorufin O-dealkylation in human and dog liver. *Biochem Pharmacol* 48: 923-936.

Pearce RE, McIntyre CJ, Madan A, Sanzgiri U, Draper AJ, Bullock PL, Cook DC, Burton LA, Latham J, Nevins C and Parkinson A. 1996. Effects of freezing, thawing, and storing human liver microsomes on cytochrome P450 activity. *Arch Biochem Biophys* 331: 145-169.

Dutton DR and Parkinson A. 1989. Reduction of 7-alkoxyresorufins by NADPH-cytochrome P450 reductase and its differential effects on their O-dealkylation by dog liver microsomal cytochrome P450. *Arch Biochem Biophys* 268: 617-629.

Huang SM, Strong JM, Zhang L. *et al.* 2008. New era in drug interaction evaluation: US Food and Drug Administration Update on CYP enzymes, transporters, and the Guidance process. *J Clin Pharmacol* 48:662-670.

Lu AYH and Levin W. 1972. Partial purification of cytochrome P-450 and cytochrome P-448 from rat liver microsomes. 1972. *Biochem Biophys Res Commun* 46:1334-1339.

Smith PK, Krohn RI, Hermanson GT, Mallia AK, Gartner FH, Provenzano MD, Fujimoto EK, Goeke NM, Olson BJ and Klenk DC. 1985. Measurement of protein using bicinchoninic acid. *Anal Biochem* 150:76-85.

Sonderfan AJ and Parkinson A. 1988. Inhibition of steroid 5 alpha-reductase and its effects on testosterone hydroxylation by rat liver microsomal cytochrome P-450. *Arch Biochem Biophys* 265:208-218.



Induction of Liver and Intestinal Cytochrome P450 (CYP) Enzymes in Male and Female Cynomolgus Monkeys

Florence Ndikum-Moffor, Michelle McBride, John Setser, C. Chengelis, Yi-Zhong Gu, Erjia Wang, Zell Woodworth, Stephanie Helmstetter, Lisa Cain and Andrew Parkinson
Presented at the 15th North American Regional ISSX Meeting, San Diego, CA, October, 2008
Abstract 194, *Drug Metabolism Reviews*, Volume 40, Supplement 3, 2008

ABSTRACT

The objective of the present study was to determine the effects of various prototypical inducers on the activity and mRNA expression of multiple CYP enzymes in monkey liver and intestine.

Male and female Cynomolgus monkeys were dosed orally (by nasogastric intubation) once daily for four days with saline (vehicle control), β-naphthoflavone (BNF, 50 mg/kg/day), phenobarbital (PB, 25 mg/kg/day), rifampin (RIF, 25 mg/kg/day), pyrazole (200 mg/kg/day) or omeprazole (OMP, 50 mg/kg/day). Microsomal samples from liver and intestine were analyzed for total cytochrome P450 content, cytochrome b₅ content, NADPH-cytochrome c reductase activity, 7-ethoxyresorufin O-dealkylation (CYP1A1/2), testosterone 16α-hydroxylation and testosterone 16β-hydroxylation (CYP2B17), 4-nitrophenol hydroxylase (CYP2E1), testosterone 2α-hydroxylation and testosterone 6β-hydroxylation (CYP3A8). CYP mRNA levels in liver were determined for CYP1A1, 1A2, 2A24, 2B6, 2C8, 2C43, 2C75, 2E1, 3A8 and 4A11 by RT-PCR. Given the amount of data collected on this project, the primary focus will be CYP mRNA data. Treatment of monkeys with BNF caused significant increases in liver CYP1A1 and CYP1A2 mRNA levels, and little or no increase (< 4-fold) in the mRNA levels of the other enzymes examined. Omeprazole caused increases in liver CYP1A1 and 1A2 mRNA levels, but as a CYP1A mRNA inducer OMP was less effective than BNF. Treatment with BNF caused an increase in CYP1A1 mRNA levels (66- and 240-fold in males and females) and CYP1A2 mRNA levels (51- and 1600-fold in males and females). Treatment with OMP caused an increase in CYP1A1 mRNA levels (2.2- and 5.8-fold in males and females) and CYP1A2 mRNA levels (14.9- and 173-fold in males and females). Similarly, BNF and OMP caused increases in liver microsomal CYP1A1/2 activity. Treatment with BNF caused significant increases in CYP1A activity in both liver and intestine. The increase in liver CYP1A activity by BNF was similar between male and female monkeys (11- and 9.4-fold, respectively) whereas induction of intestinal CYP1A activity by BNF was much lower in male than female monkeys (2.7- and 19-fold, respectively). Though to a smaller extent compared to BNF, OMP also caused increases in microsomal CYP1A1/2 activity (3.7- and 4.3-fold in the liver) and (1.5 and 4.3-fold in intestine) for male and female monkeys, respectively. Treatment with PB caused a large increase in liver CYP2B6 mRNA levels (21.6- and 52.5-fold in males and females, respectively), and caused less pronounced increases in CYP2A24 (5.6- and 7.3-fold), 2C8 (7.3- and 6.5-fold), 2C43 (6.4- and 5.7-fold), 2C75 (6.2- and 7.0-fold), and 3A8 (2.6- and 6.0-fold) mRNA levels in males and females, respectively. PB also caused an increase in liver microsomal CYP2B17 and CYP3A8 activity (3.2- and 2.2-fold, respectively). Rifampin treatment caused an increase in liver CYP2B6 (3.6- and 8.2-fold), 2C8 (8.2- and 7.5-fold), 2C43 (6.9- and 8.1-fold), 2C75 (6.7- and 8.6-fold) and 3A8 (5.9- and 5.8-fold) mRNA levels in males and females, respectively. Except for a small increase (3.5-fold in CYP2B6 and 2.7-fold in CYP4A mRNA levels), pyrazole did not cause an increase in the liver mRNA levels of any of the CYP enzymes examined. Results from this study indicate that BNF and OMP are *in vivo* inducers of CYP1A1 and CYP1A2 in monkeys, while PB and rifampin are inducers of CYP2B, 2C and CYP3A enzymes.

