



PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE
FROM DISCOVERY THROUGH CLINICAL SUPPORT

In Vitro Induction Studies: Elements of Design and Important Considerations in Data Analysis



Joanna Barbara, Ph.D.
Vice President of Scientific Operations

Overview

- Introduction
 - Enzyme induction basics
 - Relevance
- Standard study design
 - What is included and why?
- After initial data generation
- Options for next steps
- Unanswered questions
- Conclusions

Enzyme induction basics

- Drugs can cause induction (up-regulation) of drug-metabolizing enzymes
 - E.g., Cytochrome P450 (CYP) enzymes
- Primary mechanism is activation of gene transcription
 - Often by nuclear receptor activation

Enzyme	Receptor	Acronym
CYP1A2	Aryl hydrocarbon receptor	AhR
CYP2B6	Constitutive androstane receptor	CAR
CYP3A4 etc	Pregnane X receptor	PXR

- PXR activation is the most common
 - Large / flexible binding pocket
- Species differences are prevalent

Enzyme induction relevance

- Drug metabolizing enzyme induction can alter metabolic clearance
 - Of the inducer itself: auto-induction
 - Of concomitantly administered drugs
- This needs evaluating for drug-drug interaction (DDI) potential
- Efficacy of narrow therapeutic index drugs can be compromised
 - Various common antiepileptics (e.g., carbamazepine, phenobarbital, phenytoin) induce P450 enzymes and speed up clearance of comeds (e.g., warfarin), reducing exposure and so efficacy

Guidance requirements

Requirement	FDA (2017)	EMA (2013)
Test system (number of donors)	Cryopreserved or fresh human hepatocytes (other systems considered complimentary) minimum three donors	
TA concentrations	Expected or observed human plasma drug concentrations or intestinal drug concentrations (CYP3A4/5) and span therapeutic exposures and an order of magnitude above Cmax-ss unbound	50x Cmax-ss unbound or if orally dosed 1/10 th dose in 250 mL for CYP3A4
CYP emphasis	1A2, 2B6, 3A4	1A2, 2B6, 3A4
Controls	Negative: Not specified CYP specific Positive: Omeprazole (25-100 µM) Phenobarbital (500-1000 µM) Rifampin (10-50 µM)	Negative: Not required CYP specific Positive: Omeprazole (50 µM) CITCO (100 nM) Rifampin (20 µM)

FDA 2017 Guidance for Industry: In vitro metabolism- and transporter-mediated drug-drug interaction studies.

EMA 2013 Guideline on the investigation of drug interactions

A standard study

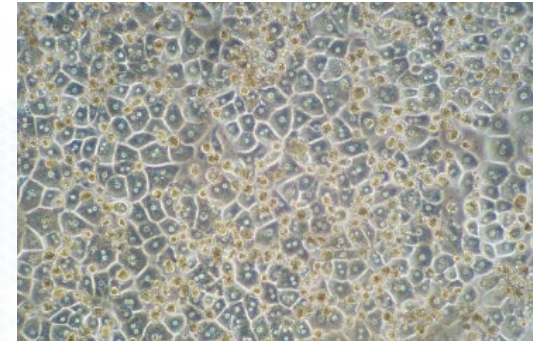
- Supporting / preliminary tests
- Cryopreserved cultured hepatocytes
 - 3 individual donors, $n = 3$
 - Lots pre-characterized for induction
- 3 day treatment at 7 concentrations (don't skimp!)
- Media drug concentration measurement
- Quantitative mRNA measurement by RT-PCR
 - CYP1A2, CYP2B6, CYP3A4)
- (Activity measurements, where applicable)
- Fold-change over vehicle control endpoint
- EC_{50}/E_{max} determination, as applicable (more later)

Supporting / preliminary assays

- 2 supporting experiments dictate test concentrations
 - Solubility testing for a high-concentration assay
 - Stock solvent solubility
 - Test system solubility (e.g., a variant of Modified Chee's Medium)
 - Cytotoxicity
 - Morphology observations
 - Lactate dehydrogenase release (LDH) assessment
- 1 more part is concurrent with the main assay
 - Spent media analysis / drug concentration analysis
 - Measuring [drug] over incubation time course on last day
 - What is it for?

Cultured hepatocytes

- Cryopreserved vs. fresh
 - Scheduling
 - Characterization
 - Continuity while the lot lasts
- Options for characterization
 - What is standard characterization?
 - Do we need to use relative induction score (RIS) characterized lots?
 - It depends. More later...



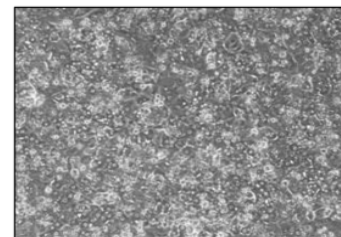
H1500.H15A+ Lot No. HC3-45

Cryopreserved Human Hepatocytes
Human, Female, Individual

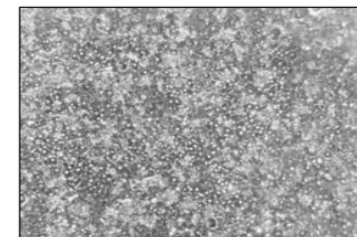
Assured Minimum Yield: 2.0×10^6 cells per vial
Viability: 83%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and OptiThaw Hepatocyte Kit.

Recommended Seeding Density: 1.4 million cells/mL



Photomicrograph (100x) of HC3-45 Day 2 of culture



Photomicrograph (100x) of HC3-45 incubation day

Induction Data

Enzyme	Inducer	mRNA Fold Induction	Marker Substrate Reaction	Enzymatic Fold Induction
CYP1A2	Omeprazole (50 μ M)	59.9	Phenacetin O-dealkylation	17.6
CYP2B6	Phenobarbital (750 μ M)	7.3	Bupropion hydroxylation	5.8
CYP3A4	Rifampin (20 μ M)	34.6	Midazolam 1'-hydroxylation	11.6

Why not immortalized cell lines?

- Acceptable per FDA 2017
- Complementary per EMA 2013
- HepaRG
 - Low CAR expression
- Fa2N4
 - Lack functional CAR
 - Low basal CYP enzyme expression
- HepG2
 - High AhR expression
 - Low basal enzyme expression
 - Little/no response to known CYP3A4 inducers

“No one cell line affords an exact reproduction of a hepatocyte and most companies felt that these cell lines are not fully understood or characterized...we do not consider immortalized hepatocyte cell lines are an inadequate replacement for primary human hepatocytes”

Target enzymes / prototypical inducers

- What are agency recommendations?

Enzyme	Receptor	FDA	EMA
CYP1A2	AhR	Omeprazole Lansoprazole	Omeprazole
CYP2B6	CAR	Phenobarbital	CITCO
CYP3A4 CYP2C8 CYP2C9 CYP2C19	PXR	Rifampin	Rifampin

- Why select phenobarbital for CYP2B6?
 - CITCO: selective for CAR, fold-change response is variable
 - Phenobarbital: consistent response, accounts for CAR/PXR crosstalk

FDA 2017 Guidance for Industry: In vitro metabolism- and transporter-mediated drug-drug interaction studies.

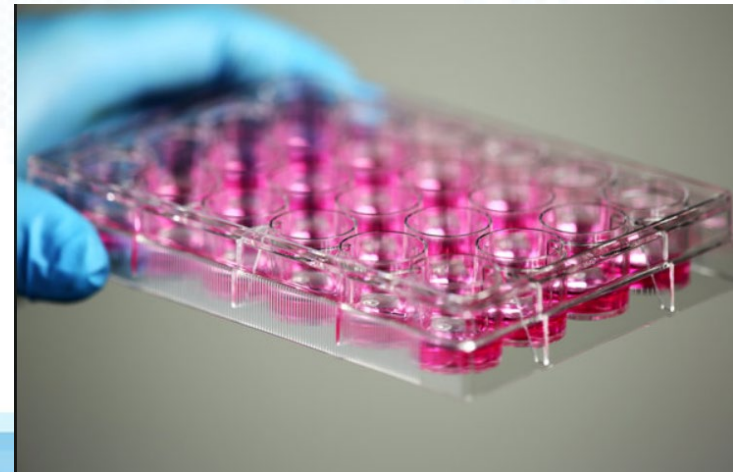
EMA 2013 Guideline on the investigation of drug interactions

Hariparsad *Drug Metab. Dispos.* 2017, 45: 1049-1059.

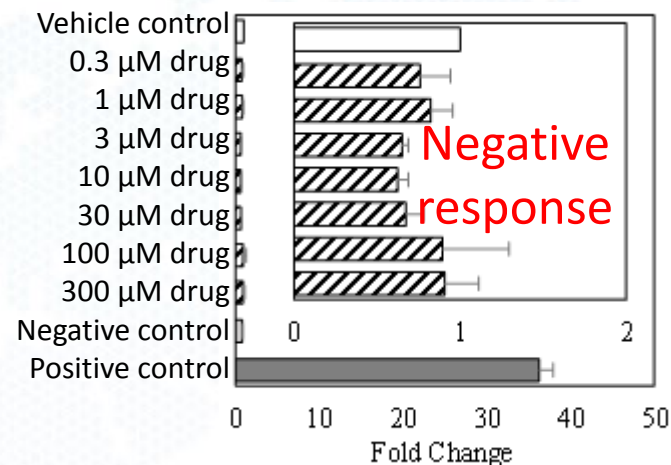
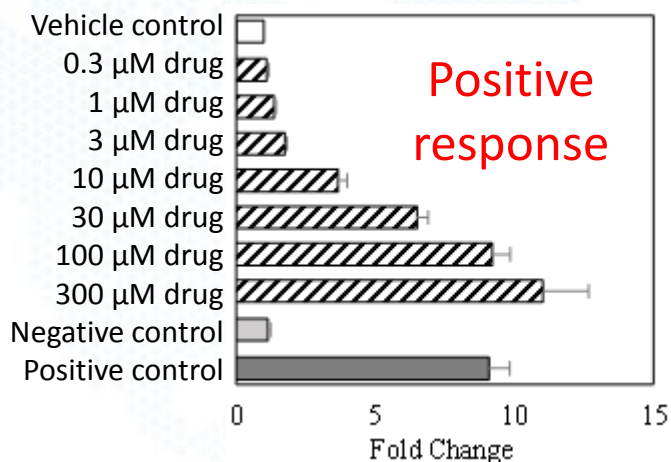
Incubation

- Sandwich-cultured hepatocytes
- 1 day adaptation
- 3 days of treatment at multiple concentrations
- Sample collection out to 24-h post last treatment
- Collect spent media for analysis
- (Perform enzyme activity incubations)
- Collect cells for mRNA harvest
- Prepare samples for analysis
- Analyze samples
 - qRT-PCR
 - LC-MS/MS

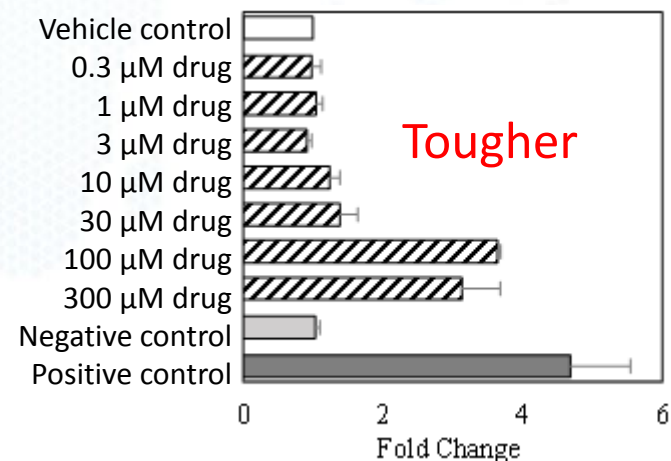
Vehicle control
0.3 μ M drug
1 μ M drug
3 μ M drug
10 μ M drug
30 μ M drug
100 μ M drug
300 μ M drug
Negative control
Positive control



First-pass data: fold-change over vehicle control

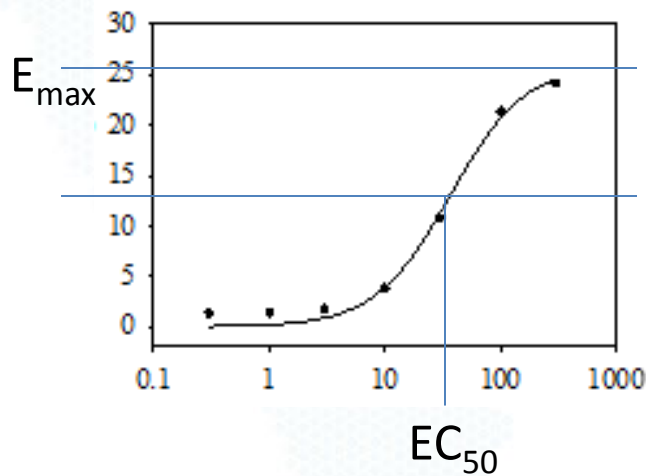


POSITIVE
 ≥ 2 -fold change over the
 vehicle control
 AND
 $\geq 20\%$ of the positive
 control response



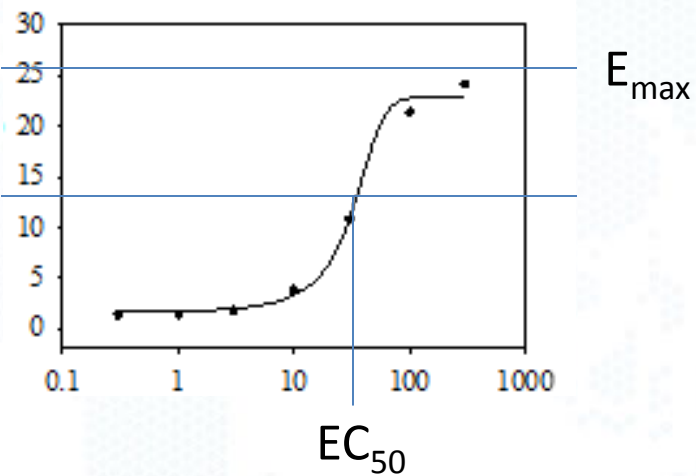
Induction constants

- EC_{50} – concentration associated with half-maximal induction
- E_{max} – maximal fold induction in vitro



Sigmoid 3-parameter

$$f = \frac{a}{1 + e^{\left(\frac{-(x - x_0)}{b}\right)}}$$



Hill 3-parameter

$$f = \frac{a \times x^b}{(x_0^b + x^b)}$$

f = fold
 $a = E_{max}$
 $x_0 = EC_{50}$
 x = drug concentration
 b = slope (Hill coefficient)

What does the data mean?

- What is a positive?
 - ≥ 2 -fold change over vehicle control and 20% of the positive control response in ANY ONE DONOR
 - FDA and EMA agree now!
- Why is one donor response a positive?
 - Interindividual variability
 - Conservative approach to safety
- What do I do about it?
 - If CYP3A4, then CYP2C evaluation warranted
 - Mathematical risk assessment is a good next step
 - More later...
 - Fourth culture? Explored in IQ IWG 2018 paper

CYP2C enzymes

- If CYP3A4 induction is observed, need to follow up with
 - CYP2C8
 - CYP2C9
 - CYP2C19
- All four enzymes can be induced through the PXR pathway
- mRNA or activity measurements?
 - mRNA fold-change highly variable for all CYP2Cs
 - mRNA rifampin fold-change often < 2
 - **Activity is a more reliable endpoint**; fold-change may still be < 2
 - Multiple possible reasons are given in IQ-IWG 2017 paper
- What about using a 'good' 2C lot?
 - Only use lots with evidence of 2C induction, but still not predictable

Data analysis for positive induction responses

- Mathematical tools for risk assessment
 - Correlation methods
 - Basic modeling
 - Static mechanistic modeling
 - Dynamic mechanistic modeling (Physiologically based pharmacokinetics - PBPK)
- Generally, need dose and $C_{\max}$ data or projections
- There are options for situations before $C_{\max}$ established

Correlation methods

Figure 3: Two Correlation Methods to Assess the Potential of an Investigational Drug to Induce Metabolizing Enzymes (Fahmi and Ripp, 2010)

Correlation Method 1: Calculate a relative induction score (RIS) using $(E_{\max} \times I_{\max,u}) / (EC_{50} + I_{\max,u})$

OR

Correlation Method 2: Calculate $I_{\max,u} / EC_{50}$ values

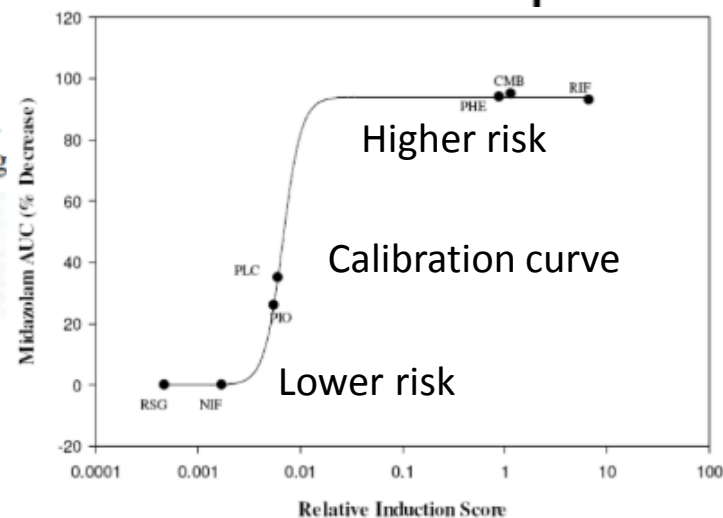
Determine the magnitude of a clinical induction effect (e.g., AUC ratio of index substrate in the presence and absence of inducers) according to a calibration curve of RIS scores or $I_{\max,u} / EC_{50}$ for a set of known inducers of the same enzyme.

$E_{\max}$ is the maximum induction effect determined in vitro.

EC_{50} is the concentration causing half-maximal effect determined in vitro.

$I_{\max,u}$ is the maximal unbound plasma concentration of the interacting drug

- EMA allows the use of AUC/F_2 for correlation
- F_2 – the [inducer] leading to 2-fold increase of the baseline levels of enzyme
- Useful if $E_{\max}/EC_{50}$ undefined



Basic modeling

Figure 4: An Equation to Calculate the Predicted Ratio of the Victim Drug's AUC in the Presence and Absence of an Inducer for Basic Models of Induction

$$R_3 = 1 / [1 + (d \times E_{\max} \times 10 \times I_{\max,u}) / (EC_{50} + (10 \times I_{\max,u}))]$$

R₃ is the predicted ratio of the victim drug's AUC in the presence and absence of an inducer for basic models of enzyme induction.

d is the scaling factor and is assumed to be 1 unless supported by prior experience with the system used.

E_{max} is the maximum induction effect determined in vitro.

I_{max,u} is the maximal unbound plasma concentration of the interacting drug.*

EC₅₀ is the concentration causing half-maximal effect determined in vitro.

*Considering uncertainties in the protein binding measurements, the unbound fraction should be set to 1% if experimentally determined to be <1%.

If $R_3 \leq 0.8$, further investigation is warranted

Static mechanistic modeling

$$AUCR = \left(\frac{1}{[A_g \times B_g \times C_g] \times (1 - F_g) + F_g} \right) \times \left(\frac{1}{[A_h \times B_h \times C_h] \times f_m + (1 - f_m)} \right)$$

A is the effect of reversible inhibitions.

B is the effect of TDI.

C is the effect of induction.

F_g is the fraction available after intestinal metabolism.

f_m is the fraction of systemic clearance of the substrate mediated by the CYP enzyme that is subject to inhibition/induction. Each value can be estimated with the following equations:

	Gut	Liver
Reversible inhibition	$A_g = \frac{1}{1 + \frac{[I]_g}{K_i}}$	$A_h = \frac{1}{1 + \frac{[I]_h}{K_i}}$
Time-dependent inhibition	$B_g = \frac{k_{deg,g}}{k_{deg,g} + \frac{[I]_g \times k_{inact}}{[I]_g + K_I}}$	$B_h = \frac{k_{deg,h}}{k_{deg,h} + \frac{[I]_h \times k_{inact}}{[I]_h + K_I}}$
Induction	$C_g = 1 + \frac{d \cdot E_{max} \cdot [I]_g}{[I]_g + EC_{50}}$	$C_h = 1 + \frac{d \cdot E_{max} \cdot [I]_h}{[I]_h + EC_{50}}$

No net effect
model now:
Why??

If AUC ratio ≤ 0.8 , further investigation is warranted

Dynamic mechanistic modeling: PBPK

- Prediction of the DDI potential as a substrate or a perpetrator

In Vitro Metabolism- and Transporter- Mediated Drug-Drug Interaction Studies Guidance for Industry

DRAFT GUIDANCE

This guidance document is being distributed for comment purposes only.

Comments and suggestions regarding this draft document should be submitted within 90 days of publication in the *Federal Register* of the notice announcing the availability of the draft guidance. Submit electronic comments to <https://www.regulations.gov>. Submit written comments to the Division of Dockets Management (HFA-305), Food and Drug Administration, 5630 Fishers Lane, rm. 1061, Rockville, MD 20852. All comments should be identified with the docket number listed in the notice of availability that publishes in the *Federal Register*.

For questions regarding this draft document, contact (CDER) Office of Clinical Pharmacology, Guidance and Policy Team at CDER_OCP_GPT@fda.hhs.gov.

Provides a
framework

U.S. Department of Health and Human Services
Food and Drug Administration
Center for Drug Evaluation and Research (CDER)

October 2017
Clinical Pharmacology

10/24/17

Physiologically Based Pharmacokinetic Analyses — Format and Content Guidance for Industry

Prescriptive
guidelines

U.S. Department of Health and Human Services
Food and Drug Administration
Center for Drug Evaluation and Research (CDER)

August 2018
Clinical Pharmacology

Physiologically Based Pharmacokinetic Analyses — Format and Content
09/05/18

Unanswered questions

- When should I place the study?
- Do I need a negative control?
- Can I use my validated method for the concentration measurements?
- More complex questions....

What if there is down-regulation?

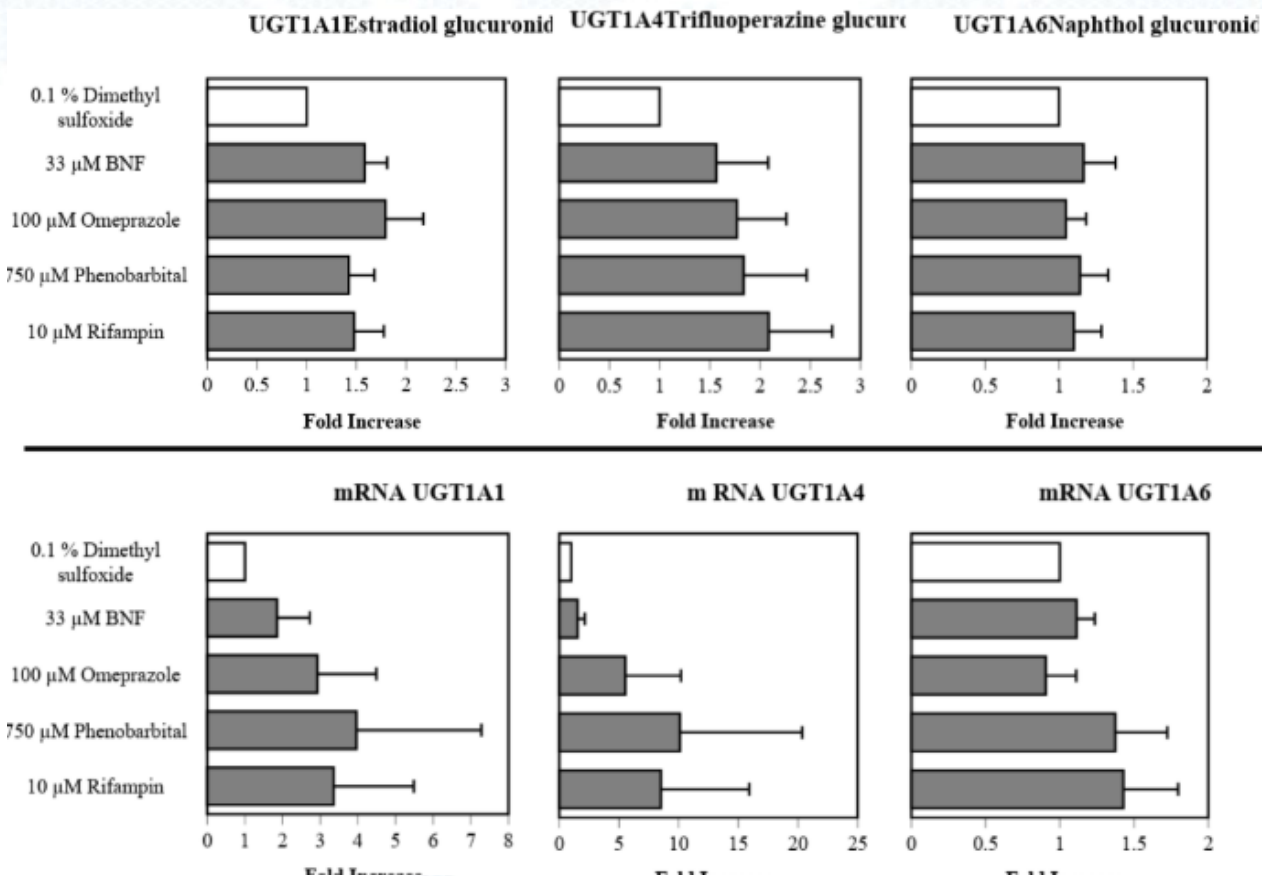
- Down-regulation does not equal clinical effects
- EMA cutoff: 50% decrease and concentration-dependence
- Evaluate relationship with
 - Mechanism of action
 - Cytotoxicity
 - Enzyme inhibition (activity measurements)
- Mechanistic studies may be needed to evaluate cause
 - May be a big undertaking
 - Scarcity of useful research in this area

IQ-IWG 2017:

Hariparsad et al. *Drug Metab. Dispos.* 2017, 45: 1049-1059.

Induction: UGT enzymes and transporters

- UGT induction usually less problematic in vivo than CYP, but...



FDA, 2017:
 “In vitro methods to evaluate induction of P-gp and other transporters are not well-established. Therefore the FDA does not currently recommend in vitro evaluation of transporter induction”

Conclusions

- Study designs should meet
 - Scientific needs
 - Regulatory needs
- A positive result (even from one donor)
 - Means some form of follow up is needed
 - Does not necessarily lead straight to the clinic
- Science is still science and some aspects are not perfect
- Make decisions carefully based on the program
- Client questions are rigorously discussed internally!

Acknowledgments

- Rebecca Campbell
- Madison Knapp
- Dr. Brian Ogilvie



www.xenotech.com

Featuring publications, posters, webinars and other useful resources.

Webinar Topic Request Form:

www.xenotech.com/scientific-resources/upcoming-webinars

XenoTech's Services & Products:

In Vitro ADMET PK & DDI

- Drug Transport
- Drug Metabolism
- Enzyme Inhibition & Induction
- Protein Binding
- Metabolite Identification
- ADME Screening
- Toxicology

In Vivo ADME/PK & Distribution

- QWBA
- Microautoradiography
- Excretion / Mass Balance
- Tissue Distribution
- Blood / Plasma & Lymphatic Partition Rate

Bioanalytical

Pharmacology

- In Vitro Ligand Binding & Radioreceptor Assays
- Immunoassays

Chemical Synthesis

- Radiolabeled Synthesis
- Metabolite Synthesis
- Peptide Synthesis

Consulting...

Cellular Products

- Hepatocytes (Cryo/Fresh, Genotyped...)
- Non-Parenchymal Cells (Kupffer Cells)

Subcellular Fractions

- Liver Microsomes
- S9 Fractions
- Cytosol
- Homogenate
- Lysosomes & Tritosomes
- Mitochondria
- Extrahepatic Fractions

Custom Products

- Various Species, Tissues & Preparations

Research Biobank

- Normal & Diseased Tissue Samples

Recombinant Enzymes

Substrates & Metabolites

Metabolite Production Kits

JCRB Cell Lines...

info@xenotechllc.com

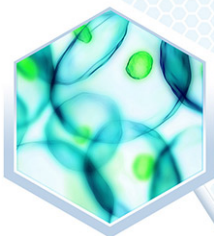


A BioIVT Company

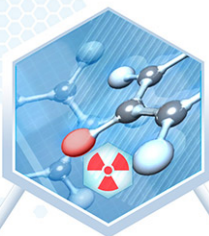
1-877-588-7530 | +913.438.7450 • WWW.XENOTECH.COM

PROVEN GLOBAL CONTRACT RESEARCH EXPERTISE
FROM DISCOVERY THROUGH CLINICAL SUPPORT

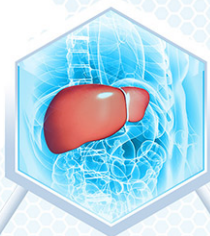
Cell & Tissue-Based Products



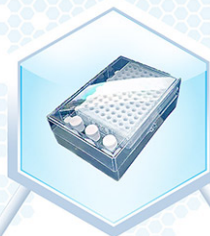
Radiolabeling



in vitro ADMET & Pharmacology



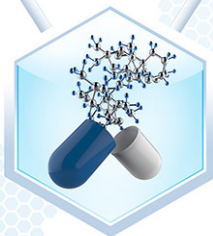
Metabolite ID & Production



Screening



API Manufacturing



in vivo ADMET & QWBA



Bioanalytical



EXPERTISE • EFFICIENCY • SUPPORT • PRECISION

Thank You!