

# Drug Metabolism: Reaction Phenotyping

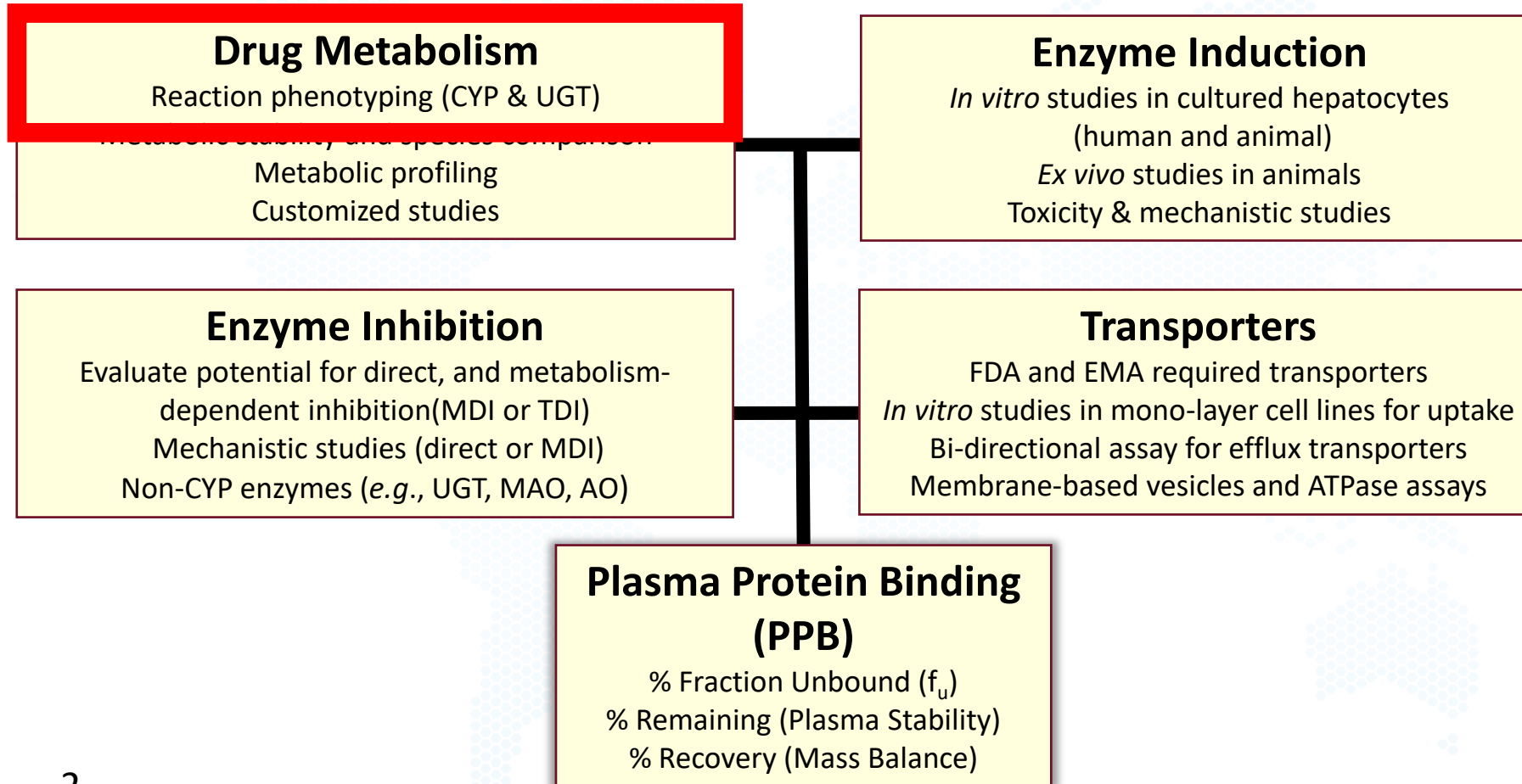
Andrew G. Taylor, Ph.D.

Manager, Technical Support for Services

XenoTech

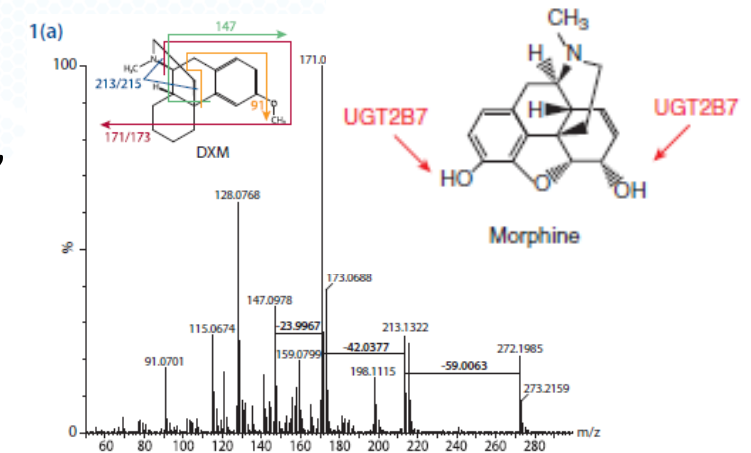
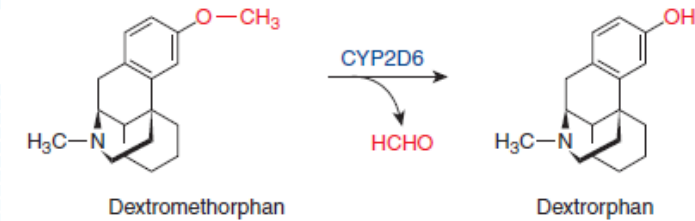


# In vitro DDI Services



# Drug Metabolism Studies

- Reaction phenotyping  
Identification of enzymes involved in the metabolism of a drug candidate
- Metabolic stability  
Comparison of metabolic stability and intrinsic clearance among drug candidates (microsomes, S9, hepatocytes)
- Metabolic profiling (AS)  
Identification of the metabolites
- Species comparison  
Provides information to select appropriate nonclinical species





## Drug Metabolism Studies

It is important to identify the enzymes that significantly contribute to the metabolism of a drug candidate and to evaluate the potential of a drug candidate to be a:

- **victim** drug (a drug which is cleared by a main route of elimination)
- **perpetrator** drug (a factor or drug which alters the clearance of a victim drug)

Test systems used depend on enzymes to be evaluated (e.g., HLM for CYP, cytosol for Phase II enzymes, S9 or hepatocytes for CYP + Phase II)

A BioIVT Company

# CYP Reaction Phenotyping: Requirements

2020 FDA

Recommendations:

Use validated analytical methods

**Triplicate determinations**

Perform at least two (rCYP & chemical inhibition) tests

CYP1A2, 2B6, 2C8, 2C9, 2C19, 2D6 and 3A.  
Other enzymes as needed

Include positive controls

In Vitro Drug Interaction Studies —  
Cytochrome P450 Enzyme- and  
Transporter-Mediated Drug Interactions  
Guidance for Industry

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

January 2020  
Clinical Pharmacology





## CYP Reaction Phenotyping: Standard Design

- **Preliminary study**
  - Time and protein
- **Optional follow-up, depending on preliminary results and/or sponsor:**
  - rCYP
  - Chemical inhibition (usually HLM)
  - Correlation analysis (rare)

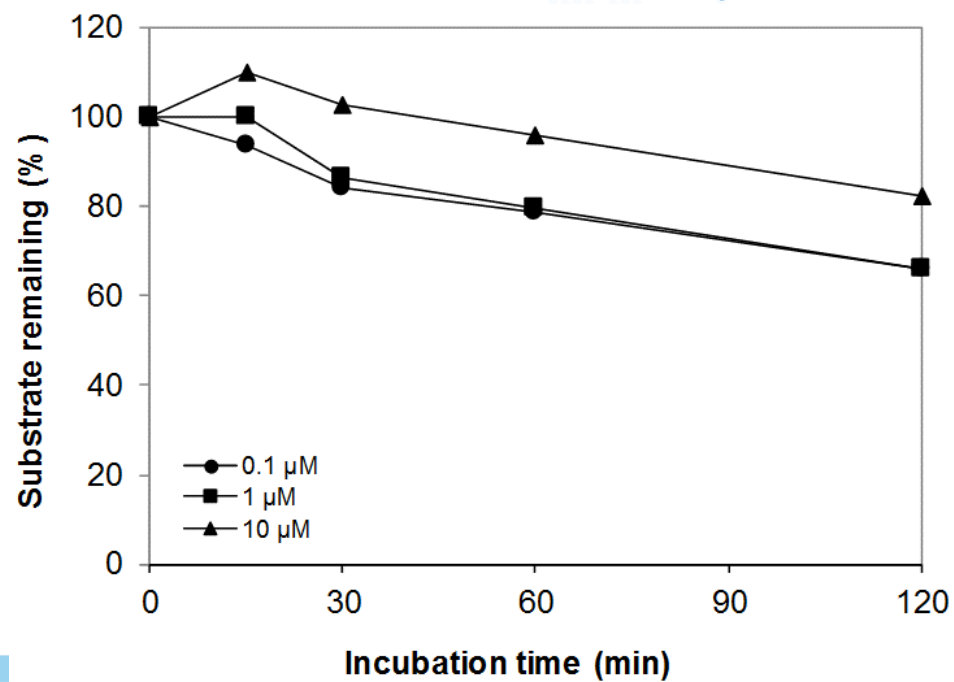


A BioIVT Company

# CYP Reaction Phenotyping: Preliminary Evaluation

## Design:

- 3 concentrations of test article
- 3 concentrations test system
- 0 + 4 incubation time points



- Establish initial rate conditions with pooled HLM
- Substrate loss (or metabolite formation) is proportional with respect to incubation time and protein concentration
- Incubation times can be long (e.g., 0.5 to 2 hr)



# Recombinant CYP (rCYP) & Chemical Inhibition

## Two complementary methods



# CYP Reaction Phenotyping: Recombinant CYP (rCYP)

- Recombinant enzymes: human enzymes expressed in bacteria or insect cells
- Enzyme expression is different than in vivo (e.g., HLM, hepatocytes)

## Design:

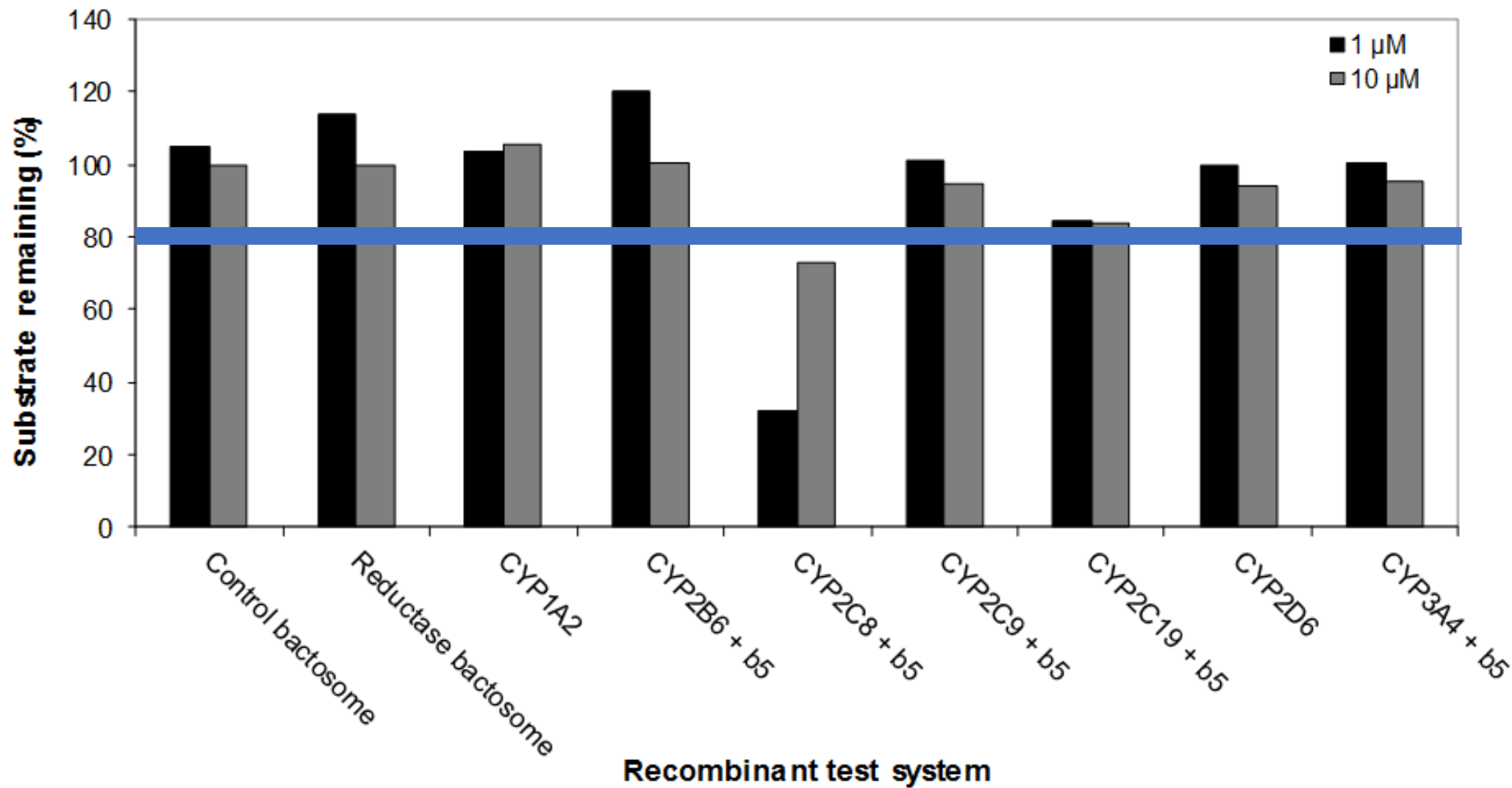
- 1 or 2 concentrations of test article
- Panel of CYP enzymes (7 standard)
- 1 concentration rCYP (time and protein samples with HLM included for comparison)
- 1 incubation time point

RECOMBINANT ENZYMES





A BioIVT Company

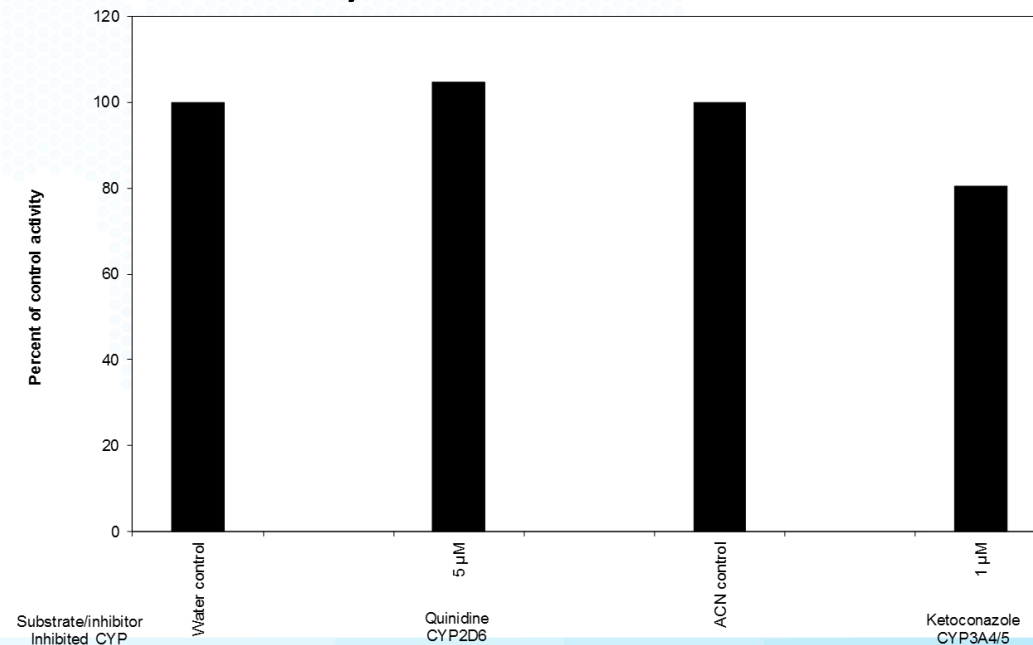


# CYP Reaction Phenotyping: Chemical Inhibition

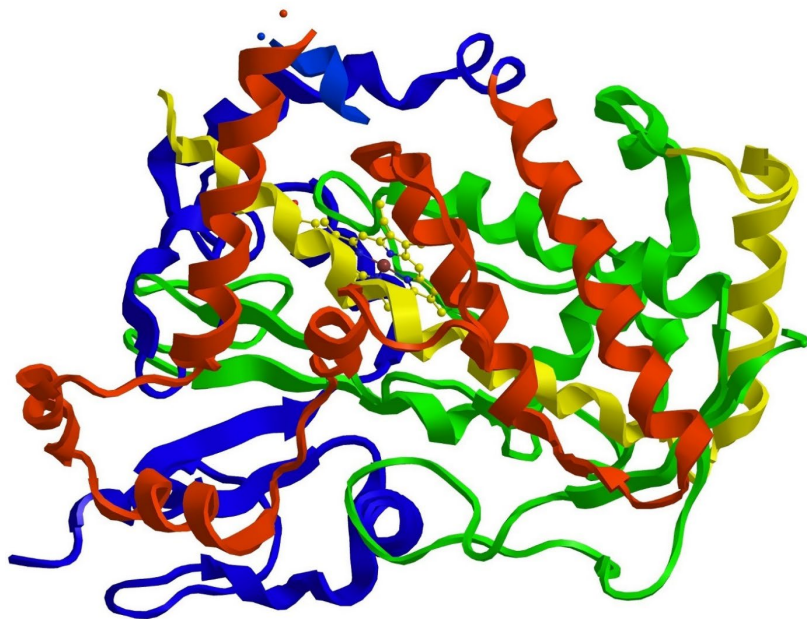
- Use chemical inhibitors specific to CYP enzymes to determine CYP enzymes responsible for clearance
- Typically use metabolism-dependent CYP inhibitors (longer lasting inhibition, better for longer incubation times)

## Design:

- 1 or 2 concentrations of test article
- Panel of CYP inhibitors
- 1 concentration HLM
- 1 incubation time point



# Reaction Phenotyping: Other Enzymes



Approach may be limited depending on enzymes evaluated

- Is recombinant test system available?
- How specific are chemical inhibitors?

For UGT enzymes:

- Recombinant human UGT enzymes
- Chemical inhibition (limited to UGT1A1, UGT1A4, UGT1A9, UGT2B10)
- Correlation analysis is not recommended due to minimal variability in UGT enzyme activity among individuals in kit.



A BioIVT Company

# XT Products (Metabolism/Reaction Phenotyping)

## Subcell Fractions:

- Microsomes, S9, and/or Cytosol
- Multiple tissues (liver and extrahepatic)
- Human, Monkey, Rat, Mouse, Dog

## Recombinant Enzymes (Distributed):

- Cypex
- Phase I and Phase II enzymes

## Primary Hepatocytes

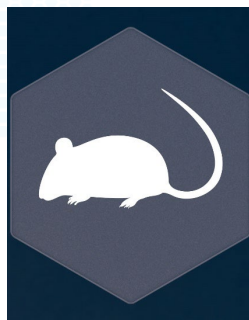
- Pooled donors
- Suspension (2-6 hours)
- Attaching (longer than 6 hours)
- Multiple species

## Reaction Phenotyping Kit (Patented)

- Well-characterized individual donor HLMs

## Support Reagents

- Rapid Start – NADPH regenerating system
  - For subcellular fractions
- Optimized hepatocyte media
  - Opti: Thaw, Plate, Incubate, and Culture
- Control metabolites (Distributed)
  - Cypex
- Technical Support
  - Recommended Protocols
  - Troubleshooting





## Thank you for watching!

Questions? Get in touch through the Contact Us tab on our website

Please contact your regional account manager if you are interested in a placing a reaction phenotyping or other drug metabolism study for your drug's development

