



# XENOTECH

A BioIVT Company

## DISCOVERY

Discovery Plus™  
Induction

*XenoTech's Discovery Plus™ Induction service provides fold change data in 10 days or less.*

XenoTech's Discovery Plus™ Induction is more than just a screening service. Our 96-well plate format allows up to 10 NCEs at 3 concentrations and positive controls in triplicate wells. Discovery Plus™ Induction is a non-GLP service that provides customers with CYP3A4 mRNA fold change data in less than 10 days. This service is best utilized for compounds that are in the early discovery / development phase. For compounds that will be submitted for regulatory agency approval, please see our definitive CYP Induction service flyer.

### *Discovery Plus™ Induction Overview*

Discovery Plus™ Induction uses XenoTech's specialized screening design to evaluate CYP3A4 fold change in mRNA. This service yields data presented in a tabular summary of fold change, compared to the vehicle for CYP3A4 (other CYPs are available). Pre-characterized cryopreserved primary hepatocytes are treated in a 96-well plate format for ~24 hours with vehicle, NCEs at 3 concentrations and positive control(s) in triplicate wells. RNA is analyzed utilizing qRT-PCR technology by the cDNA ddC<sub>i</sub> method (samples normalized to endogenous house keeping gene). The fold induction is then calculated as test compound over the vehicle control. Our Discovery Plus™ design offers the most value in the industry, providing you important data quickly, at a very attractive price point.

### Discovery Plus™ Induction Highlights

- Measures mRNA fold induction in CYP3A4 (Other CYPs may be substituted)
- Utilizes qRT-PCR techniques for mRNA measurement
- Design allows up to 10 NCEs at 3 concentrations and positive controls, performed in triplicate
- Data presented in tabular summary of fold induction
- Offered as a non-GLP service

### Benefits

- Dynamic range of fold induction data for your compound
- Rapid turnaround (10 days per 10 compounds)
- Results can aid in the design of any future induction study for your compound(s)
- Budget-friendly

### Follow-up Services Available (Additional Charge)

- Definitive enzyme induction assessment
- Consultation on results

### Additional Discovery Plus™ Services

- **Metabolic Stability**
- **Drug Inhibition**
- **Plasma Protein Binding**

Contact us to learn more about XenoTech's full suite of **Discovery Plus™** Services. Visit us online at [www.xenotech.com](http://www.xenotech.com) or call **913.438.7450**