



bioanalytical services



# XENO TECH

A BioIVT Company

Non-Regulated  
Small Molecule  
Bioanalysis

*XenoTech small molecule bioanalysis services generate drug exposure data to aid key decision making in pharmaceutical development*

Bioanalysis is a sub-discipline of analytical chemistry providing a quantitative measurement of a drug and / or its metabolite(s) in various biological matrices. XenoTech provides non-regulated small molecule bioanalysis, generating drug exposure data to aid key decision making in pharmaceutical development. Our analytical experts have supported in-house *in vitro* services for over 15 years with more than 35 validated marker assays for XenoTech products and services. They are responsible for analytical method development, qualification, validation and transfer for all *in vitro* services, as well as dose solution analysis support and metabolite profiling and characterization. Our analytical team has a combined 100 years of CRO and pharmaceutical experience spanning multiple disciplines including *in vivo* bioanalysis support for multiple clients. Let XenoTech's experts handle your bioanalytical needs, delivering consistently accurate and reliable data with rapid turnaround times.

### Small molecule non-regulated capabilities

XenoTech's knowledgeable staff can provide quality, non-GLP results in 2-4 business days. Highlights of XenoTech's bioanalytical services include:

- Protein precipitation extraction (others as required)
- Plasma, serum, blood and urine analysis
- Development of single or multiple analyte methods to support parent and metabolite(s) analysis or cassette analysis
- Method qualification (optional)
- Duplicate standard curves (bracketed samples)
- Flexible standard curve ranges up to 5000-fold (modify as needed)
- Sample analysis by LC-MS/MS
- Electronic data summary outlining analytical method and sample preparation description, calibration curve and calculated sample concentrations
- Rapid turnaround times

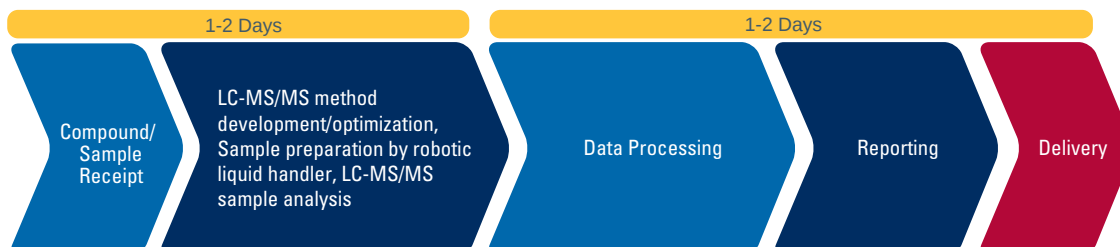
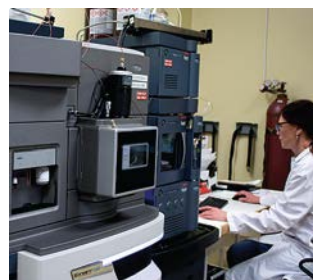


*Where Analytical and ADME Expertise Collide!*



### Equipment

- Automated liquid handlers (Tecan) for consistent sample processing
- Xevo G2-XS Qtof/Tof, Xevo TQ-S triple quadrupole, AB Sciex API 2000 and API 3000 triple quadrupole, a 4500 triple quadrupole, and API 4000 and 5500 QTrap MS/MS systems with Shimadzu Nexera or Waters Acquity UHPLC systems for high sensitivity analysis



*Timeline represents a typical non-regulated small molecule bioanalysis study. Timeline varies based on sample number.*