



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. Sekisui 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 Sekisui 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 Sekisui XenoTech's experts handle your bioanalytical needs, delivering consistently accurate and reliable data with rapid turnaround times.

Small molecule non-regulated capabilities

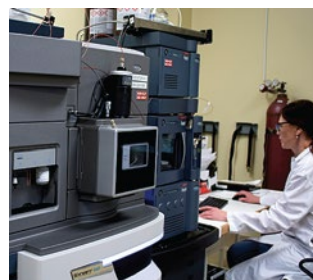
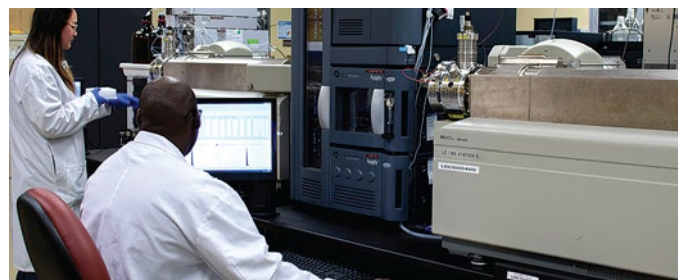
Sekisui XenoTech's knowledgeable staff can provide quality, non-GLP results in 2-4 business days. Highlights of XenoTech's bioanalytical services include:

- 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

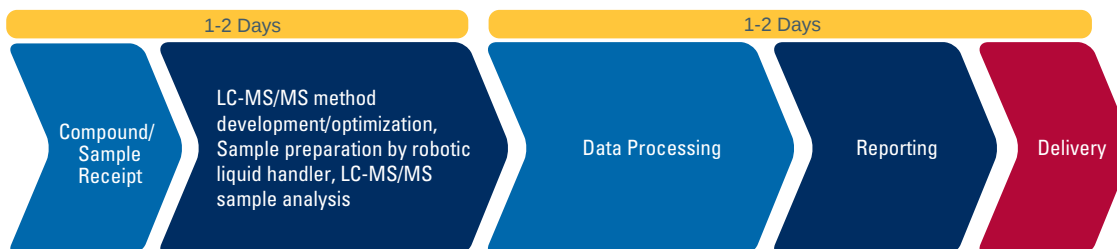


Equipment

- 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



Where Analytical and ADME Expertise Collide!



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