



Bioassay

Bioanalytical Contract Laboratory

A Reaction Biology Company

Bioassay GmbH specializes in the development, validation, and execution of biological assays in the context of pharmaceutical and chemical product development and quality control.

Overview Services Client Offerings

Potency Assay "Power House" offering complete value chain coverage

Potency Assays

- GMP potency assays of biologicals
- Focus on in-vivo and cell based potency assays
- Full value chain offered:
 - Development
 - Validation
 - Routine

Value Add

- Tailor-made Mode of Action Assays (NBEs)
- Proof of Concept
- Product/Process Development
- Biosimilar Studies
- Batch Release

Functional Bioassays

- Binding Assays
- Neutralisation AB-Assays
- ADCC and CDC
- Primary Blood Cell Assays

Value Add

- Compound Selection
- Characterization
- (Pre-) Clinical Safety

GLP & Other

- GLP Tox Safety & in-Chemico
- Pharmacokinetics
- HPLC
- Focus areas include:
 - Pharma
 - Chemicals
 - Cosmetics

Value Add

- Compound Selection
- Registration Support
- Complementary in-vitro and in-vivo services

PSL*

- Cooperation partner specialized in
 - Customized peptide synthesis
 - Peptide specific antibodies
- Production of peptides as substrates for enzymes
- Peptide modification

* <https://www.peptid.de/en>
(Acquired by Reaction Biology in 2022)

