

Preclinical & GLP Services

We execute *in vivo* and *in vitro* preclinical studies to evaluate discovery leads efficacy, toxicity and safety before entering clinical trials.



Animal Welfare

The facilities comply with the highest standards for animal welfare.



Expert team

Each study is supervised by a specialized study director.



High Transparency

We offer a full access to each stage of the study from the plan to the final report.



Certified Process

Each study is conducted in Non-GLP and GLP environments. Our QMS is ISO certified.

Discovery Leads

In vitro and *In vivo* models

- **In vitro** ADME
- **In vivo** DMPK & Tox
- Disease models
- Dose assessment

Animals: rodents

On the way to clinics

GLP regulations

- Nonclinical safety
- Toxicity studies
- Pharmacokinetics
- Histopathology & metabolism

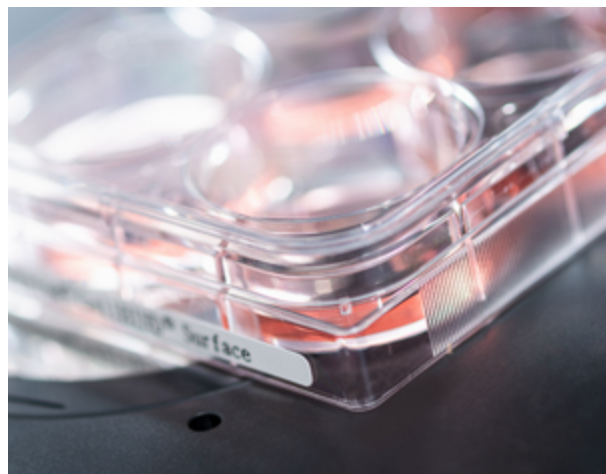
Animals: rodents + larger animals for veterinarian evaluations

Disease Models

We have various disease models to test potential treatment approach in the domains of **Neurology, oncology, nephrology and cardiometabolic.**

In vitro and *in vivo* rodent models:

- Various tumor models
- Parkinson
- Schizophrenia
- Fragile X disease
- NASH
- Diabetes
- Myocardiac infarction
- ...



Pharmacokinetics

We have **liver, lung, kidney, intestine, breast and brain models** to study how the body interacts with administered substances. Below are some of the tests we may perform:

In vitro

Pharmacology

- Migration
- Invasion
- Dose finding
- Efficacy

ADMET

- Solubility
- Permeability
- CYP450 assays
- Biocompatibility

In vivo

Pharmacology

- Efficacy
- Dose frequency
- MTD

DMPK

- Bioavailability
- Drug distribution
- Clearance

Tox & Safety

We perform tox and safety assays in **rodent and larger animals.**

In vitro

- Cyto-, Geno-, Nephro- and hepatotoxicity
- Single dose toxicity
- Toxicokinetics

In vivo

- Single dose toxicity
- Repeated dose toxicity
- Toxicokinetics
- Acute/sub-chronic and chronic toxicity
- Tolerance testing
- Sensitization testing



Learn more on www.eurogentec.com