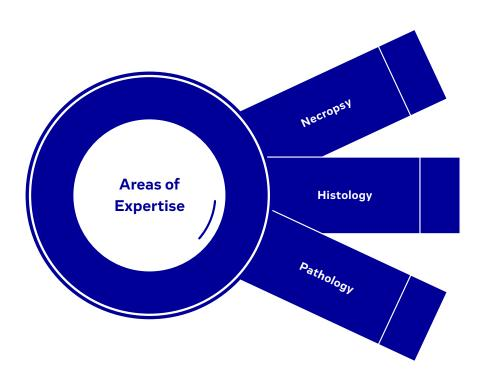


Pathology

- ▶ 25+ years' experience with a broad skillset in discovery and clinical-enabling pathology
- GLP certified with highest quality standards and an impeccable regulatory inspection history
- ▶ Unparalleled flexibility with all capabilities in-house and under one roof
- ► DACVP/DECVP, MRCPATH, DECLAM, DECVCP qualified pathologists/ clinical pathologists



Contact our experts info@evotec.com | www.evotec.com



Necropsy

- ➤ Tissue collection according to standard or customised protocols, including special procedures for immunohistochemistry, electron microscopy, neuropathology, gene expression studies, peripheral blood lymphocytes sampling for electron microscopy examination
- ► Fixation, including whole body perfusion in rodents and tissue freezing fixation
- ▶ Macroscopic and organ weight data recording
- Gross pathology reporting and interpretation
- Gross pathology photographic documentation

Histology and electron microscopy

- ► Trimming, according to standard and special procedures: enhanced brain and heart trimming
- Processing to wax and resin for light or electron microscopy
- ▶ Bone tissue decalcification
- Cryomicrotomy (frozen sectioning)
- ▶ Special histochemical stains
- ▶ Immunohistochemistry
- Grid preparation for transmission electron microscopy
- ► Peripheral lymphocytes preparation for electron microscopy
- Morphometry
- ▶ Packaging for archive of residual tissues
- ► National and international transportation and documentation

Pathology

- Pathology evaluation, interpretation, integration and reporting
- Expertise covering toxicologic pathology in drug development, research animal model validation and characterisation with focus on translation to human
- ▶ Neurovacuolation studies (Olney's test)
- International standards in terminology and diagnostic criteria addressing regulatory requirements and expectations (INHAND)
- ▶ Peer review

Pathology services applications

- Preclinical regulatory studies (acute and repeated dose studies), investigative studies (e.g. targeting attrition issues)
- ▶ Discovery research (animal model of disease: study design, validation and characterisation)
- ▶ Biomarkers identification and validation packages
- Ultra structural pathology
- Consultancy

Facilities and capabilities

- ► Rodent, non-rodent species necropsy rooms, with fully computerised working stations
- GLP-certified histology laboratory facilities and instruments
- ▶ Transmission Electron microscopy unit
- ▶ Slide scanner

Expertise

- ▶ Histopathology technicians, with extensive experience in preclinical pathology studies in GLP environments and in experimental animal models
- ► Pathologists with extensive expertise in drug development and research with competencies in risk assessment and translational research
- Clinical pathologist with expertise in biomarker setup, validation and interpretation in translational research

Regulatory compliance

Services conducted under Good Laboratory Practice (GLP) in regulatory-compliant facilities