
Fully Integrated Development: Discovery Through Clinical Proof Of Concept

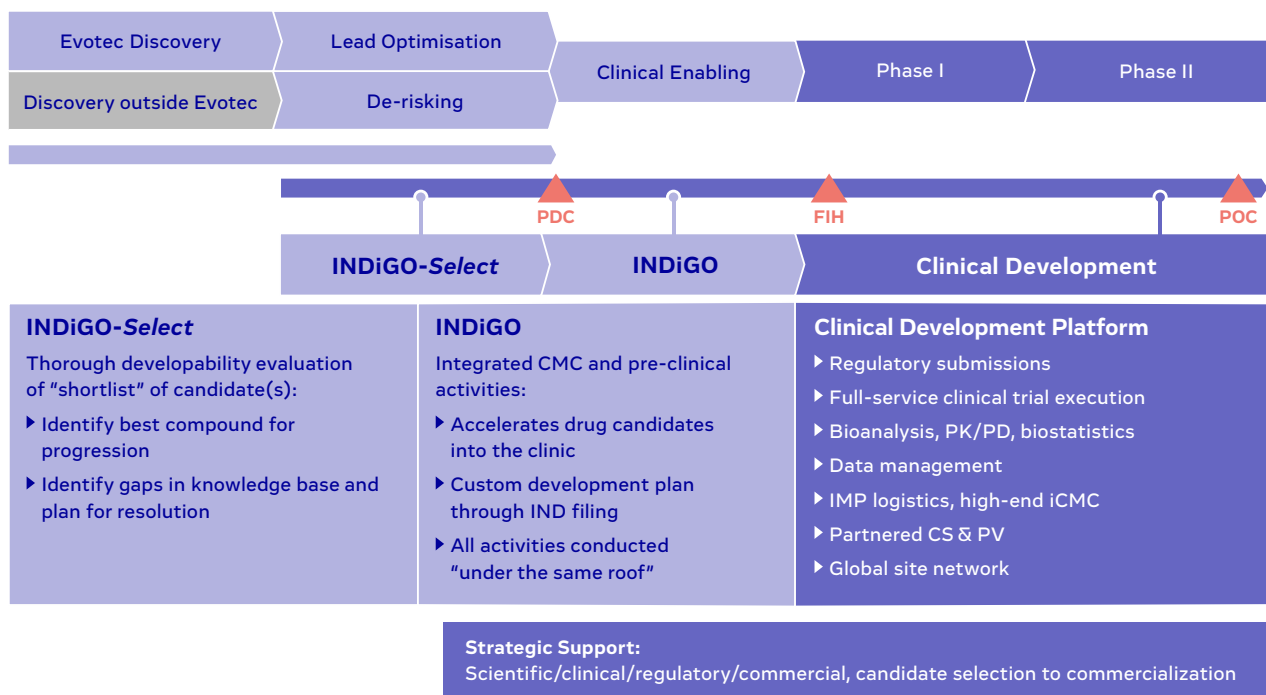
Evotec is a fully-integrated drug discovery and development partner, providing a comprehensive source of high quality scientific solutions with consultant-level expertise to our partners, which includes fully integrated research and development as well as clinical development through to Proof of Concept.

Our Clinical Development team offers:

- ▶ Deep and broad experience, derived from the perspective of the sponsor as well as the CRO, across all phases of development and multiple therapeutic areas
- ▶ Seamless integration with our pre-clinical development platform
- ▶ Industry-leading efficiency, which translates to delivery on-time and on-budget while achieving a significant reduction in time from pre-clinical to phase 1 initiation
- ▶ Regulatory and clinical experience in all global settings
- ▶ Strategic and operational support for global clinical development, regulatory, and commercial strategy, including target product profile & clinical plan development
- ▶ Strategic partnerships with established clinical sites in Italy, France, Germany, and North America - and the additional flexibility to work with any qualified site or site network our partners prefer
- ▶ Support for both healthy volunteer and patient-recruited early phase trials
- ▶ The capabilities of a large clinical CRO, with the experience and flexibility of the dedicated clinical teams found in pharmaceutical companies



Integrated Platforms Supporting All Stages Of Drug Development



While our Clinical Development team is currently focused on Phase I and Phase II trials, Evotec is actively investing in and expanding our clinical development team to a level that will support pivotal studies for marketing authorization.

Evotec Clinical Development solutions integrates seamlessly with Evotec's other core offerings

- ▶ Integrated clinical-enabling drug development
- ▶ Regulatory strategy and dossier preparation/ submission
- ▶ Full-scope clinical trial planning and execution
- ▶ Innovative translational biomarkers, ADMET and integrated CMC activities
- ▶ Clinical supplies and logistics support

The right trial at the right time can have a lasting impact on a drug's life cycle. Please contact us to discuss how we work with our partners to provide a clinical development strategy that delivers all the way through to commercialisation.

Project Team Structure

The Evotec Clinical Development leadership team is unique in its approach to trial management and communication, based on our extensive experience and coupled with areas of unmet need in the industry.

For instance, every trial for every partner, no matter how large or small, is supported by a Joint Oversight Committee to ensure quick issue identification, decision making and resolution. In addition, we employ a "Quality by Design" approach to risk management from protocol development through completion of the clinical study report.

We believe in transparency and collaboration with our partners, and chose key performance indicators that translate to meaningful representations of the highest quality deliverables.