

4 Ways De-Risking Maximizes Compound Value

Most new drugs fail because of ADME/Tox¹

With early in vitro testing, you can preemptively demonstrate low risk data and **add considerable value to your compound.**



These are facts that industry **must know and leverage** to maximize value. De-risking showcases your compound's performance by ...

1

Satisfying Regulatory Requirements



Avoid the **50% risk of IND failure** attributable to inadequate ADME



Provide **required** regulatory information

2

Avoiding Unnecessary Expenses and Approval Delays

- Each day of delay costs \$1 million or more, and over 60% of sponsors report delays of 6 months or longer.^{2,3} De-risking equates to approx. **\$230 million** of added value and that does not include the cost of additional studies and resources!
- **Prevent the losses** of time and other resources caused by clinical failures and additional studies



3

Optimizing Future Development Success



Provide insights into future in vivo study performance



Identify the **best test species** for tox studies



Hone **first in-human** dose selection



Eliminate the need for, or elevate the design of, follow-up clinical studies

4

Ensuring Patient Safety

- Reveal potentially **dangerous drug-drug interactions** before administration to humans



Investing in **high-quality ADME data** helps maximize the value of a compound. Get the return on investment you need with a custom plan from XenoTech.

Visit www.xenotech.com/contact for more information.

 **XENOTECH**
A BioIVT Company

¹ Faqi, A.S., editor. A comprehensive guide to toxicology in nonclinical drug development. Academic Press. Published 2016 November 3.

² The University of Texas at Austin. Old Drug Standards Delay New Drug Approvals. Published 2020 June 19.

³ XenoTech. Industry Survey. Conducted 2021.