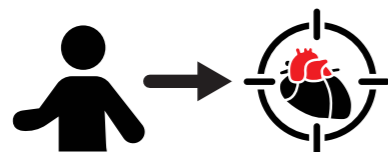


Therapeutic monoclonal antibodies

Overcoming challenges to accelerate drug development

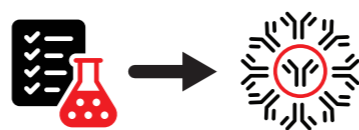
Monoclonal antibodies (mAbs) and antibody-drug conjugates (ADCs) have revolutionized the life sciences, and are crucial components of modern medicine, contributing to highly effective, personalized patient care. In less than 40 years, therapeutic mAbs have transitioned into powerful tools, transforming the lives of patients with previously untreatable diseases, including autoimmune conditions and various cancers. Hundreds of thousands of people have benefitted from these therapeutics since their entry into the market, and new mAbs and ADCs are currently in development, with many already in the final approval stages. Two Nobel Prizes in Physiology or Medicine have been awarded to scientists who made substantial contributions to mAb development, in 1984 and 2018, demonstrating their potential for improving human health.^{1,2}

Although a rapidly expanding field, the overall discovery and development of a single novel mAb is slow, with 17 years the average lag time between initial research and market launch of any drug.³ mAb development especially is still plagued by a number of challenges, both in and out of the lab.



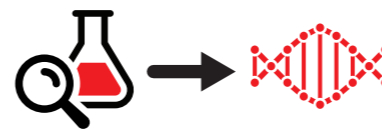
Target Selection And Validation

Identifying a promising target requires extensive and time-consuming research and a deep understanding of underlying disease biology tested in relevant disease models.¹⁰



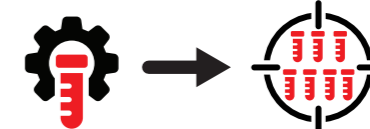
Lead Identification And Optimization

mAb binding affinity, specificity, stability, and pharmacokinetic properties must be optimized through various complex and time-consuming approaches. Candidates must also be characterized in a series of *in vivo* experiments to exclude any immunogenic glycosylation profiles.¹¹



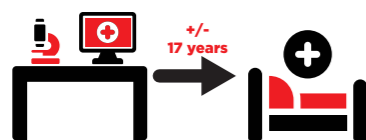
Cell Line Development

Identifying and selecting high-producing cell clones from the transfected cell population is a crucial step in mAb development, requiring quick identification of clones that exhibit desirable characteristics, such as high antibody productivity and stability.¹⁰



Process Scalability

Cell-based expression systems are costly and inefficient, and challenging to scale up. Analytical tools often have limited throughput, while downstream processes needed to remove biological contaminants can degrade the mAb, leading to product instability and loss and resulting in low yields.¹⁰

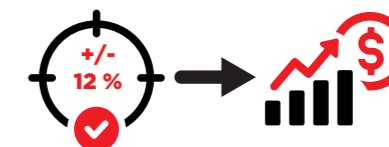


Time From Bench To Bedside

Antibody drug discovery and development is slow. Modern cell line development (e.g. clone selection) and process development (e.g. media development and feed optimization), as well as analytical solutions (e.g. clone characterization), are needed to shorten the time-to-market.

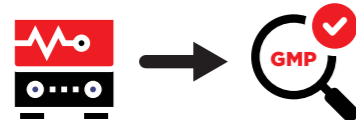
Fast-tracking mAb development

Partnering with trusted solution providers will remain a steadfast approach to combatting these challenges. Automation platforms and precision instruments can streamline mAb and ADC development - from early-stage research to analysis - helping to ensure a steady pipeline of therapeutic mAbs reach patients in need.



Success Rate

Only 12% of drugs that enter clinical trials result in an approved medicine, and each approved drug costs an estimated average of USD 2.6 billion in R&D.⁴



GMP Regulations

WHO guidance on therapeutic mAbs⁷ follows GMP regulations by the US FDA,⁸ EMA,⁹ and other regulatory agencies. Using GMP-enabling instruments from trusted providers throughout the development process can ease the transition to manufacturing.



Data Reliability In The Lab

Therapeutic mAbs require laboratory techniques and instruments that deliver accurate and reliable results across discovery, development and manufacturing, and must adhere to FDA and EMA guidelines.⁵



Instrument-To-Instrument Variability

Instrument variability directly impacts the reliability and validity of findings, and is a consideration in the ICH's analytical procedure development guidelines.⁶ A low instrument-to-instrument variability is needed to generate high quality data across multiple labs.



Cost Of Instrumentation

Laboratory automation increases speed and improves consistency of results, but high-quality instruments aren't always accessible to some organizations.



Get in touch to find out how our products can help you to overcome manual barriers, minimize downtime and reduce time-to-market.

beckman.com



© 2024 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

2024-GBL-EN-105423-v1

ICH = International Council for Harmonisation of Technical Requirements for Pharmaceuticals for Human Use, EMA = European Medicines Agency, GMP = Good Manufacturing Practices, WHO = World Health Organization, FDA = Food & Drug Administration

References

1. The Nobel Prize. The Nobel Prize in Physiology or Medicine 2018 - press release. (2018). <https://www.nobelprize.org/prizes/medicine/2018/press-release/>.
2. The Nobel Prize. The Nobel Prize in Physiology or Medicine 1984 - press release. (1984). <https://www.nobelprize.org/prizes/medicine/1984/press-release/>.
3. Morris ZS, Wooding S, Grant J. The answer is 17 years, what is the question: Understanding time lags in translational research. *J R Soc Med*. 2011;104(12):510-520. doi:10.1258/rsm.2011.10180
4. PhRMA. (2016). Modernizing Drug Discovery, Development and Approval. Accessed February 9, 2024. <https://phrma.org/en/resource-center/Topics/Research-and-Development/Modernizing-Drug-Discovery-Development-and-Approval>.
5. Alhazmi HA, Albratty M. Analytical Techniques for the Characterization and Quantification of Monoclonal Antibodies. *J Pharm*. 2023;16(2). doi:10.3390/ph16020291.
6. European Medicines Agency. (2023). ICH guideline Q14 on analytical procedure development. Accessed February 12, 2024. <https://www.ema.europa.eu/en/ich-q14-analytical-procedure-development-scientific-guideline>.
7. World Health Organization. (2022). Guidelines for the production and quality control of monoclonal antibodies and related products intended for medicinal use. Accessed February 12, 2024. <https://www.who.int/publications/m/item/guideline-for-the-safe-production-and-quality-control-of-monoclonal-antibodies>.
8. US Food & Drug Administration. (2023). Current Good Manufacturing Practice (CGMP) Regulations. Accessed February 12, 2024. <https://www.fda.gov/drugs/pharmaceutical-quality-resources/current-good-manufacturing-practice-cgmp-regulations>.
9. European Medicines Agency. Good manufacturing practice. (2023). Accessed February 12, 2024. <https://www.ema.europa.eu/en/human-regulatory-overview/research-and-development/compliance-research-and-development/good-manufacturing-practice>.
10. Sifniotis V, Cruz E, Eroglu B, Kayser V. Current Advancements in Addressing Key Challenges of Therapeutic Antibody Design, Manufacture, and Formulation. *Antibodies* 2019, Vol 8, Page 36. 2019;8(2):36. doi:10.3390/antib8020036.
11. Wolf B, Piksa M, Beley I, et al. Therapeutic antibody glycosylation impacts antigen recognition and immunogenicity. *Immunology*. 2022;166(3):380-407. doi:10.1111/imm.13481.