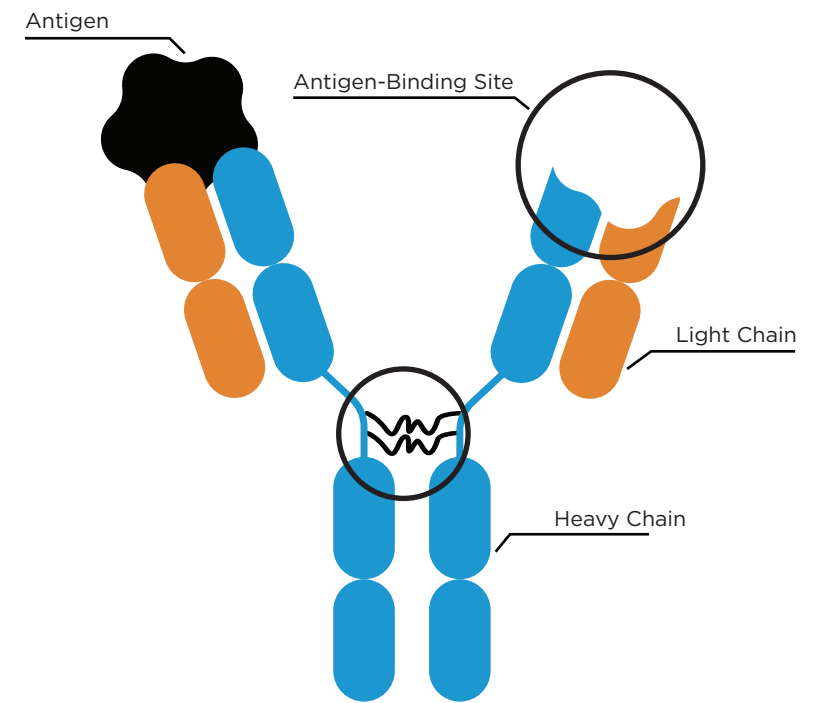


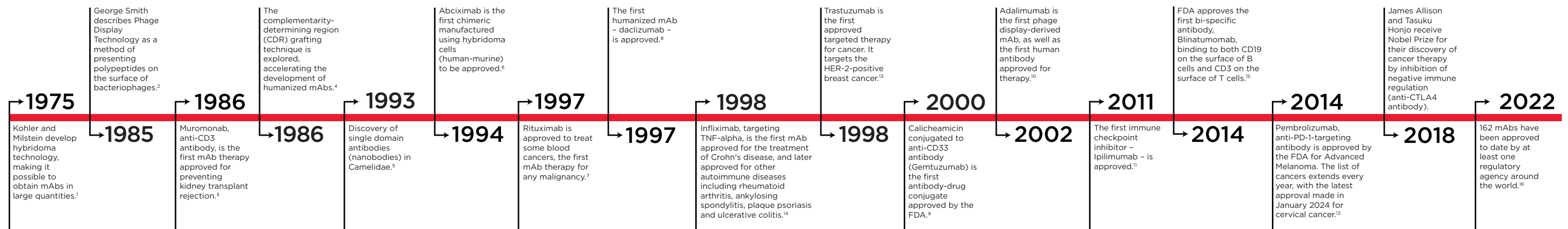
Monoclonal Antibodies: Revolutionizing Medicine

What Are Antibodies?

Monoclonal antibodies (mAbs) have revolutionized the field of therapeutics, providing highly specific and targeted treatments for a wide range of diseases. Beginning with the advent of hybridoma technology, the generation of mAbs with defined antigen specificity became possible. The subsequent advancements in antibody engineering, including the development of phage display libraries, transgenic animals, and in vitro selection techniques, expanded the repertoire of mAbs. These innovations allowed for the generation of antibodies with enhanced binding affinity, altered effector functions, and improved pharmacokinetic properties. This poster presents a chronological journey through key milestones, highlighting the scientific breakthroughs that have propelled the field forward.



Discovery Milestones



Learn more about newest trends in the discovery, development and manufacturing of mAbs. www.beckman.com/monoclonal-antibodies

References:

1. Köhler G, Milstein C. Continuous cultures of fused cells secreting antibody of predefined specificity. *Nat.* 1975;256(5517):495-497
2. Liu JKH. The history of monoclonal antibody development - Progress, remaining challenges and future innovations. *Ann Med Surg.* 2014;3(4):113
3. Smith G.P. Filamentous fusion phage: novel expression vectors that display cloned antigens on the virion surface. *Science.* 1985;228:1315-1317
4. Jones PT, Dear PH, Foote J, et al. Replacing the complementarity-determining regions in a human antibody with those from a mouse. *Nat.* 1986;321(6069):522-525
5. Hamers-Casterman C, Atarhouch T, Muyldermans S, et al. (1993) Naturally occurring antibodies devoid of light chains. *Nature* 363:446-448
6. Lu RM, Hwang YC, Liu J, et al. Development of therapeutic antibodies for the treatment of diseases. *J Biomed Sci.* 2020;27(1):1-30
7. Grillo-Lopez A, White C, Dallaire B, et al. Rituximab: the first monoclonal antibody approved for the treatment of lymphoma. *Curr Pharm Biotechnol.* 2000;1(1):1-9
8. Siegel JP. BLA approval letter: Zenapax (daclizumab BLA 97-0736). US Food and Drug Administration website. <http://www.fda.gov/downloads/Drugs/DevelopmentApprovalProcess/HowDrugsareDevelopedandApproved/ApprovalApplications/TherapeuticBiologicApplications/ucm113466.pdf>. Published December 10, 1997.
9. Bross PF, Beitz J, et al. Approval summary: gemtuzumab ozogamicin in relapsed acute myeloid leukemia. *Clin Cancer Res.* 2001 Jun;7(6):1490-6. Erratum in: *Clin Cancer Res* 2002 Jan;8(1):300.
10. Frenzel A, Schirrmann T, Hust M. Phage display-derived human antibodies in clinical development and therapy. *MAbs.* 2016;8(7):1177.
11. Lipson EJ, Drake CG. Ipilimumab: an anti-CTLA-4 antibody for metastatic melanoma. *Clin Cancer Res.* 2011 Nov 15;17(22):6958-62.
12. Przepiorka D, Ko CW, Deisseroth A, et al. FDA Approval: Blinatumomab. *Clin Cancer Res.* 2015 Sep 15;21(18):4035-9.
13. Albanell J, Baselga J. Trastuzumab, a humanized anti-HER2 monoclonal antibody, for the treatment of breast cancer. *Drugs Today (Barc).* 1999 Dec;35(12):931-46.
14. Melsheimer R, Geldhof A, Apaolaza I, Remicade® (infliximab): 20 years of contributions to science and medicine. *Biologics.* 2019 Jul 30;13:139-178.
15. Chuk MK, Chang JT, Theoret MR, et al. FDA Approval Summary: Accelerated Approval of Pembrolizumab for Second-Line Treatment of Metastatic Melanoma. *Clin Cancer Res.* 2017 Oct 1;23(19):5666-5670.
16. Lyu X, Zhao Q, Hui J, et al. The global landscape of approved antibody therapies. *Antib Ther.* 2022;5(4):233.



© 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. All other trademarks are the property of their respective owners.

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

2024-GBL-EN-105873-v1

