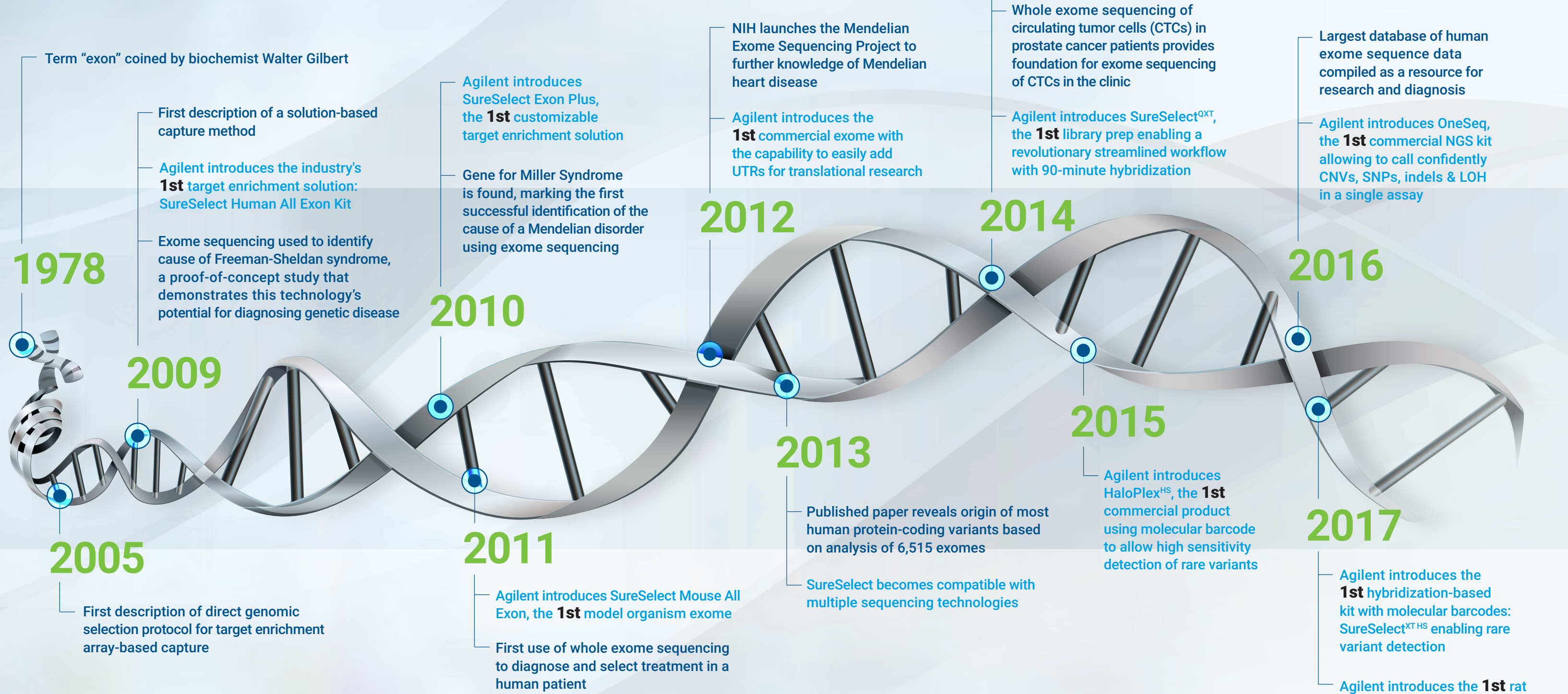


Shaping Exome Sequencing



References

1. Gilbert, W. Why genes in pieces? Nature 1978;271:501. doi: 10.1038/271501a0.
2. Stavros Basiardes; Rose Velle; Cindy Helms; Elaine R. Mardis; Anne M. Bowcock; Michael Lovett. Direct genomic selection. Nature Methods 2005. 2:63-69. doi: 10.1038/nmeth0105-63.
3. Ng SB, Turner EH, Robertson PD, Flygare SD, Bigham AW, Lee C, Shaffer T, Wong M, Bhattacharjee A, Eichler EE, Bamshad M, Nickerson DA, Shendure J. Targeted capture and massively parallel sequencing of 12 human exomes. Nature. 2009;461:272-276. doi: 10.1038/nature08250.
4. Ng SB1, Buckingham KJ, Lee C, Bigham AW, Tabor HK, Dent KM, Huff CD, Shannon PT, Jabs EW, Nickerson DA, Shendure J, Bamshad MJ. Exome sequencing identifies the cause of a mendelian disorder. Nat Genet. 2010;Jan;42(1):30-5. doi: 10.1038/ng.499.
5. Worthey E.A., Mayer A. N., Syverson G.D., Helbling D., Bonacci B.B., et al. Making a definitive diagnosis: successful clinical application of whole exome sequencing in a child with intractable inflammatory bowel disease. Genet. Med. 2011. 13: 255-262.
6. National Institute of Health: National Heart, Lung, and Blood Institute. (March 2012 to September 2015). Mendelian Exome Sequencing Project (Mendelian Exome). Retrieved from: <https://www.nhlbi.nih.gov/research/resources/genetics-genomics/mendelian>
7. Fu W1, O'Connor TD, Jun G, Kang HM, Abecasis G, Leal SM, Gabriel S, Rieder MJ, Altshuler D, Shendure J, Nickerson DA, Bamshad MJ; NHLBI Exome Sequencing Project, Akey JM. Analysis of 6,515 exomes reveals the recent origin of most human protein-coding variants. Nature. 2013 Jan 10;493(7431):216-20. doi: 10.1038/nature11690.
8. Lohr, J. et al. Whole exome sequencing of circulating tumor cells provides a window into metastatic prostate cancer. Nat Biotechnol. 2014 May;32(5): 479-484.
9. Lek, M. et al. Analysis of protein-coding genetic variation in 60,706 humans. Nature. 2016 Aug 18: 536(7616):285-91. doi: 10.1038/nature19057.