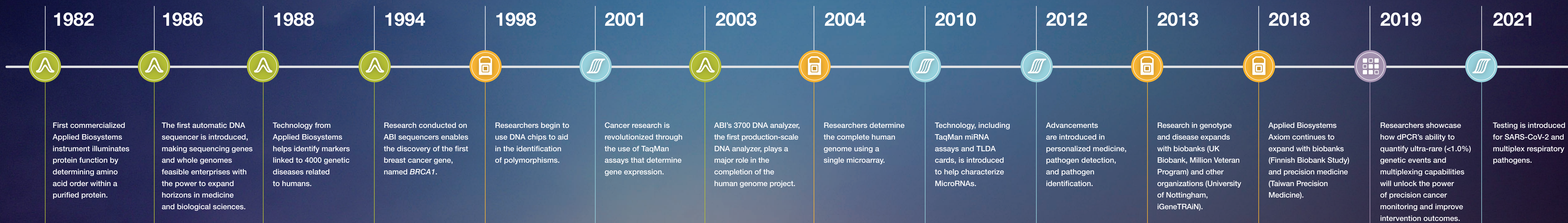




years of asking
the right questions

Biology's biggest questions. History-defining breakthroughs.

With a 40-year history of supporting breakthroughs, Applied Biosystems technologies deliver the complete genomics solutions that help scientists ask—and answer—biology's biggest questions.



= Capillary Electrophoresis (CE)
 = Microarray
 = qPCR
 = dPCR

