









Decades of DNA Discovery Now Primed for **INNOVATION**

 <ul style="list-style-type: none"> • Double helix structure of DNA discovered • First DNA successfully synthesized in test tube 	 <ul style="list-style-type: none"> • Discovery that mRNA directs protein synthesis • First restriction enzyme described 	 <ul style="list-style-type: none"> • First recombinant DNA molecules synthesized • First animal gene cloned • First genetic engineering company founded 	 <ul style="list-style-type: none"> • First transgenic mice and fruit flies • Yeast artificial chromosomes used to ferry large DNA fragments 	 <ul style="list-style-type: none"> • Human Genome Project launches • FDA approves sale of first GMO food 	 <ul style="list-style-type: none"> • Human genome sequenced • FDA approves first gene-targeted drug therapy 	 <ul style="list-style-type: none"> • First synthetic life form created • CAR-T therapy for cancer approved 	 <ul style="list-style-type: none"> • CRISPR Wins Nobel Prize
<p>1950s</p>	<p>1960s</p>	<p>1970s</p>	<p>1980s</p>	<p>1990s</p>	<p>2000s</p>	<p>2010s</p>	<p>2020s</p>
		<p>1975</p> <p>Development of Sanger method for DNA sequencing</p>	<p>1983</p> <p>Development of polymerase chain reaction (PCR) method of DNA amplification</p>	<p>1991</p> <p>Patent filed for column cleanup device and a process for the isolation of nucleic acids</p>	<p>2006</p> <p>BioBricks was established</p>	<p>2012</p> <p>Development of CRISPR/Cas9 gene editing tool</p>	<p>2021</p> <p>Launch of EMnetik 24</p>



FAST FORWARD YOUR DISCOVERIES IN 2021: Accelerate Answers and Nucleic Acid Cleanup with EMnetik 24



21.03.116.GEN © 2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service names 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. Product in development, performance characteristics have not been validated. Not intended or validated for use in diagnosis of disease or other conditions.