

Genomic Solutions Collaboration Program

Our mission today is to empower you to find answers to life's important scientific and research questions. **Our Genomic Solutions Collaboration Program is one of the ways we do that.**

Through this program, you'll have exclusive access to Beckman Coulter Life Sciences genomic technologies, data analysis tools and application scientists. It's our hope that with this collaboration, you'll be able to discover new ways to accelerate your research workflows.

How the Program Works

Step 1: Discussion	Tell us about your research goals, needs and aspirations, and we'll discuss general collaboration opportunities.
Step 2: Informal Proposal & Plan	We'll present an informal proposal and high-level action plan to you. We'll discuss and adjust where needed.
Step 3: Formal Proposal & Action Plan	We'll present an actionable proposal, generate a detailed action plan.
Step 4: Execution & Delivery	You'll conduct your research and we'll be there right alongside you with the support and genomic solutions you need. After your data is analyzed and reviewed, we'll promote your work through an application note, whitepaper or peer-reviewed paper and showcase it online and at events.

Who We've Collaborated With

We've collaborated with researchers from around the world. Here's a quick look at who they are and what they've done so far.

Collaboration Partner	About Their Collaborative Research Project
Warsaw Genomics	Built an extraction pipeline in mere weeks using the RNAdvance Viral reagent kit and Biomek i5 liquid handler.
Miami Cancer Institute	In the early days of the COVID-19 pandemic, shifted the focus of their work and completed RNA extraction validation independently using RNAdvance Viral reagents for viral research.
BridgeBio	Created a simultaneous nucleic acid extraction method, particularly relevant to gene therapies.
Henry Ford Health	Used the FormaPure XL reagent kit to generate whole transcriptome sequencing data for biomarker discovery.
Niagara Falls Memorial Medical Center	Developed an innovative collection method to preserve precious biopsy specimens, ensuring their suitability for molecular research testing.



Interested in collaborating with us?
Visit becls.co/grantly or scan the QR code to get started.