



Accelerate Answers with Beckman Coulter Life Sciences

We empower those seeking answers to life's important scientific and healthcare questions. And we're committed to accelerating answers.

Accelerating Answers

In the life sciences accelerating answers requires leading-edge solutions that you can trust. By partnering with us as an OEM supplier to develop your genomic assays, you're committing to a solution that you and your customers can rely on.

Gold Standard Reagents

Our genomic reagents are the gold standard in nucleic acid purification and clean-up technology. They've helped generate research in over 20,000 scientific publications and are suggested for use in over 200 library preparation kits, including kits from trusted sequencing companies like Illumina®, Oxford Nanopore Technologies, and Pacific Biosciences (PacBio), among others.

Our solutions are optimized to meet the stringent needs of today's genomic applications. By adding genomic reagents from Beckman Coulter Life Sciences to your assay kit, you empower your customers to maximize the recovery, consistency and speed of their entire next-generation sequencing workflow.

Help customers get the data they need

Partner with us and help your customers minimize the risk of losing important genetic data. Give your customers the gold standard and empower them to generate better results.



Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions. This protocol is for demonstration only, and is not validated by Beckman Coulter.

© 2021 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.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)
21.02.875.GEN

