



SPRI Methodology

Solid Phase Reversible Immobilization (SPRI) technology, uses paramagnetic beads to selectively bind nucleic acids by type and size

Used for

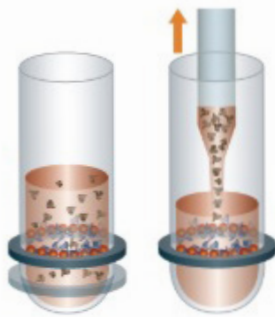
- High-performance isolation, purification and clean-up protocols
- Supporting applications such as next-generation sequencing (NGS), Sanger sequencing, qPCR, ddPCR and microarrays

Nucleic Acid Immobilization



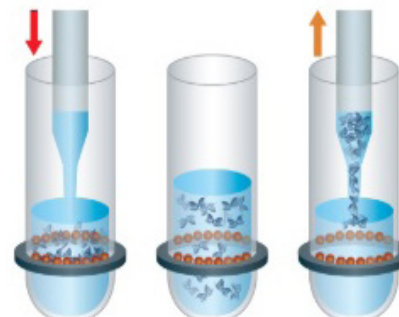
- SPRI beads are directly added to sample reactions
- Nucleic acids are selectively immobilized onto SPRI beads, leaving contaminants in solution

Contaminant Removal and Wash



- A magnetic field is used to pull the microparticles out of solution
- Contaminants are aspirated and microparticles are thoroughly washed, yielding high quality nucleic acids

Nucleic Acid Elution (Screen shot of purified DNA coming off beads bound to magnet)



- Purified nucleic acids are easily eluted from the microparticles under aqueous conditions, which provides maximum flexibility for downstream applications

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.

© 2020 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)

AAG-6504FLY02.20