



AMPure XP- Gold Standard for bead based, Next-Generation Sequencing (NGS) clean-up

AMPure XP is suggested by over 200 library prep kits, including kits from trusted sequencing companies:

- Illumina®
- Oxford Nanopore Technologies
- Pacific Biosciences
- ThermoFisher Scientific

Over 15,000 publications used AMPure XP

- Referenced in articles in Science, Nature, and PNAS

AMPure XP maximizes recovery, consistency and speed

Nucleic acid purification and clean-up are mandatory for genomic applications, such as sequencing, qPCR/ddPCR/PCR, and microarrays. Maximizing recovery, consistency, and speed, AMPure XP meets the stringent needs of today's genomic applications and minimizes the risk of losing important genetic information.

- High recovery of amplicons greater than 100bp
- Efficient removal of unincorporated dNTPs, primers, primer dimers, salts and other contaminants
- Predictable and consistent size selection

“Our findings suggest AMPure XP would be the best choice for analyses requiring very high analytical stringency.” Mikheikin, A., Olsen, A., Picco, L. et al. High-speed atomic force microscopy revealing contamination in DNA purification systems. Anal. Chem. 88:5, 2527-2523 (2016) doi: 10.1021/acs.analchem.5b04023

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

AAG-6474FLY02.20

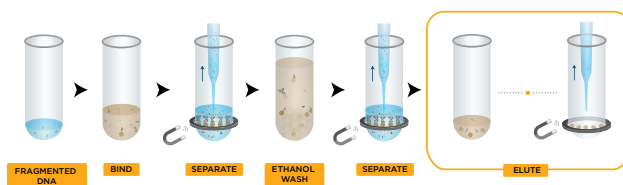


SPRIselect vs AMPure XP Feature comparison

| SPRIselect | AMPure XP |
|--|--|
| Stable at Room Temperature | Stable at 4°C |
| QC'd for size selection | QC'd for recovery |
| Suggested for use in library construction kits | Suggested for use in library construction kits |

SPRIselect and AMPure XP

- Can be substituted for each other in clean-up applications
- Use the same sample to bead ratios
- Are suggested for use by the most well-known library construction kit manufacturers:
 - Illumina®, Integrated DNA Technologies, Inc (IDT), Swift Biosciences, etc
- Have identical work flows



SPRIselect benefits

- Stability at room temperature allows:
 - Start clean-up steps without calibrating to room temperature
 - Frees up 4°C storage space
- Quality control for size selection allows:
 - Trust that lot to your size selection steps are accurate
 - Accurate and consistent size selection from lot to lot

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-6474FLY02.20

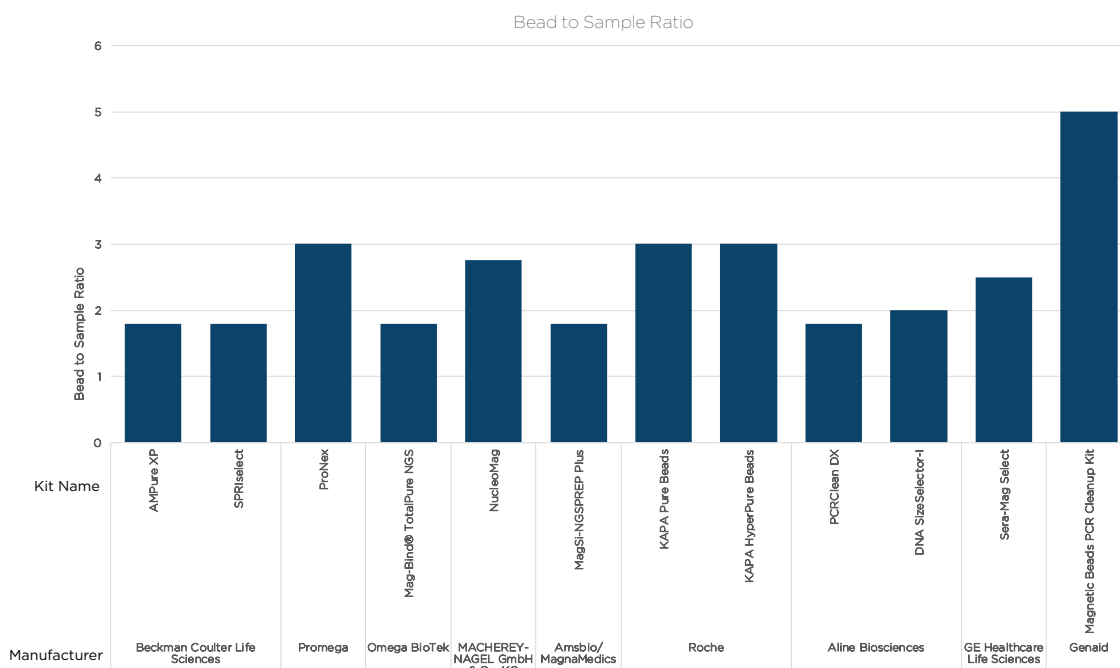


Sample to bead ratio for size selection matters

Not all bead based clean-up kits use the 1.8x ratio; the figure below shows that of 10 bead based clean-up kits only 3 suggest a 1.8x ratio for clean-up.

What switching to a kit with a different ratio could impact

- Clean-ups will require a different bead ratio that could lead to loss of fragments of interest
- Sample to bead ratios will need to be recalculated
- Lab protocols that use the standard sample to bead ratios that of AMPureXP or SPRIselect will have to be re-written



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-6474FLY02.20