



# Efficient DNA purification using SPRI technology

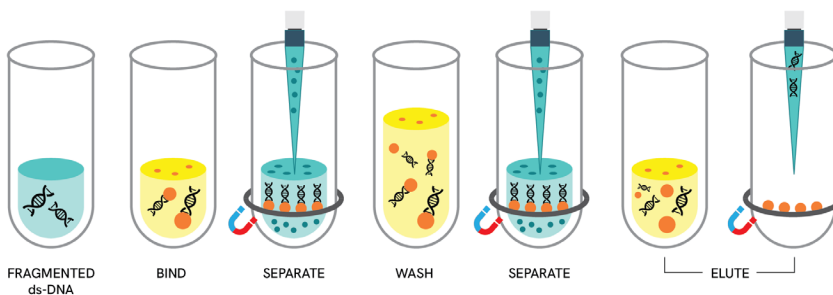
## NEXTFLEX NGS Cleanup Beads

Revvity brings you the power of SPRI in our high-performance NEXTFLEX™ NGS Cleanup Beads

- Efficiently purify nucleic acids for NGS library preparation.
- Reversible magnetic bead binding for optimal DNA recovery.
- Removes unwanted components (primers, salts, adapters).
- Precise DNA size selection capabilities.
- Compatible with manual and automated workflows.
- Drop-in replacement for most SPRI beads (no protocol changes needed).



### NEXTFLEX NGS Cleanup Beads



1. Capture DNA using magnetic beads.
2. Isolate beads from impurities.
3. Rinse magnetic beads with 70% ethanol solution to eliminate impurities.
4. Release DNA from magnetic beads.
5. Transfer eluted DNA to a fresh plate.

## Understanding SPRI technology

- SPRI technology utilizes paramagnetic beads for efficient DNA purification.
- Beads bind DNA selectively in the presence of polyethylene glycol (PEG) and salt.
- Binding is reversible, allowing for controlled DNA capture and release.
- Adjusting the bead-to-sample ratio enables size selection.
- Optimized bead composition ensures high recovery and purity.

### Size selection with bead-to-sample ratio

Fine-tune your NGS library preparation with precise size selection. By adjusting the ratio of NEXTFLEX NGS Cleanup Beads to your DNA sample, you can selectively retain fragments of the desired size range. A higher bead ratio captures smaller fragments, while a lower ratio retains larger fragments.

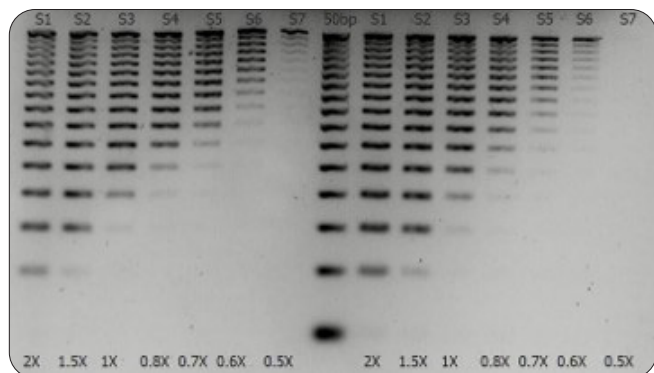


Figure 1: NEXTFLEX NGS Cleanup Beads (right) demonstrate equivalent performance to AMPure® XP beads (left) in purifying a 50 bp DNA ladder.

## NEXTFLEX NGS Cleanup Beads: performance and versatility

### Proven performance

- Achieve high DNA recovery rates.
- Ensure excellent purity for downstream NGS applications.
- Consistent and reliable results across different sample types.

### Ordering information

Part number	Size
NOVA-BEADS-1ML	1 mL
NOVA-BEADS-5ML	5 mL
NOVA-BEADS-50ML	50 mL
NOVA-BEADS-500ML	500 mL

### Storage and stability

Store at +2-8°C. Stable for at least 6 months when stored properly.

Evaluations samples available.

Contact us for more information.



Scan for detailed specifications and more.

revvity