

SPECIFICATION SHEET

NEXTFLEX® NGS Cleanup Beads: NEXTFLEX® NGS Cleanup Beads efficiently purify nucleic acids using reversible magnetic bead binding. They are ideal for NGS library preparation, removing unwanted components like primers, salts, and adapters as well as DNA size selection with consistent results. These beads are compatible with both manual and automated protocols.

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 beads at +2-8°C upon receipt. Do not freeze. The shelf life of this reagent is at least 6 months when stored properly.

Usage instructions:

- Bring beads to room temperature before use
- Resuspend beads fully before use
- Prepare 80% ethanol fresh before every use
- NEXTFLEX® NGS Cleanup Beads can be used to replace AMPure® XP beads in most NGS workflows without any protocol changes or optimizations needed (Figure 1).

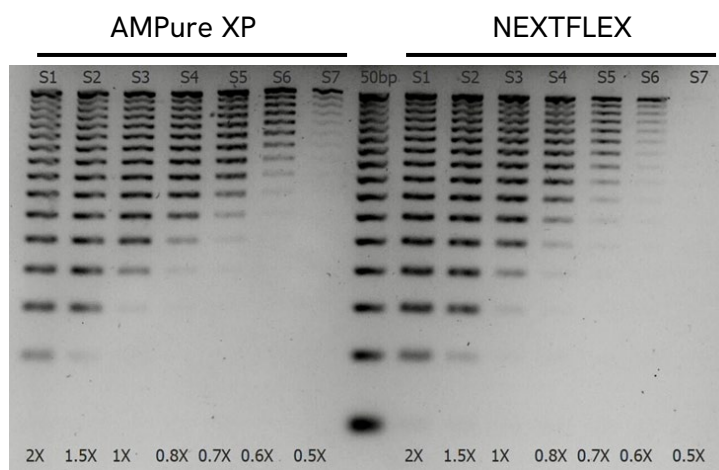


Figure 1. 625 ng of 50 bp DNA ladder (Promega) were purified with equal amount of AMPure® XP and NEXTFLEX® NGS Cleanup and run on 1% agarose gel for comparison.