

## CUTANA™ DNA Purification Beads

Catalog No	21-1407-5ML	Pack Size	5 mL
Lot No	25175001-51		

### DESCRIPTION

CUTANA™ DNA Purification Beads harness **Solid-Phase Reversible Immobilization (SPRI\*)** technology to deliver fast, high-yield DNA cleanup and precise size selection for NGS library preparation and other genomic workflows. Optimized PEG and salt concentrations drive efficient, reproducible binding of DNA to paramagnetic beads, ensuring exceptional recovery with minimal bias or carry-over contamination. The formulation **matches the bead-to-sample ratio of popular alternatives** such as AMPure XP\* and SPRIselect\*, ensuring CUTANA™ DNA Purification Beads can serve as a drop-in replacement into manual or automated cleanup workflows without any protocol re-optimization.

#### Key Advantages

- **Trusted method** – SPRI chemistry researchers rely on for DNA cleanup
- **NGS validated** – proven performance in genomic workflows including CUT&RUN, CUT&Tag, and library preparation
- **High recovery & purity** – consistently recovers DNA across a wide range of fragments and input amounts
- **Automation-compatible** – suitable for manual or automated protocols using magnetic separation
- **Drop-in AMPure XP\* alternative** – identical ratios for seamless replacement in SPRI-based cleanups
- **Room-temperature stable** – ready to use; no warmup period required

#### Validated Applications

- CUT&RUN
- CUT&Tag
- NGS Library Preparation & adapter dimer cleanup

### TECHNICAL INFORMATION

Storage	Store at 15 – 30°C (room temperature; <b>DO NOT FREEZE</b> ). Stable for 12 months from date of receipt.
Formulation	Paramagnetic bead suspension
Binding mechanism	SPRI (PEG/NaCl)
Size-selection range	Tunable via 0.8X – 1.8X bead-to-sample ratio
Handling	<b>Mix well by vortexing before use</b> to resuspend beads uniformly

### APPLICATION NOTES

Use CUTANA™ DNA Purification Beads at the following bead-to-sample ratios:

#### CUT&RUN

DNA Purification: 1.8X

#### CUT&RUN Library Preparation

Adapter Ligation Cleanup: 1.0X

PCR Cleanup: 1.0X

Adapter Dimer Removal: 1.0X

#### CUT&Tag

PCR Cleanup: 1.3X

#### Total DNA Cleanup: 1.8X

RECOMMENDED ACCESSORY PRODUCTS

Item	Cat No	Item	Cat No
CUTANA™ pAG-MNase for ChIC/CUT&RUN	15-1016	CUTANA™ pAG-Tn5 for CUT&Tag	15-1017
Magnetic Separation Rack, 0.2 mL	10-0008	Magnetic Separation Rack, 1.5 mL	10-0012
CUT&RUN 8-strip Tubes	10-0009		

CERTIFICATE OF ANALYSIS

Quality Control Metrics	<b>DNA Fragment Recovery Cleanup</b> Each lot of CUTANA™ DNA Purification Beads is subjected to a SPRI-based DNA cleanup protocol using a GeneRuler 50 bp DNA Ladder (ThermoFisher Scientific SM0371) tested at a range of bead-to-sample ratios in triplicate. A validated lot of CUTANA™ DNA Purification Beads are run in parallel. DNA Ladder only and no DNA controls are included. The resulting DNA is run on an Agilent TapeStation. Recovery of DNA fragments within 100 to 500 bp are quantified relative to DNA Ladder input control to determine percent recovery at each bead-to-sample ratio.	Pass
Approved By	Andrea Johnstone, PhD Senior Director, Product Management <i>This is a digital certificate and has not been signed by hand.</i>	
Date of Issue	July 15 2025	

This product ("Product") is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications, and/or accompanying package inserts ("Documentation") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper, and intended usage. This warranty does not extend to anyone other than the original purchase of the Product ("Buyer"). No other warranties, express or implied, are granted, including without limitation, implied warranties of merchantability, fitness for any particular purpose, or non infringement. Buyer's exclusive remedy for non-conforming Products during the warranty period is limited to replacement of or refund for the non-conforming Product(s). There is no obligation to replace Products as the result of (i) accident, disaster or event of force majeure, (ii) misuse, fault or negligence of or by Buyer, (iii) use of the Products in a manner for which they were not designed, or (iv) improper storage and handling of the Products.

\*SPRI, SPRIselect, and AMPureXP are registered trademarks of Beckman Coulter and nominative references to them are for descriptive purposes only. CUTANA™ products are not manufactured by or affiliated with Beckman Coulter.