

CELLDATA RNAsstorm™ 2.0 MagBead FFPE RNA Extraction Kit



A magnetic bead-based kit for superior RNA extraction from FFPE samples. Features proprietary technology for enhanced removal of formaldehyde-induced damage, providing RNA with higher yield and greater amplifiability.

Product Description

Formalin-fixed tissue samples are a challenge for RNA extraction, often resulting in low yields and poor performance in subsequent steps. Most existing methods rely on heat to remove crosslinks and adducts, which is only partially effective and leads to additional fragmentation of labile nucleic acids. Powered by proprietary CAT5™ catalytic technology, the RNAsstorm™ FFPE RNA Extraction Kits enhance the removal of formaldehyde-induced damage. The extracted RNA has higher yield, quality, and greater amplifiability that is more optimal for next-generation sequencing, qPCR, microarray, or other gene expression analysis.

The RNAsstorm™ 2.0 MagBead FFPE RNA Extraction Kit uses magnetic beads for RNA isolation and may be adapted for multiwell plate format. For spin column-based RNA purification, see our original [CELLDATA RNAsstorm™ 2.0 FFPE RNA Extraction Kit](#).

- Purify higher quality RNA with less fragmentation
- Extract higher yields of amplifiable RNA compared to other methods
- CAT5™ technology for chemical reversal of formaldehyde crosslinking
- Milder conditions and no harsh solvents
- Better results in downstream analysis like PCR, microarray, or next-generation sequencing
- Magnetic bead-based RNA purification adaptable for multiwell plate formats

High-Throughput and Efficient RNA Extraction from FFPE Tissues

The RNAsstorm™ 2.0 MagBead FFPE RNA Extraction Kit offers the same performance and CAT5™ technology for efficient removal of formaldehyde crosslinking found in our spin column-based [CELLDATA RNAsstorm™ 2.0 FFPE RNA Extraction Kit](#).

Advantages of CAT5™ Technology for RNA Extraction

Higher Yields of Amplifiable RNA

Significant improvements in RNA yield and quality (as measured by amount of amplifiable RNA) are seen using a prototype of the RNAsstorm™ kit featuring CAT5™ technology on FFPE samples from various tissues and with collection dates ranging from 1976 to 2015 compared to a competitor kit.

Higher Integrity RNA

See the product table below for our complete line of RNAsstorm™ and DNAsstorm™ extraction kits for FFPE tissues. We also offer the [RNAsstorm™ RNA Isolation Kit](#) for purifying RNA from fresh or frozen cells/tissue samples, and the [ExoBrite™ EV Total RNA Isolation Kit](#) for isolate RNA from extracellular vesicles.

Product attributes

Storage Conditions

See individual components for storage temperature

RNAstorm™ and DNAstorm™ FFPE Extraction Kits

Product Name	Catalog Number	Size
CELldata RNAstorm™ 2.0 FFPE RNA Extraction Kit	CD506	50 preps
CELldata DNAstorm™ 2.0 FFPE DNA Extraction Kit	CD507	50 preps
CELldata DNAstorm™/RNAstorm™ 2.0 Combination Kit	CD508	50 preps
CELldata RNAstorm™ 2.0 MagBead FFPE RNA Extraction Kit	CD510-96	96 preps
CELldata DNAstorm™ 2.0 MagBead FFPE DNA Extraction Kit	CD509-96	96 preps

References

[Download a curated list of references for CELldata DNAstorm™ and RNAstorm™ FFPE Extraction Kits](#)

This datasheet was generated on June 17, 2026 at 05:28:02 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/celldata-rnastorm-2-0-magbead-ffpe-rna-extraction-kit/>