



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## CELLDATA RNAsort™ 2.0 FFPE RNA Extraction Kit

Superior RNA extraction from FFPE (paraffin) tissue sections with 3 to 30-fold increase in yield, and 2-fold increase in DV200 (RNA longer than 200 nucleotides) compared to other methods.



### Product attributes

#### Storage Conditions

See individual components for storage temperature

## Product Description

Formalin-fixed tissue samples are a challenge for RNA extraction, often resulting in low yields of amplifiable RNA and poor performance in subsequent steps involving enzymatic manipulation, including reverse transcription and sequencing library preparation. Powered by proprietary CAT5™ technology, the RNAsort™ extraction kit enhances the removal of formaldehyde-induced damage and provides RNA with higher yield and integrity as shown by a higher RIN score and higher DV200 (% RNA >200nt). For RNA with greater amplifiability in applications such as RNA-seq, qPCR, microarray, or other gene expression analysis, the RNAsort™ 2.0 FFPE RNA Extraction kit is your best chance for success.

- CAT5™ technology for chemical reversal of formaldehyde crosslinking
- Extract higher yields of amplifiable RNA compared to other methods
- Milder conditions and no harsh solvents
- Higher quality RNA with less fragmentation
- Better results in downstream analysis like PCR, microarray, or next-generation sequencing
- Simple spin column-based workflow

**Note:** This kit is a replacement for catalog number CD501. It uses a modified deparaffinization protocol with a new Dewaxing Solution (provided in the kit). The new kit retains the same proprietary CAT5™ catalytic technology for highly efficient reversal of formaldehyde crosslinks as the original kit.

See the product table below for our complete line of RNAsort™ and DNAsort™ extraction kits for FFPE tissues. We also offer the [RNAsort™ 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.

### 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 RNAsort™ 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

# RNAstorm™ and DNAstorm™ FFPE Extraction Kits

Product Name	Catalog Number	Size
<a href="#">CELldata RNAstorm™ 2.0 FFPE RNA Extraction Kit</a>	<a href="#">CD506</a>	50 preps
<a href="#">CELldata DNAstorm™ 2.0 FFPE DNA Extraction Kit</a>	<a href="#">CD507</a>	50 preps
<a href="#">CELldata DNAstorm™/RNAstorm™ 2.0 Combination Kit</a>	<a href="#">CD508</a>	50 preps
<a href="#">CELldata RNAstorm™ 2.0 MagBead FFPE RNA Extraction Kit</a>	<a href="#">CD510-96</a>	96 preps
<a href="#">CELldata DNAstorm™ 2.0 MagBead FFPE DNA Extraction Kit</a>	<a href="#">CD509-96</a>	96 preps

## References

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

This datasheet was generated on January 19, 2026 at 07:04:53 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/celldata-rnastorm-2-ffpe-rna-extraction-kit/>