



Glowing products for science

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

CELLDATA DNASTORM™/RNAstorm™ 2.0 Combination Kit

A combination kit for dual extraction of DNA and RNA from a single FFPE sample, featuring proprietary CAT5™ technology.



Product attributes

Storage Conditions

See individual components for storage temperature

Product Description

Formalin-fixed tissue samples are a challenge for nucleic acid extraction, often resulting in low yields of amplifiable DNA/RNA and poor performance in subsequent steps involving enzymatic manipulation, including reverse transcription and sequencing library preparation. Powered by proprietary CAT5™ technology, the CELLDATA DNASTORM™/RNAstorm™ 2.0 Combination Kit enhances the removal of formaldehyde-induced damage and provides dual DNA and RNA extraction with higher yield and quality. This allows for improved amplifiability for applications such as RNA-seq, qPCR, microarray, or other gene expression analysis.

- Extract higher yields of DNA and RNA from a single FFPE tissue sample
- CAT5™ technology for chemical reversal of formaldehyde crosslinking
- Milder conditions and no harsh solvents
- Higher quality DNA/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 CD503. 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.

For extracting either RNA or DNA from FFPE tissues, please see our [CELLDATA RNAstorm™ 2.0 FFPE RNA Extraction Kit \(CD506\)](#) and [CELLDATA DNASTORM™ 2.0 FFPE DNA Extraction Kit \(CD507\)](#). We also offer the [RNAstorm™ RNA Isolation Kit](#) for purifying RNA from fresh or frozen cells/tissue samples.

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™ Extraction Kits](#)

This datasheet was generated on January 19, 2026 at 08:59:20 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/celldata-dnastorm-rnastorm-2-0-combination-kit/>