



Glowing products for science

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

## CELLDATA DNASTORM™ 2.0 MagBead FFPE DNA Extraction Kit



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

### Product Description

Formalin-fixed tissue samples are a challenge for DNA 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™ technology, the DNASTORM™ 2.0 MagBead FFPE Extraction Kit enhances the removal of formaldehyde-induced damage and provides DNA with higher yield and quality and greater amplifiability. The DNASTORM™ 2.0 MagBead FFPE DNA Extraction Kit is the best solution for next-generation sequencing and other advanced applications. This kit uses magnetic beads for DNA isolation and may be adapted for multiwell plate format. For spin column-based DNA purification, see our original [CELLDATA DNASTORM™ 2.0 FFPE DNA Extraction Kit](#).

- Purify higher quality DNA with less fragmentation
- Extract higher yields of amplifiable DNA 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 DNA purification adaptable for multiwell plate formats

The DNASTORM™ 2.0 DNA extraction kit greatly accelerates the removal of formaldehyde damage and allows recovery of higher quality nucleic acids. This results in higher yields, more intact DNA, and better results in downstream analysis like PCR, microarray, or next-generation sequencing (NGS). See the product table below for our complete line of RNASTORM™ and DNASTORM™ extraction kits for FFPE tissues.

### 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

1. Am J Surg Pathol, 43, 1720 [DOI: 10.1097/PAS.0000000000001338](https://doi.org/10.1097/PAS.0000000000001338)

This datasheet was generated on January 9, 2026 at 04:01:18 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/celldata-dnastorm-2-0-magbead-ffpe-dna-extraction-kit/>

### Product attributes

Storage Conditions

See individual components for storage temperature