CELLDATA DNAstorm™ 2.0 FFPE DNA Extraction Kit

Storage Conditions

Product attributes

See individual components for storage temperature

Call us: 800-304-5357 Email: btinfo@biotium.com

Superior DNA extraction from FFPE samples, due to the enhanced removal of formaldehyde-induced damage, providing DNA with higher yield and quality 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 CAT5TM technology, the DNAstormTM 2.0 FFPE Extraction Kit enhances the removal of formaldehyde-induced damage and provides DNA with higher yield and quality and greater amplifiability. The DNAstormTM 2.0 FFPE DNA Extraction Kit is the best solution for next-generation sequencing and other advanced applications.

- 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
- Simple workflow

Note: This kit is a replacement for catalog number CD502. It uses a modified deparaffinization protocol with a new Dewaxing Solution (provided in the kit). The new kit retains the same proprietary CAT5TM catalytic technology for highly efficient reversal of formaldehyde crosslinks as the original kit.

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
CELLDATA RNAstorm™ 2.0 FFPE RNA Extraction Kit	<u>CD506</u>	50 preps
CELLDATA DNAstorm™ 2.0 FFPE DNA Extraction Kit	<u>CD507</u>	50 preps
CELLDATA DNAstorm TM /RNAstorm TM 2.0 Combination Kit	<u>CD508</u>	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 list of Mix-n-Stain™ references.

This datasheet was generated on November 2, 2025 at 01:40:36 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/celldata-dnastorm-2-ffpe-dna-extraction-kit/