Dewaxing Solution

Dewaxing Solution provided in CELLDATA DNAstorm™ 2.0 and RNAstorm™ 2.0 FFPE extraction kits, available for separate purchase. The solution is a less hazardous alternative to xylenes for efficient removal of paraffin from FFPE tissue sections for downstream RNA or DNA extraction.



Product attributes

Storage Conditions

Store at room temperature

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

Product Description

This Dewaxing Solution (Cat. No. 99876-30ML) is supplied in our CELLDATA DNAstorm™ 2.0 and RNAstorm™ 2.0 FFPE extraction kits and is provided for separate purchase for customers who require larger volumes of the solution. For recommended procedures, please navigate to the product information sheet found on the product page for each kit linked below.

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™/RNAstorm™ 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

Learn more about our DNA and RNA extraction kits for fresh cells and FFPE tissues.

This datasheet was generated on November 6, 2025 at 09:19:33 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dewaxing-solution/