Glowing products for science

## **CELLDATA RNAstorm™ FFPE RNA Extraction Kit**

This item is discontinued.



## **Product attributes**

Storage Conditions See individual components for storage temperature

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

Kit RNAstorm™ Kit for RNA from FFPE Tissues, DNAstorm™/RNAstorm™ Combination Kit

## **Product Description**

This item is discontinued and is replaced by <u>CELLDATA RNAstorm™ 2.0 FFPE RNA Extraction Kit</u> (<u>CD506</u>).

This datasheet was generated on November 4, 2025 at 02:04:12 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/rnastorm-kit-for-isolation-of-rna-from-ffpe-tissue-samples/">https://biotium.com/product/rnastorm-kit-for-isolation-of-rna-from-ffpe-tissue-samples/</a>