## CELLDATA RNAstorm™ Fresh Cell and Tissue RNA Isolation Kit

distribution of the state of th

Storage Conditions See individual components for storage temperature

Product attributes

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

The CELLDATA RNAstorm™ Fresh Cell and Tissue RNA Isolation Kit provides a convenient and efficient method of isolating RNA from fresh or frozen cells or tissue samples.

## **Product Description**

Purify high quality total RNA from either cultured cells or tissue (fresh or fresh/frozen) in as little as 20 minutes. High yields (up to 120  $\mu$ g) can be obtained using a simple and quick column-based protocol. Contaminating DNA is removed using a DNase treatment step, and the protocol avoids toxic chemicals such as phenol or chloroform.

- Purify high quality total RNA in as little as 20 minutes
- Simple workflow yielding up to 120 µg of RNA
- DNase treatment included
- No phenol-chloroform and no ethanol precipitation
- Isolate RNA for next-gen sequencing (RNA-Seq), RT-PCR, cDNA synthesis, or microarrays

Also see the CELLDATA DNAstorm™ FFPE DNA Extraction Kit and CELLDATA RNAstorm™ FFPE RNA Extraction Kit for extraction of high-quality DNA and RNA from FFPE tissue samples.

**Higher RNA Yields** 

**High RNA Quality** 

## References

Download a curated list of references for CELLDATA DNAstorm™ and RNAstorm™ Extraction Kits

This datasheet was generated on November 2, 2025 at 06:29:03 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/rnastorm-rna-isolation-kit/">https://biotium.com/product/rnastorm-rna-isolation-kit/</a>