



Glowing products for science

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

CELLDATA RNAsort™ Fresh Cell and Tissue RNA Isolation Kit

The CELLDATA RNAsort™ 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 attributes

Storage Conditions

See individual components for storage temperature

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 µg) 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 DNAsort™ FFPE DNA Extraction Kit](#) and [CELLDATA RNAsort™ 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 DNAsort™ and RNAsort™ Extraction Kits](#)

This datasheet was generated on December 19, 2025 at 01:09:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rnasort-rna-isolation-kit/>