



Glowing products for science

CELLDATA RNAsort™ FFPE RNA Extraction Kit

This item is discontinued.



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

Product attributes

Storage Conditions	See individual components for storage temperature
Kit	RNAsort™ Kit for RNA from FFPE Tissues, DNAsort™/RNAsort™ Combination Kit

Product Description

This item is discontinued and is replaced by [CELLDATA RNAsort™ 2.0 FFPE RNA Extraction Kit \(CD506\)](#).

This datasheet was generated on January 9, 2026 at 12:05:35 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/rnasort-kit-for-isolation-of-rna-from-ffpe-tissue-samples/>