Glowing products for science

CELLDATA DNAstorm™ FFPE DNA Extraction Kit

DNA STORM Product attributes
Storage Conditions

See individual components for storage temperature

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

This item is discontinued.

Product Description

This item is discontinued and is replaced by <u>CELLDATA DNAstorm™ 2.0 FFPE DNA Extraction Kit</u> (<u>CD507</u>).

This datasheet was generated on November 11, 2025 at 10:09:25 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dnastorm-kit-for-isolation-of-dna-from-ffpe-tissue-samples/