



Glowing products for science

## CELLDATA DNASTORM™ FFPE DNA 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

### Product Description

This item is discontinued and is replaced by [CELLDATA DNASTORM™ 2.0 FFPE DNA Extraction Kit \(CD507\)](#).

This datasheet was generated on June 16, 2026 at 08:06:59 PM. 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/>