Glowing products for science

Ready-to-Use 100 bp DNA Ladder

This product is discontinued.



Product attributes

Storage Conditions

Store at -10 to -35 °C

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

Product Description

This product is discontinued. Please our replacement product 100 bp DNA Ladder Ready-to-Load (Cat. No. 31085).

This datasheet was generated on November 3, 2025 at 12:25:27 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/ready-to-use-100-bp-dna-ladder/