

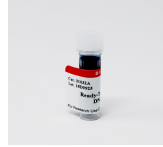


Glowing products for science

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

Ready-to-Use 100 bp DNA Ladder

This product is discontinued.



Product attributes

Storage Conditions

Store at -10 to -35 °C

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 June 5, 2026 at 07:11:49 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ready-to-use-100-bp-dna-ladder/>