

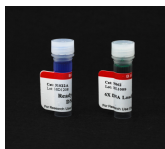


Glowing products for science

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

## Ready-to-Use 1 kb 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 [1 kb DNA Ladder, Ready-to-Load \(Cat. No. 31084\)](#).

This datasheet was generated on January 13, 2026 at 08:52:21 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/ready-to-use-1-kb-dna-ladder/>