



Glowing products for science

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

100 bp DNA Ladder, Ready-to-Load

A DNA ladder consisting of 11 DNA fragments ranging in size from 100 bp to 1500 bp, provided pre-diluted in 1X loading buffer so that it is ready to load.



Product attributes

Storage Conditions	Store at -10 to -35 °C
--------------------	------------------------

Product Description

100 bp DNA Ladder, Ready-to-Load is suitable for sizing linear double-stranded DNA fragments from 100 bp to 1500 bp. The ladder consists of 11 double stranded DNA bands. The 500 bp band has increased intensity to provide internal orientation.

- Contains 11 double-stranded DNA fragments
- Bands range from 100 bp to 1500 bp
- 500 bp band has higher intensity, for orientation
- Provided pre-diluted in 1X loading buffer, ready to load
- Ladder concentration is 20 ng/uL

Note: This product serves as a replacement for the discontinued Ready-to-Use 100 bp DNA Ladder (Cat. No. 31032)

Ready-to-Load Ladders are provided pre-diluted in 1X loading buffer at a concentration of 20 ng/uL. A separate vial of 6X loading buffer is provided to add to your DNA samples. The loading buffer provided contains density agents and two blue electrophoresis tracking dyes that run at approximately 1.5 kb and 200 bp in a 1% agarose gel. We also offer a [1 kb DNA Ladder, Ready-to-Load \(31084\)](#).

[Learn more about our products for nucleic acid gel electrophoresis workflows.](#)

This datasheet was generated on January 7, 2026 at 04:36:38 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/100-bp-dna-ladder-ready-to-load/>