1 kb DNA Ladder, Ready-to-Load

A DNA ladder consisting of 13 DNA fragments ranging in size from 250 bp to 10 kb, provided pre-diluted in 1X loading buffer so that it is ready to load.



Product attributes

Call us: 800-304-5357

Storage Conditions

Store at -10 to -35 °C

Email: btinfo@biotium.com

Product Description

1 kb DNA Ladder, Ready-to-Load is suitable for sizing linear double-stranded DNA fragments from 250 bp to 10 kb. The ladder consists of 13 double stranded DNA bands. The 1 kb and 3 kb bands have increased intensity to provide internal orientation.

- Contains 13 double-stranded DNA fragments
- Bands range from 250 bp to 10 kb
- 1 kb and 3 kb bands have 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 1 kb DNA Ladder (Cat. No. 31022).

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 100 bp DNA Ladder, Ready-to-Load (31085).

Learn more about our products for nucleic acid gel electrophoresis workflows.

This datasheet was generated on November 12, 2025 at 05:25:30 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/1-kb-dna-ladder-ready-to-load/