Glowing products for science

DNA Ladders in TE Buffer

Your choice of 100 bp DNA ladder or 1 kb DNA ladder, supplied at 100 ng/uL in TE buffer.



Product attributes

Call us: 800-304-5357

Storage Conditions

Store at -10°C to -35°C, or 2°C to 8°C

Email: btinfo@biotium.com

Product Description

DNA Ladders in TE Buffer are used for sizing linear double-stranded DNA fragments separated by gel electrophoresis. The 1 kb DNA Ladder is suitable for sizing DNA fragments from 250 bp to 10 kb, while the 100 bp DNA ladder is suitable for sizing DNA fragments from 100 bp to 1500 bp.

These DNA ladders are supplied at 100 ng/uL in TE Buffer (10 mM Tris pH 7.5, 1 mM EDTA). DNA loading buffer is not provided, but we offer <u>6X DNA Loading Buffers</u> with blue or orange tracking dyes.

The 1 kb DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 250 bp to 10 kb. The 1 kb and 3 kb bands have increased intensity to provide internal orientation. When 100 ng of 1 kb ladder is loaded, the reference bands will contain \sim 16 ng of DNA per band, while the other bands will contain \sim 6 ng of DNA per band.

The 100 bp DNA Ladder is suitable for sizing linear double-stranded DNA fragments from 100 bp to 1500 bp. The 500 bp band has increased intensity to provide internal orientation. When 100 ng of 100 bp ladder is loaded, the reference band will contain \sim 23 ng of DNA, while the other bands will contain \sim 7.7 ng of DNA per band.

Biotium also offers Ready-to-Load Ladders 1 kb (Cat. No. 31084) and 100 bp (Cat. No. 31085) that are pre-diluted in 1X loading buffer. A separate vial of 6X loading buffer is also included to add to your DNA samples.

<u>Learn more about our products for nucleic acid gel staining</u>, including safer and more sensitive <u>GelRed®</u> and <u>GelGreen®</u> nucleic acid gel stains, and <u>Go-Go™ Fast DNA Gel Running Buffer</u> for running gels 3X faster than with TAE or TBE buffer. Also see our unique <u>Gel-Bright™ Laser Diode Gel Illuminator</u> which offers exceptional sensitivity and is a safer alternative to UV transilluminators.

This datasheet was generated on November 4, 2025 at 02:49:55 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dna-ladders-in-te-buffer/