
EpiQuik 100 bp DNA Ladder

(Catalog No. R12003)

Description

EpiQuik 100 bp DNA Ladder can be used for precise visual sizing of double stranded DNA fragments from 100 to 1000 bp on agarose gels. Supplied in sufficient quantity for 100 loadings using 4-5 μ l per loading. Note: Not designed for quantifying DNA concentration in a sample.

Composition

Consists of multiple repeats of 100 bp fragments from 100 bp up to 1000 bp. (Note: This ladder needs to be mixed with loading dye of choice).

Applications

For precise sizing of double-stranded DNA fragments from 100 to 1000 bp on agarose gels.

Storage Conditions

Stable at room temperature for 12 months.

Recommendations

For best results, please load 5 μ l of the EpiQuik 100 bp DNA Ladder per well. (Do not forget to premix with appropriate loading dye.)

Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to www.epigentek.com for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

Ordering Information

Products	Size	Cat. No.
EpiQuik 100 bp DNA Ladder	500 μ l	R12003-1

This product is for research purposes only. Not intended for use in diagnostic procedures.